

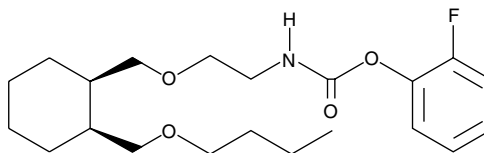
# Product Information



## (-)-AS 115

Item No. 13250

**CAS Registry No.:** 926657-43-4  
**Formal Name:** 2-fluorophenyl-(2-(((1R,2S)-2-(butoxymethyl)cyclohexyl)methoxy)ethyl)carbamate  
**MF:** C<sub>21</sub>H<sub>32</sub>FNO<sub>4</sub>  
**FW:** 381.5  
**Purity:** ≥98%  
**Stability:** ≥1 year at -20°C  
**Supplied as:** A solution in methyl acetate  
**UV/Vis.:** λ<sub>max</sub>: 205, 261, 267 nm



### Laboratory Procedures

For long term storage, we suggest that (-)-AS 115 be stored as supplied at -20°C. It should be stable for at least one year.

(-)-AS 115 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 (DMF) purged with an inert gas can be used. The solubility of (-)-AS 115 in these solvents is approximately 30 mg/ml in ethanol and approximately 12 mg/ml in DMSO and DMF.

(-)-AS 115 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of (-)-AS 115 should be diluted with the aqueous buffer of choice. (-)-AS 115 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

KIAA1363 is a 2-acetyl monoacylglycerol ether (MAGE) hydrolase that is upregulated in aggressive cancers from various tissues.<sup>1</sup> The enzyme catalyzes the hydrolysis of the 2-acetyl MAGE to MAGE and serves as a central enzyme in the PAF and LPA signaling network.<sup>2</sup> AS 115 is a potent and selective inactivator of KIAA1363, displaying an IC<sub>50</sub> value of 150 nM when tested as a racemic mixture in the invasive ovarian cancer cell line SKOV3.<sup>1</sup> Treatment of SKOV3 cells with 10 μM AS-115 for 4 hours significantly reduced the formation of MAGE, alkyl-lysophosphatidylcholine, and alkyl-lysophosphatidic acid. The activity of the individual enantiomers of AS 115, *i.e.*, (+)-AS 115 and (-)-AS 115, has not been determined.

### References

1. Jessani, N., Liu, Y., Humphrey, M., *et al.* Enzyme activity profiles of the secreted and membrane proteome that depict cancer cell invasiveness. *Proc. Natl. Acad. Sci. USA* **99**(16), 10335-10340 (2002).
2. 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).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/13250](http://www.caymanchem.com/catalog/13250)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 03/13/2015

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)