

Product Information

RNase A

Catalog Number: 99871-600UL

Unit Size: 600 uL

Concentration

Supplied at 10 mg/mL (≥ 50 U/mg (Kunitz units)) in 40 mM sodium acetate, 500 mM trehalose, and 20% glycerol.

Storage and Handling

Store at 4°C. Product is stable for at least 12 months from date of receipt when stored as recommended.

Color and Form: Clear solution

Product Description

RNase A may be useful for removing RNA from plasmid or genomic DNA preparation, recombinant protein preparations, or for use in ribonuclease protection assays. Biotium recommends this RNase A for use as a control in the RNaseReveal™ Activity Assay Kit (Cat. No. 31086). RNase A is provided as a 10 mg/mL solution and is DNase-free.

Related Products

Cat. No.	Product
31086	RNaseReveal™ Activity Assay Kit
22028	RNase-X™ Decontamination Solution
31073	AccuBlue® Broad Range RNA Quantitation Kit
31086	RNaseReveal™ Activity Assay Kit
CD507	CELLDATA DNASTORM™ 2.0 FFPE DNA Extraction Kit
31066	AccuGreen™ High Sensitivity dsDNA Quantitation Kit
31028	AccuClear® Ultra High Sensitivity dsDNA Quantitation Kit with DNA Standards
41029-5G	GelRed® Agarose LE
41003	GelRed® Nucleic Acid Gel Stain 10,000X in Water
31065	RNase-Free Calf Thymus DNA
31030-50	DNA Gel Extraction Kit

Please visit our website at www.biotium.com for information on our life science research products, including kits for RNA extraction and quantitation, stains for DNA or RNA gel electrophoresis, environmentally friendly EvaGreen® qPCR master mixes, and other products for molecular biology research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.