

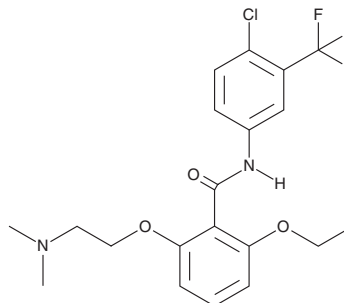
PRODUCT INFORMATION



YF-2

Item No. 31230

CAS Registry No.: 1311423-89-8
Formal Name: N-[4-chloro-3-(trifluoromethyl)phenyl]-2-[2-(dimethylamino)ethoxy]-6-ethoxy-benzamide
MF: C₂₀H₂₂ClF₃N₂O₃
FW: 430.9
Purity: ≥98%
UV/Vis.: λ_{max}: 260 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

YF-2 is supplied as a crystalline solid. A stock solution may be made by dissolving the YF-2 in the solvent of choice, which should be purged with an inert gas. YF-2 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of YF-2 in these solvents is approximately 30 mg/ml.

Description

YF-2 is a histone acetyltransferase (HAT) activator.¹ It inhibits proliferation of CCRF-CEM, Hs 578T, A549, and U251 cancer cells (IC₅₀s = 88.7-286 μM), as well as multidrug-resistant NCI/ADR-RES cells (IC₅₀ = 44.8 μM). YF-2 (5 mg/kg) increases reduced acetylated histone H3 levels in a mouse model of Alzheimer's disease induced by amyloid-β (1-42) (Aβ42; Item No. 20574). YF-2 (20 mg/kg) reduces contextual fear memory deficits in the transgenic APP/PS1 mouse model of Alzheimer's disease.

Reference

1. Feng, Y., Fa, M., Arancio, O., *et al.* Histone acetyltransferase activators and uses thereof. *Columbia University of New York*. **US20130121919A1** (2013).

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, 12/15/2022

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