

TURBO DNA-free™ Kit

TURBO™ DNase Treatment and Removal Reagents

Catalog Number AM1907

Publication Number 1907M **Revision G**

Product description

Ambion® TURBO DNA-free™ DNase Treatment and Removal Reagents are designed to remove contaminating DNA from RNA preparations, and to subsequently remove the DNase and divalent cations from the sample. The included TURBO™ DNase (patent pending) is an engineered version of wild type DNase I with 350% greater catalytic efficiency. TURBO™ DNase has a markedly higher affinity for DNA than conventional DNase I, and is thus more effective in removing trace quantities of DNA contamination. In addition, TURBO™ DNase maintains up to 50X greater activity than DNase I in solutions containing physiological salt concentrations. The TURBO™ DNase provided in the kit is overexpressed in an animal-free system, and is then extensively purified in a bovine-free process and tested. It is guaranteed to lack any contaminating RNase activity. The kit also includes an optimized DNase reaction buffer that contains a small molecule enhancer that extends the activity of the TURBO™ DNase enzyme by 100-fold or more. Using TURBO DNA-free™, contaminating DNA is digested to levels below the limit of detection by routine PCR (Figure 1). The DNase is then removed rapidly and easily using a novel method which does not require phenol/chloroform extraction, alcohol precipitation, heating, or the addition of EDTA (see Table 1 on page 2). TURBO DNA-free™ treated RNA is suitable for endpoint or real-time RT-PCR, microarray analysis, RPAs, Northern, and all other RNA analysis methods.

In addition to removing the DNase enzyme, DNase Inactivation Reagent also removes divalent cations, such as magnesium and calcium, which can catalyze RNA degradation when RNA is heated with the sample (Figure 2 on page 3).

Figure 1 TURBO DNA-free™ reduces genomic DNA contamination by greater than 5 million fold. Equal amounts of mouse spleen total RNA (purified using the RNAqueous® Kit) were either treated with 7.8 U of TURBO DNA-free™ in a 130 µL reaction for 20 min at 37°C, or were left untreated. The digestions were stopped by adding 22 µL DNase Inactivation Reagent. 5 µL (1 µg RNA) was amplified in a one step 25-µL RT-PCR using a TaqMan® primer probe set for mouse GAPDH. Treated and untreated samples were reverse transcribed with Life Technologies MessageSensor™ RT Kit. RT-minus samples were subjected to PCR to control for DNA contamination. Results are shown using a linear scale so that the amplification plot for the TURBO™ DNase-treated, RT-minus sample is visible.

The fold-removal (5.4×10^6 fold) of genomic DNA was calculated as follows: The C_T value from the untreated RNA in the RT-minus reaction is the level of gDNA contamination. The fold-removal was determined by subtracting the RT-minus reaction C_T value for the treated RNA sample, 39.5 (the other duplicate's signal was undetectable) from the C_T value of the untreated sample, 17.13, and raising the 17.13 as the exponent with a base of 2.

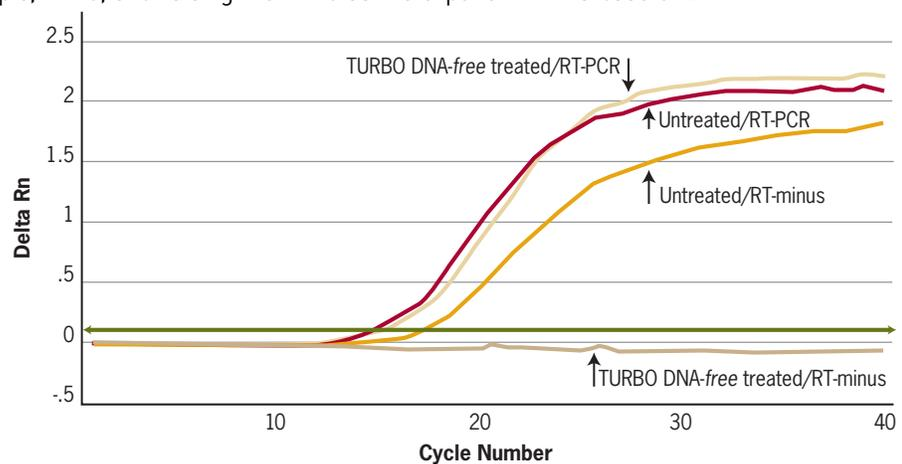
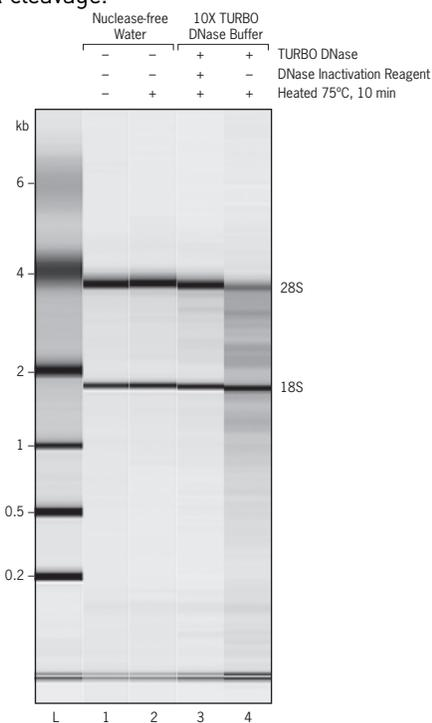


Table 1 Treatment of RNA with TURBO DNA-free™ maintains target sensitivity in real-time RT-PCR. Total RNA from HeLa S3 cells was treated with the TURBO DNA-free™ Kit following the standard protocol. 5 µL of the treated RNA was then reverse transcribed using the MessageSensor™ RT Kit, and the resulting cDNA was amplified by real-time RT-PCR using primer and probe sets for either human β-actin or CDC-2 with TaqMan® detection.

	C_T for β-actin (duplicates)	
RNA-treatment	100 pg RNA	1 pg RNA
none	24.78 / 24.67	31.83 / 31.53
TURBO DNA-free™ treated	24.50 / 24.62	30.89 / 30.88

	C_T for CDC-2 (duplicates)	
none	28.88 / 28.24	34.41 / 35.50
TURBO DNA-free™ treated	27.71 / 28.10	34.04 / 33.99

Figure 2 Removal of divalent cations by DNase inactivation reagents. HeLa-S3 total RNA (100 ng), in 50 μ L 1 \times TURBO DNase Buffer or in nuclease-free water, was treated with components from the TURBO DNA-free™ Kit as indicated. Samples were heated for 10 min at 75°C (Lanes 2, 3, & 5), or 3 min at 90°C (Lane 4), to determine if divalent cations from the TURBO DNase Buffer remained in solution, and degraded the RNA. 1 μ L of each sample was analyzed on an RNA LabChip® using the Agilent® 2100 Bioanalyzer™ Instrument. Note that RNA was degraded in the sample that contained TURBO DNase Buffer, but was not treated with the DNase Inactivation Reagent (Lane 5); this degradation is due to the presence of divalent ions that induce heat-mediated RNA cleavage.



Procedure overview

For the detailed procedure, see “TURBO DNA-free™ procedure” on page 5.

Add DNase Digestion Reagents

1. Add 0.1 volume 10X TURBO DNase Buffer and 1 μ L TURBO DNase to the RNA, and mix gently (page 5)

Incubate

2. Incubate at 37°C for 20–30 min (page 6)

Add DNase Inactivation Reagent

3. Add resuspended DNase Inactivation Reagent (typically 0.1 volume) and mix well. (page 6)

Incubate and mix

4. Incubate 5 min at room temperature, mixing occasionally. (page 6)

Centrifuge and transfer RNA

5. Centrifuge at 10,000 \times g for 1.5 min and transfer the RNA to a fresh tube (page 6)



How much RNA can be treated with TURBO DNA-free™ reagents?

This protocol is designed to remove trace to moderate amounts of contaminating DNA (up to 50 μ g DNA/mL RNA) from purified RNA to a level that is mathematically insignificant by RT-PCR. No RNA isolation method can extract RNA that is completely free from DNA contamination; in fact, RNA isolated from some tissues, such as spleen, kidney, or thymus, often contain relatively high levels of DNA. Other potential sources of DNA contamination include carryover of the interface during organic extractions, and overloaded glass-fiber filters during RNA purification.

TURBO DNA-free™ components and storage

Reagents are provided for 50 TURBO DNA-free™ treatments (up to 100 µL each).

Amount	Component	Storage
120 µL	TURBO DNase (2 Units/µL)	-20°C
600 µL	10X TURBO DNase Buffer	
600 µL	DNase Inactivation Reagent	
1.75 mL	Nuclease-free Water	any temperature†

† Store Nuclease-free Water at -20°C, 4°C, or room temperature.

Store the TURBO DNA-free™ Kit at -20°C in a non-frost-free freezer for long-term storage. For convenience, the 10X TURBO DNase Buffer and the DNase Inactivation Reagent can be stored at 4°C for up to 1 week.

TURBO DNA-free™ procedure

Procedure notes

- We recommend conducting reactions in 0.5 mL tubes to facilitate removal of the supernatant after treatment with the DNase Inactivation Reagent.
- TURBO DNA-free™ reactions can be conducted in 96-well plates. We recommend using V-bottom plates because their shape makes it easier to remove the RNA from the pelleted DNase Inactivation Reagent at the end of the procedure.
- The recommended reaction size is 10–100 µL. A typical reaction is 50 µL.

Procedure

1. Add 0.1 volume 10X TURBO DNase Buffer and 1 µL TURBO DNase to the RNA, and mix gently

There are separate DNase digestion conditions depending on the amount of contaminating DNA and the nucleic acid concentration of the sample.

- **Routine DNase treatment:** ≤200 µg nucleic acid per mL
- **Rigorous DNase treatment:** >200 µg nucleic acid per mL or RNA that is severely contaminated with DNA (i.e. >2 µg DNA/50 µL)

Routine DNase treatment: Use 1 µL TURBO™ DNase (2 U) for up to 10 µg of RNA in a 50 µL reaction. These reaction conditions will remove up to 2 µg of genomic DNA from total RNA in a 50 µL reaction volume.

Rigorous DNase treatment: If the RNA contains more than 200 µg of nucleic acid per mL, dilute the sample to 10 µg nucleic acid/50 µL before adding the TURBO DNase Buffer and TURBO™ DNase. It is also helpful to add only half of the TURBO™ DNase to the reaction initially, incubate for 30 min, then add the remainder of the enzyme and incubate for another 30 min.

If the sample cannot be diluted, simply increase the amount of TURBO DNase to 2–3 µL (4–6 U). It may be possible to successfully remove contaminating DNA from samples containing up to 500 µg/mL nucleic acid in a 10–100 µL TURBO DNA-free™ reaction. However, the efficacy of treating highly concentrated nucleic acid samples depends on the absolute level of DNA contamination, and residual DNA may or may not be detectable by PCR after 35–40 cycles.

2. Incubate at 37°C for 20–30 min

If only half of the TURBO DNase was added in the previous step, incubate for 30 min, then add the rest of the TURBO DNase and incubate for 30 min more.

3. Add resuspended DNase Inactivation Reagent (typically 0.1 volume) and mix well.

Always resuspend the DNase Inactivation Reagent by flicking or vortexing the tube before dispensing it.

- **For routine DNase treatment:** use 2 μ L or 0.1 volume DNase Inactivation Reagent, whichever is greater. For example, if the RNA volume is 50 μ L, and 1 μ L of TURBO DNase was used in the previous step, add 5 μ L of DNase Inactivation Reagent.
- **For rigorous DNase treatments:** where 2–3 μ L of TURBO DNase was used, add 0.2 volumes of DNase Inactivation Reagent.

IMPORTANT! Always use at least 2 μ L of DNase Inactivation Reagent, even if it is more than 0.1 volume.

Note: The DNase Inactivation Reagent may become difficult to pipette after multiple uses due to depletion of fluid from the interstitial spaces. If this happens, add a volume of Nuclease-free Water (supplied with the kit) equal to approximately 20–25% of the bed volume of the remaining DNase Inactivation Reagent, and vortex thoroughly to recreate a pipettable slurry.

4. Incubate 5 min at room temperature, mixing occasionally.

Flick the tube 2–3 times during the incubation period to redisperse the DNase Inactivation Reagent.

Note: If room temperature cools below 22–26°C, move the tubes to a heat block or oven to control the temperature. Cold environments can reduce the inactivation of the TURBO DNase, leaving residual DNase in the RNA sample.

5. Centrifuge at 10,000 \times g for 1.5 min and transfer the RNA to a fresh tube

Flick the tube 2–3 times during the incubation period to redisperse the DNase Inactivation Reagent.

- For 96-well plates, centrifuge at 2000 \times g for 5 min.

This centrifugation step pellets the DNase Inactivation Reagent. After centrifuging, carefully transfer the supernatant, which contains the RNA, into a fresh tube. Avoid introducing the DNase Inactivation Reagent into solutions that may be used for downstream enzymatic reactions, because it can sequester divalent cations and change the buffer conditions.

Troubleshooting

Observation	Possible cause	Recommended action
No RT-PCR product is detectable from RNA treated with TURBO DNA-free™	DNase Inactivation Reagent could inhibit RT-PCR	In step 5, remove the RNA solution from the pelleted DNase Inactivation Reagent carefully to avoid transferring it to the tube of RNA. You may have to leave a small amount of RNA behind to accomplish this. If you accidentally touch the pellet while removing the RNA, recentrifuge to pack the DNase Inactivation Reagent.
	TURBO DNA-free™ treated RNA should comprise only ~20% of an RT-PCR reaction volume.	For RT-PCR, we recommend that TURBO DNA-free™ treated RNA makes up ~20%, and no more than 40%, of the final RT-PCR volume. Otherwise, components from the TURBO DNase Buffer and the DNase Inactivation Reagent could interfere with the reaction. If necessary, RT-PCR volumes can be increased to 50 µL or more to accommodate your RNA without exceeding the 20–40% limit.
	RNA used in RT-PCR should be treated only once with TURBO DNA-free™	The salt in TURBO DNA-free™ reactions is carefully balanced for optimal TURBO DNase activity. Subjecting RNA to a second TURBO DNA-free™ treatment will introduce additional salts that could interfere with downstream enzymatic reactions.
RNA is degraded upon heating to >60°C	RNA samples that contain divalent cations, such as magnesium or calcium, will degrade when heated to temperatures above 60°C.	To ensure that divalent cations are removed in step 4, redisperse the DNase Inactivation Reagent by mixing the reaction 2–3 times during the incubation period.
The RNA absorbance spectrum has an unusual profile after treatment with TURBO DNA-free™.	If the concentration of RNA in the sample is less than about 50 ng/µL, you may notice significant absorbance at ~230 nm. A_{260}/A_{280} ratios may also be slightly lower than normal when the RNA concentration is ≤ 25 ng/µL. These differences in the absorbance profile are caused by the enhancer in the TURBO DNase Buffer. Exhaustive comparisons with both treated and untreated RNA samples indicate that the enhancer has no effect on accurate RNA quantification unless the RNA concentration is below 10 ng/µL.	

Quality control

Functional testing

The activity of the TURBO™ DNase is tested functionally in a unit activity assay. One unit is defined as the amount of enzyme required to completely degrade 1 µg DNA in 10 min at 37°C. Results are analyzed by agarose gel electrophoresis. The DNase Inactivation Reagent is tested for its ability to remove both TURBO™ DNase and TURBO™ DNase Buffer components. Results are analyzed by agarose gel electrophoresis and the Agilent 2100 bioanalyzer, respectively.

Nuclease testing

Relevant kit components are tested in the following nuclease assays:

RNase activity

A sample is incubated with labeled RNA and analyzed by PAGE.

Nonspecific endonuclease activity

A sample is incubated with supercoiled plasmid DNA and analyzed by agarose gel electrophoresis.

Exonuclease activity

A sample is incubated with labeled double-stranded DNA, followed by PAGE analysis.

Protease testing

A sample is incubated with protease substrate and analyzed by fluorescence.

Appendix A Safety

Chemical safety



WARNING! GENERAL CHEMICAL HANDLING. To minimize hazards, ensure laboratory personnel read and practice the general safety guidelines for chemical usage, storage, and waste provided below, and consult the relevant SDS for specific precautions and instructions:

- Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. To obtain SDSs, see the “Documentation and Support” section in this document.
 - Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing).
 - Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood).
 - Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer's cleanup procedures as recommended in the SDS.
 - Handle chemical wastes in a fume hood.
 - Ensure use of primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary container. Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)
 - After emptying a waste container, seal it with the cap provided.
 - Characterize (by analysis if necessary) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
 - Ensure that the waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.
 - **IMPORTANT!** Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.
-

Documentation and support

Obtaining SDSs Safety Data Sheets (SDSs) are available from www.lifetechnologies.com/support.

Note: For the SDSs of chemicals not distributed by Life Technologies, contact the chemical manufacturer.

Obtaining support For the latest services and support information for all locations, go to:

www.lifetechnologies.com/support

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)
- 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

Limited product warranty

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



For Research Use Only. Not for use in diagnostic procedures.

The information in this guide is subject to change without notice.

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.

NOTICE TO PURCHASER: LIMITED USE LABEL LICENSE: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

TRADEMARKS

The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners. TaqMan® is a registered trademark of Roche Molecular Systems, Inc., used under permission and license. LabChip is a registered trademark of Caliper Life Sciences Inc. Bioanalyzer is a trademark of Agilent Technologies, Inc.

© 2012 Life Technologies Corporation. All rights reserved.

Headquarters

5791 Van Allen Way | Carlsbad, CA 92008 USA

Phone +1 760 603 7200 | Toll Free in USA 800 955 6288

For support visit lifetechnologies.com/support or email techsupport@lifetech.com

lifetechnologies.com

05 October 2012

