

PRODUCT INFORMATION



KT474

Item No. 41053

CAS Registry No.: 2432994-31-3

Formal Name: N-[3-(difluoromethyl)-1-[*trans*-4-[[4-[[3-[1-(2,6-dioxo-3-piperidinyl)-2,3-dihydro-3-methyl-2-oxo-1H-benzimidazol-4-yl]-2-propyn-1-yl]oxy]-1-piperidinyl]methyl]cyclohexyl]-1H-pyrazol-4-yl]-5-(1R,4R)-2-oxa-5-azabicyclo[2.2.1]hept-5-yl-pyrazolo[1,5-a]pyrimidine-3-carboxamide

Synonym: SAR444656

MF: C₄₄H₄₉F₂N₁₁O₆

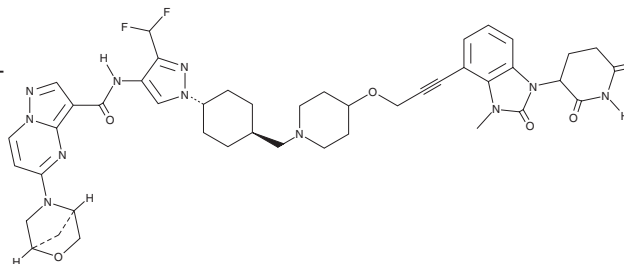
FW: 865.9

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

KT474 is supplied as a solid. A stock solution may be made by dissolving the KT474 in the solvent of choice, which should be purged with an inert gas. KT474 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

KT474 is a proteolysis-targeting chimera (PROTAC) that drives the degradation of IL-1 receptor-associated kinase 4 (IRAK4).¹ It induces the degradation of IRAK4 in primary mouse and rat splenocytes and isolated dog, monkey, and human peripheral blood mononuclear cells (PBMCs) with half-maximal degradation concentration (DC₅₀) values of 4.2, 2.1, 9.2, 2.9, and 1.5 nM, respectively. KT474 is selective for IRAK4 over a panel of 468 kinases at 1 μM. It inhibits IL-8 release induced by R-848 (Item No. 14806) in isolated human PBMCs (IC₅₀ = 3 nM). *In vivo*, KT474 (30 and 100 mg/kg, p.o.) decreases ear thickness in a mouse model of psoriasis induced by imiquimod (Item No. 14956).

Reference

1. Zheng, Z., Ji, N., Campbell, V., *et al.* Discovery of KT-474—a potent, selective, and orally bioavailable IRAK4 degrader for the treatment of autoimmune diseases. *J. Med. Chem.* **67**(20), 18022-18037 (2024).

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, 04/10/2025

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