

Molecular Grade Water

Catalog No: 37501, 37601

Format: 500 ml, 1 L

Background: Active Motif's ultra-pure molecular grade water is specifically treated to remove contaminants such as nucleases (DNases and RNases), proteins, salts, organic compounds, and microorganisms. Water is not only a universal solvent but also an essential reagent in molecular biology due to its ability to dissolve a wide range of polar molecules and salts. By providing a clean, stable medium, molecular grade water allows for accurate and reproducible results in applications such as PCR, sequencing, and enzymatic assays, where even trace impurities could compromise experimental outcomes.

Formula: H₂O

Formula Weight: 18.02 g/mol

Quality Control: This product is 0.2 µm filtered. It is designated as a Molecular Biology Reagent and is suitable for molecular biology applications. This product has been analyzed for the absence of nucleases and proteases, and has undergone bioburden analysis. It has not been treated with diethyl pyrocarbonate (DEPC).

Precautions and Disclaimer:

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

It is not recommended to autoclave the bottles in which this product is packaged.

Storage: Some products may be shipped at room temperature. This will not affect their stability or performance.

Guarantee: This product is guaranteed for 24 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

Once opened, it is recommended to use within 6 months to maintain optimal purity.

Test	Specification	Result
Appearance (Color)	Colorless	Colorless
Appearance (Form)	Liquid	Liquid
Extraneous Activities	Tested	None Detected
Dnases, Rnases, Proteases	Tested	None Detected
Bioburden aerobic / anaerobic count	<10 total CFU	Passed
Conductivity (at 25°C)	≤ 5.0 µS/cm	Passed