

PRODUCT INFORMATION

[Gly^{8,36},Glu²²]-GLP-1 (7-37) (human, bovine, guinea pig, mouse, rat) (acetate)

Item No. 40298

Formal Name:	L-histidylglycyl-L- α -glutamylglycyl-L-threonyl-L-phenylalanyl-L-threonyl-L-seryl-L- α -aspartyl-L-valyl-L-seryl-L-seryl-L-tyrosyl-L-leucyl-L- α -glutamyl-L- α -glutamyl-L-glutamyl-L-alanyl-L-alanyl-L-lysyl-L- α -glutamyl-L-phenylalanyl-L-isoleucyl-L-alanyl-L-tryptophyl-L-leucyl-L-valyl-L-lysylglycylglycyl-glycine, acetate	H—His—Gly—Glu—Gly—Thr—Phe—Thr—Ser—Asp—Val—Ser—Ser—Tyr—Leu—Glu—Glu—Gln—Ala—Ala—Lys—Glu—Phe—Ile—Ala—Trp—Leu—Val—Lys—Gly—Gly—Gly—OH • XCH ₃ CO ₂ H
Synonyms:	[Gly ^{8,36} ,Glu ²²]-Glucagon-like Peptide 1 (7-37),	
Peptide Sequence:	HGEGTFTSDVSSYLEEQAAKEFIAWLKGGG-OH	
MF:	C ₁₄₉ H ₂₂₁ N ₃₇ O ₄₉ • XC ₂ H ₄ O ₂	
FW:	3,314.6	
Purity:	≥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

[Gly^{8,36},Glu²²]-GLP-1 (7-37) (human, bovine, guinea pig, mouse, rat) (acetate) is supplied as a solid. A stock solution may be made by dissolving the [Gly^{8,36},Glu²²]-GLP-1 (7-37) (human, bovine, guinea pig, mouse, rat) (acetate) in the solvent of choice, which should be purged with an inert gas. [Gly^{8,36},Glu²²]-GLP-1 (7-37) (human, bovine, guinea pig, mouse, rat) (acetate) is soluble in methanol. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

[Gly^{8,36},Glu²²]-GLP-1 (7-37) is a derivative of GLP-1 containing alanine-to-glycine, glycine-to-glutamate, and arginine-to-glycine substitutions at positions 8, 22, and 36, respectively, and a peptide fragment of the GLP-1 receptor (GLP-1R) agonist dulaglutide.¹

Reference

1. Glaesner, W., Vick, A.M., Millican, R., *et al.* Engineering and characterization of the long-acting glucagon-like peptide-1 analogue LY2189265, an Fc fusion protein. *Diabetes Metab. Res. Rev.* **26(4)**, 287-296 (2010).

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, 04/11/2024

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