

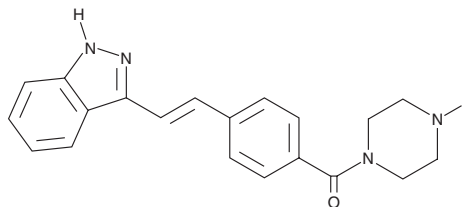
PRODUCT INFORMATION



KW 2449

Item No. 22207

CAS Registry No.: 1000669-72-6
Formal Name: [4-[2-(1H-indazol-3-yl)ethenyl]phenyl]-1-piperazinyl-methanone
MF: C₂₀H₂₀N₄O
FW: 332.4
Purity: ≥98%
UV/Vis.: λ_{max}: 266, 330 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

KW 2449 is supplied as a crystalline solid. A stock solution may be made by dissolving the KW 2449 in the solvent of choice, which should be purged with an inert gas. KW 2449 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of KW 2449 in these solvents is approximately 30 mg/ml.

KW 2449 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, KW 2449 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. KW 2449 has a solubility of approximately 0.17 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

KW 2449 is a multi-kinase inhibitor of FMS-related tyrosine kinase 3 (FLT3), Abelson tyrosine-protein kinase 1 (ABL), ABL-T315I, and Aurora kinase (IC₅₀s = 6.6, 14, 4, and 48 nM, respectively).¹ It has growth inhibitory activity against leukemia cells expressing FLT3 with activating mutations (GI₅₀s = 11-46 nM) and suppresses phosphorylation of FLT3 and STAT5 in MOLM-13 cells in a dose-dependent manner *in vitro*. KW 2499 (100 nM) inhibits colony formation of human primary acute myeloid leukemia (AML) cells with wild-type or activated mutant FLT3. It also induces MOLM-13 xenograft regression in a dose-dependent manner *in vivo*.

Reference

1. Shiotsu, Y., Kiyoi, H., Ishikawa, Y., et al. KW-2449, a novel multikinase inhibitor, suppresses the growth of leukemia cells with FLT3 mutations or T315I-mutated BCR/ABL translocation. *Blood* **114**(8), 1607-1617 (2009).

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, 01/03/2023

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