

PRODUCT INFORMATION



Hexarelin

Item No. 30716

CAS Registry No.: 140703-51-1

Formal Name: L-histidyl-2-methyl-D-tryptophyl-L-alanyl-L-tryptophyl-D-phenylalanyl-L-lysineamide

Synonyms: Examorelin, H-2-Me-WAWFK, His-2-Me-Trp-Ala-Trp-Phe-Lys-NH₂, His-(D)-Mrp-Ala-Trp-(D)-Phe-Lys-NH₂

MF: C₄₇H₅₈N₁₂O₆

FW: 887.0

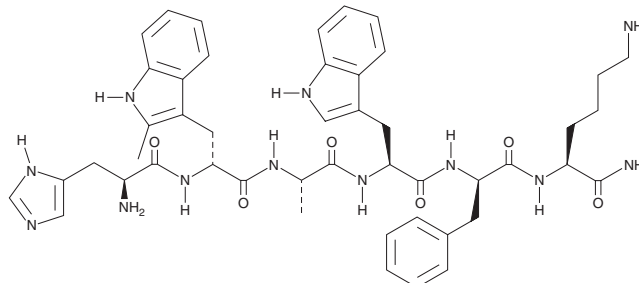
Purity: ≥98%

UV/Vis.: λ_{max}: 221 nm

Supplied as: A crystalline 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

Hexarelin is supplied as a crystalline solid. A stock solution may be made by dissolving the hexarelin in the solvent of choice, which should be purged with an inert gas. Hexarelin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of hexarelin in ethanol is approximately 5 mg/ml and approximately 30 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 hexarelin can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of hexarelin in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Hexarelin is a synthetic hexapeptide agonist of the growth hormone secretagogue receptor (GHS-R).^{1,2} It selectively binds to GHS-R over CD36 (IC₅₀s = 15.9 and 2,080 nM, respectively).¹ Hexarelin (300 µg/kg, s.c.) increases plasma growth hormone levels in neonatal rats.² It also improves left ventricular function and reduces fibrosis in a mouse model of myocardial ischemia induced by transient ligation of the left descending coronary artery (LAD) when administered at a dose of 0.3 mg/kg per day.³

References

1. Proulx, C., Picard, É., Boeglin, D., *et al.* Azapeptide analogues of the growth hormone releasing peptide 6 as cluster of differentiation 36 receptor ligands with reduced affinity for the growth hormone secretagogue receptor 1a. *J. Med. Chem.* **55**(14), 502-511 (2012).
2. Guerlavais, V., Boeglin, D., Mousseaux, D., *et al.* New active series of growth hormone secretagogues. *J. Med. Chem.* **46**(7), 1191-1203 (2003).
3. McDonald, H., Peart, J., Kurniawan, N.D., *et al.* Hexarelin targets neuroinflammatory pathways to preserve cardiac morphology and function in a mouse model of myocardial ischemia-reperfusion. *Biomed. Pharmacother.* **127**, 110165 (2020).

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, 11/15/2022

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