

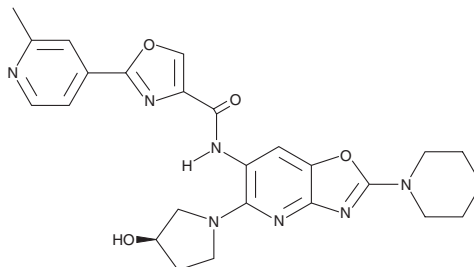
PRODUCT INFORMATION



Emavusertib

Item No. 36845

CAS Registry No.: 1801344-14-8
Formal Name: N-[5-[(3R)-3-hydroxy-1-pyrrolidinyl]-2-(4-morpholinyl)oxazolo[4,5-b]pyridin-6-yl]-2-(2-methyl-4-pyridinyl)-4-oxazolecarboxamide
Synonym: CA-4948
MF: C₂₄H₂₅N₇O₅
FW: 491.5
Purity: ≥98%
UV/Vis.: λ_{max}: 260, 345 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

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

Description

Emavusertib is an inhibitor of IL-1 receptor-associated kinase 4 (IRAK4; IC₅₀ = 31.7 nM).¹ It is selective for IRAK4 over FMS-related tyrosine kinase 3 (FLT3), cyclin-dependent kinase 2 (Cdk2), Aurora A kinase, glycogen synthase kinase 3β (GSK3β), and muscle-associated receptor tyrosine kinase (MUSK) at 1 μM. Emavusertib inhibits IL-6 release mediated by toll-like receptor 2 (TLR2), TLR5, or IL-1 receptor (IL-1R) in isolated human whole blood (IC₅₀s = 989, 696, and 1,375 nM, respectively). *In vivo*, emavusertib (25-200 mg/kg) reduces tumor volume in an OCI-Ly3 diffuse large B cell lymphoma (DLBCL) mouse xenograft model.

Reference

1. Gummadi, V.R., Boruah, A., Ainan, B.R., *et al.* Discovery of CA-4948, an orally bioavailable IRAK4 inhibitor for treatment of hematologic malignancies. *ACS Med. Chem. Lett.* **11(12)**, 2374-2381 (2020).

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, 12/21/2022

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