

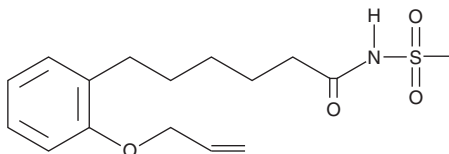
# PRODUCT INFORMATION



## MS-PPOH

Item No. 75770

CAS Registry No.: 206052-02-0  
Formal Name: N-(methylsulfonyl)-2-(2-propynyloxy)-benzenehexanamide  
MF:  $C_{16}H_{21}NO_4S$   
FW: 323.4  
Purity:  $\geq 98\%$   
UV/Vis.:  $\lambda_{max}$ : 218, 271 nm  
Supplied as: A crystalline solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

MS-PPOH is supplied as a crystalline solid. A stock solution may be made by dissolving the MS-PPOH in the solvent of choice, which should be purged with an inert gas. MS-PPOH is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of MS-PPOH in ethanol is approximately 25 mg/ml and approximately 30 mg/ml in DMSO and DMF.

MS-PPOH is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MS-PPOH should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. MS-PPOH has a solubility of approximately 0.30 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

MS-PPOH is a selective inhibitor of the epoxigenation reactions catalyzed by specific CYP450 isozymes.<sup>1</sup> MS-PPOH inhibits the formation of arachidonate 11,12-epoxides by CYP4A2 and CYP4A3 enzymes with an IC50 value of 13  $\mu M$ , but has no effect on the formation of 20-HETE, the  $\omega$ -hydroxylation product of CYP4A1.<sup>2</sup>

### References

1. Imig, J.D., Falck, J.R., and Inscho, E.W. Contribution of cytochrome P450 epoxigenase and hydroxylase pathways to afferent arteriolar autoregulatory responsiveness. *Br. J. Pharmacol.* **127**(6), 1399-1405 (1999).
2. Wang, M.H., Brand-Schieber, E., Zand, B.A., et al. Cytochrome P450-derived arachidonic acid metabolism in the rat kidney: Characterization of selective inhibitors. *J. Pharmacol. Exp. Ther.* **284**(3), 966-973 (1998).

#### 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, 04/08/2024

#### 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