

**Polyethyleneglycol 2000 - 50 % w/v**

PEG 2000

Cat. No.	Amount
CSS-245	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:**  $\text{H}(\text{OCH}_2\text{CH}_2)_n\text{OH}$ **CAS#:** 25322-68-3**EC number:** 203-473-3**Form:** liquid**Concentration:** 50 % w/v