

**Sodium Acetate - 1 M**

any pH value

useful buffer range: pH 3.6 - 5.6

Cat. No.	Amount
CSS-519	100 ml

Description:

Crystallization Stock Solution, 1M Sodium Acetate adjusted with Acetic Acid, sterile filtered

For in vitro use only!**Shipping:** shipped at ambient temperature**Storage Conditions:** store at 4 °C**Shelf Life:** 12 months**Molecular Formula:** CH₃COONa**Molecular Weight:** 82.03 g/mol**CAS#:** 127-09-3**EC number:** 204-823-8**Form:** liquid**Concentration:** 1 M**pH:** arbitrary in 0.1 increments