

# PRODUCT INFORMATION



## GLP-1 (7-13) (human, mouse, rat, bovine) (trifluoroacetate salt)

Item No. 41255

**Formal Name:** L-histidyl-L-alanyl-L- $\alpha$ -glutamylglycyl-L-threonyl-L-phenylalanyl-L-threonine, trifluoroacetate salt

**Synonyms:** Glucagon-like Peptide 1 (7-13), His-Ala-Glu-Gly-Thr-Phe-Thr-OH

**Peptide Sequence:** HAEGTFT-OH

**MF:**  $C_{33}H_{47}N_9O_{12} \cdot XCF_3COOH$

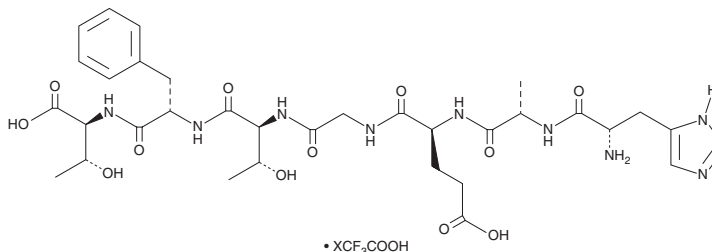
**FW:** 761.8

**Purity:**  $\geq 98\%$

**Supplied as:** A solid

**Storage:**  $-20^\circ\text{C}$

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



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

### Laboratory Procedures

GLP-1 (7-13) (human, mouse, rat, bovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the GLP-1 (7-13) (human, mouse, rat, bovine) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. GLP-1 (7-13) (human, mouse, rat, bovine) (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in DMSO.

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 (7-13) (human, mouse, rat, bovine) (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. GLP-1 (7-13) (human, mouse, rat, bovine) (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

Glucagon-like peptide 1 (GLP-1) (7-13) is a peptide fragment of the endogenous incretin hormone GLP-1 (Item Nos. 24460 | 24755).<sup>1</sup> It contains several amino acids involved in binding and activation of the GLP-1 receptor (GLP-1R).<sup>2</sup>

### References

1. Manandhar, B. and Ahn, J.-M. Glucagon-like peptide-1 (GLP-1) analogs: Recent advances, new possibilities, and therapeutic implications. *J. Med. Chem.* **58**(3), 1020-1037 (2015).
2. Adelhorst, K., Hedegaard, B.B., Knudsen, L.B., et al. Structure-activity studies of glucagon-like peptide-1. *J. Biol. Chem.* **69**(9), 6275-6278 (1994).

#### 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, 09/19/2024

#### 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