

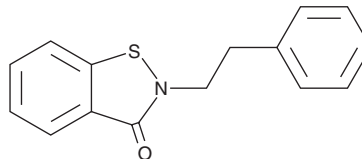
PRODUCT INFORMATION



CAY10762

Item No. 30576

CAS Registry No.: 2514-37-6
Formal Name: 2-(2-phenylethyl)-1,2-benzisothiazol-3(2H)-one
MF: C₁₅H₁₃NOS
FW: 255.3
Purity: ≥95%
UV/Vis.: λ_{max}: 227, 247 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

CAY10762 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10762 in the solvent of choice, which should be purged with an inert gas. CAY10762 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10762 in these solvents is approximately 1, 10, and 15 mg/ml, respectively.

CAY10762 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10762 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10762 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

CAY10762 is an inhibitor of monoacylglycerol lipase (MAGL; IC₅₀ = 34.1 nM).¹ It reduces hydrogen peroxide-induced lactate dehydrogenase (LDH) release from Neuro2a cells when used at a concentration of 1 μM. CAY10762 (10 mg/kg) increases levels of 2-arachidonoyl glycerol (2-AG; Item No. 62160) in mouse brain.

Reference

1. Castelli, R., Scalvini, L., Vacondio, F., *et al.* Benzisothiazolinone derivatives as potent allosteric monoacylglycerol lipase inhibitors that functionally mimic sulfenylation of regulatory cysteines. *J. Med. Chem.* **63**(3), 1261-1280 (2020).

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, 03/21/2024

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