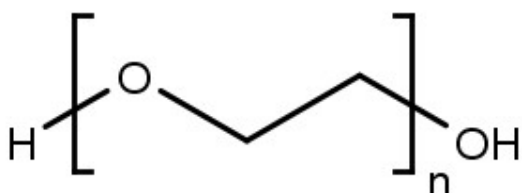


**Polyethyleneglycol 600 - 50 % w/v**

PEG 600

Cat. No.	Amount
CSS-240	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:** H(OCH₂CH₂)_nOH**CAS#:** 25322-68-3**EC number:** 203-473-3**Form:** liquid**Concentration:** 50 % w/v