

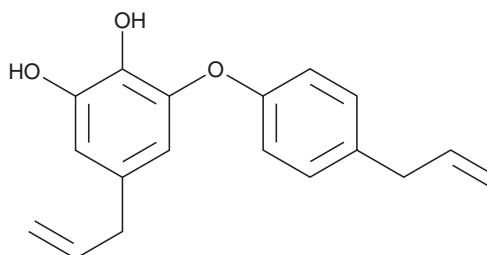
# PRODUCT INFORMATION



## Obovatol

Item No. 27553

**CAS Registry No.:** 83864-78-2  
**Formal Name:** 5-(2-propen-1-yl)-3-[4-(2-propen-1-yl)phenoxy]-1,2-benzenediol  
**Synonym:** NSC 364150  
**MF:** C<sub>18</sub>H<sub>18</sub>O<sub>3</sub>  
**FW:** 282.3  
**Purity:** ≥98%  
**Supplied as:** A neat oil  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Synthetic



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

### Laboratory Procedures

Obovatol is supplied as a neat oil. A stock solution may be made by dissolving the obovatol in the solvent of choice, which should be purged with an inert gas. Obovatol is soluble in DMSO.

### Description

Obovatol is a lignan polyphenol that has been found in *M. officinalis* and has diverse biological activities.<sup>1-3</sup> It inhibits epinephrine- or arachidonate-induced aggregation of isolated rat platelets (IC<sub>50</sub>s = 59 and 54 μM, respectively).<sup>3</sup> Obovatol (50 μM) induces cell cycle arrest at the S phase and apoptosis in SW620 colorectal cancer cells.<sup>2</sup> It reduces tumor volume and weight in a SW620 mouse xenograft model when administered at a dose of 6 mg/kg per day. Obovatol (1 mg/kg per day for three months) decreases hippocampal and cortical amyloid-β (Aβ) 1-42 (Aβ42) fibril formation and reduces latency and distance to escape in the Morris water maze in a Tg2576 mouse model of Alzheimer's disease.<sup>1</sup>

### References

1. Choi, D.-Y., Lee, J.W., Peng, J., *et al.* Obovatol improves cognitive functions in animal models for Alzheimer's disease. *J. Neurochem.* **120(6)**, 1048-1059 (2012).
2. Lee, S.-K., Kim, H.-N., Kang, Y.-R., *et al.* Obovatol inhibits colorectal cancer growth by inhibiting tumor cell proliferation and inducing apoptosis. *Bioorg. Med. Chem.* **16(18)**, 8397-8402 (2008).
3. Pyo, M.K., Lee, Y., and Yun-Choi, H.S. Anti-platelet effect of the constituents isolated from the barks and fruits of *Magnolia obovata*. *Arch. Pharm. Res.* **25(3)**, 325-328 (2002).

#### 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, 01/29/2024

#### 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