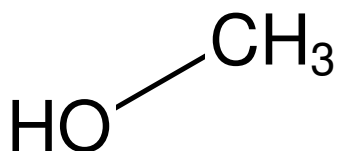


**Methanol - 50 % w/v**

Methyl alcohol

| Cat. No. | Amount    |
|----------|-----------|
| CSS-224  | 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:** CH<sub>3</sub>OH**Molecular Weight:** 32.04 g/mol**CAS#:** 67-56-1**EC number:** 200-659-6**Form:** liquid**Concentration:** 50 % w/v**Description:**  
Crystallization Stock Solution