

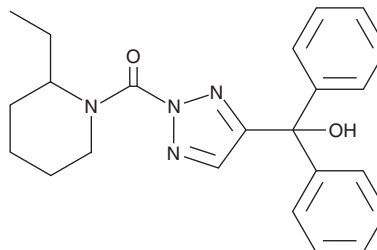
PRODUCT INFORMATION



ML-226

Item No. 25681

CAS Registry No.: 2055172-43-3
Formal Name: (2-ethyl-1-piperidinyl)
[4-(hydroxydiphenylmethyl)-2H-
1,2,3-triazol-2-yl]-methanone
MF: C₂₃H₂₆N₄O₂
FW: 390.5
Purity: ≥98%
UV/Vis.: λ_{max}: 239 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

ML-226 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-226 in the solvent of choice. ML-226 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of ML-226 in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

ML-226 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ML-226 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ML-226 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ML-226 is an inhibitor of α/β hydrolase domain-containing protein 11 (ABHD11) that inhibits ABHD11 *in vitro* and *in situ* (IC₅₀s = 15 and 0.68 nM, respectively).¹ It is 100-fold selective for ABHD11 over approximately 20 other serine hydrolases. ML-226 covalently carbamoylates the active site serine of ABHD11. It can be used as an anti-probe for ML-211 (Item No. 17630), an inhibitor of the protein palmitoyl thioesterases lysophospholipase 1 (LYPLA1) and LYPLA2.²

References

1. Adibekian, A., Hsu, K.-L., Speers, A.E., *et al.* Optimization and characterization of a triazole urea inhibitor for alpha/beta hydrolase domain-containing protein 11 (ABHD11): Anti-probe for LYPLA1/LYPLA2 dual inhibitor ML211. *Probe Reports from the NIH Molecular Libraries Program* (2010).
2. Adibekian, A., Martin, B.R., Speers, A.E., *et al.* Optimization and characterization of a triazole urea dual inhibitor for lysophospholipase 1 (LYPLA1) and lysophospholipase 2 (LYPLA2). *Probe Reports from the NIH Molecular Libraries Program* **1 R01 CA132630**, 1-42 (2013).

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, 10/16/2018

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