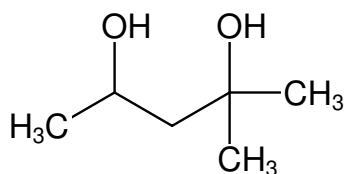


**MPD - 100 % v/v**

(+/-)-2-Methyl-2,4-Pentandiol
Hexylene glycol

Cat. No.	Amount
CSS-117	100 ml

Description:
Crystallization Stock Solution



For *in vitro* use only!

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: 12 months

Molecular Formula: CH₃CH(OH)CH₂C(CH₃)₂OH

Molecular Weight: 118.17 g/mol

CAS#: 107-41-5

EC number: 203-489-0

Form: clear solution

Concentration: 100 % v/v