

**Tris-HCl Buffer 1 M pH 8.5**

| Cat. No. | Amount |
|------------|--------|
| BU-124S-85 | 100 ml |
| BU-124L-85 | 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: 8.5 ±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 filtered, filter pore size 0.22 µm
- DNase / RNase free