

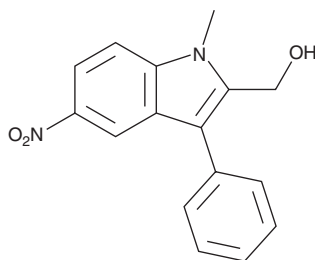
PRODUCT INFORMATION



ML-354

Item No. 19120

CAS Registry No.: 89159-60-4
Formal Name: 1-methyl-5-nitro-3-phenyl-1H-indole-2-methanol
Synonym: VU0099704
MF: C₁₆H₁₄N₂O₃
FW: 282.3
Purity: ≥95%
UV/Vis.: λ_{max}: 263, 280 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

ML-354 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-354 in the solvent of choice. ML-354 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-354 in ethanol is approximately 2 mg/ml and approximately 20 mg/ml in DMSO and DMF.

ML-354 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ML-354 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ML-354 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ML-354 is an indole-based, proteinase-activated receptor 4 (PAR4) antagonist (IC₅₀ = 140 nM) that exhibits 71-fold selectivity for PAR4 over PAR1 (IC₅₀ = 10 μM) in isolated human platelets.¹

Reference

1. Wen, W., Young, S.E., Duvernay, M.T., *et al.* Substituted indoles as selective protease activated receptor 4 (PAR-4) antagonists: Discovery and SAR of ML354. *Bioorg. Med. Chem. Lett.* **24**, 4708-4713 (2014).

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, 12/19/2022

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