



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

RNase A

A 10 mg/mL RNase solution, DNase-free. Useful for removing RNA from plasmid or genomic DNA preparation, recombinant protein preparations, or for use in ribonuclease protection assays.



Product attributes

Storage Conditions

Store at 2 to 8 °C

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 (≥50 U/mg (Kunitz units)) solution and is DNase-free.

Features

- 10 mg/mL RNase solution, DNase-free
- Removes RNA from DNA or protein preparations
- Suitable for ribonuclease protection assays
- For use with [RNaseReveal™ Activity Assay Kit \(Cat. No. 31086\)](#)

[Learn more about Biotium's unique products for RNA workflows, including RNA extraction kits, quantitation kits, and RNA stains for gel electrophoresis.](#)

This datasheet was generated on January 7, 2026 at 02:08:56 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/rnase/>