

PRODUCT INFORMATION



CPPA-TPP

Item No. 40439

Formal Name: (2-((4-cyano-4-((phenylcarbonothioyl)thio)pentanoyl)oxy)ethyl)triphenylphosphonium, monobromide

Synonyms: CPPA-Triphenylphosphonium, 4-Cyano-4-(phenylcarbonothioylthio)pentanoic Acid-TPP, 4-Cyano-4-(phenylcarbonothioylthio)pentanoic-Triphenylphosphonium

MF: $C_{33}H_{31}NO_2PS_2 \cdot Br$

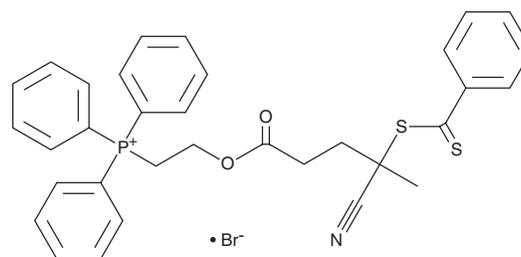
FW: 648.6

Purity: $\geq 75\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

CPPA-TPP is supplied as a solid. A stock solution may be made by dissolving the CPPA-TPP in the solvent of choice, which should be purged with an inert gas. CPPA-TPP is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

CPPA-TPP is a chain transfer agent conjugated to the mitochondria-targeting moiety triphenylphosphonium (TPP).¹ It has been used in the synthesis of reactive oxygen species-sensitive mitochondrial targeting polyprodrugs containing the topoisomerase I inhibitor camptothecin (Item No. 11694).

Reference

1. Zhang, W., Hu, X., Shen, Q., *et al.* Mitochondria-specific drug release and reactive oxygen species burst induced by polyprodrug nanoreactors can enhance chemotherapy. *Nat. Commun.* **10(1)**, 1704 (2019).

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, 10/22/2025

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