

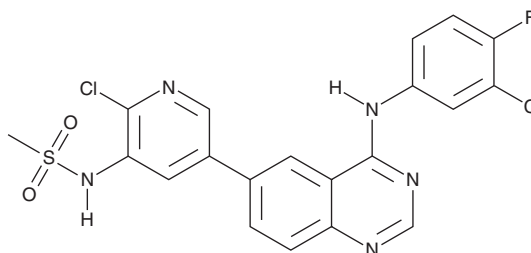
PRODUCT INFORMATION



MTX-211

Item No. 45142

CAS Registry No.: 1952236-05-3
Formal Name: N-[2-chloro-5-[4-[(3-chloro-4-fluorophenyl)amino]-6-quinazoliny]-3-pyridinyl]-methanesulfonamide
MF: C₂₀H₁₄Cl₂FN₅O₂S
FW: 478.3
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

MTX-211 is supplied as a solid. A stock solution may be made by dissolving the MTX-211 in the solvent of choice, which should be purged with an inert gas. MTX-211 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile:water (1:1).

Description

MTX-211 is a dual inhibitor of PI3K α and EGFR (IC₅₀s = <100 nM for both).¹ It reduces viability of 5637, EJ, and UM-UC-3 bladder cancer cells (IC₅₀s = 1.075, 1.514, and 2.189 μ M, respectively), an effect that can be partially inhibited by glutathione (GSH).² MTX-211 also induces apoptosis, increases the levels of the Nrf2 binding protein Keap1, and reduces the levels of Nrf2 and GSH in 5637 and EJ cells. It reduces tumor growth in a 5637 mouse xenograft model when administered at a dose of 10 mg/kg.

References

1. Whitehead, C.E. and Leopold, J.S. Small molecule inhibitors of EGFR and PI3K. *The Regents of the University of Michigan*. **WO 2016/100347 A2** (2016).
2. Hu, B., Chen, R., Jiang, M., *et al.* MTX-211 inhibits GSH synthesis through kea1/NRF2/GCLM axis and exerts antitumor effects in bladder cancer. *Int. J. Mol. Sci.* **24**, 7608 (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, 05/12/2026

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