



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
<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 January 2, 2026 at 09:05:25 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/1x-tae-1l-buffer-powder-packets/>