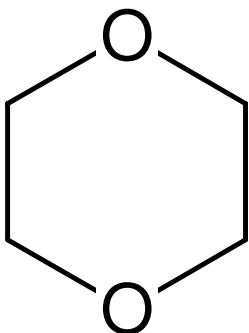


**1,4-Dioxane - 50 % w/v**

Diethylene oxide

Cat. No.	Amount
CSS-106	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:** C<sub>4</sub>H<sub>8</sub>O<sub>2</sub>**Molecular Weight:** 88.11 g/mol**CAS#:** 123-91-1**EC number:** 204-661-8**Form:** liquid**Concentration:** 50 % w/v