

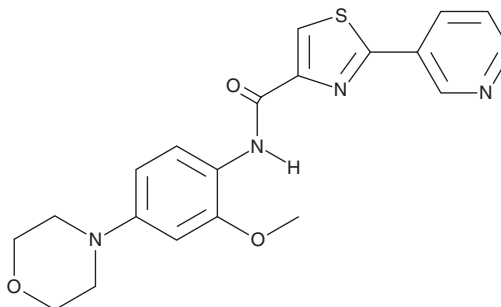
PRODUCT INFORMATION



IRAK Inhibitor 6

Item No. 33267

CAS Registry No.: 1042672-97-8
Formal Name: N-[2-methoxy-4-(4-morpholinyl)phenyl]-2-(3-pyridinyl)-4-thiazolecarboxamide
Synonyms: IL-1 Receptor-associated Kinase 4 Inhibitor 6, Interleukin-1 Receptor-associated Kinase 4 Inhibitor 6
MF: C₂₀H₂₀N₄O₃S
FW: 396.5
Purity: ≥98%
UV/Vis.: λ_{max}: 294 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.

Laboratory Procedures

IRAK inhibitor 6 is supplied as a crystalline solid. A stock solution may be made by dissolving the IRAK inhibitor 6 in the solvent of choice, which should be purged with an inert gas. IRAK inhibitor 6 is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

IRAK inhibitor 6 is an inhibitor of IL-1 receptor-associated kinase 4 (IRAK4; IC₅₀ = 160 nM).¹

Reference

1. Buckley, G.M., Gowers, L., Higuero, A.P., et al. IRAK-4 inhibitors. Part 1: A series of amides. *Bioorg. Med. Chem. Lett.* **18(11)**, 3211-3214 (2008).

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, 09/27/2024

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