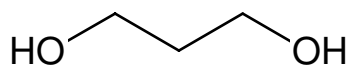


**1,3-Propanediol - 50 % v/v**

Trimethylene glycol

Cat. No.	Amount
CSS-104	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(CH₂)₃OH**Molecular Weight:** 76.06 g/mol**CAS#:** 504-63-2**EC number:** 207-997-3**Form:** liquid**Concentration:** 50 % v/v