

# PRODUCT INFORMATION



## Exendin-4 (3-39) amide (trifluoroacetate salt)

Item No. 36912

**Formal Name:** L- $\alpha$ -glutamylglycyl-L-threonyl-L-phenylalanyl-L-threonyl-L-seryl-L- $\alpha$ -aspartyl-L-leucyl-L-seryl-L-lysyl-L-glutamyl-L-methionyl-L- $\alpha$ -glutamyl-L- $\alpha$ -glutamyl-L- $\alpha$ -glutamyl-L-alanyl-L-valyl-L-arginyl-L-leucyl-L-phenylalanyl-L-isoleucyl-L- $\alpha$ -glutamyl-L-tryptophyl-L-leucyl-L-lysyl-L-asparagylglycylglycyl-L-prolyl-L-seryl-L-serylglycyl-L-alanyl-L-prolyl-L-prolyl-L-prolyl-L-serinamide, trifluoroacetate salt

**Synonym:** Exendin (3-39)

**Peptide Sequence:** EGTFTSDLKQMEEEAVRLFIEWLKNGGPSSG  
APPPS-NH<sub>2</sub>

**MF:** C<sub>176</sub>H<sub>272</sub>N<sub>46</sub>O<sub>58</sub>S • XCF<sub>3</sub>COOH

**FW:** 3,992.4

**Purity:**  $\geq$ 90%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:**  $\geq$ 4 years

H—Glu—Gly—Thr—Phe—Thr—Ser—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—NH<sub>2</sub>

• XCF<sub>3</sub>COOH

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

### Laboratory Procedures

Exendin-4 (3-39) amide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the exendin-4 (3-39) amide (trifluoroacetate salt) in the solvent of choice. Exendin-4 (3-39) amide (trifluoroacetate salt) is soluble in organic solvents such as methanol, which should be purged with an inert gas. Exendin-4 (3-39) amide (trifluoroacetate salt) is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Exendin-4 (3-39) amide is an agonist of glucagon-like peptide 1 receptor (GLP-1R) and an active metabolite of exendin-4 (48-86) amide (Item No. 11096).<sup>1,2</sup> It is formed from exendin-4 (48-86) amide in rat liver homogenates.<sup>2</sup> Exendin-4 (3-39) amide induces cAMP accumulation in HEK293 cells expressing human GLP-1R (EC<sub>50</sub> = 9 nM).<sup>1</sup>

### References

1. Patterson, J.T., Ottaway, N., Gelfanov, V.M., *et al.* A novel human-based receptor antagonist of sustained action reveals body weight control by endogenous GLP-1. *ACS Chem. Biol.* **6**(2), 135-145 (2011).
2. Liao, S., Liang, Y., Zhang, Z., *et al.* *In vitro* metabolic stability of exendin-4: Pharmacokinetics and identification of cleavage products. *PLoS One* **10**(2), e0116805 (2015).

#### 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, 12/20/2023

#### 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