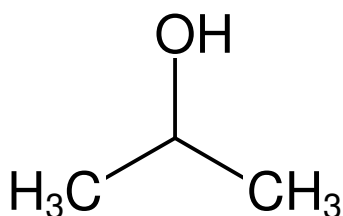


**2-Propanol - 100 % v/v**Iso-Propanol
Isopropyl alcohol

Cat. No.	Amount
CSS-119	4 x 25 ml

**For *in vitro* use only!****Description:**
Crystallization Stock Solution**Shipping:** shipped at ambient temperature**Storage Conditions:** store at ambient temperature**Shelf Life:** 12 months**Molecular Formula:** (CH₃)₂CHOH**Molecular Weight:** 60.10 g/mol**CAS#:** 67-63-0**EC number:** 200-661-7**Form:** liquid**Concentration:** 100 % v/v