

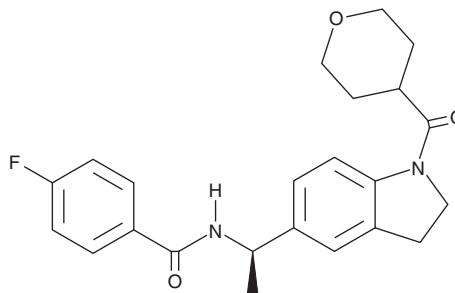
PRODUCT INFORMATION



LY3381916

Item No. 33178

CAS Registry No.: 2166616-75-5
Formal Name: N-[(1R)-1-[2,3-dihydro-1-[(tetrahydro-2H-pyran-4-yl)carbonyl]-1H-indol-5-yl]ethyl]-4-fluoro-benzamide
Synonym: IDO1-IN-5
MF: C₂₃H₂₅FN₂O₃
FW: 396.5
Purity: ≥98%
UV/Vis.: λ_{max}: 260 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

LY3381916 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY3381916 in the solvent of choice, which should be purged with an inert gas. LY3381916 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of LY3381916 in these solvents is approximately 30 mg/ml.

Description

LY3381916 is an inhibitor of indoleamine 2,3-dioxygenase 1 (IDO1).¹ It binds to apo-IDO1, which lacks heme, but does not inhibit mature heme-bound IDO1.

Reference

1. Dorsey, F.C., Benhadji, K.A., Sams, L.L., *et al.* Abstract 5245: Identification and characterization of the IDO1 inhibitor LY3381916. *Cancer Res.* **78(13 Supplement)**, 5245 (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, 11/14/2022

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