



Glowing products for science

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

1X TAE (1L) Buffer Powder Packets

TAE (Tris-acetate-EDTA) buffer powder packet for preparing 1 L of 1X TAE buffer for nucleic acid gel electrophoresis. Box includes 50 packets.



Product attributes

Product Description

TAE buffer (Tris-acetate-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. 1X TAE (1L) Buffer Powder Packets may be used as a running buffer for [Precast GelRed® Agarose TAE Gels](#).

- Convenient buffer packet for preparing 1 L of 1X TAE buffer
- Ready-to-use, with no pH adjustments needed
- Stable at room temperature
- 1X buffer will be 40 mM Tris-acetate, 1 mM EDTA, pH ~8

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 June 22, 2026 at 06:11:03 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/1x-tae-1l-buffer-powder-packets/>