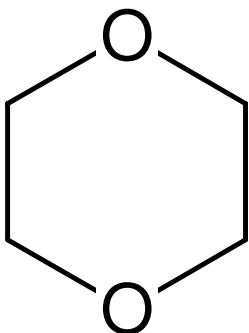


**1,4-Dioxane - 50 % v/v**

Diethylene oxide

Cat. No.	Amount
CSS-107	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: C₄H₈O₂

Molecular Weight: 88.11 g/mol

CAS#: 123-91-1

EC number: 204-661-8

Form: liquid

Concentration: 50 % v/v