

PRODUCT INFORMATION



[D-Lys³]-GHRP-6 (trifluoroacetate salt)

Item No. 36862

Formal Name: (R)-6-amino-2-((R)-2-((S)-2-amino-3-(1H-imidazol-5-yl)propanamido)-3-(1H-indol-3-yl)propanamido)-N-((S)-1-(((R)-1-(((S)-1,6-diamino-1-oxohexan-2-yl)amino)-1-oxo-3-phenylpropan-2-yl)amino)-3-(1H-indol-3-yl)-1-oxopropan-2-yl)hexanamide, trifluoroacetate salt

Synonyms: [D-Lys³]-Growth Hormone Releasing Peptide 6, His-D-Trp-D-Lys-Trp-D-Phe-Lys-NH₂

MF: C₄₉H₆₃N₁₃O₆ • CF₃COOH

FW: 930.1

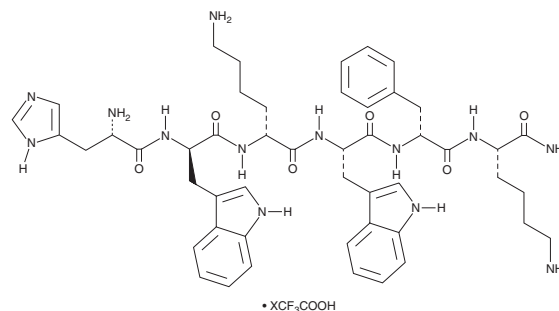
Purity: ≥95%

UV/Vis.: λ_{max}: 218 nm

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

[D-Lys³]-GHRP-6 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the [D-Lys³]-GHRP-6 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. [D-Lys³]-GHRP-6 (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of [D-Lys³]-GHRP-6 (trifluoroacetate salt) in ethanol is approximately 1 mg/ml and approximately 2 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of [D-Lys³]-GHRP-6 (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of [D-Lys³]-GHRP-6 (trifluoroacetate salt) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

[D-Lys³]-GHRP-6 is an antagonist of the growth hormone secretagogue receptor 1a (GHS-R1a; K_i = 4.25 nM).¹ It inhibits ghrelin-induced cAMP accumulation in human aortic smooth muscle cells when used at a concentration of 20 nM. [D-Lys³]-GHRP-6 (12 mg/kg) decreases food intake in Zucker FA/FA lean and fa/fa obese rats.²

References

1. Rossi, F., Castelli, A., Bianco, M.J., *et al.* Ghrelin inhibits contraction and proliferation of human aortic smooth muscle cells by cAMP/PKA pathway activation. *Atherosclerosis* **203**(1), 97-104 (2008).
2. Beck, B., Richy, S., and Sticker-Krongrad, A. Feeding response to ghrelin agonist and antagonist in lean and obese Zucker rats. *Life Sci.* **76**(4), 473-478 (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, 12/21/2022

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