

---

## C646, p300/CBP Inhibitor

(Catalog No. M31000)

---

### Alternative Names

(E)-4-(4-((5-(4,5-Dimethyl-2-nitrophenyl)furan-2-yl)-methylene)-3-methyl-5-oxo-4,5-dihydro-1H-pyrazol-1-yl)benzoic acid

### Description

Cell-permeable. A reversible, p300/CBP (CREB-binding protein) histone acetyltransferase (HAT) inhibitor ( $K_i = 400$  nM). Specific inhibition to p300 (86%) compared to N-acetyltransferase, PCAF, GCN5, Rtt109, Sas or MOZ histone acetyltransferases (<10%). Suppresses histone H3 and H4 acetylation in mouse fibroblast cell lines.

### CAS Number

328968-36-1

### Chemical Formula

$C_{24}H_{19}N_3O_6$

### Molecular Weight

445.4

### Size

1 mg

### Purity

≥98% by NMR

### Solubility

DMSO (25 mM)

### Storage Temp

-20°C

### Ordering Information:

Products	Size	Cat. No.
C646, p300/CBP Inhibitor	1 mg	M31000 -1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*