

1X PBS (2L) Buffer Powder Packets

PBS (phosphate-buffered saline) buffer powder packet for preparing 2 L of 1X PBS buffer. Box includes 50 packets.



Product attributes

Storage Conditions	Store at room temperature
Shelf life	Guaranteed for at least 12 months from date of receipt when stored as recommended

Product Description

PBS buffer (phosphate-buffered saline) is widely used buffer in biology and biochemical research. 1X PBS (2L) Buffer Powder Packets are a ready-to-use format for easy and reliable preparation of PBS buffer. Simply dissolve the contents in 2 L dH₂O to prepare 2 L of 1X PBS.

- Convenient buffer packet for preparing 2 L of 1X PBS buffer
- Ready-to-use, with no pH adjustments needed
- Stable at room temperature
- 1X buffer will be 137 mM NaCl, 2.7 mM KCl, 10 mM phosphate, pH ~7.3

Procedure

Prepare 1X PBS buffer by dissolving the contents of the buffer packet in 2 L of dH₂O. 10X PBS may also be prepared by dissolving the contents in 200 mL of dH₂O. The final composition of the 1X buffer will be 137 mM NaCl, 2.7 mM KCl, 10 mM phosphate, pH ~7.3 (do not adjust pH). PBS buffer may be stored at room temperature indefinitely as long as the solution remains clear.

Buffer Powder Packets

Product Name	Catalog Number	Size
1X TAE (1L) Buffer Powder Packets	22031	50 packets
1X TBE (1L) Buffer Powder Packets	22032	50 packets
1X PBS (2L) Buffer Powder Packets	22033	50 packets

Browse Biotium's wide range of [buffers, background reducers and other accessory products for research](#).

This datasheet was generated on June 5, 2026 at 02:54:08 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/1x-pbs-2l-buffer-powder-packets/>