

C646

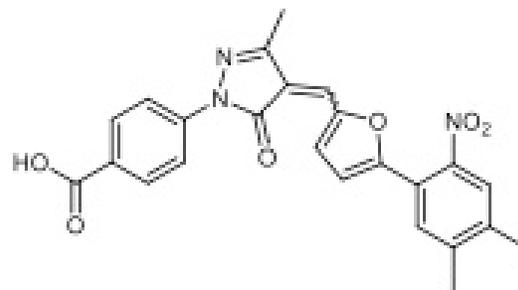
Catalog No: 14052, 14053**Format:** 25 mg, 5 mg**Chemical Properties:**

MW = 445.4

C₂₄H₁₉N₃O₆

CAS 328968-36-1

Physical Properties: Orange powder

Names: C-646, 4-[4-[[5-(4,5-Dimethyl-2-nitrophenyl)-2-furanyl]methylene]-4,5-dihydro-3-methyl-5-oxo-1H-pyrazol-1-yl]benzoic acid**Pharmacology:** Potent and selective inhibitor of the histone acetyltransferase p300/CBP (K_i = 400 nM)(ref 1). Induces apoptosis in prostate cancer cells (ref 2). Enhances fear extinction memory and synaptic plasticity in mice (ref 3). Promotes tau deacetylation reducing levels of pathogenic p-tau (ref 4). Cell permeable.**Solubilization:** May be dissolved in DMSO (10 mg/ml)**Fluorescent Properties:** N/A**Quality Control:**

>98% (TLC); NMR (Conforms)

References:

1. EM Bowers *et al. Chem. Biol.* 2010, 17:471
2. FR Santer *et al. Mol. Cancer Ther.* 2011, 10:1644
3. R Marek *et al. J. Neurosci.* 2011, 31:7486
4. S-W Min *et al. Neuron* 2010, 67:953

Storage and Guarantee: Store desiccated as supplied at 4°C for up to 2 years. Store solutions at -20°C for up to 2 months. This product is guaranteed for 6 months from date of arrival.