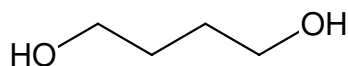


**1,4-Butanediol - 50 % v/v**

1,4-Butylene glycol

Cat. No.	Amount
CSS-386	100 ml

**For *in vitro* use only!****Description:**  
Crystallization Stock Solution**Shipping:** shipped at ambient temperature**Storage Conditions:** store at 4 °C**Shelf Life:** 12 months**Molecular Formula:** C<sub>4</sub>H<sub>10</sub>O<sub>2</sub>**Molecular Weight:** 90.12 g/mol**CAS#:** 110-63-4**EC number:** 203-786-5**Form:** liquid**Concentration:** 50 % v/v