

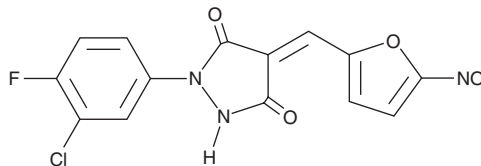
PRODUCT INFORMATION



PYZD-4409

Item No. 44131

CAS Registry No.: 423148-78-1
Formal Name: 1-(3-chloro-4-fluorophenyl)-4-[(5-nitro-2-furanyl)methylene]-3,5-pyrazolidinedione
MF: C₁₄H₇ClFN₃O₅
FW: 351.7
Purity: ≥98%
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

PYZD-4409 is supplied as a solid. A stock solution may be made by dissolving the PYZD-4409 in the solvent of choice, which should be purged with an inert gas. PYZD-4409 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

PYZD-4409 is an inhibitor of ubiquitin-activating enzyme (UAE; IC₅₀ = 20 μM in a cell-free assay).¹ It decreases levels of ubiquitinated cell division cycle 34 (Cdc34) in K562 chronic myeloid leukemia (CML) cells when used at a concentration of 50 μM. PYZD-4409 increases levels of cyclin D3 and p53 in HCT116 colorectal cancer cells. It selectively reduces colony formation in primary patient-derived acute myeloid leukemia (AML) cells over primary human hematopoietic cells when used at a concentration of 10 μM. PYZD-4409 (10 mg/kg) decreases tumor weight in an MDAY-D2 murine lymphoma model.

Reference

1. Xu, G.W., Ali, M., Wood, T.E., *et al.* The ubiquitin-activating enzyme E1 as a therapeutic target for the treatment of leukemia and multiple myeloma. *Blood* **115(11)**, 2251-2259 (2021).

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, 09/02/2025

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