



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## 10X Phosphate-Buffered Saline (PBS)

This item is discontinued.



### Product attributes

Storage Conditions
Store at room temperature

### Product Description

This item is discontinued. See our [1X PBS \(2L\) Buffer Powder Packets](#) as a replacement.

This datasheet was generated on June 20, 2026 at 04:43:34 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/10x-phosphate-buffered-saline-pbs/>