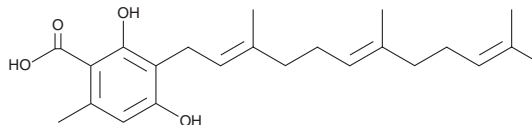


PRODUCT INFORMATION



Grifolic Acid Item No. 17398

CAS Registry No.: 80557-12-6
Formal Name: 2,4-dihydroxy-6-methyl-3-
[(2E,6E)-3,7,11-trimethyl-2,6,10-
dodecatrien-1-yl]-benzoic acid
MF: C₂₃H₃₂O₄
FW: 372.5
Purity: ≥95%
UV/Vis.: λ_{max}: 221, 268, 303 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

Grifolic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the grifolic acid in the solvent of choice. Grifolic acid is soluble in organic solvents such as DMSO, which should be purged with an inert gas. The solubility of grifolic acid in this solvent is approximately 100 mg/ml.

Description

GPR120 (free fatty acid receptor 4/FFAR4) is a G protein-coupled receptor expressed in intestine, adipocytes, and pro-inflammatory macrophages that is activated by long-chain free fatty acids. It has been linked to the ability of fatty acids to produce anti-inflammatory effects and to acutely potentiate insulin secretion. Grifolic acid is a selective partial agonist of GPR120 that induces ERK activation and intracellular calcium responses in mouse enteroendocrine STC-1 cells that express GPR120 endogenously, but not those expressing GPR40.¹

Reference

1. Sun, Q., Hirasawa, A., Hara, T., *et al.* Structure-activity relationships of GPR120 agonists based on a docking simulation. *Mol. Pharm.* **78**, 804-810 (2010).

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, 11/17/2022

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