

## DNA & RNA Cleaner

<b>Catalogue number</b>	<b>Presentation</b>
MB46201	250 mL
MB46202	500 mL (Refill)

### Features

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

### Description

The DNA & RNA Cleaner by NZYtech offers a comprehensive solution for combating DNA contamination and eradicating unwanted DNA and RNA molecules from various 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. Our advanced formula is specifically designed to target and eliminate nucleic acid contaminants, providing you with a highly effective approach to ensure the integrity of your molecular biology techniques, resulting in trustworthy and reproducible outcomes.

The DNA & RNA Cleaner is meticulously designed to meet the demanding requirements of a wide range of molecular biology techniques, including highly sensitive qPCR assays. It is compatible with PCR, qPCR, RT-qPCR, LAMP, RT-LAMP, and other DNA-centric methods commonly employed in research and diagnostic laboratories. The risk of genetic material contamination during PCR amplification, a crucial technique in DNA-based detection workflows, is well-recognized. Even a single undesired DNA molecule (contaminant) can undergo exponential amplification, resulting in millions of identical copies during the PCR process. DNA & RNA Cleaner effectively tackles this challenge, ensuring optimal decontamination and allowing you to focus on the core aspects of your work with enhanced efficiency and productivity.

Extensive testing and validation have been conducted on the DNA & RNA 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 DNA & RNA 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 DNA & RNA Cleaner for 1 minute.
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 DNA & RNA 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.