

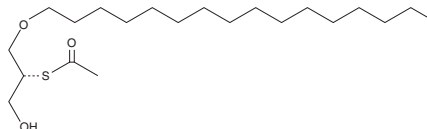
PRODUCT INFORMATION



2-thio-Acetyl MAGE

Item No. 10009651

CAS Registry No.: 112014-15-0
Formal Name: S-[2-(hexadecyloxy)-1-(hydroxymethyl)ethyl]ester, ethanethioic acid
MF: C₂₁H₄₂O₃S
FW: 374.6
Purity: ≥98%
Supplied as: A solution in methyl acetate
Storage: -80°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

2-thioacetyl MAGE is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 2-thioacetyl MAGE in these solvents is approximately 10 mg/ml.

2-thioacetyl MAGE is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, first evaporate the methyl acetate under a gentle stream of nitrogen and then immediately add the aqueous buffer of choice. 2-thioacetyl MAGE has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

KIAA1363 is a recently characterized 2-acetyl monoacylglycerol ether hydrolase that is critical to the survival and proliferation of many cancer cell lines.¹ 2-thioacetyl MAGE is a colorimetric substrate for KIAA1363. It can readily be adapted to perform activity/inhibition studies of this new enzyme using standard DTNB detection of the free thiol as a readout for enzymatic activity.¹

Reference

1. Chiang, K.P., Niessen, S., Saghatelian, A., *et al.* An enzyme that regulates ether lipid signaling pathways in cancer annotated by multidimensional profiling. *Chemistry & Biology* **13**, 1041-1050 (2006).

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, 10/10/2023

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