

**Magnesium Acetate Tetrahydrate - 1 M**

Cat. No.	Amount
CSS-210	100 ml

Description:

Crystallization Stock Solution, sterile filtered

For *in vitro* use only!**Shipping:** shipped at ambient temperature**Storage Conditions:** store at 4 °C**Shelf Life:** 12 months**Molecular Formula:** $(\text{CH}_3\text{COO})_2\text{Mg} \cdot 4\text{H}_2\text{O}$ **Molecular Weight:** 214.45 g/mol**CAS#:** 16674-78-5**EC number:** 205-554-9**Form:** liquid**Concentration:** 1 M