

PRODUCT INFORMATION



GLPG1837

Item No. 9003143

CAS Registry No.: 1654725-02-6
Formal Name: N-[3-(aminocarbonyl)-4,7-dihydro-5,5,7,7-tetramethyl-5H-thieno[2,3-c]pyran-2-yl]-1H-pyrazole-3-carboxamide

Synonym: ABBV-974

MF: C₁₆H₂₀N₄O₃S

FW: 348.4

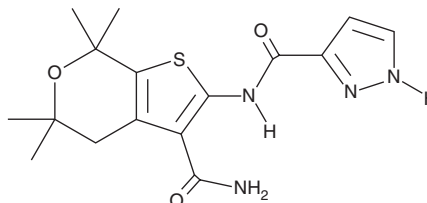
Purity: ≥98%

UV/Vis.: λ_{max}: 226, 326 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

GLPG1837 is supplied as a crystalline solid. A stock solution may be made by dissolving the GLPG1837 in the solvent of choice, which should be purged with an inert gas. GLPG1837 is soluble in DMSO.

Description

GLPG1837 is a potentiator of cystic fibrosis transmembrane conductance regulator (CFTR).^{1,2} It increases forskolin-induced opening of wild-type and mutant CFTRs expressing the F508 deletion (F508del) and G551D missense mutations with EC₅₀ values of 88, 3, and 339 nM, respectively, in a reporter assay. GLPG1837 increases forskolin-induced conductivity of patient-derived bronchial epithelial cells expressing both F508del- and G551D-mutant CFTRs (EC₅₀ = 159 nM in a transepithelial clamp circuit (TECC) assay).¹

References

1. Gees, M., Musch, S., Van der Plas, S., *et al.* Identification and characterization of novel CFTR potentiators. *Front. Pharmacol.* **9:1221**, (2018).
2. Van der Plas, S., Kelgtermans, H., De Munck, T., *et al.* Discovery of N (3-carbamoyl-5,5,7,7-tetramethyl-5,7-dihydro 4H thieno[2,3 c]pyran-2-yl) 1H pyrazole-5-carboxamide (GLPG1837), a novel potentiator which can open class III mutant cystic fibrosis transmembrane conductance regulator (CFTR) channels to a high extent. *J. Med. Chem.* **61(4)**, 1425-1435 (2018).

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, 10/31/2019

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