

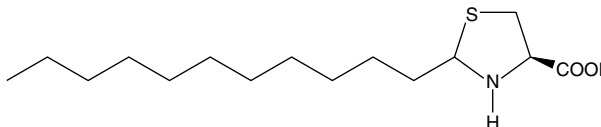
# Product Information



## CAY10444

Catalog No. 10005033

**CAS Registry No.:** 298186-80-8  
**Formal Name:** 2-undecyl-thiazolidine-4-carboxylic acid  
**Synonym:** BML-241  
**MF:** C<sub>15</sub>H<sub>29</sub>NO<sub>2</sub>S  
**FW:** 287.5  
**Purity:** ≥98%  
**Stability:** ≥1 year at -20°C  
**Supplied as:** A crystalline solid



### Laboratory Procedures

For long term storage, we suggest that CAY10444 be stored as supplied at -20°C. It should be stable for at least one year.

CAY10444 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10444 in an organic solvent purged with an inert gas. CAY10444 is soluble in dimethyl formamide (DMF). The solubility of CAY10444 in DMF is approximately 0.5 mg/ml.

CAY10444 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10444 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10444 has a solubility of approximately 0.15 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

A family of related G-protein coupled receptors that bind sphingosine-1 phosphate (S1P) as a high-affinity ligand have recently been cloned.<sup>1,2</sup> S1P<sub>3</sub>/EDG-3 is a member of this family that is widely expressed in many tissues, signaling *via* the ERK-1/2 and PLC/IP<sub>3</sub> pathways. CAY10444 is a selective antagonist of S1P binding to the S1P<sub>3</sub> receptor, blocking the calcium increase in HeLa cells by about 40% when present at 10 μM.<sup>3</sup>

### References

1. Kluk, M.J. and Hla, T. Signaling of sphingosine-1-phosphate *via* the S1P/EDG-family of G-protein-coupled receptors. *Biochim. Biophys. Acta* **1582**, 72-80 (2002).
2. Chun, J., Goetzl, E.J., Hla, T., *et al.* International union of pharmacology. XXXIV. Lysophospholipid receptor nomenclature. *Pharmacol. Rev.* **54**, 265-269 (2002).
3. Koide, Y., Hasegawa, T., Takahashi, A., *et al.* Development of novel EDG3 antagonists using a 3D database search and their structure-activity relationships. *J. Med. Chem.* **45**, 4629-4638 (2002).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/10005033](http://www.caymanchem.com/catalog/10005033)

## 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)

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

#### MATERIAL 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 Material 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, 12/05/2011