

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

## PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: TBE buffer, 5X

**CATALOG #** 41006

**DESCRIPTION:** TBE buffer is used for polyacrylamide and agarose gel electrophoresis. It is suited for

DNA and RNA separation on longer run, higher voltage and amperage gel. A 1X working solution is obtained by adding 1 part of the 5X TBE buffer to 4 parts of

deionized water

**1X COMPOSITION:** 89mM Tris base, 89mM Boric acid and 2mM EDTA

**STORAGE:** Store at room temperature. Gently heat the solution if a precipitate is formed.

FIRST AID: Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on

clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing,

move individual to fresh air and seek medical advice immediately.

**Disclaimer:** Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.