

Product Information

10X Phosphate-Buffered Saline (PBS)

Catalog Number: 22020

Unit Size: 4L

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

10X Phosphate-Buffered Saline (PBS) is a concentrated buffer solution designed to be used at a final concentration of 1X. In the 1X working concentration the buffer contains 137 mM NaCl, 2.7 mM KCl, 10 mM Na₂HPO₄, 1.8 mM KH₂PO₄, at a pH of 7.4.

This molecular biology grade 10X PBS is suitable for a variety of molecular biology or cell biology applications. It has been tested to ensure it is RNase-, DNase- and protease-free. It is not designed for sterile cell culture, as it is not sterile-filtered.

Related Products

Catalog number	Product
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
22003	Mini Cell Scrapers
23001, 23002	EverBrite™ Mounting Medium
23003, 23004	EverBrite™ Hardset Mounting Medium
23005	CoverGrip™ Coverslip Sealant
23007	TrueBlack™ Lipofuscin Autofluorescence Quencher, 20X
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.