

Heparin

Heparin sodium salt from porcine intestinal

Cat.-No.	Amount
CPP-A04S	100 mg
CPP-C04L	500 mg

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

Substance not yet fully tested. (EU)
Avoid contact and inhalation.
Target organ(s): Blood.
For R&D use only. Not for drug, household or other use.

Product Information

CAS Number: 9041-08-1

Form

White to off-white powder

Solubility

soluble in water (50 mg/ml)

Description

If microscopically indicated remove cargo from outside of cell membrane (proteoglycane scaffold) with heparin (0.1%), which is dissolved in PBS or glycine buffer.

Because of its polarity and large molecular size Heparin poorly crosses cell membranes. Heparin has been traditionally used as an anticoagulant.

More detailed protocols are given in the **general manual (4.2.2)**.