

**Ammonium Sulfate - 4 M**

Cat. No.	Amount
CSS-143	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:** $(\text{NH}_4)_2\text{SO}_4$ **Molecular Weight:** 132.14 g/mol**CAS#:** 7783-20-2**EC number:** 231-984-1**Form:** liquid**Concentration:** 4 M