

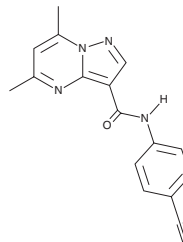
PRODUCT INFORMATION



ML-198

Item No. 37755

CAS Registry No.: 1380716-06-2
Formal Name: N-(4-ethynylphenyl)-5,7-dimethyl-pyrazolo[1,5-a]pyrimidine-3-carboxamide
Synonym: NCGC 758
MF: C₁₇H₁₄N₄O
FW: 290.3
Purity: ≥98%
UV/Vis.: λ_{max}: 215, 244, 293 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-198 is supplied as a solid. A stock solution may be made by dissolving the ML-198 in the solvent of choice, which should be purged with an inert gas. ML-198 is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

Description

ML-198 is a chemical chaperone of glucocerebrosidase.¹ It activates hydrolysis of the fluorescent glucocerebrosidase substrate 4-methylumbelliferyl-β-D-glucopyranoside (4-MUG; Item No. 20948) in spleen homogenate isolated from patients with Gaucher disease with the homozygous N370S mutant of glucocerebrosidase (glucocerebrosidase^{N370S}; AC₅₀ = 0.36 μM). ML-198 (5 μM) increases the translocation of glucocerebrosidase^{N370S} to lysosomes in dermal fibroblasts isolated from patients with Gaucher disease.

Reference

1. Patnaik, S., Zheng, W., Choi, J.H., *et al.* Discovery, structure-activity relationship, and biological evaluation of noninhibitory small molecule chaperones of glucocerebrosidase. *J. Med. Chem.* **55**(12), 5734-5748 (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, 04/20/2023

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