

**Hepes Buffer 1 M pH 7.5**

Cat. No.	Amount
BU-106-75	100 ml

For *in vitro* use only!

Shipping: shipped at ambient temperature

Storage Conditions: store at 4 °C

Shelf Life: 12 months

Molecular Formula: C₈H₁₈N₂O₄S

Molecular Weight: 238.30 g/mol

CAS#: 7365-45-9

Form: liquid

Concentration: 1 M

pH: 7.5 ±0.1 (22 °C)

Description:

Hepes Buffer is one of the best general purpose buffers available for biological research. It has been used to advantage in tissue culture, oxidative phosphorylation and protein synthesis with cell-free bacterial systems. Hepes is a recommended buffer for the glutamate binding assay, prevents binding to non-receptor materials.

Quality:

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