# **PRODUCT** INFORMATION



## (4-Carboxybutyl-d<sub>4</sub>)triphenylphosphonium (bromide)

Item No. 31254

CAS Registry No.:	42932-63-8	
Formal Name:	(4-carboxybutyl-2,2,3,3-d <sub>4</sub> )triphenyl-	
	phosphonium, monobromide	
Synonyms:	TPP-d <sub>4</sub> , 5-Triphenylphosphoniovaleric Acid-d <sub>4</sub>	
MF:	$C_{23}H_{20}D_4O_2P \bullet Br$	
FW:	447.3	
Chemical Purity:	≥98% ((4-Carboxybutyl)triphenylphosphonium)	
Deuterium		
Incorporation:	≥99% deuterated forms (d <sub>1</sub> -d <sub>4</sub> ); ≤1% d <sub>0</sub>	
Supplied as:	A solid	$\setminus$ /
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-Carboxybutyl-d<sub>4</sub>)triphenylphosphonium (bromide) is intended for use as an internal standard for the quantification of (4-carboxybutyl)triphenylphosphonium by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

(4-Carboxybutyl-d<sub>4</sub>)triphenylphosphonium (bromide) is supplied as a solid. A stock solution may be made by dissolving the (4-carboxybutyl- $d_{d}$ )triphenylphosphonium (bromide) in the solvent of choice, which should be purged with an inert gas. (4-Carboxybutyl- $d_{4}$ )triphenylphosphonium (bromide) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (4-carboxybutyl-d<sub>4</sub>)triphenylphosphonium (bromide) in ethanol is approximately 1 mg/ml and approximately 5 mg/ml in DMSO and DMF.

#### Description

(4-Carboxybutyl)triphenylphosphonium is a Wittig reagent.<sup>1</sup> It has been used in the synthesis of prostaglandin  $F_{2a}$  (PGF<sub>2a</sub>; Item No. 16010) and the monounsaturated fatty acid 7-methyl-6-octadecenoic acid.<sup>1,2</sup> (4-Carboxybutyl) triphenylphosphonium has also been used to enhance mitochondrial localization of albumin nanoparticles loaded with the anticancer agent docetaxel (Item No. 11637).<sup>3</sup>

#### References

- 1. Stork, G., Sher, P.M., and Chen, H.-L. Radical cyclization-trapping in the synthesis of natural products. A simple, stereocontrolled route to prostaglandin F<sub>2n</sub>. J. Am. Chem. Soc. 108(20), 6384-6385 (1986).
- 2. Carballeira, N.M. Recent developments in the total synthesis of bioactive marine fatty acids. Studies in Natural Products Chemistry. Atta-ur-Rahman, editor, 1st edition, Elsevier (2002).
- 3. Battogtokh, G., Gotov, O., Kang, J.H., et al. Triphenylphosphine-docetaxel conjugate-incorporated albumin nanoparticles for cancer treatment. Nanomedicine (Lond.) 13(3), 325-338 (2018).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFFTY 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

uyer 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, 11/04/2022

### 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