

PRODUCT INFORMATION

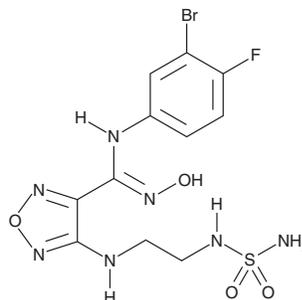


Epacadostat

Item No. 19875

CAS Registry No.: 1204669-58-8
Formal Name: [C(Z)]-4-[[2-[(aminosulfonyl)amino]ethyl]amino]-N-(3-bromo-4-fluorophenyl)-N'-hydroxy-1,2,5-oxadiazole-3-carboximidamide

Synonym: INCB024360
MF: C₁₁H₁₃BrFN₇O₄S
FW: 438.2
Purity: ≥98%
UV/Vis.: λ_{max}: 290 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

Epacadostat is supplied as a crystalline solid. A stock solution may be made by dissolving the epacadostat in the solvent of choice, which should be purged with an inert gas. Epacadostat is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of epacadostat in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Epacadostat is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, epacadostat should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Epacadostat has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Epacadostat is an inhibitor of indoleamine 2,3-dioxygenase 1 (IDO1) with an IC₅₀ value of 7.1 nM in HeLa cells that demonstrates little activity against the related enzymes IDO2 or tryptophan 2,3-dioxygenase (TDO).¹ It has been shown to promote T and natural killer-cell growth, to increase IFN-γ production, and to reduce conversion to regulatory T (Treg)-like cells in a coculture system of human allogeneic lymphocytes with either dendritic cells or tumor cells.¹ Epacadostat can also inhibit tumor growth in tumor-bearing mice in a lymphocyte-dependent manner.¹

Reference

1. Liu, X., Shin, N., Koblisch, H.K., *et al.* Selective inhibition of IDO1 effectively regulates mediators of antitumor immunity. *Blood* **115**(17), 3520-3530 (2010).

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, 10/26/2022

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