

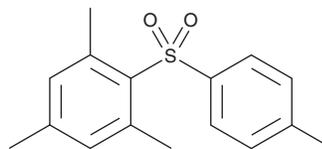
PRODUCT INFORMATION



ESI-05

Item No. 29702

CAS Registry No.: 5184-64-5
Formal Name: 1,3,5-trimethyl-2-[(4-methylphenyl)sulfonyl]-benzene
Synonym: NSC 116966
MF: C₁₆H₁₈O₂S
FW: 274.4
Purity: ≥95%
UV/Vis.: λ_{max}: 243 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

ESI-05 is supplied as a crystalline solid. A stock solution may be made by dissolving the ESI-05 in the solvent of choice, which should be purged with an inert gas. ESI-05 is soluble in DMSO.

Description

ESI-05 is an inhibitor of exchange protein activated by cAMP 2 (Epac2) that inhibits cAMP-induced Epac2 guanine nucleotide exchange factor (GEF) activity with an IC₅₀ value of 0.43 μM in a cell-free assay.¹ It is selective for Epac2 over Epac1 and PKA at 25 μM. ESI-05 (1, 5, 10, and 25 μM) reduces activation of the GTPase RAP1 induced by the Epac-selective cAMP analog 007-AM in HEK293 cells expressing Epac2 but has no effect in HEK293 cells expressing Epac1.

Reference

1. Tsalkova, T., Mei, F.C., Li, S., *et al.* Isoform-specific antagonists of exchange proteins directly activated by cAMP. *Proc. Natl. Acad. Sci. USA* **109**(45), 18613-18618 (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, 12/09/2022

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