

---

## Protease Inhibitor Cocktail

(Catalog No. R-1101)

---

### Description

Protease Inhibitor Cocktail (PIC) is a mixture of several protease inhibitors with broad inhibitory specificity. This product is in an easy-to-use solution form and can be directly employed for inhibition of a broad range of proteases.

### Concentration/Volume

1000 X, 1 ml

### Storage Temp & Conditions

Store aliquots at 4°C. Avoid multiple freeze/thaw cycles.

### Application Notes

Used for the inhibition of proteolytic activity in cell culture and protein isolation.

### Ordering Information

Products	Size	Cat. No.
Protease Inhibitor Cocktail	1 ml	R-1101-1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*