

LY171883

Catalog #: 27211

Lot #: 120605

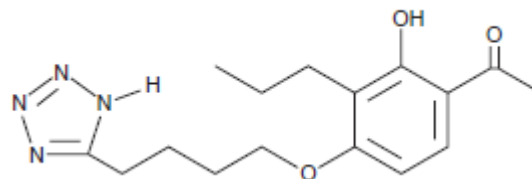
Size: 10 mg

Structure:

CAS Registry #: 88107-10-2

Purity: >98%

Chemical Formula: C₁₆H₂₂N₄O₃



Molecular Weight: 318.4

Description: LY171883 is a PDE inhibitor and a selective, potent inhibitor of the leukotriene D₄ receptor. Also, LY171883 binds to the PPAR γ nuclear receptor at 50-100 μ M and induces adipogenesis in cultured NIH3T3 fibroblasts.

Appearance: A crystalline solid

Solubility: Soluble in ethanol, DMSO and DMF, purged with an inert gas. Solubility in ethanol, DMSO and DMF is approximately 25 mg/ml. For maximum solubility in aqueous buffers, dissolve directly in 0.5 M Na₂CO₃ (15 mg/ml) then dilute with PBS (pH 7.2) to the desired concentration or pH. Do not store aqueous solutions for more than one day.

Biological Activity: LY171883 inhibits PDEs from human polymorphonuclear leukocytes with an IC₅₀ of 22.6 μ M, and from guinea pig tissues with an IC₅₀ range from 6.9 – 209 μ M.

Storage/Stability: Store at or below -20°C for up to two years.

Quality Control: The purity was determined by HPLC analysis.

References:

1. Fleisch, J.H., *et al.*, *J. Pharmacol. Exp. Ther.* 1985; **233**: 148-157.
2. Forman, B.M., *et al.*, *Cell.* 1995; **83**: 803-812.
3. Tontonoz, P., *et al.*, *Cell.* 1994; **79**: 1147-1156.