

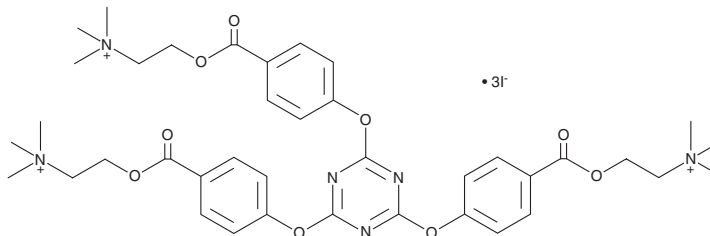
PRODUCT INFORMATION



TAE-1

Item No. 20464

CAS Registry No.: 1414469-59-2
Formal Name: 2,2',2'-[1,3,5-triazine-2,4,6-triyltris(oxy-4,1-phenylenecarbonyloxy)]tris[N,N,N-trimethyl-ethanaminium, triiodide
MF: C₃₉H₅₁N₆O₉ • 3I
FW: 1,128.6
Purity: ≥95%
UV/Vis.: λ_{max}: 224 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

TAE-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the TAE-1 in the solvent of choice, which should be purged with an inert gas. TAE-1 is soluble in organic solvents such as acetonitrile, DMSO, and dimethyl formamide (DMF). The solubility of TAE-1 in acetonitrile is approximately 5 mg/ml and approximately 20 mg/ml in DMSO and DMF. TAE-1 is also slightly soluble in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of TAE-1 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of TAE-1 in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

TAE-1 is a sym-triazine with diverse biological activities.¹ It inhibits *in vitro* amyloid-β (1-42) (Aβ₄₂; Item No. 20574) fibrilization when used at a concentration of 20 μM. TAE-1 inhibits acetylcholinesterase (AChE; IC₅₀ = 0.465 μM for the human erythrocyte enzyme). It also increases neuronal cellular process length and branching and synaptophysin protein levels in differentiated human SH-SY5Y neuronal cells.

Reference

1. Veloso, A.J., Chow, A.M., Dhar, D., *et al.* Biological activity of sym-triazines with acetylcholine-like substitutions as multitarget modulators of Alzheimer's disease. *ACS Chem. Neurosci.* **4(6)**, 924-929 (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, 12/16/2022

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