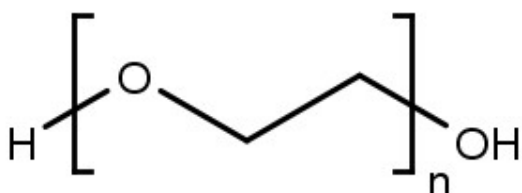


**Polyethyleneglycol 400 - 100 % v/v**

PEG 400

| Cat. No. | Amount |
|----------|--------|
| CSS-252 | 100 ml |

Description:
Crystallization Stock Solution**For general laboratory use.****Please centrifuge briefly before opening (volume ≤ 2 ml).****Shipping:** shipped at ambient temperature**Storage Conditions:** store at ambient temperature**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:** 100 % v/v