

CTPB

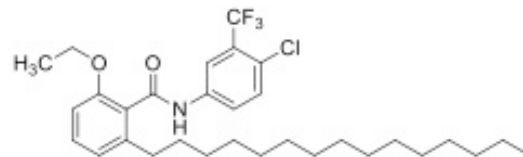
Catalog No: 14064, 14065**Format:** 25 mg, 5 mg**Chemical Properties:**

MW = 554.1

 $C_{31}H_{43}ClF_3NO_2$

CAS 586976-24-1

Physical Properties: Off white powder

Names: CTPB, N-(4-chloro-3-trifluoromethyl-phenyl)-2-ethoxy-6-pentadecyl-benzamide**Pharmacology:** A selective activator of p300 HAT (histone acetyltransferase). Selective over PCAF (p300 CPB-associated factor) (ref 1, 2).**Solubilization:** May be dissolved in DMSO (100 mg/ml); Ethanol (50 mg/ml)**Fluorescent Properties:** N/A**Quality Control:**98% (TLC; 1:10 EtOAc:Hexane; R_f = 0.26); NMR (Conforms)

References:

1. K Balasubramanyam *et al. J. Biol. Chem.* 2003, 278:19134
2. K Mantelingu *et al. J. Phys. Chem.* 2007, 111:4527

Storage and Guarantee: Store desiccated as supplied at -20°C for up to 3 years. Store solutions at -20°C for up to 3 months. This product is guaranteed for 6 months from date of arrival.