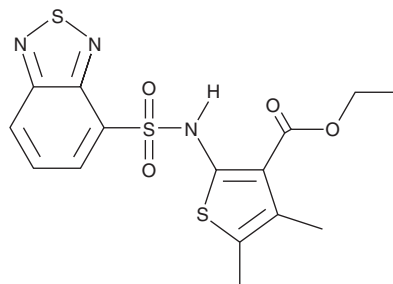


PRODUCT INFORMATION

Calcium Influx Inducer Compound 634

Item No. 41986

CAS Registry No.: 882291-37-4
Formal Name: 2-[(2,1,3-benzothiadiazol-4-ylsulfonyl)amino]-4,5-dimethyl-3-thiophenecarboxylic acid, ethyl ester
MF: C₁₅H₁₅N₃O₄S₃
FW: 397.5
Purity: ≥95%
Supplied as: A 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

Calcium influx inducer compound 634 is supplied as a solid. A stock solution may be made by dissolving the calcium influx inducer compound 634 in the solvent of choice, which should be purged with an inert gas. Calcium influx inducer compound 634 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Calcium influx inducer compound 634 is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Calcium influx inducer compound 634 is an inducer of calcium influx.¹ It enhances the release of extracellular vesicles (EVs) from mouse bone marrow-derived dendritic cells (mBMDCs) when used at a concentration of 10 μM. Calcium influx inducer compound 634 (10 μM) also increases the levels of CD86 and CD80 on the surface of mBMDCs, an effect that can be blocked by the store-operated calcium entry (SOCE) inhibitor BTP2 (YM-58483; Item No. 13246), and increases CD86 and CD80 levels on mBMDC-derived EVs.

Reference

1. Sako, Y., Sato-Kaneko, F., Shukla, N.M., *et al.* Identification of a novel small molecule that enhances the release of extracellular vesicles with immunostimulatory potency via induction of calcium influx. *ACS Chem. Biol.* **18**(4), 982-993 (2023).

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, 09/19/2024

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