

**Tris-EDTA Buffer pH 7.6 - 100x conc.**

Cat. No.	Amount
BU-121S-100	100 ml
BU-121L-100	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: 100x

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