

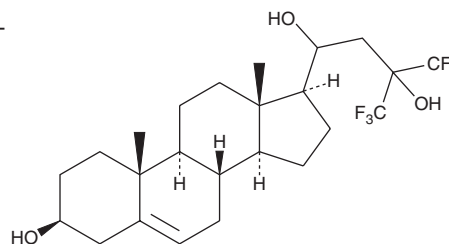
PRODUCT INFORMATION



NSC 12

Item No. 20117

CAS Registry No.: 102586-30-1
Formal Name: (3 β)-24,24,24-trifluoro-23-(trifluoromethyl)-21-norchol-5-ene-3,20,23-triol
Synonym: NSC 172285
MF: C₂₄H₃₄F₆O₃
FW: 484.5
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

NSC 12 is supplied as a crystalline solid. A stock solution may be made by dissolving the NSC 12 in the solvent of choice, which should be purged with an inert gas. NSC 12 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of NSC 12 in these solvents is approximately 20, 0.1, and 2 mg/ml, respectively.

NSC 12 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NSC 12 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. NSC 12 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

NSC 12 is an extracellular trap for fibroblast growth factor 2 (FGF2) that binds FGF2 ($K_d = 51 \mu\text{M}$) and interferes with its interaction with FGFR1, without affecting the ability of FGF2 to bind heparin or heparin sulfate proteoglycans (HSPGs).¹ NSC 12 also binds several other, but not all, FGF isoforms with K_d values ranging between 16 and 120 μM , preventing them from forming HSPG/FGF/FGFR ternary complexes.¹ NSC 12 inhibits FGF-dependent angiogenesis and tumor cell proliferation *in vitro* and reduces tumor growth in mice.¹

Reference

1. Ronca, R., Giacomini, A., Di Salle, E., *et al.* Long-pentraxin 3 derivative as a small-molecule FGF trap for cancer therapy. *Cancer Cell* **28**(2), 225-238 (2015).

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, 10/17/2022

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