

**Calcium Chloride Dihydrate - 5 M**

sterile filtered

Cat. No.	Amount
CSS-155	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:** CaCl₂ * 2H₂O**Molecular Weight:** 147.01 g/mol**CAS#:** 10035-04-8**EC number:** 233-140-8**Form:** liquid**Concentration:** 5 M