



Glowing products for science

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

## RNase-X™ Decontamination Solution

A decontamination solution that effectively removes RNases and other contamination from surfaces and instruments for RNA research.



### Product attributes

## Product Description

Working with RNA samples can be very difficult due to the ubiquitous presence of RNase contamination on benchtops, pipettors, and other surfaces. RNases are notoriously stable and resistant to many heat and chemical treatments. Therefore it is crucial to clean work surfaces and implements with a cleanser that is proven to eliminate RNases.

RNase-X™ Decontamination Solution protects precious RNA samples by effectively eliminating RNase contamination on working surfaces such as benchtops, pipets, tip boxes, and instruments. It is provided in a convenient spray bottle: simply spray the surface, wipe, and rinse with RNase-free water or 70% ethanol.

- Effectively eliminates RNase contamination on work surfaces
- Ideal for cleaning benchtops, pipettors, and instruments that must be RNase-free
- Provided in a convenient and ready-to-use spray bottle
- Excellent performance and more cost-effective than RNaseZap™

### Other RNA Research Solutions

For more RNA research tools and solutions, please view our [AccuBlue® Broad Range RNA Quantitation Kit](#), [EMBER500™ RNA Prestain Loading Dye](#), and our [RNA extraction kits for FFPE tissues or fresh tissues and cells](#).

This datasheet was generated on January 7, 2026 at 05:29:39 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/rnase-x-decontamination-solution/>