

## CP 80633

**Catalog #:** 27301

**Lot #:** 120628

**Size:** 10 mg

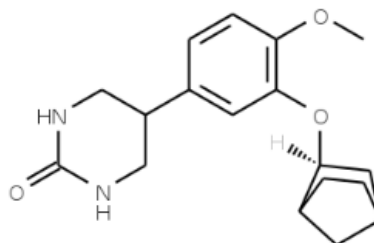
**Structure:**

**CAS Registry #:** 135637-46-6

**Purity:** ≥95%

**Chemical Formula:** C<sub>18</sub>H<sub>24</sub>N<sub>2</sub>O<sub>3</sub>

**Molecular Weight:** 316.39



**Description:** CP 80633 is a selective inhibitor of PDE4, which does not display significant isozyme selectivity. It inhibits hydrolysis of cAMP in isolated monocytes, eosinophils, human peripheral blood and T-cells. It has been shown to exhibit anti-inflammatory and bronchodilatory activity *in vivo*.

**Appearance:** A crystalline solid

**Solubility:** Soluble in 100 mM DMSO or 100 mM Ethanol.

**Biological Activity:** CP-80633 inhibits PDE4 isozymes (human lung IC<sub>50</sub> = 1.27 μM) in the absence of effects on PDE1, PDE2, PDE3 and PDE5 isozymes (IC<sub>50</sub> > 100 μM).

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

**Quality Control:** The purity was determined by HPLC analysis.

### Reference(s):

1. Turner, CR., *et al.*, *J. Pharmacol. Exp. Ther.* 1996 Sept;**278**(3):1349-55.
2. Cohan, VL., *et al.*, *J. Pharmacol. Exp. Ther.* 1996 Sept;**278**(3):1356-61.
3. Cheng, JB., *et al.*, *J. Pharmacol. Exp. Ther.* 1997 Feb;**280**(2):621-6.