

PRODUCT INFORMATION



[Leu¹⁵]-Gastrin I amide (human) (trifluoroacetate salt)

Item No. 31535

Formal Name:	5-oxo-L-prolylglycyl-L-prolyl-L-tryptophyl-L-leucyl-L- α -glutamyl-L- α -glutamyl-L- α -glutamyl-L- α -glutamyl-L- α -glutamyl-L-alanyl-L-tyrosylglycyl-L-tryptophyl-L-leucyl-L- α -aspartyl-L-phenylalaninamide, trifluoroacetate salt	H—Glp—Gly—Pro—Trp—Leu—Glu—Glu—Glu—Glu—Glu—
Synonyms:	15-Leu-Gastrin-17, GPWLEEEEEAYGWLDF-NH ₂	Ala—Tyr—Gly—Trp—Leu—Asp—Phe—NH ₂
MF:	C ₉₈ H ₁₂₆ N ₂₀ O ₃₁ • XCF ₃ COOH	• XCF ₃ COOH
FW:	2,080.2	
Purity:	≥95%	
Supplied as:	A solid	
Storage:	-20°C	
Stability:	≥4 years	

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

Laboratory Procedures

[Leu¹⁵]-Gastrin I amide (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the [Leu¹⁵]-gastrin I amide (human) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. [Leu¹⁵]-Gastrin I amide (human) (trifluoroacetate salt) is soluble in 0.1 M NaOH.

Description

[Leu¹⁵]-Gastrin I amide is a synthetic peptide fragment of the peptide hormone gastrin I (1-17) (Item No. 24457) that contains a substitution of leucine for methionine at position 15.^{1,2}

References

1. Wunsch, E. Synthesis of pure polypeptide hormones with full biological activity. *The chemistry of polypeptides*. Katsoyannis, P.G., editor, 1st edition, Plenum Press (1973).
2. Morley, J.S. and Smith, J.M. Polypeptides. Part IX. Variations of the methionyl position in the C-terminal tetrapeptide amide sequence of the gastrins. *J. Chem. Soc. Perkin 1* 6, 726-733 (1968).

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, 10/07/2022

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