

PRODUCT INFORMATION



Exendin-3 (9-39) amide

Item No. 46369

CAS Registry No.: 133514-43-9

Formal Name: L- α -aspartyl-L-leucyl-L-seryl-L-lysyl-L-glutaminyl-L-methionyl-L- α -glutamyl-L- α -glutamyl-L- α -glutamyl-L-alanyl-L-valyl-L-arginyl-L-leucyl-L-phenylalanyl-L-isoleucyl-L- α -glutamyl-L-tryptophyl-L-leucyl-L-lysyl-L-asparaginylglycylglycyl-L-prolyl-L-seryl-L-serylglycyl-L-alanyl-L-prolyl-L-prolyl-L-prolyl-L-serinamide

H—Asp—Leu—Ser—Lys—Gln—Met—Glu—Glu—Glu—Ala—

Val—Arg—Leu—Phe—Ile—Glu—Trp—Leu—Lys—Asn—

Gly—Gly—Pro—Ser—Ser—Gly—Ala—Pro—Pro—Pro—

Ser—NH₂

Synonyms: Avexitide, Exendin (9-39) amide, Exendin (9-39)NH₂

Peptide Sequence: DLSKQMEEEAVRLFIEWLKNGGPSSGAPPPS-NH₂

MF: C₁₄₉H₂₃₄N₄₀O₄₇S

FW: 3,369.8

Purity: $\geq$ 98%

Supplied as: A solid

Storage: -20°C

Stability: $\geq$ 4 years

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

Laboratory Procedures

Exendin-3 (9-39) amide is supplied as a solid. A stock solution may be made by dissolving the exendin-3 (9-39) amide in the solvent of choice, which should be purged with an inert gas. Exendin-3 (9-39) amide is sparingly soluble (1-10 mg/ml) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Exendin-3 (9-39) amide is an antagonist of the glucagon-like peptide 1 receptor (GLP-1R) and a truncated form of exendin-4 amide (Item No. 11096).¹ It inhibits cAMP production induced by GLP-1 or glucose-dependent insulinotropic polypeptide (GIP) in Chinese hamster lung fibroblasts expressing GLP-1R or the GIP receptor (GIPR) at 10 nM. Exendin-3 (9-39) amide (25 nmol/kg) reduces plasma insulin levels in fed or fasted *ob/ob* mice.

Reference

1. Gault, V.A., O'Harte, F.P., Harriott, P., *et al.* Effects of the novel (Pro³)GIP antagonist and exendin(9-39) amide on GIP- and GLP-1-induced cyclic AMP generation, insulin secretion and postprandial insulin release in obese diabetic (*ob/ob*) mice: Evidence that GIP is the major physiological incretin. *Diabetologia* 46(2), 222-230 (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, 07/22/2026

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