

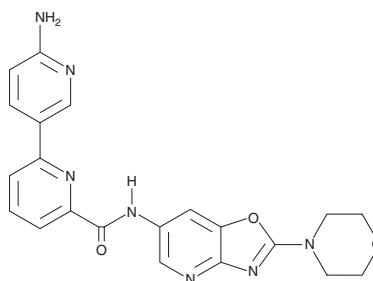
PRODUCT INFORMATION



CA-4948

Item No. 27598

CAS Registry No.: 1801343-74-7
Formal Name: 6'-amino-N-[2-(4-morpholinyl)oxazolo[4,5-b]pyridin-6-yl]-[2,3'-bipyridine]-6-carboxamide
MF: C₂₁H₁₉N₇O₃
FW: 417.4
Purity: ≥98%
UV/Vis.: λ_{max}: 315 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.

Description

CA-4948 is an orally bioavailable inhibitor of interleukin-1 receptor associated kinase-4 (IRAK-4).¹ It has antitumor activity in ABC-type diffuse large B cell lymphoma patient-derived xenograft (PDX) mouse models.

Reference

1. Booher, R.N., Samson, M.E., Xu, G.-X., *et al.* Abstract 1168: Efficacy of the IRAK4 inhibitor CA-4948 in patient-derived xenograft models of diffuse large B cell lymphoma. *Cancer Res.* **77(13 Supp)**, 1168 (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, 12/19/2022

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