



Glowing products for science

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

## 1X TBE (1L) Buffer Powder Packets

TBE (Tris-borate-EDTA) buffer powder packet for preparing 1L of 1X TBE buffer for nucleic acid gel electrophoresis. 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

TBE buffer (Tris-borate-EDTA) is widely used for electrophoresis of nucleic acids in agarose and polyacrylamide gels. The buffer may be used as running buffer for gel electrophoresis, or when casting agarose gels.

- Convenient buffer packet for preparing 1 L of 1X TBE buffer
- Ready-to-use, with no pH adjustments needed
- Stable at room temperature
- 1X buffer will be 89 mM Tris-borate, 2 mM EDTA, pH ~8

### Procedure

Prepare 1X TBE buffer by dissolving the contents of the buffer packet in 1 L of dH<sub>2</sub>O. The final composition of the 1X buffer will be 89 mM Tris, 89 mM boric acid, 2 mM EDTA, pH ~8 (do not adjust pH). 1X TBE buffer may be stored at room temperature indefinitely as long as the solution remains clear.

**Note:** We recommend using the same buffer for both casting and running gels for electrophoresis.

## 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

Learn more about Biotium's safer and more sensitive [GelRed® and GelGreen® nucleic acid gel stains](#), as well as other unique products for [DNA/RNA gel electrophoresis](#).

This datasheet was generated on February 4, 2026 at 03:45:15 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/1x-tbe-1l-buffer-powder-packets/>