

# PRODUCT INFORMATION

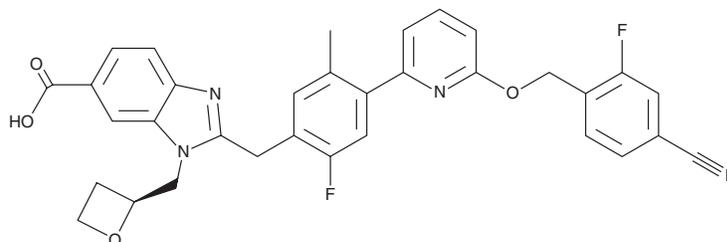


## Naperiglipron

Item No. 45151

**CAS Registry No.:** 2572566-11-9  
**Formal Name:** 2-[[4-[6-[(4-cyano-2-fluorophenyl)methoxy]-2-pyridinyl]-2-fluoro-5-methylphenyl)methyl]-1-[(2S)-2-oxetanylmethyl]-1H-benzimidazole-6-carboxylic acid

**Synonym:** LY3549492  
**MF:** C<sub>33</sub>H<sub>26</sub>F<sub>2</sub>N<sub>4</sub>O<sub>4</sub>  
**FW:** 580.6  
**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

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

### Description

Naperiglipron is a glucagon-like peptide 1 receptor (GLP-1R) agonist.<sup>1</sup> It induces cAMP production in HEK293 cells expressing human GLP-1R (EC<sub>50</sub> = 1.14 nM). *In vivo*, naperiglipron reduces blood glucose levels in human GLP1R knock-in mice (ED<sub>50</sub> = 0.07 mg/kg).

### Reference

1. Coates, D.A., Fielgs, T., Ho, J.D., *et al.* Glucagon-like peptide1 receptor agonists. *Eli Lilly and Company, US2020/0407347A1* (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, 02/27/2026

#### 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