

# PRODUCT INFORMATION

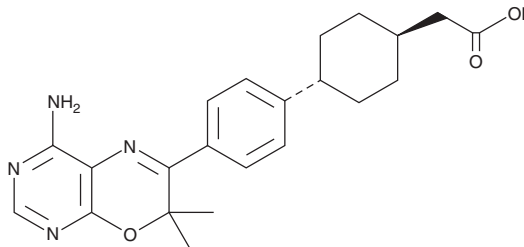


**T-863**

Item No. 25807

**CAS Registry No.:** 701232-20-4  
**Formal Name:** *trans*-4-[4-(4-amino-7,7-dimethyl-7H-pyrimido[4,5-b][1,4]oxazin-6-yl)phenyl]-cyclohexaneacetic acid

**MF:** C<sub>22</sub>H<sub>26</sub>N<sub>4</sub>O<sub>3</sub>  
**FW:** 394.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 238, 335 nm  
**Supplied as:** A crystalline 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

T-863 is supplied as a crystalline solid. A stock solution may be made by dissolving the T-863 in the solvent of choice, which should be purged with an inert gas. T-863 is slightly soluble in DMSO and dimethyl formamide.

## Description

T-863 is a potent inhibitor of diacylglycerol acetyltransferase-1 (DGAT-1; IC<sub>50</sub> = 15 nM).<sup>1</sup> It is selective for DGAT-1 over DGAT-2 and acyl-coenzyme A:cholesterol acyltransferase-1 (ACAT-1; IC<sub>50</sub>s = >10 and 12 μM, respectively). T-863 inhibits triacylglycerol synthesis in HuTu 80 cells (IC<sub>50</sub> = 4 nM). *In vivo*, T-863 (20 mg/kg) reduces plasma triacylglycerol levels, the ratio of triacylglycerols to diacylglycerols, and body weight in a mouse model of diet-induced obesity.

## Reference

1. Birch, A.M., Birtles, S., Buckett, L.K., *et al.* Discovery of a potent, selective, and orally efficacious pyrimidinooxazinyl bicyclooctaneacetic acid diacylglycerol acyltransferase-1 inhibitor. *J. Med. Chem.* **52**(6), 1558-1568 (2009).

### 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, 11/04/2022

## 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