

## 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

|                           |   |
|---------------------------|---|
| <b>Storage Conditions</b> | Store at room temperature   |
| <b>Shelf life</b>         | 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 dH<sub>2</sub>O 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 dH<sub>2</sub>O. 10X PBS may also be prepared by dissolving the contents in 200 mL of dH<sub>2</sub>O. 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       |
|---|-----------------------|------------|
| <a href="#">1X TAE (1L) Buffer Powder Packets</a> | <a href="#">22031</a> | 50 packets |
| <a href="#">1X TBE (1L) Buffer Powder Packets</a> | <a href="#">22032</a> | 50 packets |
| <a href="#">1X PBS (2L) Buffer Powder Packets</a> | <a href="#">22033</a> | 50 packets |

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

This datasheet was generated on January 15, 2025 at 05:33:11 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/1x-pbs-2l-buffer-powder-packets/>