

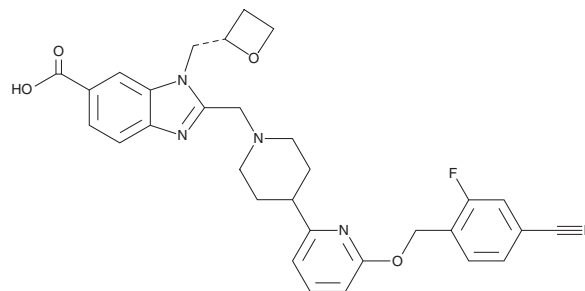
PRODUCT INFORMATION



PF-06882961

Item No. 30657

CAS Registry No.: 2230198-02-2
Formal Name: 2-[[4-[6-[(4-cyano-2-fluorophenyl)methoxy]-2-pyridinyl]-1-piperidinyl)methyl]-1-[(2S)-2-oxetanylmethyl]-1H-benzimidazole-6-carboxylic acid
Synonym: Danuglipron
MF: C₃₁H₃₀FN₅O₄
FW: 555.6
Purity: ≥95%
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

PF-06882961 is supplied as a solid. A stock solution may be made by dissolving the PF-06882961 in the solvent of choice, which should be purged with an inert gas. PF-06882961 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

PF-06882961 is an agonist of the glucagon-like peptide 1 receptor (GLP-1R).¹ It induces cAMP accumulation in CHO cells expressing the human receptor (EC₅₀ = 0.71 nM). Oral administration of PF-06882961 enhances insulin secretion in a glucose tolerance test, and subcutaneous administration decreases food intake in cynomolgus monkeys.

Reference

1. Griffith, D.A., Edmonds, D.J., Fortin, J.P., *et al.* A small-molecule oral agonist of the human glucagon-like peptide-1 receptor. *J. Med. Chem.* **65**(12), 8208-8226 (2022).

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, 07/22/2026

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