

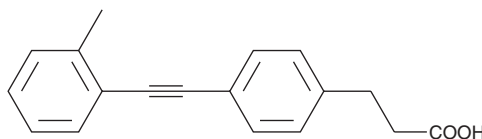
# PRODUCT INFORMATION



**CAY10587**

Item No. 13143

**CAS Registry No.:** 1082058-99-8  
**Formal Name:** 4-[2-(2-methylphenyl)ethynyl]-benzenepropanoic acid  
**MF:** C<sub>18</sub>H<sub>16</sub>O<sub>2</sub>  
**FW:** 264.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 286, 295, 305 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

## Laboratory Procedures

CAY10587 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10587 in the solvent of choice, which should be purged with an inert gas. CAY10587 is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

## Description

CAY10587 is a selective FFAR1 (GPR40) agonist (EC<sub>50</sub> = 32 nM) that does not exhibit activity on the related FFARs: FFAR2 (GPR43) or FFAR3 (GPR41).<sup>1</sup> At 100 nM, CAY10587 increases glucose-stimulated insulin secretion in the rat insulin-secreting cell line INS-1E.<sup>1</sup>

## Reference

- Christiansen, E., Urban, C., Merten, N., *et al.* Discovery of potent and selective agonists for the free fatty acid receptor 1 (FFA1/GPR40), a potential target for the treatment of type II diabetes. *J. Med. Chem.* **51**(22), 7061-7064 (2008).

### 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, 03/23/2022

## 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](http://WWW.CAYMANCHEM.COM)