

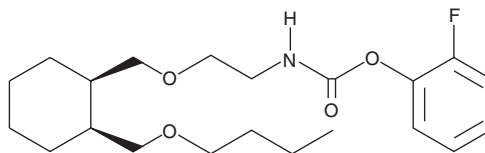
# PRODUCT INFORMATION



## (-)-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%  
**UV/Vis.:** λ<sub>max</sub>: 261,267 nm  
**Supplied as:** A solution in methyl acetate  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

(-)-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 ethanol is approximately 30 mg/ml 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 0.25 mg/ml/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.

### Description

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

- 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).
- 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**(10), 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, 10/13/2022

#### 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