

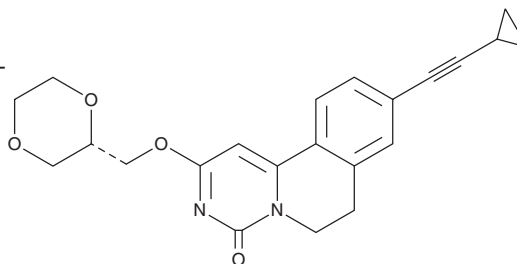
# PRODUCT INFORMATION



## GLPG1205

Item No. 33185

**CAS Registry No.:** 1445847-37-9  
**Formal Name:** 9-(2-cyclopropylethynyl)-2-[(2S)-1,4-dioxan-2-ylmethoxy]-6,7-dihydro-4H-pyrimido[6,1-a]isoquinolin-4-one  
**MF:** C<sub>22</sub>H<sub>22</sub>N<sub>2</sub>O<sub>4</sub>  
**FW:** 378.4  
**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

GLPG1205 is supplied as a solid. A stock solution may be made by dissolving the GLPG1205 in the solvent of choice, which should be purged with an inert gas. GLPG1205 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

### Description

GLPG1205 is a G protein-coupled receptor 84 (GPR84) antagonist.<sup>1</sup> It inhibits GTPγS binding induced by 3,3'-diindolylmethane (DIM) in cells overexpressing human GPR84 (IC<sub>50</sub> = 54 nM). GLPG1205 inhibits embelin-induced chemotaxis of isolated human neutrophils (IC<sub>50</sub> = 11 nM). *In vivo*, GLPG1205 (3 mg/kg) reduces body weight loss and rectal bleeding and improves stool consistency in a rat model of colitis induced by dextran sulfate sodium (DSS; Item No. 23250).

### Reference

1. Labéguère, F., Dupont, S., Alvey, L., *et al.* Discovery of 9-cyclopropylethynyl-2-[(S)-1-[1,4]dioxan-2-ylmethoxy]-6,7-dihydropyrimido[6,1-a]isoquinolin-4-one (GLPG1205), a unique GPR84 negative allosteric modulator undergoing evaluation in a phase II clinical trial. *J. Med. Chem.* **63(22)**, 13526-13545 (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, 11/13/2024

#### 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