

Certificate of Analysis

RQ1 RNase-Free DNase (Cat.# M6101):

Contents

Part No.	Name	Size (units)
M610A	RQ1 RNase-Free DNase	1,000
M199A	Stop Solution	1ml
M198A	RQ1 DNase 10X Reaction Buffer	1ml

Storage Conditions: Store at -20°C . Avoid exposure to frequent temperature changes. See the expiration date on the Product Information Label.

Description: RQ1 (RNA-Qualified) RNase-Free DNase is a DNase I (endonuclease) that degrades both double-stranded and single-stranded DNA, producing 3'-OH oligonucleotides (1). (RQ1 RNase-Free DNase may be used in applications where maintaining RNA integrity is critical.) This DNase is suited for applications such as nick translation (2), production of random fragments (3), cleavage of genomic DNA for footprinting (3), removal of DNA template after in vitro transcription (4) and removal of DNA from RNA samples prior to applications such as RT-PCR (5).

Unit Definition: One unit of RQ1 RNase-Free DNase is defined as the amount required to completely degrade $1\mu\text{g}$ of lambda DNA in 10 minutes at 37°C in $50\mu\text{l}$ of a buffer containing 40mM Tris-HCl (pH 7.9), 10mM NaCl, 6mM MgCl_2 and 10mM CaCl_2 . See the unit concentration on the Product Information Label.

Biological Source: Bovine pancreas.

Requirement: Ca^{2+} and Mg^{2+} or Mn^{2+} (6).

Inhibitors: EGTA; EDTA (6); salt concentrations $>100\text{mM}$ will reduce DNase activity.

RQ1 DNase 10X Reaction Buffer (M198A): 400mM Tris-HCl (pH 8.0), 100mM MgSO_4 and 10mM CaCl_2 .

Stop Solution (M199A): 20mM EGTA (pH 8.0).

Enzyme Storage Buffer: RQ1 DNase is supplied in 10mM HEPES (pH 7.5), 50% glycerol (v/v), 10mM CaCl_2 and 10mM MgCl_2 .

Usage Notes:

1. This DNase solution does not contain an RNase inhibitor. Observe caution in handling the product to avoid contaminating it with RNase.
2. Under different buffer conditions the amount of DNase required to completely digest a given amount of DNA may need to be empirically determined. For example, salt concentrations $>100\text{mM}$ will reduce DNase activity.
3. To inactivate, heat for 10 minutes at 65°C in the presence of Stop Solution.

Quality Control Assays

Contaminant Activity

RNase Assay: To test for RNase activity, 50ng of $[^3\text{H}]\text{RNA}$ is incubated with 5 units of RQ1 RNase-Free DNase in Transcription Optimized 1X Buffer (Cat.# P1181, diluted fivefold) for 1 hour at 37°C , and the release of radiolabeled nucleotides is monitored by scintillation counting of TCA-soluble material. The specification is $<3\%$ release.

Part# 9PIM610
Revised 12/16



AF9PIM610 1216M610



Promega

Promega Corporation

2800 Woods Hollow Road	
Madison, WI 53711-5399	USA
Telephone	608-274-4330
Toll Free	800-356-9526
Fax	608-277-2516
Internet	www.promega.com

PRODUCT USE LIMITATIONS, WARRANTY, DISCLAIMER

Promega manufactures products for a number of intended uses. Please refer to the product label for the intended use statements for specific products. Promega products contain chemicals which may be harmful if misused. Due care should be exercised with all Promega products to prevent direct human contact.

Each Promega product is shipped with documentation stating specifications and other technical information. Promega products are warranted to meet or exceed the stated specifications. Promega's sole obligation and the customer's sole remedy is limited to replacement of products free of charge in the event products fail to perform as warranted. Promega makes no other warranty of any kind whatsoever, and SPECIFICALLY DISCLAIMS AND EXCLUDES ALL OTHER WARRANTIES OF ANY KIND OR NATURE WHATSOEVER, DIRECTLY OR INDIRECTLY, EXPRESS OR IMPLIED, INCLUDING, WITHOUT LIMITATION, AS TO THE SUITABILITY, PRODUCTIVITY, DURABILITY, FITNESS FOR A PARTICULAR PURPOSE OR USE, MERCHANTABILITY, CONDITION, OR ANY OTHER MATTER WITH RESPECT TO PROMEGA PRODUCTS. In no event shall Promega be liable for claims for any other damages, whether direct, incidental, foreseeable, consequential, or special (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including negligence) or strict liability arising in connection with the sale or the failure of Promega products to perform in accordance with the stated specifications.

Riboprobe is a registered trademark of Promega Corporation.

© 1996–2016 Promega Corporation. All Rights Reserved.

All specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.

Signed by:

R. Wheeler, Quality Assurance

Part# 9PIM610
Printed in USA. Revised 12/16

I. DNase Treatment of RNA Samples Prior to RT-PCR

1. Set up the DNase digestion reaction as follows:

RNA in water or TE buffer	1–8µl
RQ1 RNase-Free DNase 10X Reaction Buffer	1µl
RQ1 RNase-Free DNase	1µ/µg RNA
Nuclease-free water to a final volume of	10µl

Note: Use 1 unit of RQ1 RNase-Free DNase per microgram of RNA. For smaller amounts of RNA, use 1 unit of RQ1 RNase-Free DNase per reaction.

2. Incubate at 37°C for 30 minutes.

Note: If analyzing RNA samples by gel electrophoresis, perform a phenol:chloroform extraction and ethanol precipitation before loading the samples on the gel because salts in the RQ1 DNase Reaction Buffer and Stop Solution may cause aberrant migration or smearing of RNA on gels. Steps 3 and 4 may be omitted if a phenol:chloroform extraction is performed.

3. Add 1µl of RQ1 DNase Stop Solution to terminate the reaction.
4. Incubate at 65°C for 10 minutes to inactivate the DNase.
5. Add all, or a portion of, the treated RNA to the RT-PCR. See the *Access RT-PCR System Technical Bulletin #TB220* (5).

Note: The RQ1 RNase-Free DNase digestion contains a final concentration of 10mM MgSO₄. When adding DNase-treated RNA to an RT-PCR, carryover of magnesium must be considered. For example, adding 1µl of treated RNA to a 50µl RT-PCR will raise the magnesium concentration by 0.2mM, and adding 5µl of treated RNA will raise the magnesium concentration by 1mM. The requirement for magnesium may be different in the RQ1 DNase digestion step and the amplification reaction.

- RQ1 DNase activity increases as Mg²⁺ concentration increases up to 5–10mM. At a concentration of 1mM Mg²⁺, RQ1 DNase is expected to be at least fourfold less active than at the optimal Mg²⁺ concentration.
- For some templates, the yield from the amplification reaction is highly dependent on Mg²⁺ concentration, and the optimal Mg²⁺ concentration may be as low as 1mM.

If an increased Mg²⁺ concentration is not tolerable in the amplification reaction, the following alternatives may be used.

- The RQ1 RNase-Free DNase 10X Reaction Buffer may be diluted 1:10 with 400mM Tris (pH 8.0), 10mM CaCl₂ prior to DNase digestion. (Note that under these conditions, the RQ1 DNase will be approximately fourfold less active than under standard reaction conditions.)
- An alternative DNase reaction buffer may be used (such as the RT or PCR reaction buffer) if that buffer contains at least 1mM Mg²⁺.
- The RNA sample may be diluted in water prior to RT-PCR, allowing dilution of the MgSO₄ to a concentration that is compatible with this application.
- The RNA may be purified with a standard phenol:chloroform extraction followed by an ethanol precipitation.

II. Other Applications

RQ1 RNase-Free DNase may be used in a number of other applications where maintaining RNA integrity is important. These include in vitro transcription, nick translation and DNase I footprinting.

A. In Vitro Transcription

To remove template DNA, RQ1 RNase-Free DNase may be added directly to the transcription reaction. Refer to the *Riboprobe® in vitro Transcription Systems Technical Manual #TM016* (4) for specific protocol information.

B. Nick Translation

For protocol information on the use of RQ1 RNase-Free DNase for this application, refer to the *Protocols and Application Guide* (2).

C. Production of Random Fragments

For protocol information on the use of DNase I for random fragmentation of DNA, see reference 7.

Note: Under different buffer conditions, the amount of RQ1 RNase-Free DNase required to completely digest a given amount of DNA must be empirically determined. For example, salt concentrations >100mM will result in reduced DNase activity. Ca²⁺ and Mg²⁺ are essential for RQ1 DNase activity.

III. References

1. Moore, S. (1981) Pancreatic DNase In: *The Enzymes*, Volume 14A, P.D. Boyer, Ed., Academic Press, New York, 281.
2. *Protocols and Applications Guide* (1996) Promega Corporation.
3. Cobianchi, F. and Wilson S.H. (1987) *Meth. Enzymol.* **152**, 94–110.
4. *Riboprobe® in vitro Transcription Systems Technical Manual #TM016*, Promega Corporation.
5. *Access RT-PCR System Technical Bulletin #TB220*, Promega Corporation.
6. Ausubel, F.M. (1994) *Current Protocols in Molecular Biology*, John Wiley and Sons, New York, 3.12.
7. Sambrook, J., Fritsch, E.F. and Maniatis, T. (1989) *Molecular Cloning: A Laboratory Manual* 2nd ed., Cold Spring Harbor Laboratory, Cold Spring Harbor, New York.