

# PRODUCT INFORMATION



## GLPG3970

Item No. 45730

**CAS Registry No.:** 2403733-82-2  
**Formal Name:** 3,4-dihydro-8-methoxy-6-[7-[2-(4-morpholinyl)ethoxy]imidazo[1,2-a]pyridin-3-yl]-2-(2,2,2-trifluoroethyl)-1(2H)-isoquinolinone

**MF:** C<sub>25</sub>H<sub>27</sub>F<sub>3</sub>N<sub>4</sub>O<sub>4</sub>

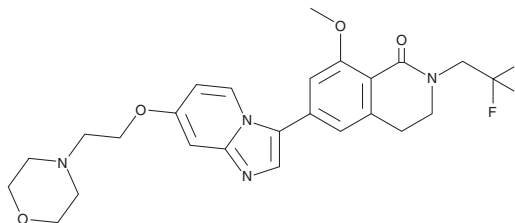
**FW:** 504.5

**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

GLPG3970 is supplied as a solid. A stock solution may be made by dissolving the GLPG3970 in the solvent of choice, which should be purged with an inert gas. GLPG3970 is soluble (≥10 mg/ml) in ethanol and DMSO.

### Description

GLPG3970 is an inhibitor of salt-inducible kinase 1 (SIK1), SIK2, SIK3, and receptor-interacting serine/threonine kinase 2 (RIPK2; IC<sub>50</sub>s = 282.8, 7.8, 3.8, and 78.4 nM, respectively).<sup>1</sup> It is selective for SIK1, SIK2, SIK3, and RIPK2 over a panel of 371 other kinases at 1 μM. GLPG3970 reduces the levels of TNF-α and IL-12 and increases the level of IL-10 in LPS-stimulated primary human monocytes and whole blood. It decreases disease severity in a mouse model of colitis induced by dextran sodium sulfate (DSS) when administered at a dose of 30 mg/kg.

### Reference

1. Peixoto, C., Joncour, A., Temal-Laib, T., *et al.* Discovery of clinical candidate GLPG3970: A potent and selective dual SIK2/SIK3 inhibitor for the treatment of autoimmune and inflammatory diseases. *J. Med. Chem.* **67**(7), 5233-5258 (2024).

#### 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, 05/28/2026

#### 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](http://WWW.CAYMANCHEM.COM)