

Catestatin (human)

Human Chromogranin A [352-372]). Wild type catestatin, initially described as a potent catecholamine release inhibitory peptide. Cst located in the skin is active against Gram-negative and Gram-positive bacteria as well as against fungi and yeast. Also proposed as a modulator or regulator of cardiac function and as an anti-hypertensive agent.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-CP6420-0001	1mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Chromogranin A (352-372)

Appearance Solid.

Formula $C_{104}H_{164}N_{32}O_{27}S_1$

Formulation Lyophilized.

MW 2326.7

Purity ≥98% (HPLC)

Sequence H-Ser-Ser-Met-Lys-Leu-Ser-Phe-Arg-Ala-Arg-Ala-Tyr-Gly-Phe-Arg-Gly-Pro-Gly-Pro-Gln-Leu-OH (human Chromogranin A[352-372])

Solubility Soluble in Water.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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