

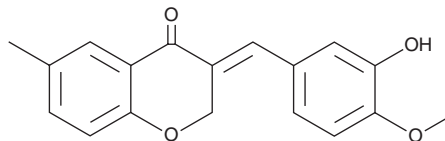
PRODUCT INFORMATION



IMS 2186

Item No. 31157

CAS Registry No.: 1031206-36-6
Formal Name: (3E)-2,3-dihydro-3-[(3-hydroxy-4-methoxyphenyl)methylene]-6-methyl-4H-1-benzopyran-4-one
MF: $C_{18}H_{16}O_4$
FW: 296.3
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

IMS 2186 is supplied as a solid. A stock solution may be made by dissolving the IMS 2186 in the solvent of choice, which should be purged with an inert gas. IMS 2186 is soluble in organic solvents such as ethanol and DMSO. The solubility of IMS 2186 in these solvents is approximately 1 and 5 mg/ml, respectively.

Description

IMS 2186 is an anti-choroidal neovascularization agent.¹ It inhibits proliferation of NIH3T3 mouse and IMR-90 human fibroblasts ($IC_{50}s = 0.5\text{--}3\text{ }\mu\text{M}$), as well as in a panel of eight human cancer cell lines, including ovarian, lung, colon, and pancreatic cancer cells ($IC_{50}s = 0.3\text{--}3\text{ }\mu\text{M}$). IMS 2186 ($10\text{ }\mu\text{M}$) decreases migration of THP-1 macrophages in wound healing and chamber assays. Subcutaneous administration of IMS 2186 ($20\text{ }\mu\text{g/animal}$) inhibits VEGF-induced blood vessel formation in a MatrigelTM plug assay in mice. It reduces ocular leakage, as well as cellular proliferation and infiltration in a rat model of laser-induced choroidal neovascularization when administered intravitreally at a dose of $100\text{ }\mu\text{g/eye}$.

Reference

1. Falkenstein, I.A., Cheng, L., Wong-Staal, F., *et al.* Toxicity and intraocular properties of a novel long-acting anti-proliferative and anti-angiogenic compound IMS2186. *Curr. Eye Res.* **33**(7), 599-609 (2008).

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, 11/04/2022

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