

PRODUCT INFORMATION



GSA 10

Item No. 16164

CAS Registry No.: 300833-95-8
Formal Name: 4-[[[(1-hexyl-1,2-dihydro-4-hydroxy-2-oxo-3-quinolinyl) carbonyl]amino]-benzoic acid, propyl ester

MF: C₂₆H₃₀N₂O₅

FW: 450.5

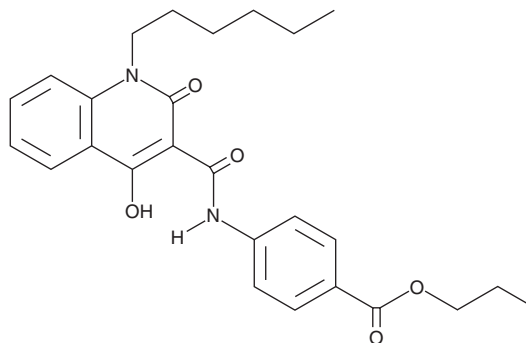
Purity: ≥95%

UV/Vis.: λ_{max}: 224, 237, 259, 314 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

GSA 10 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSA 10 in the solvent of choice, which should be purged with an inert gas. GSA 10 is soluble in the organic solvent dimethyl formamide at a concentration of approximately 1 mg/ml.

Description

GSA 10 is a quinolinecarboxamide derivative that acts as an agonist of SMO.¹ By acting at SMO, GSA 10 has been shown to promote the differentiation of multipotent mesenchymal progenitor cells into osteoblasts (EC₅₀ = 1.2 μM).¹ Most notably, GSA 10 does not recognize the classic cycloamine binding site as do other SMO agonists and as such has been used to characterize a novel SMO active site.¹

Reference

1. Gorojankina, T., Hoch, L., Faure, H., *et al.* Discovery, molecular and pharmacological characterization of GSA-10, a novel small-molecule positive modulator of Smoothed. *Mol. Pharmacol.* **83**(5), 1020-1029 (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, 06/29/2020

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