

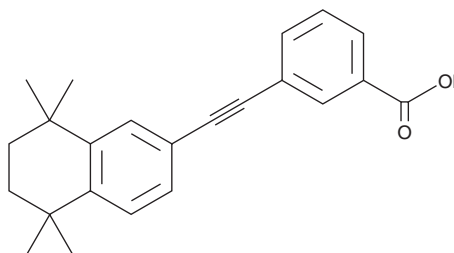
PRODUCT INFORMATION



EC 19

Item No. 44117

CAS Registry No.: 1010694-08-2
Formal Name: 3-[2-(5,6,7,8-tetrahydro-5,5,8,8-tetramethyl-2-naphthalenyl)ethynyl]-benzoic acid
MF: $C_{23}H_{24}O_2$
FW: 332.4
Purity: $\geq 98\%$
Supplied as: A 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

EC 19 is supplied as a solid. A stock solution may be made by dissolving the EC 19 in the solvent of choice, which should be purged with an inert gas. EC 19 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

EC 19 is a synthetic retinoid.¹ It induces epithelial differentiation in TERA2.cl.SP12 pluripotent stem cells when used at a concentration of 10 μM . EC 19 also decreases the proliferation of HepG2 hepatocellular carcinoma, Caco-2 colon carcinoma, and MCF-7 breast cancer cells (IC_{50}s = 42.2, 10.8, and 9.7 μM , respectively).² It inhibits migration and invasion and induces cell cycle arrest at the G_2/M and G_0/G_1 phases in Caco-2 cells when used at a concentration of 42.2 μM . EC 19 reduces total antioxidant levels in Caco-2 cells.

References

1. Christie, V.B., Barnard, J.H., Batsanov, A.S., *et al.* Synthesis and evaluation of synthetic retinoid derivatives as inducers of stem cell differentiation. *Org. Biomol. Chem.* **6**(19), 3497-3507 (2008).
2. Abdelaal, M.R., Soror, S.H., Elnagar, M.R., *et al.* Revealing the potential application of EC-synthetic retinoid analogues in anticancer therapy. *Molecules* **26**(2), 506 (2021).

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, 09/16/2025

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