

PRODUCT INFORMATION



Albiglutide Fragment (sodium salt)

Item No. 42542

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 -glutamylglycyl-L-glutamyl-L-alanyl-L-alanyl-L-lysyl-L- α -glutamyl-L-phenylalanyl-L-isoleucyl-L-alanyl-L-tryptophyl-L-leucyl-L-valyl-L-lysylglycyl-L-argininamide, sodium salt

Synonyms: GLP-1 Gly⁸, GLP-1 8-glycine, [Gly⁸]-GLP-1 (7-36) amide

Peptide Sequence: HGEFTSDVSSYLEGQAAKEFIAWLVKGR-NH₂

MF: C₁₄₈H₂₂₄N₄₀O₄₅ • XNa

FW: 3,283.6

Purity: $\geq 95\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

H—His—Gly—Glu—Gly—Thr—Phe—Thr—Ser—Asp—Val—

Ser—Ser—Tyr—Leu—Glu—Gly—Gln—Ala—Ala—Lys—

Glu—Phe—Ile—Ala—Trp—Leu—Val—Lys—Gly—Arg—NH₂

• XNa

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

Laboratory Procedures

Albiglutide fragment (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the albiglutide fragment (sodium salt) in the solvent of choice, which should be purged with an inert gas. Albiglutide fragment (sodium salt) is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Albiglutide fragment is a peptide fragment of albiglutide, a fusion peptide with antidiabetic activity that contains two copies of [Gly⁸]-GLP-1 (7-36), which contains an alanine-to-glycine substitution that confers resistance to hydrolysis by dipeptidyl peptidase 4 (DPP-4), fused to albumin.¹ Albiglutide fragment comprises one copy of the DPP-4-resistant [Gly⁸]-GLP-1 (7-36) sequence.

Reference

1. Rendell, M.S. Albiglutide: A unique GLP-1 receptor agonist. *Expert Opin. Biol. Ther.* **16**(12), 1557-1569 (2016).

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, 01/29/2025

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