

CTPB

p300 HAT activator

Potent activator of p300 HAT (histone acetyltransferase), but not of PCAF (p300/CBP-associated factor) HAT activities.

Citations: 1

[View Online »](#)

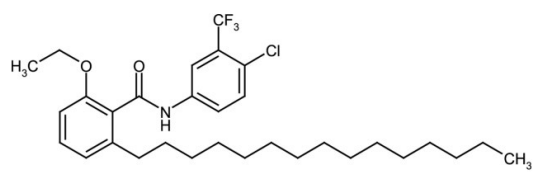
Ordering Information

[Order Online »](#)

ALX-420-033-M001	1mg
ALX-420-033-M005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-(4-Chloro-3-trifluoromethyl-phenyl)-2-ethoxy-6-pentadecyl-benzamide
Appearance	White to off-white crystalline solid.
CAS	586976-24-1
Couple Target	HAT
Couple Type	Activator
Formula	$C_{31}H_{43}ClF_3NO_2$
MW	554.1
Purity	≥97% (TLC)
Solubility	Soluble in DMSO.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com