

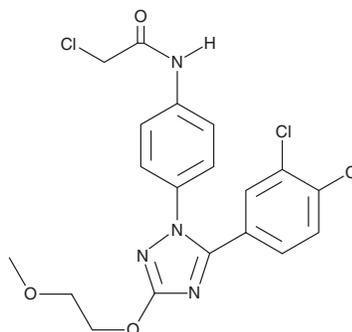
PRODUCT INFORMATION



MI 2

Item No. 23437

CAS Registry No.: 1047953-91-2
Formal Name: 2-chloro-N-[4-[5-(3,4-dichlorophenyl)-3-(2-methoxyethoxy)-1H-1,2,4-triazol-1-yl]phenyl]-acetamide
MF: C₁₉H₁₇Cl₃N₄O₃
FW: 455.7
Purity: ≥98%
UV/Vis.: λ_{max}: 253 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

MI 2 is supplied as a crystalline solid. A stock solution may be made by dissolving the MI 2 in the solvent of choice, which should be purged with an inert gas. MI 2 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of MI 2 in ethanol is approximately 16 mg/ml and approximately 33 mg/ml in DMSO and DMF.

MI 2 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MI 2 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. MI 2 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

MI 2 is an inhibitor of the paracaspase MALT1 with an IC₅₀ value of 5.84 μM in a fluorescence assay.¹ It inhibits the growth of MALT1-dependent activated B cell-like diffuse large B cell lymphoma (ABC-DLBCL) cell lines (GI₅₀s = 0.2, 0.5, 0.4, and 0.4 μM for HBL-1, TMD8, OCI-Ly3, and OCI-Ly10 cells, respectively). MI 2 inhibits cleavage of the MALT1 target protein CYLD in a dose-dependent manner in HBL-1 cells but does not inhibit caspase-3, caspase-8, and caspase-9, which are structurally similar to MALT1. MI 2 (25 mg/kg, i.v.) reduces tumor size in TMD8 and HBL-1 mouse xenograft models.

Reference

1. Fontan, L., Yang, C., Kabaleeswaran, V., et al. MALT1 small molecule inhibitors specifically suppress ABC-DLBCL *in vitro* and *in vivo*. *Cancer Cell*. **22(6)**, 812-824 (2012).

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, 11/15/2022

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