

# Ni-NTA Purification System

For purification of polyhistidine-containing recombinant proteins

**Catalog Numbers** K950-01, K953-01, R901-01, R901-10, R901-15

**Document Part Number** 25-0496

**Publication Number** MAN0000270

**Revision** A.0



# Contents

Kit Contents and Storage.....	4
Accessory Products .....	6
<b>Introduction .....</b>	<b>7</b>
Overview .....	7
<b>Methods .....</b>	<b>8</b>
Preparing Cell Lysates .....	8
Purification Procedure – Native Conditions.....	13
Purification Procedure – Denaturing Conditions .....	17
Purification Procedure – Hybrid Conditions.....	19
Troubleshooting.....	21
<b>Appendix.....</b>	<b>23</b>
Additional Protocols .....	23
<i>In Situ</i> Digestion .....	24
Recipes .....	25
Frequently Asked Questions.....	28
Technical Support.....	29
Purchaser Notification.....	29
References.....	30

## Kit Contents and Storage

### Types of Products

This manual is supplied with the following products:

Kit Name	Catalog No.
Ni-NTA Purification System	K950-01
Ni-NTA Purification System with Antibody, <i>with Anti-His(C-term)-HRP Antibody</i>	K953-01
Ni-NTA Agarose (10 mL)	R901-01
Ni-NTA Agarose (25 mL)	R901-15
Ni-NTA Agarose (100 mL)	R901-10

### System Components

The Ni-NTA Purification System components are listed in the following table and include enough resin, reagents, and columns for six purifications.

Component	Composition	Quantity
Ni-NTA Agarose	50% slurry in 30% ethanol	10 mL
5X Native Purification Buffer	250 mM NaH <sub>2</sub> PO <sub>4</sub> , pH 8.0 2.5 M NaCl	1 × 125 mL bottle
Guanidinium Lysis Buffer	6 M Guanidine HCl 20 mM sodium phosphate, pH 7.8 500 mM NaCl	1 × 60 mL bottle
Denaturing Binding Buffer	8 M Urea 20 mM sodium phosphate, pH 7.8 500 mM NaCl	2 × 125 mL bottles
Denaturing Wash Buffer	8 M Urea 20 mM sodium phosphate, pH 6.0 500 mM NaCl	2 × 125 mL bottles
Denaturing Elution Buffer	8 M Urea 20 mM NaH <sub>2</sub> PO <sub>4</sub> , pH 4.0 500 mM NaCl	1 × 60 mL bottle
Imidazole	3 M Imidazole 20 mM sodium phosphate, pH 6.0 500 mM NaCl	1 × 8 mL bottle
Purification columns	10 mL columns	6 each

## Kit Contents and Storage, Continued

---

### Ni-NTA Purification System with Antibody

The Ni-NTA Purification System with Antibody includes resin, reagents, and columns as described for the Ni-NTA Purification System and 50  $\mu$ L of the appropriate purified mouse monoclonal antibody. Sufficient reagents are included to perform six purifications and 25 Western blots with the antibody.

For more details on the antibody specificity, subclass, and protocols for using the antibody, refer to the antibody manual supplied with the system.

---

### Storage

Store Ni-NTA Agarose at 4°C. Store buffers and columns at room temperature.

Store the antibody at 4°C. Avoid repeated freezing and thawing of the antibody as it may result in loss of activity.

The product is guaranteed for 6 months when stored properly.

---

### Note

All native purification buffers are prepared from the 5X Native Purification Buffer and the 3 M Imidazole, as described on page 13.

The Denaturing Wash Buffer pH 5.3 is prepared from the Denaturing Wash Buffer (pH 6.0), as described on page 17.

---

### Resin and Column Specifications

Ni-NTA Agarose is precharged with Ni<sup>2+</sup> ions and appears blue in color. It is provided as a 50% slurry in 30% ethanol.

Ni-NTA Agarose and purification columns have the following specifications:

- Binding capacity of Ni-NTA Agarose: 5–10 mg of protein per mL of resin
  - Average bead size: 45–165 microns
  - Pore size of purification columns: 30–35 microns
  - Recommended flow rate: 0.5 mL/min
  - Maximum linear flow rate: 700 cm/h
  - Maximum pressure: 2.8 psi (0.2 bar)
  - Column material: Polypropylene
  - pH stability (long term): 3–13
  - pH stability (short term): 2–14
-

## Accessory Products

---

### Additional Products

The following products are also available for order from Life Technologies:

Product	Quantity	Catalog No.
ProBond™ Nickel-Chelating Resin	50 mL	R801-01
	150 mL	R801-15
Polypropylene columns (empty)	50 each	R640-50
Ni-NTA Agarose	10 mL	R901-01
	25 mL	R901-15
	100 mL	R901-10
ProBond™ Purification System	6 purifications	K850-01
Anti- <i>myc</i> Antibody	50 µL	R950-25
Anti-V5 Antibody	50 µL	R960-25
Anti-Xpress™ Antibody	50 µL	R910-25
Anti-His(C-term) Antibody	50 µL	R930-25
InVision™ His-tag In-gel Stain	500 mL	LC6030
InVision™ His-tag In-gel Staining Kit	1 kit	LC6033

---

### Pre-Cast Gels and Pre-made Buffers

A large variety of pre-cast gels for SDS-PAGE and pre-made buffers for your convenience are available from Life Technologies. For details, visit our website at [www.lifetechnologies.com](http://www.lifetechnologies.com) or contact Technical Support (page 29).

---

# Introduction

## Overview

---

### Introduction

The Ni-NTA Purification System is designed for purification of 6xHis-tagged recombinant proteins expressed in bacteria, insect, and mammalian cells. The system is designed around the high affinity and selectivity of Ni-NTA Agarose for recombinant fusion proteins that are tagged with six tandem histidine residues.

The Ni-NTA Purification System is a complete system that includes purification buffers and resin for purifying proteins under native, denaturing, or hybrid conditions. The resulting proteins are ready for use in many target applications.

This manual is designed to provide generic protocols that can be adapted for your particular proteins. The optimal purification parameters will vary with each protein being purified.

---

### Ni-NTA Resin

Ni-NTA Agarose is used for purification of recombinant proteins expressed in bacteria, insect, and mammalian cells from any 6xHis-tagged vector. The resin exhibits high affinity and selectivity for 6xHis-tagged recombinant fusion proteins.

Proteins can be purified under native, denaturing, or hybrid conditions using the Ni-NTA Agarose. Proteins bound to the resin are eluted with low pH buffer or by competition with imidazole or histidine. The resulting proteins are ready for use in target applications.

---

### Note

The protocols provided in this manual are generic, and may not result in 100% pure protein. These protocols should be optimized based on the binding characteristics of your particular proteins.

---

### Binding Characteristics

Ni-NTA Agarose uses nitrilotriacetic acid (NTA), a tetradentate chelating ligand, in a highly cross-linked 6% agarose matrix. NTA binds Ni<sup>2+</sup> ions by four coordination sites.

---

### Native Versus Denaturing Conditions

The decision to purify 6xHis-tagged proteins under native or denaturing conditions depends on the solubility of the protein and the need to retain biological activity for downstream applications.

- Use **native conditions** if your protein is soluble (in the supernatant after lysis) and you want to preserve protein activity.
  - Use **denaturing conditions** if the protein is insoluble (in the pellet after lysis) or if your downstream application does not depend on protein activity.
  - Use **hybrid protocol** if your protein is insoluble but you want to preserve protein activity. Prepare the lysate and columns under denaturing conditions and then use native buffers during the wash and elution steps to refold the protein. Note that this protocol may not restore activity for all proteins. See page 20.
-

# Methods

## Preparing Cell Lysates

---

### Introduction

Instructions for preparing lysates from bacteria, insect, and mammalian cells using native or denaturing conditions are described in the following sections.

---

### Materials Needed

You will need the following items:

- Native Binding Buffer (recipe is on page 14) for preparing lysates under native conditions
  - Sonicator
  - *(Optional)* 10 µg/mL RNase and 5 µg/mL DNase I
  - Guanidinium Lysis Buffer (supplied with the system) for preparing lysates under denaturing conditions
  - 18-gauge needle
  - Centrifuge
  - Sterile, distilled water
  - SDS-PAGE sample buffer
  - Lysozyme for preparing bacterial cell lysates
  - Bestatin or leupeptin, for preparing mammalian cell lysates
- 

### Processing Higher Amount of Starting Material

Instructions for preparing lysates from specific amount of starting material (bacteria, insect, and mammalian cells) and purification using 2 mL resin under native or denaturing conditions are described in this manual.

If you wish to purify your protein of interest from higher amounts of starting material, you may need to optimize the lysis protocol and purification conditions (amount of resin used for binding). The optimization depends on the expected yield of your protein and amount of resin to use for purification.

Perform a pilot experiment to optimize the purification conditions and then based on the pilot experiment results, scale up accordingly.

---

## Preparing Cell Lysates, Continued

---

### Preparing Bacterial Cell Lysate – Native Conditions

Use the following procedure to prepare bacterial cell lysate under native conditions. Scale up or down as necessary.

1. Harvest cells from a 50 mL culture by centrifugation (e.g., 5000 rpm for 5 minutes in a Sorvall SS-34 rotor). Resuspend the cells in 8 mL of Native Binding Buffer (recipe on page 14).
2. Add 8 mg lysozyme and incubate on ice for 30 minutes.
3. Using a sonicator equipped with a microtip, sonicate the solution on ice using six 10-second bursts at high intensity with a 10-second cooling period between each burst.

Alternatively, sonicate the solution on ice using two or three 10-second bursts at medium intensity, then flash freeze the lysate in liquid nitrogen or a methanol dry ice slurry. Quickly thaw the lysate at 37°C and perform two more rapid sonicate-freeze-thaw cycles.

4. (*Optional*) If the lysate is very viscous, add RNase A (10 µg/mL) and DNase I (5 µg/mL) and incubate on ice for 10–15 minutes. Alternatively, draw the lysate through an 18-gauge syringe needle several times.
5. Centrifuge the lysate at  $3000 \times g$  for 15 minutes to pellet the cellular debris. Transfer the supernatant to a fresh tube.

**Note:** Some 6xHis-tagged protein may remain insoluble in the pellet, and can be recovered by preparing a denatured lysate (page 8) followed by the denaturing purification protocol (page 18). To recover this insoluble protein while preserving its biological activity, you can prepare the denatured lysate and then follow the hybrid protocol on page 20. Note that the hybrid protocol may not restore activity in all cases, and should be tested with your particular protein.

6. Remove 5 µL of the lysate for SDS-PAGE analysis. Store the remaining lysate on ice or freeze at –20°C. When ready to use, proceed to the protocol on page 16.
-

## Preparing Cell Lysates, Continued

---

### Preparing Bacterial Cell Lysate – Denaturing Conditions

Use the following procedure to prepare bacterial cell lysate under denaturing conditions:

1. Equilibrate the Guanidinium Lysis Buffer, pH 7.8 (supplied with the system or see page 26 for recipe) to 37°C.
2. Harvest cells from a 50 mL culture by centrifugation (e.g., 5000 rpm for 5 minutes in a Sorvall SS-34 rotor).
3. Resuspend the cell pellet in 8 mL of Guanidinium Lysis Buffer from Step 1.
4. Slowly rock the cells for 5–10 minutes at room temperature to ensure thorough cell lysis.
5. Sonicate the cell lysate on ice with three 5-second pulses at high intensity.
6. Centrifuge the lysate at  $3000 \times g$  for 15 minutes to pellet the cellular debris. Transfer the supernatant to a fresh tube.
7. Remove 5  $\mu$ L of the lysate for SDS-PAGE analysis. Store the remaining lysate on ice or at  $-20^{\circ}\text{C}$ . When ready to use, proceed to the denaturing protocol on page 17 or hybrid protocol on page 19.

**Note:** To perform SDS-PAGE with samples in Guanidinium Lysis Buffer, you need to dilute the samples, dialyze the samples, or perform TCA precipitation prior to SDS-PAGE to prevent the precipitation of SDS.

---

### Harvesting Insect Cells

For detailed protocols dealing with insect cell expression, consult the manual for your particular system. The following lysate protocols are for baculovirus-infected cells and are intended to be highly generic. They should be optimized for your cell lines.

For baculovirus-infected insect cells, when the time point of maximal expression has been determined, large scale protein expression can be carried out. Generally, the large-scale expression is performed in 1 liter flasks seeded with cells at a density of  $2 \times 10^6$  cells/mL in a total volume of 500 mL and infected with high titer viral stock at an MOI of 10 pfu/cell. At the point of maximal expression, harvest cells in 50 mL aliquots. Pellet the cells by centrifugation and store at  $-70^{\circ}\text{C}$  until needed. Proceed to preparing cell lysates using native or denaturing conditions as described on page 11.

---

## Preparing Cell Lysates, Continued

---

### Preparing Insect Cell Lysate – Native Condition

1. Prepare 8 mL Native Binding Buffer (recipe on page 14) containing leupeptin (a protease inhibitor) at a concentration of 0.5  $\mu\text{g}/\text{mL}$ .
  2. After harvesting the cells (page 10), resuspend the cell pellet in 8 mL Native Binding Buffer containing 0.5  $\mu\text{g}/\text{mL}$  Leupeptin.
  3. Lyse the cells by two freeze-thaw cycles using a liquid nitrogen or dry ice/ethanol bath and a 42°C water bath.
  4. Shear DNA by passing the preparation through an 18-gauge needle four times.
  5. Centrifuge the lysate at  $3000 \times g$  for 15 minutes to pellet the cellular debris. Transfer the supernatant to a fresh tube.
  6. Remove 5  $\mu\text{L}$  of the lysate for SDS-PAGE analysis. Store remaining lysate on ice or freeze at  $-20^\circ\text{C}$ . When ready to use, proceed to the protocol on page 13.
- 

### Preparing Insect Cell Lysate – Denaturing Condition

1. After harvesting insect cells (page 10), resuspend the cell pellet in 8 mL Guanidinium Lysis Buffer (supplied with the system or see page 26 for recipe).
  2. Pass the preparation through an 18-gauge needle four times.
  3. Centrifuge the lysate at  $3000 \times g$  for 15 minutes to pellet the cellular debris. Transfer the supernatant to a fresh tube.
  4. Remove 5  $\mu\text{L}$  of the lysate for SDS-PAGE analysis. Store remaining lysate on ice or freeze at  $-20^\circ\text{C}$ . When ready to use, proceed to the denaturing protocol on page 17 or hybrid protocol on page 20.  
**Note:** To perform SDS-PAGE with samples in Guanidinium Lysis Buffer, you need to dilute the samples, dialyze the samples, or perform TCA precipitation prior to SDS-PAGE to prevent the precipitation of SDS.
-

## Preparing Cell Lysates, Continued

---

### Preparing Mammalian Cell Lysate – Native Conditions

For detailed protocols dealing with mammalian expression, consult the manual for your particular system. The following protocols are intended to be highly generic, and should be optimized for your cell lines.

To produce recombinant protein, you need between  $5 \times 10^6$  and  $1 \times 10^7$  cells. Seed cells and grow in the appropriate medium until they are 80–90% confluent. Harvest the cells by trypsinization. You can freeze the cell pellet in liquid nitrogen and store at  $-70^\circ\text{C}$  until use.

1. Resuspend the cell pellet in 8 mL Native Binding Buffer (recipe on page 14). The addition of protease inhibitors such as bestatin and leupeptin may be necessary depending on the cell line and expressed protein.
  2. Lyse the cells by two freeze-thaw cycles using a liquid nitrogen or dry ice/ethanol bath and a  $42^\circ\text{C}$  water bath.
  3. Shear the DNA by passing the preparation through an 18-gauge needle four times.
  4. Centrifuge the lysate at  $3000 \times g$  for 15 minutes to pellet the cellular debris. Transfer the supernatant to a fresh tube.
  5. Remove 5  $\mu\text{L}$  of the lysate for SDS-PAGE analysis. Store the remaining lysate on ice or freeze at  $-20^\circ\text{C}$ . When ready to use, proceed to the protocol on page 13.
- 

### Preparing Mammalian Cell Lysates – Denaturing Conditions

For detailed protocols dealing with mammalian expression, consult the manual for your particular system. The following protocols are intended to be highly generic, and should be optimized for your cell lines.

To produce recombinant protein, you need between  $5 \times 10^6$  and  $1 \times 10^7$  cells. Seed cells and grow in the appropriate medium until they are 80–90% confluent. Harvest the cells by trypsinization. You can freeze the cell pellet in liquid nitrogen and store at  $-70^\circ\text{C}$  until use.

1. Resuspend the cell pellet in 8 mL Guanidinium Lysis Buffer (supplied with the system or see page 26 for recipe).
2. Shear the DNA by passing the preparation through an 18-gauge needle four times.
3. Centrifuge the lysate at  $3000 \times g$  for 15 minutes to pellet the cellular debris. Transfer the supernatant to a fresh tube.
4. Remove 5  $\mu\text{L}$  of the lysate for SDS-PAGE analysis. Store the remaining lysate on ice or freeze at  $-20^\circ\text{C}$  until use. When ready to use, proceed to the denaturing protocol on page 17 or hybrid protocol on page 20.

**Note:** To perform SDS-PAGE with samples in Guanidinium Lysis Buffer, you need to dilute the samples, dialyze the samples, or perform TCA precipitation prior to SDS-PAGE to prevent the precipitation of SDS.

---

# Purification Procedure – Native Conditions

---

## Introduction

In the following procedure, use the prepared Native Binding, Wash, and Elution Buffers, columns, and cell lysate prepared under native conditions for purification. Be sure to check the pH of your buffers before starting.

---

## Buffers for Native Purification

All buffers for purification under native conditions are prepared from the 5X Native Purification Buffer supplied with the system. Dilute and adjust the pH of the 5X Native Purification Buffer to create 1X Native Purification Buffer (see page 14). From this, you can create the following buffers:

- Native Binding Buffer
- Native Wash Buffer
- Native Elution Buffer

The recipes described in this section will create sufficient buffers to perform one native purification using one kit-supplied purification column. Scale up accordingly.

If you are preparing your own buffers, see page 25 for recipe.

---

## Materials Needed

You will need the following items:

- 5X Native Purification Buffer (supplied with the system or see page 25 for recipe)
  - 3 M Imidazole (supplied with the system or see page 25 for recipe)
  - NaOH
  - HCl
  - Sterile distilled water
  - Prepared Ni-NTA columns with native buffers (page 15)
  - Lysate prepared under native conditions (page 8)
- 

## Imidazole Concentration in Native Buffers

Imidazole is included in the Native Wash and Elution buffers to minimize the binding of untagged, contaminating proteins and increase the purity of the target protein with fewer wash steps. Note that, if your level of contaminating proteins is high, you may add imidazole to the Native Binding Buffer.

If your protein does not bind well under these conditions, you can experiment with lowering or eliminating the imidazole in the buffers and increasing the number of wash and elution steps.

---

## Purification Procedure – Native Conditions, Continued

---

### 1X Native Purification Buffer

To prepare 100 mL 1X Native Purification Buffer, combine:

- 80 mL of sterile distilled water
- 20 mL of 5X Native Purification Buffer (supplied with the system or see page 25 for recipe)

Mix well and adjust pH to 8.0 with NaOH or HCl.

---

### Native Binding Buffer

#### Without Imidazole

Use 30 mL of the 1X Native Purification Buffer (see above for recipe) for use as the Native Binding Buffer (used for column preparation, cell lysis, and binding).

#### With Imidazole (Optional):

You can prepare the Native Binding Buffer with imidazole to reduce the binding of contaminating proteins. (Note that some His-tagged proteins may not bind under these conditions.)

To prepare 30 mL Native Binding Buffer with 10 mM imidazole, combine:

- 30 mL of 1X Native Purification Buffer
- 100  $\mu$ L of 3 M Imidazole, pH 6.0

Mix well and adjust pH to 8.0 with NaOH or HCl.

---

### Native Wash Buffer

To prepare 50 mL Native Wash Buffer with 20 mM imidazole, combine:

- 50 mL of 1X Native Purification Buffer
- 335  $\mu$ L of 3 M Imidazole, pH 6.0

Mix well and adjust pH to 8.0 with NaOH or HCl.

---

### Native Elution Buffer

To prepare 15 mL Native Elution Buffer with 250 mM imidazole, combine:

- 13.75 mL of 1X Native Purification Buffer
- 1.25 mL of 3 M Imidazole, pH 6.0

Mix well and adjust pH to 8.0 with NaOH or HCl.

---

## Purification Procedure – Native Conditions, Continued

---

### Note

Do not use strong reducing agents such as DTT with Ni-NTA Agarose columns. DTT reduces the nickel ions in the resin. In addition, do not use strong chelating agents such as EDTA or EGTA in the loading buffers or wash buffers, as these will strip the nickel from the columns.

Be sure to check the pH of your buffers before starting.

---

### Preparing Ni-NTA Column

When preparing a column as described below, make sure that the snap-off cap at the bottom of the column remains **intact**. To prepare a column:

1. Resuspend the Ni-NTA Agarose in its bottle by inverting and gently tapping the bottle repeatedly.
  2. Pipet or pour 1.5 mL of the resin into a 10-mL Purification Column. Allow the resin to settle completely by gravity (5–10 minutes) or gently pellet it by low-speed centrifugation (1 minute at  $800 \times g$ ). Gently aspirate the supernatant.
  3. Add 6 mL sterile, distilled water and resuspend the resin by alternately inverting and gently tapping the column.
  4. Allow the resin to settle using gravity or centrifugation as described in Step 2, and gently aspirate the supernatant.
  5. For purification under **Native Conditions**, add 6 mL Native Binding Buffer (recipe on previous page).
  6. Resuspend the resin by alternately inverting and gently tapping the column.
  7. Allow the resin to settle using gravity or centrifugation as described in Step 2, and gently aspirate the supernatant.
  8. Repeat Steps 5 through 7.
- 

### Storing Prepared Columns

To store a column containing resin, add 0.02% azide or 20% ethanol as a preservative and cap or parafilm the column. Store at room temperature.

---

## Purification Procedure – Native Conditions, Continued

---

### Purification Under Native Conditions

Using the native buffers, columns and cell lysate, follow the procedure below to purify proteins under native conditions:

1. Add 8 mL lysate prepared under native conditions to a prepared Purification Column (page 15).
2. Bind for 30–60 minutes using gentle agitation to keep the resin suspended in the lysate solution.
3. Settle the resin by gravity or low-speed centrifugation ( $800 \times g$ ), and carefully aspirate the supernatant. Save supernatant at  $4^{\circ}\text{C}$  for SDS-PAGE analysis.
4. Wash with 8 mL Native Wash Buffer (see page 14). Settle the resin by gravity or low-speed centrifugation ( $800 \times g$ ), and carefully aspirate the supernatant. Save supernatant at  $4^{\circ}\text{C}$  for SDS-PAGE analysis.
5. Repeat Step 4 three more times.
6. Clamp the column in a vertical position and snap off the cap on the lower end. Elute the protein with 8–12 mL Native Elution Buffer (see page 14). Collect 1 mL fractions and analyze with SDS-PAGE.

**Note:** Store the eluted fractions at  $4^{\circ}\text{C}$ . If  $-20^{\circ}\text{C}$  storage is required, add glycerol to the fractions. For long term storage, add protease inhibitors to the fractions.

If you wish to reuse the resin to purify the same recombinant protein, wash the resin with 0.5 M NaOH for 30 minutes and equilibrate the resin in a suitable binding buffer. If you need to recharge the resin, see page 23.

---

# Purification Procedure – Denaturing Conditions

---

## Introduction

Instructions to perform purification using denaturing conditions with prepared denaturing Buffers, columns, and cell lysate are described below. Be sure to check the pH of your buffers before starting.

---

## Materials Needed

You will need the following items:

- Denaturing Binding Buffer (supplied with the system or see page 26 for recipe)
  - Denaturing Wash Buffer, pH 6.0 (supplied with the system or see page 27 for recipe) and Denaturing Wash Buffer, pH 5.3 (see recipe below)
  - Denaturing Elution Buffer (supplied with the system or see page 27 for recipe)
  - Prepared Ni-NTA Agarose with denaturing buffers (below)
  - Lysate prepared under denaturing conditions (page 8)
- 

## Note

Be sure to check the pH of your buffers before starting. Note that the denaturing buffers containing urea will become more basic over time.

---

## Preparing the Denaturing Wash Buffer pH 5.3

Using a 10 mL aliquot of the kit-supplied Denaturing Wash Buffer (pH 6.0), adjust the pH to 5.3 using HCl. Use this for the Denaturing Wash Buffer pH 5.3 in Step 5 next page.

---

## Preparing Ni-NTA Column

When preparing a column as described below, make sure that the snap-off cap at the bottom of the column remains **intact**.

If you are reusing the Ni-NTA Agarose, see page 23 for recharging protocol.

To prepare a column:

1. Resuspend the Ni-NTA Agarose in its bottle by inverting and gently tapping the bottle repeatedly.
  2. Pipet or pour 2 mL of the resin into a 10-mL Purification Column supplied with the kit. Allow the resin to settle completely by gravity (5–10 minutes) or gently pellet it by low-speed centrifugation (1 minute at  $800 \times g$ ). Gently aspirate the supernatant.
  3. Add 6 mL of sterile, distilled water and resuspend the resin by alternately inverting and gently tapping the column.
  4. Allow the resin to settle using gravity or centrifugation as described in Step 2, and gently aspirate the supernatant.
  5. For purification under **Denaturing Conditions**, add 6 mL of Denaturing Binding Buffer.
  6. Resuspend the resin by alternately inverting and gently tapping the column.
  7. Allow the resin to settle using gravity or centrifugation as described in Step 2, and gently aspirate the supernatant. Repeat Steps 5 through 7.
-

## Purification Procedure – Denaturing Conditions, Continued

---

### Purification Under Denaturing Conditions

Using the denaturing buffers, columns, and cell lysate, follow the provided procedure to purify proteins under denaturing conditions:

1. Add 8 mL lysate to a prepared Purification Column.
2. Bind for 15–30 minutes at room temperature using gentle agitation (e.g., using a rotating wheel) to keep the resin suspended in the lysate solution. Settle the resin by gravity or low-speed centrifugation ( $800 \times g$ ), and carefully aspirate the supernatant.
3. Wash the column with 4 mL Denaturing Binding Buffer by resuspending the resin and rocking for two minutes. Settle the resin by gravity or low-speed centrifugation ( $800 \times g$ ), and carefully aspirate the supernatant. Save supernatant at  $4^\circ\text{C}$  for SDS-PAGE analysis. Repeat this step one more time.
4. Wash the column with 4 mL Denaturing Wash Buffer (pH 6.0) by resuspending the resin and rocking for two minutes. Settle the resin by gravity or low-speed centrifugation ( $800 \times g$ ), and carefully aspirate the supernatant. Save supernatant at  $4^\circ\text{C}$  for SDS-PAGE analysis. Repeat this step one more time.
5. Wash the column with 4 mL Denaturing Wash Buffer pH 5.3 (see previous page) by resuspending the resin and rocking for two minutes. Settle the resin by gravity or low-speed centrifugation ( $800 \times g$ ), and carefully aspirate the supernatant. Save supernatant at  $4^\circ\text{C}$  for SDS-PAGE analysis. Repeat this step once more for a total of two washes with Denaturing Wash Buffer pH 5.3.
6. Clamp the column in a vertical position and snap off the cap on the lower end. Elute the protein by adding 5 mL Denaturing Elution Buffer. Collect 1 mL fractions and monitor the elution by taking  $\text{OD}_{280}$  readings of the fractions. Pool fractions that contain the peak absorbance and dialyze against 10 mM Tris, pH 8.0, 0.1% Triton X-100 overnight at  $4^\circ\text{C}$  to remove the urea. Concentrate the dialyzed material by any standard method (i.e., using 10,000 MW cut-off, low-protein binding centrifugal instruments or vacuum concentration instruments).

If you wish to reuse the resin to purify the same recombinant protein, wash the resin with 0.5 M NaOH for 30 minutes and equilibrate the resin in a suitable binding buffer. If you need to recharge the resin, see page 23.

---

# Purification Procedure – Hybrid Conditions

---

## Introduction

For certain insoluble proteins, the following protocol can be used to restore protein activity following cell lysis and binding under denaturing conditions. Note that this procedure will not work for all proteins, and should be tested using your particular recombinant proteins.

---

## Note

Be sure to check the pH of your buffers before starting. Note that the denaturing buffers containing urea will become more basic over time.

---

## Materials Needed

You will need the following items:

- Denaturing Binding Buffer (supplied with the system or see page 26 for recipe)
  - Denaturing Wash Buffer, pH 6.0 (supplied with the system or see page 27 for recipe)
  - Native Wash Buffer (page 14 for recipe)
  - Native Elution Buffer (page 14 for a recipe)
  - Prepared Ni-NTA Agarose Columns under denaturing conditions (page 17)
  - Lysate prepared under denaturing conditions (page 8)
- 

## Ni-NTA Columns

Prepare the Ni-NTA columns using Denaturing Binding Buffer as described on page 17.

---

## Purification Procedure – Hybrid Conditions, Continued

---

### Purification Under Hybrid Conditions

Using the denaturing buffers and columns and cell lysate prepared under denaturing conditions, follow the provided purification procedure to purify and renature target proteins:

1. Add 8 mL lysate to a prepared Purification Column.
2. Bind for 15–30 minutes at room temperature using gentle agitation (e.g., on a rotating wheel) to keep the resin suspended in the lysate solution. Settle the resin by gravity or low-speed centrifugation ( $800 \times g$ ) and carefully aspirate the supernatant.
3. Wash the column with 4 mL Denaturing Binding Buffer by resuspending the resin and rocking for two minutes. Settle the resin by gravity or low-speed centrifugation ( $800 \times g$ ) and carefully aspirate the supernatant. Save supernatant at  $4^\circ\text{C}$  for SDS-PAGE analysis. Repeat this step one more time.
4. Wash the column with 4 mL Denaturing Wash Buffer (pH 6.0) by resuspending the resin and rocking for two minutes. Settle the resin by gravity or low-speed centrifugation ( $800 \times g$ ) and carefully aspirate the supernatant. Save supernatant at  $4^\circ\text{C}$  for SDS-PAGE analysis. Repeat this step one more time.
5. Wash the column with 8 mL Native Wash Buffer (see page 14) by resuspending the resin and rocking for two minutes. Settle the resin by gravity or low-speed centrifugation ( $800 \times g$ ) and carefully aspirate the supernatant. Save supernatant at  $4^\circ\text{C}$  for SDS-PAGE analysis. Repeat this step three more times for a total of four native washes.
6. Clamp the column in a vertical position and snap off the cap on the lower end. Elute the protein with 8–12 mL Native Elution Buffer (see page 14). Collect 1 mL fractions and analyze with SDS-PAGE.

If you wish to reuse the resin to purify the same recombinant protein, wash the resin with 0.5 M NaOH for 30 minutes and equilibrate the resin in a suitable binding buffer. If you need to recharge the resin, see page 23.

---

# Troubleshooting

## Introduction

Review the information below to troubleshoot your experiments with the Ni-NTA Purification System.

For troubleshooting problems with antibody detection, see the antibody manual supplied with the system.

Problem	Probable Cause	Possible Solution
No recombinant protein recovered following elution.	Nothing bound because of protein "folding" .	Try denaturing conditions.
	Expression levels too low.	Optimize expression levels using the guidelines in your expression manual.
	Protein washed out by too stringent washing.	Raise pH of wash buffer in high-stringency wash step. Wash less extensively in high-stringency wash step.
	Not enough sample loaded.	Increase amount of sample loaded or lysate used.
	Recombinant protein has very high affinity for Ni-NTA Agarose.	Increase stringency of elution by decreasing the pH or increasing the imidazole concentration. To preserve activity, use EDTA or EGTA (10–100 mM) to strip resin of nickel ions and elute protein.
	Protein degraded.	Perform all purification steps at 4°C. Check to make sure that the 6xHis-tag is not cleaved during processing or purification. Include protease inhibitors during cell lysis.
Some recombinant protein is in the flow through and wash fractions.	Protein overload.	Load less protein on the column or use more resin for purification.
Good recombinant-protein recovery but contaminated with non-recombinant proteins.	Wash conditions not stringent enough.	Lower pH of wash buffer in high-stringency wash step. Wash more extensively.
	Other His-rich proteins in sample.	Consider an additional high stringency wash at a lower pH (i.e., between pH 6 and pH 4) before the elution step. Further purify the eluate on a new Ni-NTA Agarose column after dialysis of the eluate against the binding buffer and equilibrating the column with binding buffer. Perform second purification over another type of column.
	Recombinant protein has low affinity for resin; comes off in wash with many contaminating proteins.	Try denaturing conditions. Try an imidazole step gradient elution. Try a pH gradient with decreasing pH.

## Troubleshooting, Continued

Problem	Probable Cause	Possible Solution
Low recombinant protein recovery and contaminated with non-recombinant proteins.	Recombinant protein not binding tightly to resin.	Try denaturing conditions. Try "reverse-chromatography": bind lysate, including recombinant protein; allow recombinant protein to come off in low stringency washes; collect these fractions; re-do chromatography on saved fractions on new or stripped and recharged column. Works for native purification only.
	Expression levels too low.	Consider an additional high stringency wash at a lower pH (i.e., between pH 6 and pH 4) before elution step.
Column turns reddish brown.	DTT is present in buffers.	Use $\beta$ -mercaptoethanol as a reducing agent.
Column turns white.	Chelating agents present in buffer that strip the nickel ions from the column.	Recharge the column as described on page 23.
Protein precipitates during binding.	Temperature is too low.	Perform purification at room temp.
	Protein forms aggregates.	Add solubilization reagents such as 0.1% Triton X-100 or Tween-20 or stabilizers such as $Mg^{2+}$ . These may be necessary in all buffers to maintain protein solubility. Run column in drip mode to prevent protein from dropping out of solution.

# Appendix

## Additional Protocols

---

### Cleavage of the Fusion Peptide

If your recombinant fusion protein contains the recognition sequence for enterokinase (EnterokinaseMax™ enzyme) or AcTEV™ Protease between the 6xHis-tag and the protein, you may cleave the 6xHis-tag from the fusion protein using the specific protease. You may cleave the tag after obtaining the purified recombinant fusion protein or while the protein is bound to the nickel-chelating resin (see *In Situ Digestion*).

**EnterokinaseMax™** (EKMax™ enzyme) is a recombinant preparation of the catalytic subunit of enterokinase. This enzyme recognizes -Asp-Asp-Asp-Asp-Lys- and cleaves after the lysine. It has high specific activity, leading to more efficient cleavage, and requires less enzyme.

Description	Cat. No.
EKMax™ Enterokinase, 250 units	E180-01
EKMax™ Enterokinase, 1000 units	E180-02

**AcTEV™ Protease** is an enhanced form of Tobacco Etch Virus (TEV) protease that is highly site-specific, active, and more stable than native TEV protease. AcTEV™ Protease recognizes the seven-amino-acid sequence Glu-Asn-Leu-Tyr-Phe-Gln-Gly- and cleaves between Gln and Gly with high specificity.

Description	Cat. No.
AcTEV™ Protease, 1000 units	12575-015
AcTEV™ Protease, 10,000 units	12575-023

---

### Recharging Ni-NTA Resin

Ni-NTA resin can be used for up to three or four purifications of the same protein without recharging. Wash the resin with 0.5 M NaOH for 30 minutes and equilibrate the resin with the appropriate binding buffer, if you are reusing the resin.

We recommend not recharging the resin more than three times and only reusing it for purification of the same recombinant protein. If the resin turns white due to the loss of nickel ions from the column, recharge the resin.

To recharge 2 mL of resin in a purification column:

1. Wash the column two times with 8 mL 50 mM EDTA to strip away the chelated nickel ions.
  2. Wash the column two times with 8 mL 0.5 M NaOH.
  3. Wash the column two times with 8 mL sterile, distilled water.
  4. Recharge the column with two washes of 8 mL NiCl<sub>2</sub> hexahydrate at a concentration of 5 mg/mL prepared in sterile, distilled water.
  5. Wash the column two times with 8 mL distilled water.
  1. Add 0.02% azide or 20% ethanol as a preservative and cap or apply a parafilm to the column. Store at room temperature.
-

# In Situ Digestion

---

## Introduction

EKMax™ enzyme can be used to cleave Xpress™ fusion proteins while they are bound to the Ni-NTA resin. The leader peptide will remain bound to the resin while the cleaved native protein is collected in the flow-through fraction. This in situ digestion simplifies purification by removing the leader peptide and undigested fusion protein from the native protein.

It is necessary to exchange the column buffer with 1X EKMax™ buffer prior to digestion since EKMax™ enzyme is known to be inhibited by >2 M urea, >20 mM β-mercaptoethanol (β-ME), >0.1% SDS, >50 mM Imidazole, and pH values below 6 and above 9.

---

## Materials Needed

You will need the following items:

- 5X Native Binding Buffer **without** NaCl (see page 25 for recipe)
  - 1X Native Binding Buffer **without** NaCl
  - 10X EKMax™ Reaction Buffer (supplied with EKMax™ Enzyme)
  - EKMax™ Enterokinase
- 

## In Situ Digestion Protocol

Following step 5 of Purification Under Native Conditions (page 16).

1. Wash the resin 2 times with 2X bed volumes of 1X Native Binding Buffer **without** NaCl and save the supernatant. (This is very important to do as NaCl inhibits EKMax™.)
  2. After the last wash, securely cap the bottom of the column. Add 450 μL 1X Native Binding Buffer without NaCl and 50 μL 10X EKMax™ Reaction Buffer. (EKMax™ enzyme requires CaCl<sub>2</sub> and Tween-20 for maximum activity.)
  3. Add twice the units of EKMax™ enzyme required to cleave the fusion protein in solution.  
**Note:** See EKMax™ User Guide for pilot reaction protocol to determine the amount of EKMax™ enzyme required.
  4. Seal the column and rock or rotate overnight at room temperature or 4°C.
  5. Settle the resin by gravity or low-speed centrifugation (800 × g), and carefully aspirate the supernatant. Save supernatant at 4°C for SDS-PAGE analysis. This fraction contains your cleaved protein.
  6. Resuspend the resin in 2X bed volume of Native Wash Buffer (with NaCl) and rock or rotate for 2 minutes. (Inclusion of NaCl will reduce non-specific interactions between the cleaved protein and the resin.)
  7. Settle the resin by gravity or low-speed centrifugation (800 × g), and carefully aspirate the supernatant. Repeat Steps 6 and 7 twice for a total of 3 washes. Store these fractions on ice. These fractions contain your cleaved fusion protein.
  8. Resuspend the resin in 2X bed volume Native Elution Buffer containing 250 mM Imidazole and rock or rotate for 2 minutes.
  9. Settle the resin by gravity or low-speed centrifugation (800 × g), and carefully aspirate the supernatant. Store the fractions on ice.
  10. (Steps 8 and 9 elute the Xpress™ fusion partner from the resin in order to evaluate the extent of digestion and recovery of your cleaved protein, and to enable re-use of the resin.)
-

# Recipes

---

## Buffer Stock Solutions (10X)

To prepare the buffer solutions described below, you need to prepare sodium phosphate stock solutions:

### Stock Solution A (10X)

200 mM sodium phosphate, monobasic ( $\text{NaH}_2\text{PO}_4$ )

5 M NaCl

Dissolve 27.6 g sodium phosphate, monobasic ( $\text{NaH}_2\text{PO}_4$ ) and 292.9 g NaCl in 900 mL of deionized water. Mix well and adjust the volume to 1 L with deionized water. Store solution at room temperature.

### Stock Solution B (10X)

200 mM sodium phosphate, dibasic ( $\text{Na}_2\text{HPO}_4$ )

5 M NaCl

Dissolve 28.4 g sodium phosphate, dibasic ( $\text{Na}_2\text{HPO}_4$ ) and 292.9 g of NaCl in 900 mL of deionized water. Mix well and adjust the volume to 1 L with deionized water. Store solution at room temperature.

---

## 5X Native Purification Buffer

250 mM  $\text{NaH}_2\text{PO}_4$ , pH 8.0

2.5 M NaCl

Prepare 200 mL solution as follows:

1. To 180 mL deionized water, add

Sodium phosphate, monobasic	7 g
NaCl	29.2 g
2. Mix well and adjust the pH with NaOH to pH 8.0.
3. Bring the final volume to 200 mL with water.
4. Store buffer at room temperature.

**Note:** Omit the NaCl to prepare Native Purification Buffer without NaCl.

---

## 3 M Imidazole pH 6.0

3 M Imidazole

500 mM NaCl

20 mM Sodium Phosphate Buffer, pH 6.0

Prepare 100 mL solution as follows:

2. To 80 mL deionized water, add

Imidazole	20.6 g
Stock Solution A (10X)	8.77 mL
Stock Solution B (10X)	1.23 mL
  3. Mix well and adjust the pH to 6.0 with concentrated HCl or NaOH as necessary.
  4. Bring the final volume to 100 mL with water. If the solution forms a precipitate, heat solution until the precipitate dissolves.
  5. Store buffer at room temperature.
-

## Recipes, Continued

---

### Guanidinium Lysis Buffer

6 M Guanidine Hydrochloride  
20 mM Sodium Phosphate, pH 7.8  
500 mM NaCl

Prepare 100 mL solution as follows:

1. To 60 mL deionized water, add

Stock Solution A (10X)	0.58 mL
Stock Solution B (10X)	9.42 mL
Guanidine Hydrochloride	57.3 g
  2. Stir the solution until completely dissolved. Adjust the pH to 7.8 using 1 N NaOH or 1 N HCl.
  3. Bring the volume to 100 mL and filter sterilize the buffer using a 0.45  $\mu$ m filter (autoclaving the solution will alter the pH of the buffer).
  4. Store buffer at room temperature.
- 

### Denaturing Binding Buffer

8 M Urea  
20 mM Sodium Phosphate pH 7.8  
500 mM NaCl

Prepare 100 mL solution as follows:

1. To 60 mL deionized water, add

Stock Solution A (10X)	0.58 mL
Stock Solution B (10X)	9.42 mL
Urea	48.1g
  2. Stir the solution with gentle heating (50–60°C, do not overheat) until completely dissolved. When cooled to room temperature, adjust the pH to 7.8 using 1 N NaOH or 1 N HCl.
  3. Bring the volume to 100 mL and filter sterilize the buffer using a 0.45  $\mu$ m filter (autoclaving the solution will alter the pH of the buffer).
  4. Store buffer at room temperature.
-

## Recipes, Continued

---

### Denaturing Wash Buffer

8 M Urea  
20 mM Sodium Phosphate, pH 6.0  
500 mM NaCl

Prepare 100 mL solution as follows:

1. To 60 mL deionized water, add

Stock Solution A (10X)	7.38 mL
Stock Solution B (10X)	2.62 mL
Urea	48.1g
  2. Stir the solution with gentle heating (50–60°C, do not overheat) until completely dissolved. Adjust the pH to 6.0 using 1 N NaOH or 1 N HCl.
  3. Bring the volume to 100 mL and filter sterilize the buffer using a 0.45 µm filter (autoclaving the solution will alter the pH of the buffer).
  4. Store buffer at room temperature.
- 

### Denaturing Elution Buffer

8 M Urea  
20 mM Sodium Phosphate, pH 4.0  
500 mM NaCl

Prepare 100 mL as follows:

1. To 60 mL deionized water, add

Stock Solution A (10X)	10 mL
Urea	48.1g
  2. Stir the solution with gentle heating (50–60°C, do not overheat) until completely dissolved. Adjust the pH to 4.0 using 1 N NaOH or 1 N HCl.
  3. Bring the volume to 100 mL and filter sterilize the buffer using a 0.45 µm filter (autoclaving the solution will alter the pH of the buffer).
  4. Store buffer at room temperature.
-

## Frequently Asked Questions

---

**For denatured conditions, why is Guanidinium used for lysis of cells?**

We have found that guanidinium works better for cell lysis than urea; however, urea works well for the remaining steps.

**Can proteins bind to the resin at a pH lower than 7.8?**

The optimal binding range is pH 7.2–7.8. However, we have performed purifications with columns equilibrated to pH 6.0. Some proteins bind well under these conditions and will remain bound to the column following a pH 6.0 wash.

**Can glycine be used instead of sodium phosphate in the purification system binding buffers?**

No, because glycine is a competitive ligand for nickel. People have successfully used:

- Tris-HCl
- Tris-Phosphate
- Tris-Acetate
- Sodium Acetate
- Sodium Borate
- MES-NaOH
- Pipes-HCl
- HEPES

**Can I use the resin to purify a protein with fewer than six histidine residues?**

We have not tried to purify proteins with less than six histidines. However, if several histidines are near each other, you may be able to attach the protein to the resin well enough for purification.

**Is there a cell lysis procedure that will liberate microsome-bound proteins for subsequent purification using Ni-NTA?**

If solubility is a problem, you can include up to 0.2% Sarkosyl in the 6 M Guanidinium Lysis Buffer—this should solubilize everything and may still be compatible with purification on the Ni-NTA columns. In general, anionic detergents are incompatible with nickel chelating columns, but up to 0.2% Sarkosyl has been used in some cases.

**What are recommended elution conditions for His-tagged proteins that are unstable at a pH < 7.0?**

You can elute with a stepped imidazole gradient at a neutral pH (pH 7.0–7.5). Use 10 mM imidazole, then 50 mM, 75 mM and so on until the protein elutes. Note that more contaminating proteins that would have been washed off at pH 6.0 will remain on the resin at pH 7.0.

**What is the importance of NaCl in the binding buffer?**

Ni-NTA resin has a net positive charge, and 500 mM NaCl is used to prevent the nonspecific binding of negatively charged proteins.

---

# Technical Support

---

**Obtaining Support** For the latest services and support information for all locations, go to [www.lifetechnologies.com](http://www.lifetechnologies.com).

At the website, you can:

- Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
  - Search through frequently asked questions (FAQs)
  - Submit a question directly to Technical Support ([techsupport@lifetech.com](mailto:techsupport@lifetech.com))
  - Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents
  - Obtain information about customer training
  - Download software updates and patches
- 

## **Safety Data Sheets (SDS)**

Safety Data Sheets (SDSs) are available at [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support).

---

## **Certificate of Analysis**

The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support) and search for the Certificate of Analysis by product lot number, which is printed on the box.

---

## **Limited Product Warranty**

Life Technologies and/or its affiliate(s) warrant their products as set forth in the Life Technologies General Terms and Conditions of Sale found on the Life Technologies website at [www.lifetechnologies.com/termsandconditions](http://www.lifetechnologies.com/termsandconditions). If you have any questions, please contact Life Technologies at [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support).

---

# Purchaser Notification

---

## **Limited Use Label License No. 102: Ni-NTA Resin**

The Ni-NTA Agarose contained in this product is manufactured by QIAGEN® under a license from Hoffmann-LaRoche Inc., Nutley, NJ and/or Hoffmann-LaRoche Ltd., Basel, Switzerland and is provided only for use in research. Information about licenses for commercial use is available from QIAGEN GmbH, Max-Volmer-Strasse 4, D-40724 Hilden, Germany.

---

## References

---

- Ausubel, F.M., Brent, R., Kingston, R.E., Moore, D.D., Seidman, J.G., Smith, J.A. and Struhl, K. (1994) *Current Protocols in Molecular Biology Vol. 1*. John Wiley and Sons, New York.
- Blochlinger, K. and Diggelmann, H. (1984) *Mol. Cell Biol.* **4**: 2929.
- Frost, E. and Williams, J. (1978) *Virology* **91**: 39.
- Goeddel, D.V. ed. (1991) "Expression in Mammalian Cells." *Methods in Enzymology Vol. 185*. Academic Press, San Diego, California.
- Graham, F.L. and van der Ebb, A.J. (1973) *Virology* **52**: 456.
- Kawasaki Y., *et al.* (2000) *Science* 289(5482):1194–7.
- Laemmli, U.K. (1970) *Nature* **227**: 680-685.
- Lopata, M.A., Cleveland, D.W. and Sollner-Webb, B. (1984) *Nuc. Acids Res.* **12**: 5707.
- Lowry, O. H., Rosebrough, N. J., Farr, A. L., and Randall, R. J. (1951) *J. Biol. Chem.*, **193**: 265-275.
- Lowy, D.R., Rands, E. and Scolnick, E.M. (1978) *J. Virology* **26**: 291.
- Lewis, W.H., *et al.* (1980) *Somat. Cell Genet.* **6**: 333.
- Maniatis, T., Fritsch, E.F. and Sambrook, M.D. (1989) *Molecular Cloning: A Laboratory Manual*. Cold Spring Harbor Laboratory, Cold Spring Harbor, New York.
- Wigler, M. *et al.*, (1977) *Cell* **11**: 223.
- Zhou *et al.*, (1990). *Biotechniques*, **8(2)**: 172.

---

©2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

**DISCLAIMER:** LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.



