

PRODUCT INFORMATION



UNC-CA2-103

Item No. 41241

Formal Name: (Z)-N-(3-((4-methyl-1H-imidazol-5-yl)methylene)-2-oxoindolin-5-yl)benzamide

MF: C₂₀H₁₆N₄O₂

FW: 344.4

Purity: ≥98%

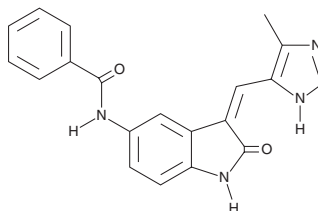
Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Special Conditions: Light sensitive; sensitive to protic solvents

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.



Laboratory Procedures

UNC-CA2-103 is supplied as a solid. A stock solution may be made by dissolving the UNC-CA2-103 in the solvent of choice, which should be purged with an inert gas. UNC-CA2-103 is soluble (≥10 mg/ml) in DMSO.

Description

UNC-CA2-103 is an inhibitor of tousel-like kinase 2 (TLK2; IC₅₀ = 18 nM).¹ It is selective for TLK2 over TLK1, tropomyosin-related kinase C (TrkC), and FMS-related tyrosine kinase 3 (FLT3) at 500 nM and a panel of 468 kinases at 1 μM, but it does inhibit 11 kinases by greater than 90% in the same panel.

Reference

1. Asquith, C.R.M., East, M.P., Laitinen, T., *et al.* Discovery and optimization of narrow spectrum inhibitors of Tousel like kinase 2 (TLK2) using quantitative structure activity relationships. *BioRxiv* **271(6)**, 116357 (2023).

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, 11/08/2024

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