

PRODUCT INFORMATION



MM-11253

Item No. 21800

CAS Registry No.: 345952-44-5
Formal Name: 6-[2-(5,6,7,8-tetrahydro-5,5,8,8-tetra methyl-2-naphthalenyl)-1,3-dithiolan-2-yl]-2-naphthalenecarboxylic acid

MF: C₂₈H₃₀O₂S₂

FW: 462.7

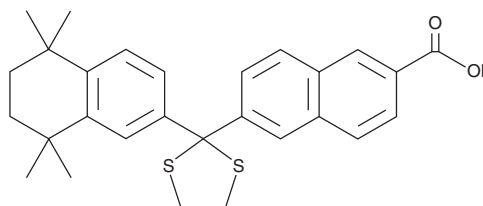
Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Stability: Store in desiccating conditions



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

Laboratory Procedures

MM-11253 is supplied as a solid. A stock solution may be made by dissolving the MM-11253 in the solvent of choice. MM-11253 is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 100 mM.

Description

MM-11253 is a synthetic, conformationally restricted retinoid that acts as a selective antagonist of retinoic acid receptor γ (RAR γ).¹ It dose-dependently blocks the ability of RAR γ -selective agonists, but not retinoic acid (all-*trans* retinoic acid; Item No. 11017), to inhibit the growth of oral squamous cell carcinoma cells *in vitro*.¹

Reference

1. Le, Q.T., Dawson, M.I., Soprano, D.R., *et al.* Modulation of retinoic acid receptor function alters the growth inhibitory response of oral SCC cells to retinoids. *Oncogene* **19(11)**, 1457-1465 (2000).

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, 12/22/2022

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