

PRODUCT INFORMATION



Oxyntomodulin (human, mouse, rat) (trifluoroacetate salt)

Item No. 24758

Formal Name: L-histidyl-L-seryl-L-glutaminyglycyl-L-threonyl-L-phenylalanyl-L-threonyl-L-seryl-L- α -aspartyl-L-tyrosyl-L-seryl-L-lysyl-L-tyrosyl-L-leucyl-L- α -aspartyl-L-seryl-L-arginyl-L-arginyl-L-alanyl-L-glutaminy-L- α -aspartyl-L-phenylalanyl-L-valyl-L-glutaminy-L-tryptophyl-L-leucyl-L-methionyl-L-asparaginy-L-threonyl-L-lysyl-L-arginyl-L-asparaginy-L-arginyl-L-asparaginy-L-asparaginy-L-isoleucyl-L-alanine, trifluoroacetate salt

H—His—Ser—Gln—Gly—Thr—Phe—Thr—Ser—Asp—Tyr—
Ser—Lys—Tyr—Leu—Asp—Ser—Arg—Arg—Ala—Gln—
Asp—Phe—Val—Gln—Trp—Leu—Met—Asn—Thr—Lys—
Arg—Asn—Arg—Asn—Asn—Ile—Ala—OH

Synonyms: Glicentin (33-69), Proglucagon (33-69)

MF: $C_{192}H_{295}N_{61}O_{60}S \cdot XCF_3COOH$

FW: 4,449.8

Purity: $\geq 95\%$

Supplied as: A lyophilized powder

Storage: $-20^{\circ}C$

Stability: ≥ 4 years

$\cdot XCF_3COOH$

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

Laboratory Procedures

Oxyntomodulin (human, mouse, rat) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the oxyntomodulin (human, mouse, rat) (trifluoroacetate salt) in water. The solubility of oxyntomodulin (human, mouse, rat) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Oxyntomodulin is a peptide hormone involved in regulation of food intake, energy expenditure, and glucose metabolism.¹ It is produced in gut endocrine L-cells *via* post-translational processing of proglucagon by prohormone convertase 1/3 following nutrient uptake. Oxyntomodulin is an agonist of the glucagon-like peptide-1 receptor (GLP-1R) and the glucagon receptor (GCGR; $IC_{50}s = 4.3$ and 355 nM for rat GCGR and GLP-1R, respectively).² It is selective for GCGR and GLP-1R over GLP-2R and glucose-dependent insulinotropic peptide receptor (GIPR), as it stimulates cAMP formation in BHK cells expressing rat GCGR and GLP-1R when used at concentrations of 10 or 100 nM, but has no effect on cells expressing rat GLP-2R or GIPR at either concentration.³ Oxyntomodulin also stimulates cAMP formation in CHO cells expressing mouse GLP-1R and GCGR ($EC_{50}s = 2.5$ and 6.2 nM, respectively).⁴ *In vivo*, oxyntomodulin reduces refeeding in 24-hour fasted rats when injected intracerebroventricularly or into the paraventricular nucleus (PVN) of the hypothalamus at doses of 3 and 1 nmol, respectively.⁵ Oxyntomodulin (50 nmol/kg twice per day for 7 days) reduces cumulative daily food intake and rate of body weight gain in rats.⁶

References

1. Pocai, A. *Mol. Metab.* **3**(3), 241-251 (2013).
2. Price, S.L., Minnion, J.S., and Bloom, S.R. *Curr. Ther. Res. Clin. Exp.* **77**, 111-115 (2015).
3. Baggio, L.L., Huang, Q., Brown, T.J., et al. *Gastroenterology* **127**(2), 546-558 (2004).
4. Pocai, A., Carrington, P.E., Adams, J.R., et al. *Diabetes* **58**(10), 2258-2266 (2009).
5. Dakin, C.L., Gunn, I., Small, C.J., et al. *Endocrinology* **142**(10), 4244-4250 (2001).
6. Dakin, C.L., Small, C.J., Batterham, R.L., et al. *Endocrinology* **145**(6), 2687-2695 (2004).

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, 12/06/2022

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