

PRODUCT INFORMATION



BI-2545

Item No. 24285

CAS Registry No.: 2162961-71-7

Formal Name: 3,5-bis(trifluoromethyl)benzyl
(1R,5S,6s)-6-((1H-benzo[d][1,2,3]
triazole-6-carboxamido)methyl)-
3-azabicyclo[3.1.0]hexane-3-
carboxylate

MF: C₂₃H₁₉F₆N₅O₃

FW: 527.4

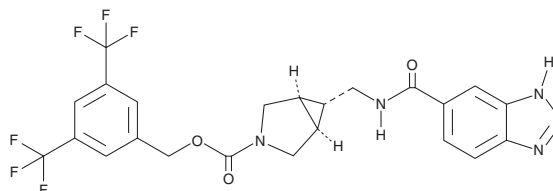
Purity: ≥98%

UV/Vis.: λ_{max}: 213 nm

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

BI-2545 is supplied as a solid. A stock solution may be made by dissolving the BI-2545 in the solvent of choice, which should be purged with an inert gas. BI-2545 is soluble in DMSO.

Description

BI-2545 is an inhibitor of autotaxin (IC₅₀s = 2.2 and 3.4 for the human and rat enzyme, respectively).¹ It is selective for autotaxin over a panel of 48 enzymes, transporters, ion channels, and receptors at 10 μM. BI-2545 (10 mg/kg) reduces plasma lysophosphatidic acid (LPA) levels in rats.

Reference

1. Kuttruff, C.A., Ferrara, M., Bretschneider, T., *et al.* Discovery of BI-2545: A novel autotaxin inhibitor that significantly reduces LPA levels *in vivo*. *ACS Med. Chem. Lett.* **8**(12), 1252-1257 (2017).

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, 12/17/2022

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