

# 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.



#### Call us: 800-304-5357 Email: techsupport@biotium.com

#### 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**

<b>Product Name</b>	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	<u>22033</u>	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 October 30, 2025 at 11:53:28 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/1x-tbe-1I-buffer-powder-packets/">https://biotium.com/product/1x-tbe-1I-buffer-powder-packets/</a>