

**Tris-HCl Buffer 1 M pH 7.5**

Cat. No.	Amount
BU-125S	100 ml
BU-125L	1 l

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at 4 °C

Shelf Life: 12 months

Molecular Formula: C₄H₁₁NO₃·HCl

Molecular Weight: 157.60 g/mol

CAS#: 1185-53-1

Form: liquid

Concentration: 1 M

pH: 7.5 ±0.1 (22°C)

Description:

Tris-HCl Buffer is often included in PCR reaction buffer. It is used for cell culture in vitro, enzyme assays and some electrophoretic applications at physiological pH.

Quality:

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