

Polyethyleneglycol 550 Monomethylether

PEG 550 MME, 50 % w/v

Cat. No.	Amount
CSS-237	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% 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: 9004-74-4

Storage: 4°C

Expiry: 12 months