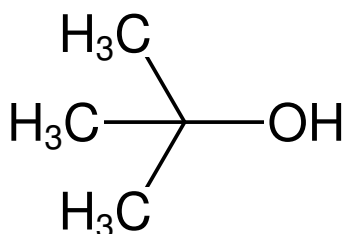


**tert-Butanol - 50 % w/v**2-Methyl-2-propanol
Trimethyl carbinol

Cat. No.	Amount
CSS-310	4 x 25 ml

**For *in vitro* use only!****Shipping:** shipped at ambient temperature**Storage Conditions:** store at 4 °C**Shelf Life:** 12 months**Molecular Formula:** (CH₃)₃COH**Molecular Weight:** 74.12 g/mol**CAS#:** 75-65-0**EC number:** 200-889-7**Form:** liquid**Concentration:** 50 % w/v**Description:**
Crystallization Stock Solution