

BAY-60-7550

Catalog #: 27215

Lot #: 120607

Size: 5 mg

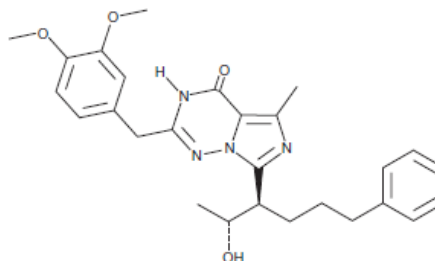
Structure:

CAS Registry #: 439083-90-6

Purity: ≥95%

Chemical Formula: C₂₇H₃₂N₄O₄

Molecular Weight: 476.6



Description: BAY-60-7550 is a potent inhibitor of PDE2. It is more selective for PDE2 than PDE1 by 50-fold. Similarly, it is 100-fold more selective for PDE4B, PDE5B, PDE7B, PDE8A, PDE9A, PDE10A, and PDE11A.

Appearance: A crystalline solid

Solubility: Soluble in ethanol, DMSO and DMF, purged with an inert gas. Solubility in ethanol, DMSO and DMF is ~10 mg/ml. Do not store aqueous solutions for more than one day.

Biological Activity: BAY-60-7550 inhibits PDE2 with an IC₅₀ of 2.0 nM in bovine, and 4.7 nM in humans.

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

Quality Control: The purity was determined by HPLC analysis.

References:

1. Boess, F.G., *et al.*, *Neuropharmacology*. 2004; **47**: 1081-1092.
2. Masood, A., *et al.*, *J. Pharmacol. Exp. Ther.* 2008; **326(2)**: 369-379.
3. Rutten, K., *et al.*, *Neuropsychopharmacology*. 2009; **34**: 1914-1925.