TBE Buffer, 5X

A 5X concentrated buffer solution in deionized water. TBE is used for polyacrylamide and agarose gel electrophoresis.



Product attributes

Storage Conditions

Store at room temperature

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

Product Description

5X TBE (Tris-Borate-EDTA) is a concentrated buffer solution in deionized water. TBE is used for polyacrylamide and agarose gel electrophoresis. 1X solution: 89 mM Tris, 89 mM boric acid, 2 mM EDTA.

- 4L Cubitainer® with detachable spigot for easy dispensing
- Non-glug design for a continuous, uninterrupted stream
- Opening can be raised or depressed for pouring

Biotium also offers ready-to-use <u>1X TBE (1L) Buffer Powder Packets</u> for quickly preparing 1L of 1X TBE buffer.

Learn about our <u>Go-GoTM Fast DNA Gel Running Buffer</u>, a novel low ionic strength running buffer that allows you to run DNA agarose gels at higher voltage to get results up to 3X faster than with TAE or TBE buffer, streamlining routine DNA gel analysis.

Cubitainer is a registered trademark of Hedwin Corporation.

This datasheet was generated on November 12, 2025 at 03:51:32 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/tbe-buffer-5x/