

Protease Inhibitor Cocktail

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-KI103-0001	1Set
----------------	------

Manuals, SDS & CofA

[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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)