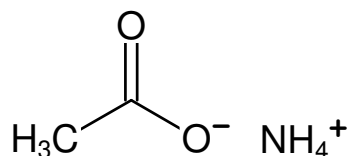


**Ammonium Acetate - 5 M**

sterile filtered

Cat. No.	Amount
CSS-129	100 ml

Description:
Crystallization Stock Solution**For *in vitro* use only!****Shipping:** shipped at ambient temperature**Storage Conditions:** store at 4 °C**Shelf Life:** 12 months**Molecular Formula:** CH₃CO₂NH₄**Molecular Weight:** 77.08 g/mol**CAS#:** 631-61-8**EC number:** 211-162-9**Form:** liquid**Concentration:** 5 M