

Product Information

Water, Ultrapure Molecular Biology Grade

Catalog Number: 41024-4L

Unit Size: 4L Cubitainer® with spigot

Storage and Handling

Store at room temperature. Product is stable for at least 2 years from date of receipt when stored as recommended.

Product Description

Water, Ultrapure Molecular Biology Grade is suitable for use in molecular biology applications that require high quality water. Type I Milli-Q (18 megOhm) water is subjected to ultrafiltration to eliminate RNase, DNase, protease, pyrogens/endotoxins (< 0.001 EU/mL) and bacteria (< 1 cfu/mL). Each lot is quality tested for the absence of RNase, DNase and protease contamination. The product is not certified as sterile, and is therefore not recommended for cell culture. Water, Ultrapure Molecular Biology Grade is free of toxic chemical additives, such as diethyl pyrocarbonate (DEPC).

Applications

- Buffer preparation
- Protein and nucleic acid electrophoresis
- DNA and RNA preparation
- Southern and northern blotting
- PCR, DNA sequencing, cloning, and more

Related Products

Catalog number	Product
41003	GelRed™ Nucleic Acid Gel Stain
41005	GelGreen™ Nucleic Acid Gel Stain
31042	Forget-Me-Not™ EvaGreen® qPCR Master Mix
31028	AccuClear Ultra High Sensitivity dsDNA Quantitation Kit with 7 DNA Standards
21003-1L	One-Step Blue® Protein Gel Stain
22020	10X Phosphate-Buffered Saline (PBS)
41006	TBE Buffer, 5X
22014	Bovine Serum Albumin (BSA) 30% solution
22013	Bovine Serum Albumin (BSA) Fraction V
22012	Dry Milk Powder
22011	Fish Gelatin Powder
90082	DMSO, anhydrous
22015	Fixation Buffer
22016	Permeabilization Buffer
22017	Permeabilization and Blocking Buffer (5X)
23006	Flow Cytometry Fixation/Permeabilization Kit
23001	EverBrite™ Mounting Medium
23002	EverBrite™ Mounting Medium with DAPI
23005	CoverGrip™ Coverslip Sealant
22002	Tween-20
99821	1X Passive Lysis Buffer 2.0 for Single-Tube Luciferase Kit
99912	5X Passive Lysis Buffer for Renilla Luciferase Kit
99923	5X Firefly Luciferase Lysis Buffer

Please visit our website at www.biotium.com for information on our life science research products, including fluorescent CF™ dye primary and secondary antibodies, environmentally friendly EvaGreen® qPCR master mixes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

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