

**TBE Buffer - 10x conc.**

| Cat. No. | Amount |
|----------|--------|
| BU-120 | 1 l |

For in vitro use only!

Shipping: shipped at ambient temperature

Storage Conditions: store at 4 °C

Additional Storage Conditions: do not store below 4 °C to avoid precipitation of the buffer

Shelf Life: 12 months

Form: liquid

Concentration: 10x

Description:

A 10 x concentration of TBE (Tris-Borate-EDTA) Buffer contains:

108 g/l Tris

55 g/l Boric Acid

7 g/l EDTA

TBE Buffer is optimized for DNA electrophoresis techniques using both acrylamide and agarose. The TBE Buffer has a greater buffering capacity and high ionic strength.

Quality:

- Molecular biology grade
- sterile
- DNase / RNase free

