

# PRODUCT INFORMATION

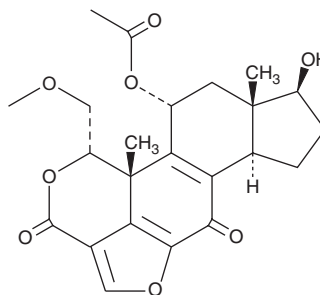


## 17 $\beta$ -hydroxy Wortmannin

Item No. 13812

**CAS Registry No.:** 58053-83-1  
**Formal Name:** (1S,6bR,9S,9aS,11R,11bR)-11-(acetyloxy)-6b,7,8,9,9a,10,11,11b-octahydro-9-hydroxy-1-(methoxymethyl)-9a,11b-dimethyl-3H-furo[4,3,2-de]indeno[4,5-h]-2-benzopyran-3,6(1H)-dione

**MF:** C<sub>23</sub>H<sub>26</sub>O<sub>8</sub>  
**FW:** 430.4  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 259, 292 nm  
**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

17 $\beta$ -hydroxy Wortmannin is supplied as a crystalline solid. A stock solution may be made by dissolving the 17 $\beta$ -hydroxy wortmannin in the solvent of choice, which should be purged with an inert gas. 17 $\beta$ -hydroxy Wortmannin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 17 $\beta$ -hydroxy wortmannin in these solvents is approximately 0.15, 2.5, and 3 mg/ml, respectively.

17 $\beta$ -hydroxy Wortmannin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 17 $\beta$ -hydroxy wortmannin should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 17 $\beta$ -hydroxy Wortmannin has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

17 $\beta$ -hydroxy Wortmannin is an analog of wortmannin. It irreversibly binds phosphoinositide 3-kinase (PI3K) and potently blocks fMLP-stimulated respiratory burst in neutrophils (IC<sub>50</sub> = 5 nM).<sup>1</sup> 17 $\beta$ -hydroxy Wortmannin inhibits recombinant PI3K and mTOR (IC<sub>50</sub> = 2.7 and 193 nM, respectively) and prevents the growth of LNCap prostate cancer cells (IC<sub>50</sub> = 1.46  $\mu$ M).<sup>2</sup> The 17-hydroxyl group has been used to further modify this compound, e.g., by pegylation and conjugation with rapamycin.<sup>3,4</sup>

### References

1. Thelen, M., Wymann, M.P., and Langen, H. Wortmannin binds specifically to 1-phosphatidylinositol 3-kinase while inhibiting guanine nucleotide-binding protein-coupled receptor signaling in neutrophil leukocytes. *Proc. Natl. Acad. Sci. USA* **91**(11), 4960-4964 (1994).
2. Zask, A., Kaplan, J., Toral-Barza, L., et al. Synthesis and structure-activity relationships of ring-opened 17-hydroxywortmannins: Potent phosphoinositide 3-kinase inhibitors with improved properties and anticancer efficacy. *J. Med. Chem.* **51**(5), 1319-1323 (2008).
3. Zhu, T., Gu, J., Yu, K., et al. Pegylated wortmannin and 17-hydroxywortmannin conjugates as phosphoinositide 3-kinase inhibitors active in human tumor xenograft models. *J. Med. Chem.* **49**(4), 1373-1378 (2006).
4. Ayral-Kaloustian, S., Gu, J., Lucas, J., et al. Hybrid inhibitors of phosphatidylinositol 3-kinase (PI3K) and the mammalian target of rapamycin (mTOR): Design, synthesis, and superior antitumor activity of novel wortmannin-rapamycin conjugates. *J. Med. Chem.* **53**(1), 452-459 (2010).

#### 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, 03/28/2024

#### 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