

# PRODUCT INFORMATION



**Ki16425**

Item No. 10012659

**CAS Registry No.:** 355025-24-0  
**Formal Name:** 3-[[[4-[4-[[[1-(2-chlorophenyl)ethoxy]carbonyl]amino]-3-methyl-5-isoxazolyl]phenyl]methyl]thio]-propanoic acid

**MF:** C<sub>23</sub>H<sub>23</sub>ClN<sub>2</sub>O<sub>5</sub>S

**FW:** 475.0

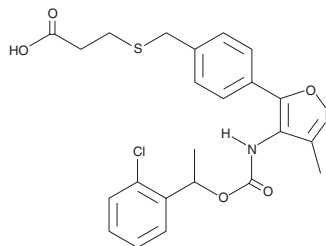
**Purity:** ≥95%

**UV/Vis.:** λ<sub>max</sub>: 276 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

Ki16425 is supplied as a crystalline solid. A stock solution may be made by dissolving the Ki16425 in an organic solvent purged with an inert gas. Ki16425 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Ki16425 in these solvents is approximately 10 mg/ml.

Ki16425 is sparingly soluble in aqueous buffers. Ki16425 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Ki16425 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

Lysophosphatidic acid (LPA) is a bioactive lipid mediator that signals through five distinct G protein-coupled receptors (LPA<sub>1-5</sub>).<sup>1</sup> Ki16425 is a LPA receptor antagonist with selectivity for LPA<sub>1</sub> and LPA<sub>3</sub>. It exhibits K<sub>i</sub> values of 0.34, 6.5, and 0.93 μM for the human LPA<sub>1</sub>, LPA<sub>2</sub>, and LPA<sub>3</sub> receptors, respectively, as determined by measuring inositol phosphate production in RH7777-transfected cells.<sup>2</sup> Ki1642, at 10 μM, significantly blocks the response of a variety of cancer cell lines to LPA-induced cell migration.<sup>3</sup>

## References

1. Choi, J.W., Lee, C.-W., and Chun, J. Biological roles of lysophospholipid receptors revealed by genetic null mice: An update. *Biochim. Biophys. Acta* **1781**(9), 531-539 (2008)
2. Ohta, H., Sato, K., Murata, N., et al. Ki16425, a subtype-selective antagonist for EDG-family lysophosphatidic acid receptors. *Mol. Pharmacol.* **64**(4), 994-1005 (2003).
3. Yamada, T., Sato, K., Komachi, M., et al. Lysophosphatidic acid (LPA) in malignant ascites stimulates motility of human pancreatic cancer cells through LPA. *J. Biol. Chem.* **279**(8), 6595-6605 (2004).

### 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, 12/13/2022

## 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