

**Sodium Potassium Phosphate pH 7.5 - 1 M**

Cat. No.	Amount
CSS-351	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:** Na₂HPO₄ (di-Sodium hydrogen Phosphate);
KH₂PO₄ (Potassium dihydrogen Phosphate)**Molecular Weight:** 141.96 g/mol (di-Sodium hydrogen Phosphate);
136.09 g/mol (Potassium dihydrogen Phosphate)**CAS#:** 7558-79-4 (di-Sodium hydrogen Phosphate), 7778-77-0
(Potassium dihydrogen Phosphate)**EC number:** 231-448-7 (di-Sodium hydrogen Phosphate), 231-913-4
(Potassium dihydrogen Phosphate)**Form:** liquid**Concentration:** 1 M**pH:** 7.5, 1M Potassium dihydrogen Phosphate adjusted with 1M
di-Sodium hydrogen Phosphate