

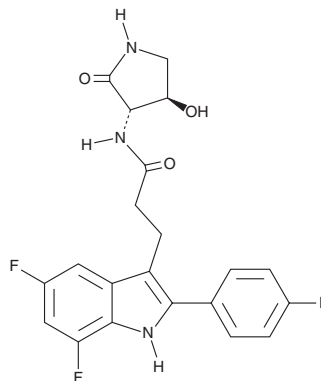
PRODUCT INFORMATION



Inaxaplin

Item No. 40852

CAS Registry No.: 2446816-88-0
Formal Name: 5,7-difluoro-2-(4-fluorophenyl)-N-[(3S,4R)-4-hydroxy-2-oxo-3-pyrrolidinyl]-1H-indole-3-propanamide
Synonym: VX-147
MF: C₂₁H₁₈F₃N₃O₃
FW: 417.4
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

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

Description

Inaxaplin is an apolipoprotein L1 (ApoL1) inhibitor.¹ It inhibits cytotoxicity induced by overexpression of the *APOL1* variants G1 and G2 in cells when used at a concentration of 3 μM. Inaxaplin also reduces ApoL1 G1-induced inhibition of global protein synthesis in podocytes isolated from *APOL1* transgenic mice. *In vivo*, inaxaplin (10 mg/kg) reduces podocyte injury and proteinuria in *APOL1* transgenic mice.

Reference

1. Datta, S., Antonio, B.M., Zahler, N.H., *et al.* *APOL1*-mediated monovalent cation transport contributes to *APOL1*-mediated podocytopathy in kidney disease. *J. Clin. Invest.* **134**(5), (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, 10/08/2024

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