

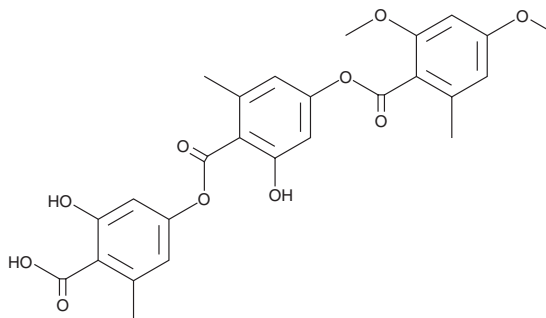
# PRODUCT INFORMATION



## Amidepsine D

Item No. 17775

**CAS Registry No.:** 79786-34-8  
**Formal Name:** 4-[(2,4-dimethoxy-6-methylbenzoyl)oxy]-2-hydroxy-6-methyl-benzoic acid, 4-carboxy-3-hydroxy-5-methylphenyl ester  
**MF:** C<sub>26</sub>H<sub>24</sub>O<sub>10</sub>  
**FW:** 496.5  
**Purity:** ≥95%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Fungi/*Humicola* sp. FO-2942



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

### Laboratory Procedures

Amidepsine D is supplied as a crystalline solid. A stock solution may be made by dissolving the amidepsine D in the solvent of choice, which should be purged with an inert gas. Amidepsine D is soluble in the organic solvent DMSO.

### Description

Diacylglycerol acyltransferases (DGATs) catalyze the final step in triglyceride synthesis. DGAT-1 deficient mice demonstrate a phenotype that is protective against the development of diet-induced obesity or insulin resistance, implicating the actions of this enzyme in the development of such metabolic disorders. Amidepsine D is a fungal metabolite isolated from the culture broth of *Humicola* sp. FO-2942 that inhibits DGAT activity (IC<sub>50</sub> = 17.5 μM in rat liver microsomes).<sup>1</sup> It inhibits triacylglycerol formation in Raji cells with an IC<sub>50</sub> value of 2.8 μM.<sup>1</sup>

### Reference

1. Tomoda, H., Ito, M., Tabata, N., *et al.* Amidepsines, inhibitors of diacylglycerol acyltransferase produced by *Humicola* sp. FO-2942. I. Production, isolation and biological properties. *J. Antibiot. (Tokyo)* **48(9)**, 937-941 (1995).

#### 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, 11/10/2022

#### 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