

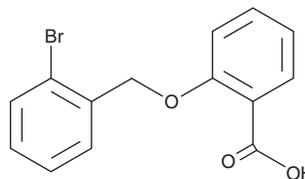
PRODUCT INFORMATION



ML-097

Item No. 15174

CAS Registry No.: 743456-83-9
Formal Name: 2-[(2-bromophenyl)methoxy]-benzoic acid
Synonym: CID-2160985
MF: C₁₄H₁₁BrO₃
FW: 307.1
Purity: ≥95%
UV/Vis.: λ_{max}: 292 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-097 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-097 in the solvent of choice, which should be purged with an inert gas. ML-097 is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

Description

ML-097 is a small molecule pan activator of Ras-related GTPases, activating cell division cycle 42, Ras, and Rac1 with EC₅₀ values of 102, 109, and 151 nM, respectively.¹

References

1. Surviladze, Z., Ursu, O., Miscisocia, F., et al. UNMCMD Probe Report: Three small molecule pan activator families of Ras-related GTPases, in *Probe Reports from the NIH Molecular Libraries Program*, 1 (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, 12/13/2022

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