

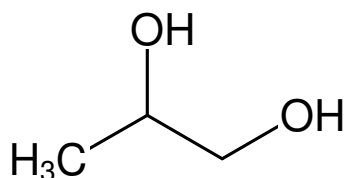
**Propylene Glycol - 50 % v/v**

1,2-Propanediol

Cat. No.	Amount
CSS-280	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:** CH<sub>3</sub>CH(OH)CH<sub>2</sub>OH**Molecular Weight:** 76.09 g/mol**CAS#:** 57-55-6**EC number:** 200-338-0**Form:** liquid**Concentration:** 50 % v/v