

**Magnesium Sulfate Solution 1 M**

| Cat. No.  | Amount |
|-----------|--------|
| BU-111-1M | 100 ml |

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at 4 °C

**Shelf Life:** 12 months

**Molecular Formula:**  $\text{MgSO}_4 \cdot 7 \text{H}_2\text{O}$

**Molecular Weight:** 246.48 g/mol

**CAS#:** 10034-99-8

**Form:** liquid

**Concentration:** 1 M

**Description:**

Magnesium Sulfate is used as a source of magnesium ions and is an essential Co-factor in many enzymes (DNase).

**Quality:**

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