

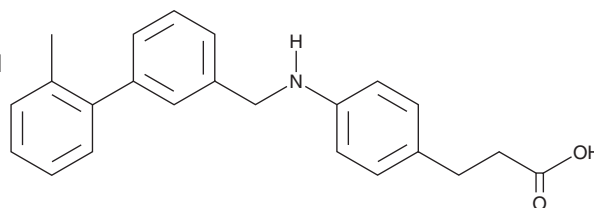
# PRODUCT INFORMATION



**TUG-469**

Item No. 44272

**CAS Registry No.:** 1236109-67-3  
**Formal Name:** 4-[[[2'-methyl[1,1'-biphenyl]-3-yl)methyl]amino]-benzenepropanoic acid  
**MF:** C<sub>23</sub>H<sub>23</sub>NO<sub>2</sub>  
**FW:** 345.4  
**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

TUG-469 is supplied as a solid. A stock solution may be made by dissolving the TUG-469 in the solvent of choice, which should be purged with an inert gas. TUG-469 is slightly soluble (0.1-1 mg/ml) in ethanol and slightly soluble (0.1-1 mg/ml) in DMSO.

## Description

TUG-469 is an agonist of free fatty acid receptor 1 (FFAR1/GPR40).<sup>1</sup> It induces calcium mobilization in a reporter assay using cells expressing human FFAR1/GPR40 (EC<sub>50</sub> = 18.6 nM). TUG-469 (10 μM) induces insulin secretion in INS-1E insulinoma cells. It inhibits increases in blood glucose levels in a glucose tolerance test (GTT) in mice when administered at a dose of 5 mg/kg.<sup>2</sup>

## References

- Christiansen, E., Due-Hansen, M.E., Urban, C., *et al.* Structure-activity study of dihydrocinnamic acids and discovery of the potent FFA1 (GPR40) agonist TUG-469. *ACS Med. Chem. Lett.* **1(7)**, 345-349 (2010).
- Urban, C., Hamacher, A., Partke, H.J., *et al.* In vitro and mouse in vivo characterization of the potent free fatty acid 1 receptor agonist TUG-469. *Naunyn Schmiedebergs Arch Pharmacol.* **386(12)**, 1021-1030 (2013).

### 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, 09/18/2025

## 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