

## Polyethyleneglycol 550 Monomethylether

PEG 550 MME, 50 % v/v

Cat. No.	Amount
CSS-236	100 ml

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

**JBScreen Single Stocks - PEG 550 MME** is a ready-made solution.

The single stock is provided as a 50% v/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:** 9004-74-4

**Storage:** 4°C

**Expiry:** 12 months