

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 dH₂O. 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
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

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 January 15, 2025 at 05:43:52 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/1x-tbe-1l-buffer-powder-packets/>