

**TAE Buffer - 10x conc.**

| Cat. No. | Amount |
|-----------|--------|
| BU-119-10 | 1 l |

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at 4 °C

Shelf Life: 12 months

Form: liquid

Concentration: 10x

pH: 8.2 - 8.4 (22 °C)

Description:

A 10 x concentration of TAE (Tris-Acetate-EDTA) Buffer contains:

0.4 M Tris

0.01 M EDTA

0.2 M Acetic Acid

TAE is an extensively used buffer for agarose gel electrophoresis applications requiring high resolution and separation of high molecular weight, double-stranded DNA. TAE Buffer is more compatible with in-gel manipulations and band recovery procedures than TBE Buffer.

Quality:

- Molecular biology grade
- sterile filtered, filter pore size 0.22 µm
- DNase / RNase free