

PRODUCT INFORMATION



Adipokinetic Hormone

Item No. 43823

Formal Name: N-((5S,8S,11S,14S,17S,20S)-5-((1H-indol-3-yl)methyl)-1-amino-14-benzyl-11,17-bis((R)-1-hydroxyethyl)-8-(hydroxymethyl)-22-methyl-1,4,7,10,13,16,19-heptaaxo-3,6,9,12,15,18-hexaazatricosan-20-yl)-5-oxopyrrolidine-2-carboxamide

Synonyms: AKH, Pyr-Leu-Thr-Phe-Thr-Ser-Trp-Gly-NH₂ Pyr—Leu—Thr—Phe—Thr—Ser—Trp—Gly—NH₂

Peptide Sequence: XLTFTSWG-NH₂ (X = pyroglutamyl)

MF: C₄₄H₆₀N₁₀O₁₂

FW: 921.0

Purity: ≥98%

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

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

Adipokinetic hormone is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Adipokinetic hormone is a peptide derivative of adipokinetic hormones, which are insect neuropeptides that regulate metabolism.¹

Reference

1. Van der Horst, D.J. Insect adipokinetic hormones: Release and integration of flight energy metabolism. *Comp. Biochem. Physiol. B Biochem. Mol. Biol.* **136(2)**, 217-226 (2003).

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, 09/08/2025

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