

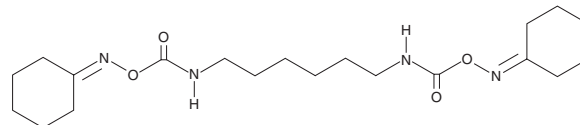
# PRODUCT INFORMATION



**RHC-80267**

Item No. 16432

**CAS Registry No.:** 83654-05-1  
**Formal Name:** 1,1'-[O,O']-[1,6-hexanediylbis(iminocarbonyl)]dioxime  
**Synonym:** U-57908  
**MF:** C<sub>20</sub>H<sub>34</sub>N<sub>4</sub>O<sub>4</sub>  
**FW:** 394.5  
**Purity:** ≥95%  
**Supplied as:** A crystalline 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

RHC-80267 is supplied as a crystalline solid. A stock solution may be made by dissolving the RHC-80267 in the solvent of choice, which should be purged with an inert gas. RHC-80267 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of RHC-80267 in ethanol and DMSO is approximately 5 mg/ml and approximately 15 mg/ml in DMF.

RHC-80267 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, RHC-80267 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. RHC-80267 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

RHC-80267 is an inhibitor of diacylglycerol lipase (DAGL; IC<sub>50</sub> = 4 μM for the canine platelet enzyme).<sup>1</sup> It also inhibits the serine hydrolases KIAA1363, fatty acid amide hydrolase (FAAH), BAT5, phospholipase A<sub>2</sub> (PLA<sub>2</sub>), hormone-sensitive lipase (HSL), and acetylcholinesterase (AChE) by greater than 60% when used at a concentration of 50 μM.<sup>2</sup> RHC-80267 (0.1-10 μM) potentiates ACh-induced relaxation of isolated rat mesenteric arteries.<sup>3</sup>

## References

1. Sutherland, C.A. and Amin, D. Relative activities of rat and dog platelet phospholipase A<sub>2</sub> and diglyceride lipase. Selective inhibition of diglyceride lipase by RHC 80267. *J. Biol. Chem.* **257(23)**, 14006-14010 (2014).
2. Hoover, H.S., Blankman, J.L., Niessen, S., *et al.* Selectivity of inhibitors of endocannabinoid biosynthesis evaluated by activity-based protein profiling. *Bioorg. Med. Chem. Lett.* **18**, 5838-5841 (2008).
3. Ghisdal, P., Vandenberg, G., Hamaide, M.-C., *et al.* The diacylglycerol lipase inhibitor RHC-80267 potentiates the relaxation to acetylcholine in rat mesenteric artery by anti-cholinesterase action. *Eur. J. Pharmacol.* **517(1-2)**, 97-102 (2005).

### 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, 12/19/2022

## 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