

ezDNase

Catalog Number 11766051

Pub. No. MAN0015899 Rev. B



WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](https://www.thermofisher.com/support).

Product description

Invitrogen™ ezDNase™ Enzyme is a recombinant DNase for the fast removal of contaminating genomic DNA from RNA preparations. It eliminates double-stranded DNA specifically by cleaving phosphodiester bonds to yield 2–8 bp oligonucleotides with 5'-phosphate and 3'-hydroxyl termini.

The high specificity of ezDNase™ Enzyme for double-stranded DNA enables efficient and fast genomic DNA removal without reduction in the quality or quantity of RNA or single-stranded DNA present in the reaction such as cDNA and primers. ezDNase™ Enzyme is heat-labile and so can be easily deactivated by heat treatment at moderate temperature (55°C). These features make ezDNase™ Enzyme a superior choice for genomic DNA removal prior to reverse transcription reactions.

Contents and storage

Contents	Amount	Storage
ezDNase™ Enzyme	50 µL	Store at –20°C
10X ezDNase™ Enzyme Buffer	100 µL	
Nuclease-free Water	1.25 mL	

Source

The enzyme is purified from a *Pichia pastoris* strain expressing the recombinant DNase.

Unit definition

One unit increases the absorbance at 260 nm by 0.001 OD per min at 25°C and pH 5.0 with large molecular weight DNA as the substrate according to the assay method of Kunitz.

Storage buffer

The enzyme is supplied in 20 mM Tris-HCl (pH 7.5 at 25°C), 2 mM MgCl₂, 10 mM NaCl, 0.01% (v/v) Tween 20, 50% (v/v) glycerol.

Inhibition and inactivation

- The enzyme is inhibited by metal chelators, transition metals (e.g., Zn) in millimolar concentrations, reducing agents (DTT and β-mercaptoethanol), SDS (even at concentrations <0.1%), and high salt concentration.
- Incubate at 55°C for 5 minutes in the presence of 10 mM DTT to inactivate the enzyme.

Procedural guidelines

- Use RNase-free tubes, gloves and filter tips.
- If preparing several parallel reactions, use a master mix of water, buffer and ezDNase™ Enzyme to minimize the possibility of pipetting errors.

Two-step gDNA digestion and RT PCR protocol

The following protocol allows gDNA digestion and cDNA synthesis to be combined in a two-step procedure within a single reaction tube. Ensure that the reaction volume does not exceed the allowable volume limits of the RT reaction.

1. Mix the following components in a RNase-free tube.

Component	Volume
10X ezDNase™ Enzyme Buffer	1 µL
ezDNase™ Enzyme	1 µL
Template RNA	up to 8 µL
Nuclease-free Water	to 10 µL
Total volume	10 µL

2. Gently mix the samples then centrifuge the tube.
3. Incubate the sample for 2 minutes at 37°C.
4. (*Optional*) If the RNA sample is to be used for RT-PCR of fragments ≥3 kb, incubate the sample at 55°C for 5 minutes in the presence of 10 mM DTT to inactivate the enzyme.
5. Chill the tube on ice to bring the sample to room temperature, then briefly centrifuge and place the tube on ice.
6. Add reverse transcription reagents directly to the same tube and proceed with first strand cDNA synthesis protocol according to the manufacturer's instructions for your cDNA synthesis kit.

The product of the first strand cDNA synthesis can be used directly in PCR and qPCR or stored at -20°C for up to one week. For longer storage, -70°C is recommended. Avoid freeze/thaw cycles of cDNA.

Troubleshooting

Observation	Possible cause	Recommended action
PCR product is observed in the qPCR negative control reaction	The gDNA contamination was incompletely digested in the RNA sample.	Increase 37°C incubation step up to 5 minutes.
	Trace amounts of contaminants (SDS, EDTA, guanidine salts, phosphate, pyrophosphate, polyamines, spermidine) from the RNA purification process are inhibiting the DNase.	Re-precipitate the RNA with ethanol, wash the pellet with 75% ethanol then dissolve in nuclease-free water.

Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have questions, contact Life Technologies at www.thermofisher.com/support.

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2016-2025 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.