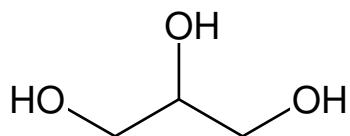


**Glycerol - 100 % v/v**Glycerin
1,2,3-Propanetriol

Cat. No.	Amount
CSS-188	100 ml

Description:
Crystallization Stock Solution**For *in vitro* use only!****Shipping:** shipped at ambient temperature**Storage Conditions:** store at ambient temperature**Shelf Life:** 12 months**Molecular Formula:** HOCH₂CH(OH)CH₂OH**Molecular Weight:** 92.09 g/mol**CAS#:** 56-81-5**EC number:** 200-289-5**Form:** liquid**Concentration:** 100 % v/v