

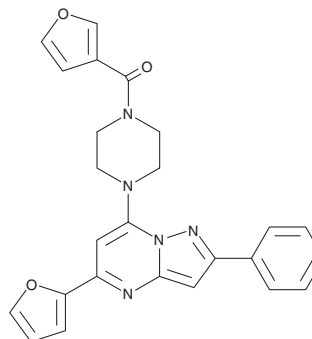
PRODUCT INFORMATION



ML-230

Item No. 32552

CAS Registry No.: 1776055-05-0
Formal Name: 3-furanyl[4-[5-(2-furanyl)-2-phenylpyrazolo[1,5-a]pyrimidin-7-yl]-1-piperazinyl]-methanone
Synonym: CID-44640177
MF: C₂₅H₂₁N₅O₃
FW: 439.5
Purity: ≥98%
UV/Vis.: λ_{max}: 285, 296 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

ML-230 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-230 in the solvent of choice, which should be purged with an inert gas. ML-230 is soluble in DMSO.

Description

ML-230 is an inhibitor of ATP-binding cassette family G member 2 (ABCG2).¹ It selectively inhibits JC-1 efflux in Ig-MXP3 cells overexpressing ABCG2 over Jurkat cells overexpressing ABCB1 (IC₅₀s = 0.13 and 4.65 μM, respectively). ML-230 enhances inhibition of IGROV-1/T8 ovarian cancer cell colony formation induced by topotecan (Item No. 14129).² Intratumoral injection of ML-230 (100 nM) and topotecan reduces tumor size in an IGROV-1/T8 mouse xenograft model.

References

1. Strouse, J.J., Ivnitski-Steele, I., Khawaja, H.M., *et al.* A selective ATP-binding cassette subfamily G member 2 efflux inhibitor revealed via high-throughput flow cytometry. *J. Biomol. Screen.* **18(1)**, 26-38 (2013).
2. Ricci, J.W., Lovato, D.M., Severns, V., *et al.* Novel ABCG2 antagonists reverse topotecan-mediated chemotherapeutic resistance in ovarian carcinoma xenografts. *Mol. Cancer Ther.* **15(12)**, 2853-2862 (2016).

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, 04/13/2023

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