

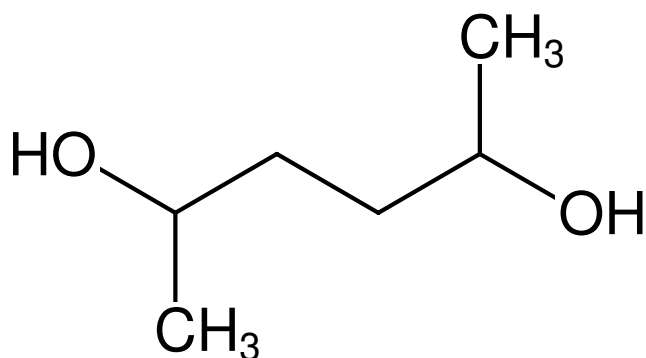
**2,5-Hexanediol - 8 % v/v**

2,5-Hexylene glycol

Cat. No.	Amount
CSS-419	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:** C₆H₁₄O₂**Molecular Weight:** 118.17 g/mol**CAS#:** 2935-44-6**EC number:** 220-910-3**Form:** liquid**Concentration:** 8 % v/v