

# PRODUCT INFORMATION



## GLP-1 (1-32) (bullfrog) (trifluoroacetate salt)

Item No. 40718

<b>Formal Name:</b>	L-histidyl-L-alanyl-L- $\alpha$ -aspartylglycyl-L-threonyl-L-phenylalanyl-L-threonyl-L-seryl-L- $\alpha$ -aspartyl-L-methionyl-L-seryl-L-seryl-L-tyrosyl-L-leucyl-L- $\alpha$ -glutamyl-L- $\alpha$ -glutamyl-L-lysyl-L-alanyl-L-alanyl-L-lysyl-L- $\alpha$ -glutamyl-L-phenylalanyl-L-valyl-L- $\alpha$ -aspartyl-L-tryptophyl-L-leucyl-L-isoleucyl-L-lysylglycyl-L-arginyl-prolyl-L-lysineamide, trifluoroacetate salt	H — His — Ala — Asp — Gly — Thr — Phe — Thr — Ser — Asp — Met — Ser — Ser — Tyr — Leu — Glu — Glu — Lys — Ala — Ala — Lys — Glu — Phe — Val — Asp — Trp — Leu — Ile — Lys — Gly — Arg —
<b>Synonym:</b>	Glucagon-like Peptide 1 (1-32)	
<b>Peptide Sequence:</b>	HADGTFSDMSSYLEEKAAKEFVDWLKGRPK-NH <sub>2</sub>	Pro — Lys — NH <sub>2</sub>
<b>MF:</b>	C <sub>164</sub> H <sub>250</sub> N <sub>42</sub> O <sub>51</sub> S • XCF <sub>3</sub> COOH	
<b>FW:</b>	3,658.1	• XCF <sub>3</sub> COOH
<b>Purity:</b>	≥98%	
<b>Supplied as:</b>	A solid	
<b>Storage:</b>	-20°C	
<b>Stability:</b>	≥4 years	

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

### Laboratory Procedures

Glucagon-like peptide 1 (GLP-1) (1-32) (bullfrog) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the GLP-1 (1-32) (bullfrog) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. GLP-1 (1-32) (bullfrog) (trifluoroacetate salt) is soluble (≥10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

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 GLP-1 (1-32) (bullfrog) (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. GLP-1 (1-32) (bullfrog) (trifluoroacetate salt) is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

GLP-1 (1-32) is an endogenous incretin hormone.<sup>1</sup> It increases intracellular cAMP levels using HEK293 cells expressing human GLP-1 receptor (GLP-1R; EC<sub>50</sub> = 0.051 nM). *In vivo*, GLP-1 (1-32) (30 mg/kg) reduces blood glucose levels in an oral glucose tolerance test (OGTT) in fasted mice.

### Reference

1. Jiang, N., Su, D., Chen, D., *et al.* Discovery of a novel glucagon-like peptide-1 (GLP-1) analogue from bullfrog and investigation of its potential for designing GLP-1-based multiagonists. *J. Med. Chem.* **67**(1), 180-198 (2024).

#### 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, 06/03/2024

#### 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