

o-Phenantroline **1,10-Phenanthroline Monohydrate**

Cat.-No.	Amount
CPP-A06S	50 mg
CPP-A06L	250 mg

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

Protect from light.

Toxic if swallowed. Very toxic to aquatic organisms; may cause long-term adverse effects in the aquatic environment. Wear suitable gloves. In case of accident or if you feel unwell, seek medical advice immediately. This material and its container must be disposed of as hazardous waste. Avoid release to the environment.

Product Information

CAS Number: 5144-89-8

Molecular Formula

$C_{12}H_8N_2 \cdot H_2O$

Molecular Weight

198.2 Da

Solubility

DMSO (100 mg/ml) or water (3 mg/ml)

Description

Membrane bound and into the medium secreted proteases may inactivate CPPs indicating a need for using protease inhibitors.

o-Phenantroline acts as a general metalloprotease inhibitor.

Add the inhibitor to the serum free incubation medium to achieve a concentration of 1 mM.

For more information see **general manual (4.2.2)**.