

DMSO

Dimethylsulfoxid $\geq 99,5\%$

Cat.-No.	Amount
CPP-A01S	25 ml
CPP-A01L	125 ml

For *in vitro* use only
Quality guaranteed for 12 months
Store at 4°C

Product Information

CAS Number: 67-68-5

Molecular Formula

C₂H₆OS

Molecular Weight

78.13 Da

Boiling point

189 °C

Melting point

18.5 °C

Flash point

87 °C

Density

1.1 g/ml

Solubility

soluble in water

Description

The addition of DMSO may enhance penetration dependent on the membrane composition of cells. We recommend the use of 10% DMSO in serum free transduction medium.

For more information see **general manual (4.2.1)**.