

Cystatin C (human), (purified)

Protease inhibitor

Cystatin C is a major member of a large family of endogenous cysteine protease inhibitors found throughout the body. It inhibits proteases such as cathepsin B, H, DPPI, and papain, and is inactivated by neutrophil elastase. It is involved in cancer, rheumatoid arthritis, and other inflammatory diseases, and serves as a marker for renal function. Recently it was shown to bind and inhibit amyloid fibril formation, indicating therapeutic potential for Alzheimer's disease.

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE479-0100	100µg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Although this is in highly purified form, and the source individual tested negative for all diseases tested, handle this product with all due precautions.
Handling	Aliquot product into individual tubes and freeze at -20 or -70°C. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Cystatin-3, Neuroendocrine basic polypeptide, CysC
Application Notes	Can be used for the inhibition of cysteine protease activity or of amyloid fibril formation in vitro, in vivo, or in tissue culture.
Formulation	Lyophilized from 0.02M ammonium bicarbonate buffer.
MW	~13kDa
Purity	≥95% (SDS-PAGE)
Reconstitution	Resuspend in water and freeze at -20°C in aliquots.
Source	Isolated from human urine.
UniProt ID	P01034



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-
fr@enzolifesciences.com](mailto:info-fr@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)