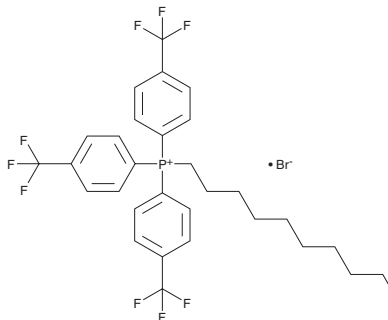


PRODUCT INFORMATION



4-CF₃-TPP-DC Item No. 33513

CAS Registry No.: 2661108-93-4
Formal Name: decyltris[4-(trifluoromethyl)phenyl]-
phosphonium, monobromide
Synonym: 4-CF₃-Triphenylphosphonium-DC
MF: C₃₁H₃₃F₉P • Br
FW: 687.5
Purity: ≥98%
UV/Vis.: λ_{max}: 225 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

4-CF₃-TPP-DC is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-CF₃-TPP-DC in the solvent of choice, which should be purged with an inert gas. 4-CF₃-TPP-DC is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 4-CF₃-TPP-DC in ethanol and DMF is approximately 5 mg/ml and approximately 10 mg/ml in DMSO.

Description

4-CF₃-TPP-DC is an inert mitochondriotropic carrier.¹ It does not induce mitochondrial depolarization or cell death in C2C12 cells.

Reference

1. Kulkarni, C.A., Fink, B.D., Gibbs, B.E., *et al.* A novel triphenylphosphonium carrier to target mitochondria without uncoupling oxidative phosphorylation. *J. Med. Chem.* **64(1)**, 662-676 (2021).

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, 10/02/2025

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