

00-4303 1 Kit
Avidin/Biotin Blocking Kit
Lot No.: See product label
Exp. Date: See product label

ENDOGENOUS AVIDIN/BIOTIN BLOCKING KIT

INVITROGEN HISTOCHEMICAL REAGENTS

INTENDED USE

General purpose reagent. For laboratory use.

FORM:

The following reagents are supplied with the kit:

- A. One dropper bottle of Avidin solution (18 mL).
- B. One dropper bottle of d-Biotin solution (18 mL).

Both reagents contain 0.05% sodium azide (NaN₃).

APPLICATION:

This kit is designed to eliminate background caused by endogenous avidin or biotin-like substances.

RECOMMENDED GUIDELINES FOR USE:

These blocking steps should be performed before the primary antibody incubation step.

1. Apply 2 drops (50 µL/drop) of reagent A to the tissue.
2. Incubate for 10 minutes at room temperature.
3. Rinse extensively with phosphate buffered saline (PBS).
4. Apply 2 drops of reagent B to the tissue.
5. Incubate for 10 minutes at room temperature.
6. Rinse extensively with PBS.

The endogenous avidin/biotin activity is now blocked and the tissue section is ready for immunohistochemical staining.

STORAGE:

Storage at 2-8°C.

Authorized Representative for IVDD 98/79/EC: Invitrogen Ltd.
Inchinnan Business Park
3 Fountain Drive
Paisley
PA4 9RF
UK

Explanation of symbols

Symbol	Description	Symbol	Description
	Catalogue Number		Batch code
	Research Use Only		<i>In vitro</i> diagnostic medical device
	Use by		Temperature limitation
	Manufacturer		European Community authorised representative
	Without, does not contain		With, contains
	Protect from light		Consult accompanying documents
	Directs the user to consult instructions for use (IFU), accompanying the product.		

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288

E-mail: techsupport@invitrogen.com

PI004303

(Rev 08/10) DCC-10-1984

© 2007 Invitrogen Corporation. All right reserved. These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses.

00-4303 1 kit
Avidina/Biotina
Equipo de bloqueo

EQUIPO DE BLOQUEO DE LA AVIDINA/BIOTINA ENDÓGENA
REACTIVOS HISTOQUÍMICOS INVITROGEN

USO PREVISTO

Reactivos de uso general. Para uso en laboratorio.

FORMA:

Los siguientes reactivos son proporcionados con el equipo:

- A. Un frasco gotero de solución de avidina (18 ml).
- B. Un frasco gotero de solución de biotina (18 ml).

Ambos reactivos contienen azida de sodio (NaN_3) al 0,05%.

APLICACIÓN:

Este equipo está diseñado para eliminar la base provocada por sustancias endógenas del tipo de la avidina o la biotina.

RECOMENDACIONES DE USO:

Estos pasos de bloqueo deben ser realizados antes de proceder al paso de incubación del anticuerpo primario.

1. Aplicar 2 gotas (50 μl /gota) de reactivo A al tejido.
2. Incubar durante 10 minutos a temperatura ambiente.
3. Enjuagar muy bien con solución salina tampón de fostato (PBS).
4. Aplicar 2 gotas de reactivo B al tejido.
5. Incubar durante 10 minutos a temperatura ambiente.
6. Enjuagar muy bien con PBS.

La actividad endógena de la avidina/biotina está bloqueada y el corte tisular queda listo para la inmunotinción.

ALMACENAMIENTO:

Guardar a 2-8 °C.

Representante autorizado de IVDD 98/79/EC: Invitrogen Ltd.
Inchinnan Business Park
3 Fountain Drive
Paisley
PA4 9RF
UK

00-4303 1 kit
Kit de blocage
de l'avidine/biotine

**KIT DE BLOCAGE DE L'AVIDINE/BIOTINE ENDOGÈNE
RÉACTIFS HISTOCHIMIQUES DE INVITROGEN**

L'UTILISATION PREVUE

General Purpose réactif. Pour une utilisation en laboratoire.

FORME :

Les réactifs suivants sont fournis avec le kit :

- A. Un flacon compte-gouttes de solution d'avidine (18 ml).
- B. Un flacon compte-gouttes de solution de d-biotine (18 ml).

Les deux réactifs contiennent de l'azoture de sodium (NaN_3) à 0,05 %.

APPLICATION :

Ce kit a été conçu pour éliminer tout mouvement propre causé par des substances ressemblant à des substances d'avidine ou de biotine endogène.

DIRECTIVES D'UTILISATION RECOMMANDÉES :

Ces étapes de blocage doivent être effectuées avant l'étape primaire d'incubation des anticorps.

1. Déposer 2 gouttes (50 µl/goutte) de réactif A sur le tissu.
2. Incuber pendant 10 minutes à température ambiante.
3. Rincer abondamment avec du soluté tampon de phosphate (PBS).
4. Déposer 2 gouttes de réactif B sur le tissu.
5. Incuber pendant 10 minutes à température ambiante.
6. Rincer abondamment avec du PBS.

L'activité d'avidine/biotine endogène est bloquée et la section de tissu est prête pour la coloration immunohistochimique.

STOCKAGE :

Conserver entre 2 et 8 °C.

Représentant Autorisé pour IVDD 98/79/EC:

Invitrogen Ltd.
Inchinnan Business Park
3 Fountain Drive
Paisley
PA4 9RF
UK

00-4303 1 kit
Kit per il blocco
dell'avidina/biotina

**KIT PER IL BLOCCO DELLA AVIDINA/BIOTINA ENDOGENA
REAGENTI ISTOCIMICI INVITROGEN**

DESTINAZIONE D'USO

Reagente di uso generale. Per uso di laboratorio.

FORMA:

Nel kit sono acclusi i seguenti reagenti:

- A. Una boccetta contagocce (18 ml) di soluzione di avidina.
- B. Una boccetta contagocce (18 ml) di soluzione di d-biotina.

Entrambi i reagenti contengono azoturo di sodio allo 0,05% (NaN_3).

APPLICAZIONE:

Questo kit è da usarsi per l'eliminazione del fondo causato da sostanze simili all'avidina o alla biotina endogena.

RACCOMANDAZIONI PER L'USO:

Prima di eseguire la procedura di incubazione degli anticorpi primari, bisognerebbe eseguire le fasi riportate qui di seguito.

1. Applicare sul tessuto 2 gocce (50 μL /goccia) di reagente A.
2. Incubare per 10 minuti a temperatura ambiente.
3. Risciacquare accuratamente con soluzione fisiologica tamponata a base di fosfato (PBS).
4. Applicare sul tessuto 2 gocce di reagente B.
5. Incubare per 10 minuti a temperatura ambiente.
6. Risciacquare accuratamente con PBS.

L'attività dell'avidina/biotina endogena è ora bloccata e la sezione tissutale è pronta per la colorazione immunoistochimica.

CONSERVAZIONE:

Conservare a 2-8°C.

Rappresentante Autorizzato per IVDD 98/79/EC: Invitrogen Ltd.
Inchinnan Business Park
3 Fountain Drive
Paisley
PA4 9RF
UK

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288
E-mail: techsupport@invitrogen.com

PI004303

(Rev 08/10) DCC-10-1984

© 2007 Invitrogen Corporation. All right reserved. These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses.

00-4303 1 Kit
Avidin/Biotin-
Blocking-Kit

**ENDOGENES AVIDIN/BIOTIN-BLOCKING-KIT
HISTOCHEMISCHE REAGENZIEN VON INVITROGEN**

VERWENDUNGSZWECK

Allzweck-Reagens. Für das Labor.

FORMULIERUNG:

Folgende Reagenzien sind im Kit enthalten:

- A. Eine Tropfflasche mit Avidinlösung (18 ml).
- B. Eine Tropfflasche mit d-Biotinlösung (18 ml).

Beide Reagenzien enthalten 0,05% Natriumazid (NaN_3).

ANWENDUNG:

Dieses Kit ist dafür vorgesehen, den durch endogene Avidin- oder Biotin-ähnliche Substanzen verursachten Hintergrund zu beseitigen.

EMPFOHLENE RICHTLINIEN FÜR DEN GEBRAUCH:

Diese Blocking-Schritte sollten vor dem primären Antikörper-Inkubationsschritt durchgeführt werden.

1. 2 Tropfen (50 µl/Tropfen) Reagenz A auf das Gewebe auftragen.
2. Für 10 Minuten bei Raumtemperatur inkubieren.
3. Gründlich mit phosphatgepufferter Kochsalzlösung (PBS) spülen.
4. 2 Tropfen Reagenz B auf das Gewebe auftragen.
5. Für 10 Minuten bei Raumtemperatur inkubieren.
6. Gründlich mit PBS spülen.

Die endogene Avidin-/Biotin-Aktivität ist jetzt blockiert und der Gewebeschnitt ist bereit für immunohistochemische Färbung.

AUFBEWAHRUNG:

Bei 2-8°C aufbewahren.

Bevollmächtigter Repräsentant für IVDD 98/79/EC: Invitrogen Ltd.
Inchinnan Business Park
3 Fountain Drive
Paisley
PA4 9RF
UK

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288
E-mail: techsupport@invitrogen.com

PI004303

(Rev 08/10) DCC-10-1984

© 2007 Invitrogen Corporation. All right reserved. These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses.