

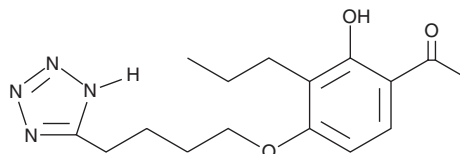
# PRODUCT INFORMATION



**LY171883**

Item No. 70710

**CAS Registry No.:** 88107-10-2  
**Formal Name:** 1-[2-hydroxy-3-propyl-4-[4-(1H-tetrazol-5-yl)butoxy]phenyl]-ethanone  
**MF:** C<sub>16</sub>H<sub>22</sub>N<sub>4</sub>O<sub>3</sub>  
**FW:** 318.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 218, 284 nm  
**Supplied as:** A crystalline solid  
**Storage:** Room temperature  
**Stability:** ≥4 years



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

## Laboratory Procedures

LY171883 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY171883 in the solvent of choice, which should be purged with an inert gas. LY171883 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of LY171883 in these solvents is approximately 25 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. For greater aqueous solubility, LY171883 can be directly dissolved in 0.5 M sodium bicarbonate (15 mg/ml) and then diluted with PBS (pH 7.2) to achieve the desired concentration or pH. We do not recommend storing the aqueous solution for more than one day.

## Description

LY171883 is a selective, potent, orally active antagonist of the leukotriene D<sub>4</sub> receptor.<sup>1</sup> Dissociation constants (K<sub>D</sub>) for LY171883 on guinea pig ileum and parenchyma are 0.07 and 0.34 μM, respectively.<sup>1</sup> LY171883 is an inhibitor of phosphodiesterase obtained from human polymorphonuclear leukocytes (IC<sub>50</sub> of 22.6 μM) and various guinea pig tissues (IC<sub>50</sub>s range from 6.9-209 μM).<sup>1</sup> At a concentration of 50-100 μM, LY171883 binds to the PPARγ nuclear receptor, inducing adipogenesis in cultured NIH3T3 fibroblasts.<sup>2,3</sup>

## References

1. Fleisch, J.H., Rinkema, L.E., Haisch, K.D., *et al.* LY171883, 1-[2-hydroxy-3-propyl-4-[4-(1H-tetrazol-5-yl)butoxy]phenyl]-ethanone, an orally active leukotriene D<sub>4</sub> antagonist. *J. Pharmacol. Exp. Ther.* **233**, 148-157 (1985).
2. Sala, A., Rossoni, G., Buccellati, C., *et al.* Formation of sulphidopeptide-leukotrienes by cell-cell interaction causes coronary vasoconstriction in isolated, cell-perfused heart of rabbit. *Br. J. Pharmacol.* **110(3)**, 1206-1212 (1993).
3. Tontonoz, P., Hu, E., Spiegelman, B.M. Stimulation of adipogenesis in fibroblasts by PPARγ2, a lipid-activated transcription factor. *Cell* **79(7)**, 1147-1156 (1994).

### 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, 11/28/2022

## 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