

## Polyethyleneglycol 3350

PEG 3350, 50%w/v

Cat. No.	Amount
CSS-249	100 ml

**Application:** Reagent for optimization of crystallization conditions for proteins.

**JBScreen Single Stocks - PEG 3350** is a ready-made solution.

The single stock is provided as a 50% w/v sterile filtered stock solution and supplied in 100 ml aliquots.

**Molecular Formula:**  $\text{CH}_3(\text{OCH}_2\text{CH}_2)_n\text{OH}$

**CAS Number:** 25322-68-3

**EC Number:** 2034733

**Storage:** 4°C

**Expiry:** 12 months