

Polyethyleneglycol 300

Polyethyleneglycol 300, 50 % v/v

Cat. No.	Amount
CSS-247	100 ml

Application: Reagent for optimization of crystallization conditions for proteins.

JBScreen Single Stocks - PEG 300 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: 25322-68-3

EC Number: 2034733

Storage: 4°C

Expiry: 12 months