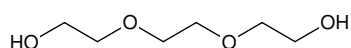


**Triethyleneglycol - 50 % w/v**

Triglycol

Cat. No.	Amount
CSS-314	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:** HO(CH₂CH₂O)₂CH₂CH₂OH**Molecular Weight:** 150.17 g/mol**CAS#:** 112-27-6**EC number:** 203-953-2**Form:** liquid**Concentration:** 50 % w/v