

PRODUCT INFORMATION



GLP-1R Agonist 2

Item No. 41319

CAS Registry No.: 2230197-64-3
Formal Name: 2-[[[(2S)-4-[6-[(4-chloro-2-fluorophenyl)methoxy]-2-pyridinyl]-2-methyl-1-piperazinyl]methyl]-1-[(2S)-2-oxetanylmethyl]-1H-benzimidazole-6-carboxylic acid

Synonyms: GLP-1 Receptor Agonist 2,
Glucagon-like Peptide 1 Receptor Agonist 2

MF: C₃₀H₃₁ClFN₅O₄

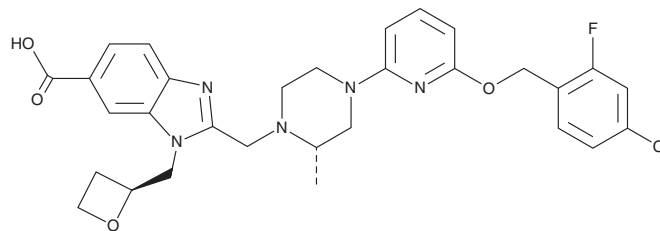
FW: 580.1

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

GLP-1R agonist 2 is supplied as a solid. A stock solution may be made by dissolving the GLP-1R agonist 2 in the solvent of choice, which should be purged with an inert gas. GLP-1R agonist 2 is sparingly soluble (1-10 mg/ml) in acetonitrile and DMSO.

Description

GLP-1R agonist 2 is a glucagon-like peptide 1 receptor agonist.¹ It induces cAMP accumulation in CHO cells expressing human GLP-1R (EC₅₀ = 0.28 nM).

Reference

1. Aspnes, G.E., Bagley, S.W., Curto, J.M., *et al.* GLP-1 receptor agonists and uses thereof. *Pfizer, Inc.* CA2988721A1 (2018).

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, 07/17/2024

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