

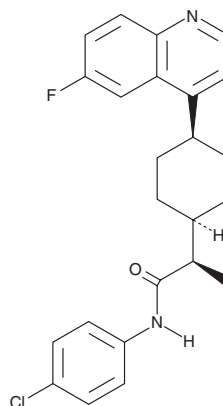
PRODUCT INFORMATION



BMS 986205

Item No. 25025

CAS Registry No.: 1923833-60-6
Formal Name: cis-(αR)-N-(4-chlorophenyl)-4-(6-fluoro-4-quinolinyl)-α-methyl-cyclohexaneacetamide
Synonym: Linrodostat
MF: C₂₄H₂₄ClFN₂O
FW: 410.9
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 251 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

BMS 986205 is supplied as a crystalline solid. A stock solution may be made by dissolving the BMS 986205 in the solvent of choice, which should be purged with an inert gas. BMS 986205 is soluble in the organic solvent DMSO at a concentration of approximately 90 mg/ml.

Description

BMS 986205 is an inhibitor of indoleamine 2,3-dioxygenase 1 (IDO1; IC₅₀ = <50 nM).¹ It inhibits kynurenine synthesis in HeLa cells and IDO1-overexpressing HEK293 cells (IC₅₀s = 1.1 and 1.7 nM, respectively).

Reference

1. Cheong, J.E., Ekkati, A., and Sun, L. A patent review of IDO1 inhibitors for cancer. *Exp. Opin. Ther. Pat.* 28(4), 317-330 (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, 10/21/2022

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