

RNase & DNase Cleaner

Catalogue number	Presentation
MB46301	250 mL
MB46302	500 mL (Refill)

Features

- Ready-to-use solution specifically designed to remove nucleases from stainless-steel surfaces, glassware, plastics, laboratory equipment, and pipettors.
- The cleaner comes in a user-friendly, lightweight bottle equipped with a trigger spray for convenient application.
- Using this cleaner simplifies the decontamination process, saving both time and effort.
- A refill option is available for added convenience.

Description

RNase & DNase Cleaner is the ultimate solution for effectively eliminating both RNase and DNase enzymes from laboratory surfaces, instruments, and apparatus. It is particularly suitable for decontaminating pipettes, gel boxes, as well as RNA or DNA preparation areas. By incorporating RNase & DNase Cleaner into your laboratory protocols, you can enhance the accuracy and reliability of your results with utmost confidence. Our advanced formula is specifically designed to target and eliminate the ubiquitous contaminant enzymes that degrade nucleic acids, providing you with a highly effective approach to ensure the integrity of your molecular biology techniques, resulting in trustworthy and reproducible results.

The RNase & DNase Cleaner is meticulously designed to meet the demanding requirements of a wide range of molecular biology techniques, including highly sensitive qPCR assays. It is particularly advantageous for labware that cannot undergo autoclaving, eliminating the need for time-consuming procedures such as DEPC treatments.

Extensive testing and validation have been conducted on the RNase & DNase Cleaner to ensure optimal performance in diverse laboratory settings. The effectiveness of our product is scientifically proven and supported by robust evidence, offering you the assurance you need for critical decontamination requirements.

Storage temperature

Store at 15 °C to 25 °C.

Protocol of usage

Use gloves as personal protective equipment when manipulating this product.

Decontamination of surfaces and laboratory devices:

1. Apply RNase & DNase Cleaner onto a paper towel.
2. Vigorously rub the surfaces, appliances, or equipment with the soaked paper towel.
3. Optionally, rinse well with DNA-free water.
4. Thoroughly dry the surface with a clean paper towel.

Note: Prior to use, it is recommended to perform a surface compatibility test.

Decontamination of pipettors:

1. Follow the manufacturer's instructions to disassemble the pipettor.
2. Remove gaskets and seals, and immerse the stem of the pipette in RNase & DNase Cleaner for 1 min.
3. Rinse thoroughly in DNA-free water.
4. Ensure the pipette is completely dry before reassembling it.

Please note that while our recommendations provide a robust decontamination solution, we encourage users to verify the suitability of RNase & DNase Cleaner for their specific purposes and applications. Do not add the cleaner to the reaction.

For life science research only. Not for use in diagnostic procedures.