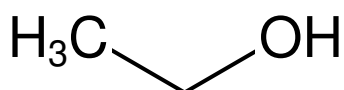


**Ethanol - 50 % v/v**

Ethylalkohol

Cat. No.	Amount
CSS-330	4 x 25 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:** CH₃CH₂OH**Molecular Weight:** 46.07 g/mol**CAS#:** 64-17-5**EC number:** 200-578-6**Form:** liquid**Concentration:** 50 % v/v