

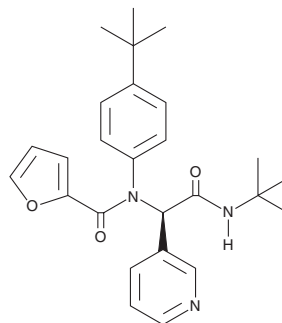
PRODUCT INFORMATION



ML-188

Item No. 37346

CAS Registry No.: 1417700-13-0
Formal Name: (αR)-N-(1,1-dimethylethyl)-α-[[4-(1,1-dimethylethyl)phenyl](2-furanylcarbonyl)amino]-3-pyridineacetamide
MF: C₂₆H₃₁N₃O₃
FW: 433.5
Purity: ≥98%
UV/Vis.: λ_{max}: 262 nm
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

ML-188 is supplied as a solid. A stock solution may be made by dissolving the ML-188 in the solvent of choice, which should be purged with an inert gas. ML-188 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of ML-188 in these solvents is approximately 30 mg/ml.

Description

ML-188 is an inhibitor of severe acute respiratory coronavirus (SARS-CoV) main protease (M^{pro}), also known as 3C-like protease (3CL^{pro}; IC₅₀ = 1.5 μM).¹

Reference

1. Jacobs, J., Grum-Tokars, V., Zhou, Y., *et al.* Discovery, synthesis, and structure-based optimization of a series of N-(tert-butyl)-2-(N-arylamido)-2-(pyridin-3-yl) acetamides (ML188) as potent noncovalent small molecule inhibitors of the severe acute respiratory syndrome coronavirus (SARS-CoV) 3CL protease. *J. Med. Chem.* **56**(2), 534-546 (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, 05/15/2023

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