

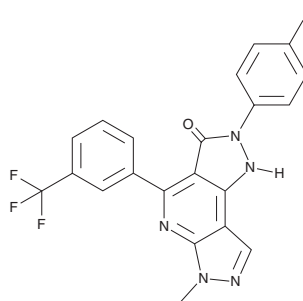
PRODUCT INFORMATION



B7-CD28 Interaction Inhibitor

Item No. 42398

CAS Registry No.: 635324-72-0
Formal Name: 2-(4-fluorophenyl)-1,6-dihydro-6-methyl-4-[3-(trifluoromethyl)phenyl]-dipyrzolo[3,4-b:3',4'-d]pyridin-3(2H)-one
MF: C₂₁H₁₃F₄N₅O
FW: 427.4
Purity: ≥95%
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

B7-CD28 interaction inhibitor is supplied as a solid. A stock solution may be made by dissolving the B7-CD28 interaction inhibitor in the solvent of choice, which should be purged with an inert gas. B7-CD28 interaction inhibitor is soluble (≥10 mg/ml) in DMSO.

Description

B7-CD28 interaction inhibitor is an inhibitor of the interaction between B7-1/CD80 and CD28 (IC₅₀ = 50 nM).¹

Reference

1. Green, N.J., Xiang, J., Chen, J., *et al.* Structure-activity studies of a series of dipyrzolo[3,4-b:3',4'-d]pyridin-3-ones binding to the immune regulatory protein B7.1. *Bioorg. Med. Chem.* **11(13)**, 2991-3013 (2003).

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, 11/27/2024

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