

Aprotinin

Pancreatic Trypsin Inhibitor, Trypsin-Kallikrein Inhibitor Competitive Inhibitor for Esterases and certain Serin Proteases

Cat.-No.	Amount
CPP-A05S	5 mg
CPP-C05L	25 mg

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

Product Information

CAS Number: 9087-70-1

Form

White crystalline powder

Purity

≥ 95%

Molecular Weight

6.5 kDa

Solubility

soluble in water (10 mg/ml)

Activity

≥ 5,500 KIU/mg

Description

Aprotinin is used for inhibiting proteolytic degradation of CPPs and peptide cargos. It can be combined with BSA and o-phenanthroline. We recommend a concentration of 50 µg per 1 ml serum free incubation medium.

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