

Nucleases & Nucleic Acid Cleaner

Catalogue number	Presentation
MB48301	250 mL
MB48302	500 mL (Refill)

Features

- All-in-one ready-to-use solution designed to remove nucleases as well as nucleic acids 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

Nucleases & Nucleic Acid Cleaner is cutting-edge solution for daily laboratory operations involving nucleic acids manipulation and highly sensitive enzymatic reactions. This all-in-one ready-to-use solution does not just remove RNase and DNase enzymes, as it also wipes out unwanted DNA and RNA guaranteeing the utmost purity of laboratory surfaces, instruments, and equipment. By incorporating this cleaner into your laboratory protocols, you can enhance the accuracy and reliability of your results with utmost confidence.

The Nucleases & Nucleic Acid Cleaner is purpose-built to meet the rigorous demands of a broad spectrum of molecular biology techniques, including the most sensitive qPCR, LAMP assays, and other DNA-centric methods commonly embraced by cutting-edge research and diagnostic laboratories. It is also particularly advantageous for labware that cannot undergo autoclaving, eliminating the need for time-consuming procedures such as DEPC treatments typically used for RNase removal.

Extensive testing and validation have been conducted on the Nucleases & Nucleic Acid 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 Nucleases & Nucleic Acid 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 Nucleases & Nucleic Acid 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 Nucleases & Nucleic Acid 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.