

PRODUCT INFORMATION



Catestatin (bovine)

Item No. 42518

CAS Registry No.: 142211-96-9

Formal Name: L-arginyl-L-seryl-L-methionyl-L-arginyl-L-leucyl-L-seryl-L-phenylalanyl-L-arginyl-L-alanyl-L-arginylglycyl-L-tyrosylglycyl-L-phenylalanyl-L-arginylglycyl-L-prolylglycyl-L-leucyl-L-glutaminy-L-leucine

Synonyms: bCgA₃₄₄₋₃₆₄, Bovine Chromogranin A₃₄₄₋₃₆₄

Peptide Sequence: RSMRLSFRARGYGFRGPGQL-OH

MF: C₁₀₇H₁₇₃N₃₇O₂₆S

FW: 2,425.8

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

H—Arg—Ser—Met—Arg—Leu—Ser—Phe—Arg—Ala—Arg—Gly—Tyr—Gly—Phe—Arg—Gly—Pro—Gly—Leu—Gln—Leu—OH

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Catestatin (bovine) is supplied as a solid. A stock solution may be made by dissolving the catestatin (bovine) in the solvent of choice, which should be purged with an inert gas. Catestatin (bovine) is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Catestatin is a peptide derivative of chromogranin A and an inhibitor of norepinephrine release.¹ It inhibits nicotine-induced secretion of norepinephrine and nicotine-induced sodium intake in PC12 rat adrenal medulla cells (IC₅₀s = ~0.2 and ~0.25 μM, respectively). Catestatin decreases mean arterial blood pressure and current-induced pressor response and increases plasma levels of epinephrine and histamine in pithed rats.²

References

1. Mahata, S.K., O'Connor, D.T., Mahata, M., *et al.* Novel autocrine feedback control of catecholamine release. A discrete chromogranin a fragment is a noncompetitive nicotinic cholinergic antagonist. *J. Clin. Invest.* **100(6)**, 1623-1633 (1997).
2. Kennedy, B.P., Mahata, S.K., O'Connor, D.T., *et al.* Mechanism of cardiovascular actions of the chromogranin A fragment catestatin in vivo. *Peptides* **19(7)**, 1241-1248 (1998).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/10/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM