

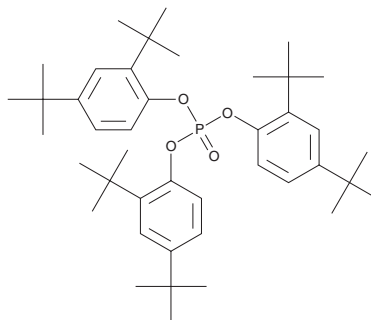
PRODUCT INFORMATION



TDTBPP

Item No. 39138

CAS Registry No.: 95906-11-9
Formal Name: 2,4-bis(1,1-dimethylethyl)-phenol,
1,1',1''-phosphate
Synonyms: AAN06119, TBPP-ox,
Tris(2,4-di-*t*-butylphenyl) Phosphate
MF: C₄₂H₆₃O₄P
FW: 662.9
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

TDTBPP is supplied as a solid. A stock solution may be made by dissolving the TDTBPP in the solvent of choice, which should be purged with an inert gas. TDTBPP is slightly soluble in methanol and chloroform.

TDTBPP is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

TDTBPP is a degradation product of the phenolic antioxidant tri(2,4-di-*t*-butylphenyl) phosphite (TDTBPPO), which has been used in plastics manufacturing.¹ It has been found in the dust from discarded electronic parts (e-waste dust), house dust, sediment, water, and air.^{1,2} TDTBPP has also been found in *V. negundo* and has anti-inflammatory activity.³ It reduces paw edema in a rat model of carrageenan-induced acute inflammation when administered at doses of 50 and 70 mg/kg.

References

1. Dorival-García, N., and Bones, J. Evaluation of solvent systems for optimized extractables studies of single use bioprocessing solutions. *J. Chromatogr. A* **1513**, 69-77 (2017).
2. Venier, M., Stubbings, W.A., Guo, J., *et al.* Tri(2,4-di-*t*-butylphenyl) phosphate: A previously unrecognized, abundant, ubiquitous pollutant in the built and natural environment. *Environ. Sci. Technol.* **52(22)**, 12997-13003 (2018).
3. Vinuchakkaravarthy, T., Kumaravel, K.P., Ravichandran, S., *et al.* Active compound from the leaves of *Vitex negundo* L. shows anti-inflammatory activity with evidence of inhibition for secretory phospholipase A₂ through molecular docking. *Bioinformation* **7(4)**, 199-206 (2011).

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, 10/25/2023

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