

10X Phosphate-Buffered Saline (PBS)

This item is discontinued.



Product attributes

Storage Conditions

Call us: 800-304-5357

Store at room temperature

Product Description

This item is discontinued. See our $\underline{\text{1X PBS (2L)}}$ Buffer Powder Packets as a replacement.

This datasheet was generated on November 22, 2025 at 09:00:45 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/10x-phosphate-buffered-saline-pbs/