## Protease Inhibitor Cocktail

Citations: 7

Ordering Information

BML-KI103-0001

1Set

View Online »

View Online »

## **Handling & Storage**

**Use/Stability** As indicated on product label or CoA when stored as recommended. The tablets are

stable at +4°C, stored dry. The stock solution is stable for one to two weeks at +4°C, or

at least 12 weeks at -20°C.

Long Term Storage +4°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Appearance Two tablets

**Application Notes**Mixture of several protease inhibitors. For the inhibition of

serine, cysteine, but not metalloproteases, and protects proteins isolated from animal tissues, plants, yeast and bacteria. 1 tablet contains protease inhibitors sufficient for

a 10 ml cell extract.

**Solubility** Soluble in aqueous buffers, or add directly to extraction

media. Alternatively, prepare 7X stock solutions in 1.5ml

water or 100mM phosphate buffer, pH 7.0.

Last modified: May 29, 2024