

**Hepes Buffer 1 M pH 7.0**

Cat. No.	Amount
BU-106-70	100 ml

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at 4 °C

**Shelf Life:** 12 months

**Molecular Formula:**  $C_8H_{18}N_2O_4S$

**Molecular Weight:** 238.30 g/mol

**CAS#:** 7365-45-9

**Form:** liquid

**Concentration:** 1 M

**pH:** 7.0  $\pm$  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  $\mu$ m
- DNase / RNase free