

# PRODUCT INFORMATION

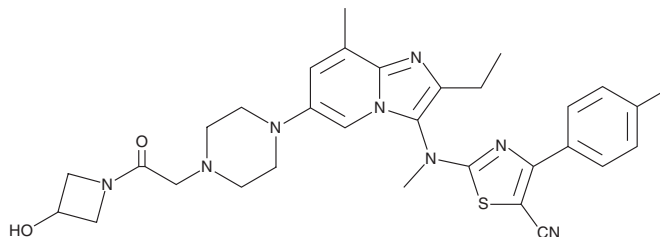


## GLPG1690

Item No. 25498

**CAS Registry No.:** 1628260-79-6  
**Formal Name:** 2-[[2-ethyl-6-[4-[2-(3-hydroxy-1-azetidiny)]-2-oxoethyl]-1-piperazinyl]-8-methylimidazo[1,2-a]pyridin-3-yl]methylamino]-4-(4-fluorophenyl)-5-thiazolecarbonitrile

**Synonym:** Ziritaxestat  
**MF:** C<sub>30</sub>H<sub>33</sub>FN<sub>8</sub>O<sub>2</sub>S  
**FW:** 588.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 252 nm  
**Supplied as:** A crystalline 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

GLPG1690 is supplied as a crystalline solid. A stock solution may be made by dissolving the GLPG1690 in the solvent of choice, which should be purged with an inert gas. GLPG1690 is soluble in the organic solvent DMSO at a concentration of approximately 80 mg/ml.

### Description

GLP1690 is a potent inhibitor of autotaxin ( $K_i = 15 \text{ nM}$ ).<sup>1</sup> It inhibits the production of lysophosphatidic acid (LPA) in mouse, rat, and human plasma ( $IC_{50}$ s = 418, 542, and 242 nM, respectively). GLPG1690 (30 mg/kg) reduces extracellular matrix deposition in the lung as well as LPA 18:2 levels in bronchoalveolar lavage fluid in a mouse model of pulmonary fibrosis induced by bleomycin (Item No. 13877).

### Reference

1. Liu, Y., Laufer, R., Patel, N.K., *et al.* Discovery of pyrazolo[1,5-a]pyrimidine TTK inhibitors: CFI-402257 is a potent, selective, bioavailable anticancer agent. *ACS Med. Chem. Lett.* **7**(7), 671-675 (2016).

#### 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, 08/11/2023

#### 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