

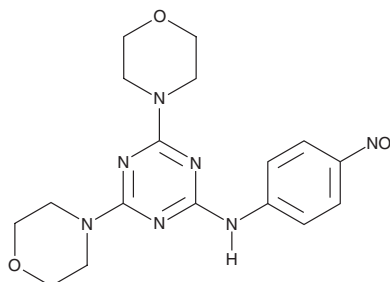
PRODUCT INFORMATION



MHY1485

Item No. 18008

CAS Registry No.: 326914-06-1
Formal Name: 4,6-di-4-morpholinyl-N-(4-nitrophenyl)-
1,3,5-triazin-2-amine
MF: C₁₇H₂₁N₇O₄
FW: 387.4
Purity: ≥95%
UV/Vis.: λ_{max}: 233, 349 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

MHY1485 is supplied as a crystalline solid. A stock solution may be made by dissolving the MHY1485 in the solvent of choice, which should be purged with an inert gas. MHY1485 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of MHY1485 in these solvents is approximately 0.5 and 5 mg/ml, respectively.

Description

MHY1485 is a cell-permeable activator of mTOR that has been shown to increase levels of cellular mTOR Ser²⁴⁴⁸ and downstream substrate 4E-BP Thr^{37/46} phosphorylation in rat liver Ac2F cells.¹ It can induce cellular LC3-II accumulation in Ac2F cells and has been shown to inhibit autophagy by suppressing the fusion between autophagosomes and lysosomes.¹

Reference

1. Choi, Y.J., Park, Y.J., Park, J.I., *et al.* Inhibitory effect of mTOR activator MHY1485 on autophagy: suppression of lysosomal fusion. *PLoS One* **7(8)**, e43418 (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, 10/28/2022

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