

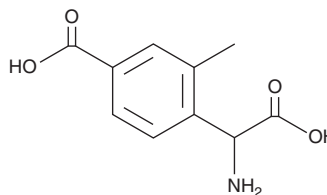
# PRODUCT INFORMATION



(±)-LY367385

Item No. 33158

CAS Registry No.: 198419-90-8  
Formal Name: α-amino-4-carboxy-2-methyl-  
benzeneacetic acid  
MF: C<sub>10</sub>H<sub>11</sub>NO<sub>4</sub>  
FW: 209.2  
Purity: ≥98%  
Supplied as: A 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

(±)-LY367385 is supplied as a solid. A stock solution may be made by dissolving the (±)-LY367385 in water. The solubility of (±)-LY367385 in water is approximately 20 mg/ml. We do not recommend storing the aqueous solution for more than one day.

## Description

(±)-LY367385 is an antagonist of metabotropic glutamate receptor 1α (mGluR1α).<sup>1</sup> It inhibits phosphoinositide (PI) hydrolysis selectively in AV-12 cells expressing mGluR1α over those expressing mGluR5a (IC<sub>50</sub>s = 19 and >300 μM, respectively).

## Reference

1. Salt, T.E., Jones, H.E., Copeland, C.S., *et al.* Function of mGlu1 receptors in the modulation of nociceptive processing in the thalamus. *Neuropharmacology* **79(100)**, 405-411 (2014).

### 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, 11/30/2022

## 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