

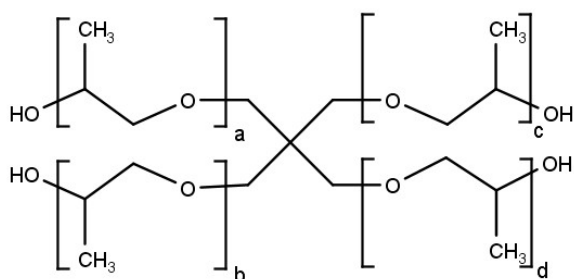


Pentaerythritol propoxylate (17/8 PO/OH) - 60 % w/v

PEP 629

Cat. No.	Amount
CSS-376	100 ml

Description:
Crystallization Stock Solution



For *in vitro* use only!

Shipping: shipped at ambient temperature

Storage Conditions: store at 4 °C

Shelf Life: 12 months

Molecular Formula: $C[CH_2[OCH_2CH(CH_3)]_nOH]_4$

CAS#: 9051-49-4

EC number: 500-030-9

Form: liquid

Concentration: 60 % w/v