

PRODUCT INFORMATION

Elsiglutide (acetate)

Item No. 41273

MF: $C_{196}H_{323}N_{53}O_{56} \cdot XC_2H_4O_2$
FW: 4,318.0
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years

H — His — Gly — Glu — Gly — Ser — Phe — Ser — Ser — Glu — Leu —
 Ser — Thr — Ile — Leu — Asp — Ala — Leu — Ala — Ala — Arg —
 Asp — Phe — Ile — Ala — Trp — Leu — Ile — Ala — Thr — Lys —
 Ile — Thr — Asp — Lys — Lys — Lys — Lys — Lys — Lys — NH₂
 • XCH_3CO_2H

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

Laboratory Procedures

Elsiglutide (acetate) is supplied as a solid. Aqueous solutions of elsiglutide (acetate) can be prepared by directly dissolving the solid in aqueous buffers. Elsiglutide (acetate) is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Elsiglutide is a peptide derivative of glucagon-like peptide 2 (GLP-2).¹ It increases jejunum and ileum crypt depth and villus height without disrupting gut microbiota diversity in rats when administered at a dose of 0.9 mg/kg four days per week. Elsiglutide decreases diarrhea severity and incidence induced by the dual inhibitor of EGFR and HER2 lapatinib (Item No. 11493) in rats. It prevents lapatinib-induced loss of body weight in rats.

Reference

- Mayo, B.J., Secombe, K.R., Wignall, A.D., *et al.* The GLP-2 analogue elsiglutide reduces diarrhoea caused by the tyrosine kinase inhibitor lapatinib in rats. *Cancer Chemother. Pharmacol.* **85(4)**, 793-803 (2020).

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, 08/21/2024

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