

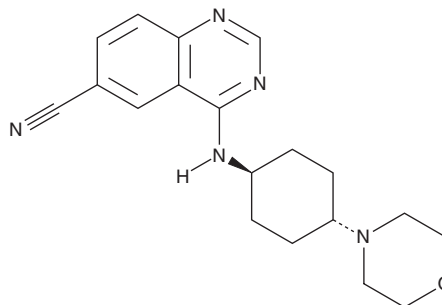
PRODUCT INFORMATION



IRAK4 Inhibitor 1

Item No. 37548

CAS Registry No.: 1820787-94-7
Formal Name: 4-[[*trans*-4-(4-morpholinyl)cyclohexyl]amino]-6-quinazolinecarbonitrile
Synonym: IRAK4-IN-1
MF: C₁₉H₂₃N₅O
FW: 337.4
Purity: ≥98%
UV/Vis.: λ_{max}: 217 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

IRAK4 Inhibitor 1 (IRAK4-IN-1) is supplied as a solid. A stock solution may be made by dissolving the IRAK4-IN-1 in the solvent of choice, which should be purged with an inert gas. IRAK4-IN-1 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of IRAK4-IN-1 in these solvents is approximately 10 and 0.5 mg/ml, respectively.

Description

IRAK4-IN-1 is an inhibitor of IL-1 receptor-associated kinase 4 (IRAK4; IC₅₀ = 7 nM).¹ It reduces TNF-α level increases induced by the toll-like receptor 7 (TLR7) agonist resiquimod (R-848; Item No. 14806) in isolated human peripheral blood mononuclear cells (PBMCs; EC₅₀ = 300 nM).

Reference

1. Smith, G.F., Altman, M.D., Andresen, B., *et al.* Identification of quinazoline based inhibitors of IRAK4 for the treatment of inflammation. *Bioorg. Med. Chem. Lett.* **27(12)**, 2721-2726 (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, 06/08/2023

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