

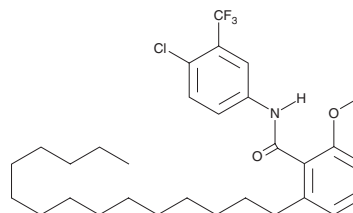
PRODUCT INFORMATION



CTPB

Item No. 19570

CAS Registry No.: 586976-24-1
Formal Name: N-[4-chloro-3-(trifluoromethyl)phenyl]-2-ethoxy-6-pentadecyl-benzamide
MF: C₃₁H₄₃ClF₃NO₂
FW: 554.1
Purity: ≥98%
UV/Vis.: 258 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Special Conditions: Keep cool and dry. Protect from light.

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

CTPB is supplied as a crystalline solid. A stock solution may be made by dissolving the CTPB in the solvent of choice, which should be purged with an inert gas. CTPB is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CTPB in these solvents is approximately 15, 20, and 30 mg/ml, respectively.

CTPB is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CTPB should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CTPB has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CTPB is an amide derivative that selectively activates the histone acetyltransferase (HAT) p300, with maximal activation at 275 μM.¹ It directly binds p300.² CTPB does not alter PCAF HAT or histone deacetylase activity.¹ It dose-dependently increases the acetylation of histones H3 and H4, with maximal effect at 200 μM.¹

References

1. Balasubramanyam, K., Swaminathan, V., Ranganathan, A., *et al.* Small molecule modulators of histone acetyltransferase p300. *J. Biol. Chem.* **278**(21), 19134-19140 (2003).
2. Devipriya, B., Parameswari, A.R., Rajalakshmi, G., *et al.* Exploring the binding affinities of p300 enzyme activators CTPB and CTB using docking method. *Indian J. Biochem. Biophys.* **47**(6), 364-369 (2010).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/29/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM