

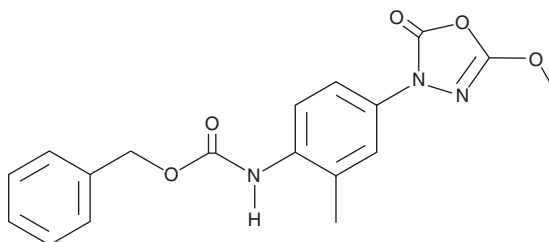
# Product Information



**CAY10499**

Catalog No. 10007875

**CAS Registry No.:** 359714-55-9  
**Formal Name:** [4-(5-methoxy-2-oxo-1,3,4-oxadiazol-3(2H)-yl)-2-methylphenyl]-carbamic acid, phenylmethyl ester  
**MF:** C<sub>18</sub>H<sub>17</sub>N<sub>3</sub>O<sub>5</sub>  
**FW:** 355.3  
**Purity:** ≥98%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid



## Laboratory Procedures

For long term storage, we suggest that CAY10499 be stored as supplied at -20°C. It should be stable for at least two years.

CAY10499 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10499 in an organic solvent purged with an inert gas. CAY10499 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10499 in ethanol is approximately 2 mg/ml and it is approximately 20 mg/ml in DMSO and DMF.

CAY10499 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10499 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. CAY10499 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Hormone sensitive lipase (HSL) catalyzes the hydrolysis of tri-, di-, and monoacylglycerols, as well as cholesterol esters, thus mobilizing fatty acids as a primary source of energy in mammals.<sup>1</sup> The enzyme is highly expressed in adipose tissue and steroidogenic tissues, and less abundantly in skeletal muscle, heart, brain, pancreatic beta cells, adrenal gland, ovaries, testes, and macrophages. Its presence in various tissues indicates the enzyme plays diverse roles including those in steroidogenesis and spermatogenesis, foam cell formation in atherosclerosis, and diabetes.<sup>1,2</sup> CAY10499 is a potent inhibitor of human HSL exhibiting an IC<sub>50</sub> of 90 nM for the recombinant enzyme.<sup>3</sup> The *in vivo* pharmacological efficacy of CAY10499 has not been reported.

## References

1. Saltiel, A.R. Another hormone-sensitive triglyceride lipase in fat cells. *Proc. Natl. Acad. Sci. USA* **97**(2), 535-537 (2000).
2. Kraemer, F.B. and Shen, W.-J. Hormone-sensitive lipase: Control of intracellular tri-(di-)acylglycerol and cholesteryl ester hydrolysis. *J. Lipid Res.* **43**, 1585-1594 (2002).
3. Ali, Y.B., Müller, G., Petry, S., *et al.* Kinetic characterization and specific inhibition of hormone sensitive lipase. *Chem. Phys. Lipids* **136**, (2005).

## Related Products

Endothelial Lipase Polyclonal Antibody - Cat. No. 100030 • Monoacylglycerol Lipase Polyclonal Antibody - Cat. No. 100035 • Hormone Sensitive Lipase Polyclonal Antibody - Cat. No. 10006371

**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, 05/06/2010

## 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

### Web

www.caymanchem.com