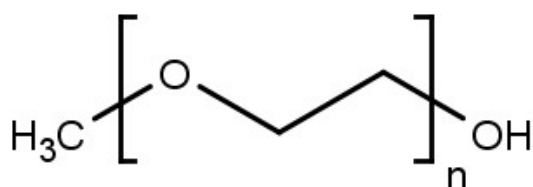


**Polyethyleneglycol 5000 Monomethylether - 50 % w/v**

PEG 5000 MME

Cat. No.	Amount
CSS-235	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>(OCH<sub>2</sub>CH<sub>2</sub>)<sub>n</sub>OH**CAS#:** 9004-74-4**Form:** liquid**Concentration:** 50 % w/v