

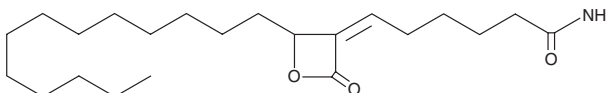
PRODUCT INFORMATION



KC01

Item No. 17364

CAS Registry No.: 1646795-59-6
Formal Name: 6Z-(2-oxo-4-tridecyl-3-oxetanylidene), hexanamide
MF: $C_{22}H_{39}NO_3$
FW: 365.6
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 213 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

KC01 is supplied as a crystalline solid. A stock solution may be made by dissolving the KC01 in the solvent of choice. KC01 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of KC01 in ethanol is approximately 16 mg/ml, and approximately 5 mg/ml in DMSO and DMF.

KC01 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, KC01 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. KC01 has a solubility of approximately 0.5 mg/ml in a 1:5 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ABHD16A (also known as HLA-B associated transcript 5 (BAT5)) is a phosphatidylserine hydrolase that regulates the formation of lysophosphatidylserines (lyso-PSs) and has been implicated in certain immunological and neurological diseases.¹ KC01 is a covalent inhibitor of ABHD16A (IC_{50} s = 90 and 520 nM for human and mouse, respectively).¹ It has been shown to deplete lyso-PS from various human cancer cell lines, a lymphoblast cell line, and mouse brain membrane lysates and to decrease LPS-induced cytokine production in macrophages.¹ This compound has been used to characterize the role of ABHD16A as a principle phosphatidylserine lipase *in vivo*.¹

Reference

1. Kamat, S.S., Camara, K., Parsons, W.H., *et al.* Immunomodulatory lysophosphatidylserines are regulated by ABHD16A and ABHD12 interplay. *Nat. Chem. Biol.* **11**, 164-175 (2015).

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, 12/02/2022

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