

Product Information

RNase-X™ Decontamination Solution

Catalog Number: 22028

Unit Size: 250 mL

Storage and Handling

Store at room temperature. Product is stable for at least 12 months from date of receipt when stored as recommended.

Always wear gloves and exercise universal laboratory safety precautions during handling. Prolonged skin contact with RNase-X™ Decontamination Solution may cause irritation. Use in a well-ventilated area and take care to avoid breathing the aerosol which may cause irritation to lungs and mucous membranes. RNase-X™ Decontamination Solution may corrode metal surfaces, so care should be taken to rinse treated surfaces well with water after use.

Product Description

RNase-X™ Decontamination Solution is a ready-to-use cleaning agent for effective removal of RNase and other contaminants from working surfaces. The solution features an optimized formulation of key ingredients for maximum removal of RNase to protect precious RNA samples during purification, quantitation, amplification, and other analytical workflows. The solution may be applied to benchtops, pipets, tip boxes, and instruments as well as the outside of reaction vessels such as microfuge tubes. RNase-X™ is provided in a convenient spray bottle and performs as well as RNaseZap™ and other common decontamination reagents.

Instructions for Use

RNase-X™ Decontamination Solution is ready-to-use. Simply spray the surface, wipe, and rinse with RNase-free ultrapure water or 70% ethanol. See below for detailed instructions on cleaning specific surfaces or items.

Cleaning work surfaces

Apply RNase-X™ Decontamination Solution directly to the work area to be cleaned and wipe thoroughly with a laboratory wipe. Rinse with ultrapure water. Wipe with 70% ethanol (optional), and then wipe dry.

Cleaning instruments

Apply RNase-X™ Decontamination Solution to a laboratory wipe and wipe all exposed surfaces of the instrument thoroughly. Rinse with ultrapure water, then wipe with 70% ethanol (optional), then wipe dry. Small parts may be cleaned by briefly soaking them in RNase-X™, followed by rinsing with ultrapure water. Wipe with 70% ethanol (optional), and then wipe dry.

Cleaning pipettes

Apply RNase-X™ Decontamination Solution to a laboratory wipe and wipe all exposed surfaces of the pipettor thoroughly. If applicable, remove ejector according to manufacturer's instructions to clean separately. Rinse thoroughly with ultrapure water. Wipe with 70% ethanol (optional), dry, and reattach ejector if needed. Alternatively, the shaft of the pipettor may be removed and briefly soaked in RNase-X™ to decontaminate, followed by rinsing and drying as described above. Do not expose inner gaskets or seals to RNase-X™.

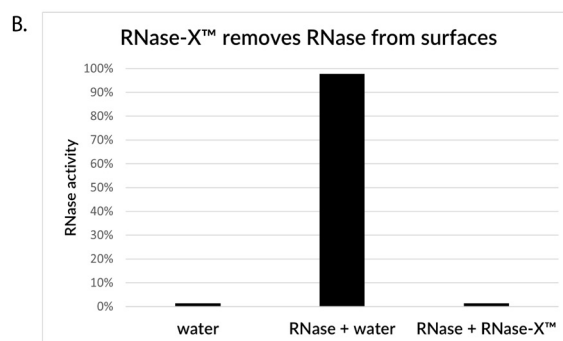
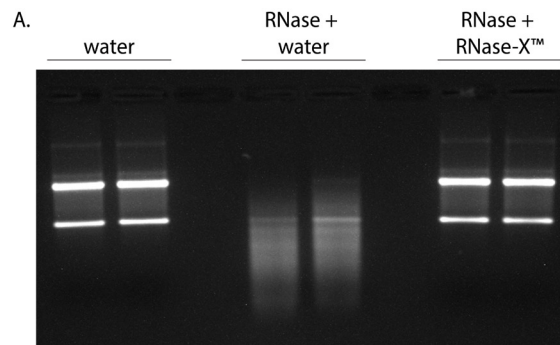


Figure 1. RNase-X™ effectively removes RNase from surfaces. To simulate RNase-contaminated surfaces, the insides of microfuge tubes were left untreated (water), or coated with RNase A, then rinsed with either water or RNase-X™, followed by water washes. A. The final water washes were incubated with total human RNA, which was then stained with EMBER500™ RNA Prestain Loading Dye and analyzed on an agarose gel. First two lanes: RNA incubated with water alone. Middle two lanes: RNA incubated with the water from RNase-treated, water-rinsed tube. Last two lanes: RNA incubated with the water from RNase-treated, RNase-X™-rinsed tube. B. Analysis of RNase activity in the final water washes using the RNaseAlert™ assay from IDT.

Related Products

Catalog number	Product
CD501	RNastorm™ Kit for Isolation of RNA from FFPE Tissue Samples
CD504	RNastorm™ RNA Isolation Kit
41032	EMBER500™ RNA Prestain Loading Dye
31073	AccuBlue® Broad Range RNA Quantitation Kit
99850	RNA Broad Range Standard, 100 ng/μL
31065	RNase-free Calf Thymus DNA, 1 mg/mL
41024	Water, Ultrapure Molecular Biology Grade, RNase-Free

Please visit our website at www.biotium.com for information on our products for RNA research and applications, including RNA extraction kits for fresh cells and FFPE tissues, RNA quantitation kits, and RNA gel stains.

RNaseZap is a trademark of Thermo Fisher Scientific; RNaseAlert is a trademark of Integrated DNA Technologies.

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