

**Tris-HCl Buffer 1 M pH 7.0**

Cat. No.	Amount
BU-124S-70	100 ml
BU-124L-70	1 l

For *in vitro* use only!

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: 7.0 ±0.1 (22 °C)

Description:

Tris-HCl Buffer is included in PCR reaction buffer, used for *in vitro* cell culture, enzyme assays and electrophoretic applications.

Quality:

- Molecular biology grade
- sterile
- DNase / RNase free