PRODUCT INFORMATION



KC02

Item No. 17365

CAS Registry No.:	1646795-60-9	
Formal Name:	6Z-[2-oxo-4-(2-phenylethyl)-3-	
	oxetanylidene], hexanamide	
MF:	C ₁₇ H ₂₁ NO ₃	
FW:	287.4	
Purity:	≥98%	
Supplied as:	A solution in ethyl acetate	0 \0
Storage:	-20°C	
Stability:	≥2 years	

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

Laboratory Procedures

KC02 is supplied as a solution in ethyl acetate. To change the solvent, simply evaporate the KC02 under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of KC02 in ethanol is 16 mg/ml, and approximately 5 mg/ml in DMSO and DMF.

KC02 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethyl acetate solution of KC02 should be diluted with the aqueous buffer of choice. KC02 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

KC02 is an inactive control probe for KC01 (Item No. 17364), the potent inhibitor of ABHD16A, a phosphatidylserine hydrolase that regulates the formation of lysophosphatidylserines in mammalian systems. KC02 demonstrates IC₅₀ values greater 10 μ M against human and mouse ABHD16A.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

uyer 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, 02/15/2024

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