

**Tris-EDTA Buffer pH 7.6 - 1x conc.**

Cat. No.	Amount
BU-121S-1	100 ml
BU-121L-1	1 l

**For *in vitro* use only!**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at 4 °C

**Shelf Life:** 12 months

**Form:** liquid

**Concentration:** 1x

**pH:** 7.6 ±0.1

**Description:**

The 1 x concentration of Tris-EDTA Buffer contains:

10 mM Tris

1 mM EDTA

Tris-EDTA Buffer is commonly used as storage buffer for DNA and RNA samples. The Buffer protects DNA and RNA from degradation.

**Quality:**

- Molecular biology grade
- sterile
- DNase / RNase free