

PRODUCT INFORMATION



Bamadutide

Item No. 41269

CAS Registry No.: 1596343-09-7

Formal Name: L-histidyl-D-seryl-L-glutaminyglycyl-L-threonyl-L-phenylalanyl-L-threonyl-L-seryl-L- α -aspartyl-L-leucyl-L-seryl-L-lysyl-L-glutaminy-L- γ -glutamyl-L-lysyl-L- α -glutamyl-L-seryl-L-lysyl-L-alanyl-L-alanyl-L-glutaminy-L- α -aspartyl-L-phenylalanyl-L-iso-leucyl-L- α -glutamyl-L-tryptophyl-L-leucyl-L-lysyl-L-alanylglycylglycyl-L-prolyl-L-seryl-L-serylglycyl-L-alanyl-L-prolyl-L-prolyl-L-prolyl-L-serinamide

Synonym: SAR425899

Peptide Sequence: HsQGTFSTDLKQXESKAAQDFI
EWLKAGGPSSGAPPPS-NH₂ (X = (S)-4-carboxy-4-hexadecanoylamino-butyl-lysine)

MF: C₂₀₀H₃₁₃N₅₁O₆₃

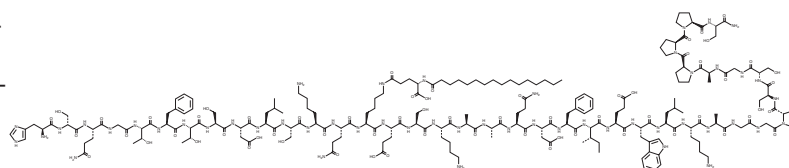
FW: 4,440.0

Purity: $\geq 98\%$

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

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

Bamadutide 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

Bamadutide is a peptide agonist of glucagon receptor (GCGR) and glucagon-like peptide 1 receptor (GLP-1R).¹ It induces cAMP accumulation in HEK293 cells expressing human GCGR or GLP-1R (EC₅₀s = 1 and 3.5 pM, respectively). Bamadutide (0.1 mg/kg) reduces food intake in female mice. It decreases blood levels of glucose and hemoglobin A1c (HbA1c) in female *db/db* mice when administered at a dose of 0.1 mg/kg per day. Bamadutide (5 or 15 μ g/kg twice per day) reduces body weight and total fat mass in male mice fed a high-fat diet.

Reference

1. Haack, T., Wagner, M., Henkel, B., *et al.* Extendin-4 derivatives as dual GLP1/glucagon agonists. *Sanofi*. US20140100156A1 (2014).

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, 07/15/2024

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