

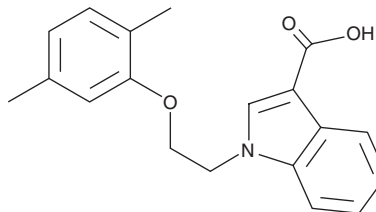
PRODUCT INFORMATION



ML-098

Item No. 15175

CAS Registry No.: 878978-76-8
Formal Name: 1-[2-(2,5-dimethylphenoxy)ethyl]-
1H-indole-3-carboxylic acid
Synonym: CID-7345532
MF: C₁₉H₁₉NO₃
FW: 309.4
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 282 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-098 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-098 in the solvent of choice, which should be purged with an inert gas. ML-098 is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

Description

The Ras family of GTPases, including the Arf, Rho, Ras, and Rab GTPase subfamilies, regulates various cellular processes ranging from membrane trafficking to the control of cell proliferation. They regulate cell physiology by switching between active and inactive conformational states by binding either GTP or GDP. Alterations in GTPase function are a known hallmark of certain cancers and genetic diseases. ML-098 is an activator of the GTP-binding protein Rab7 (EC₅₀ = 77.6 nM) that demonstrates selectivity against the related GTPases cdc42, Ras, Rab-2A, and Rac1 (EC₅₀s = 588.8, 346.7, 158.5, and 794.3 nM, respectively).¹ It has been shown to function by increasing the affinity of the GTPase for guanine nucleotides.¹

Reference

1. Surviladze, Z., Ursu, O., Miscisocia, F., *et al.* Three small molecule pan activator families of Ras-related GTPases. *Probe Reports from the NIH Molecular Libraries Program* (2009).

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, 05/16/2023

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