

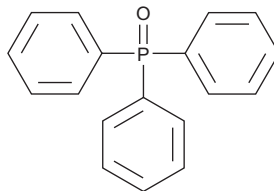
# PRODUCT INFORMATION



## Triphenylphosphine oxide

Catalog No. 9000289

**CAS Registry No.:** 791-28-6  
**Formal Name:** triphenyl-phosphine oxide  
**Synonyms:** NSC 398, TPPO  
**MF:** C<sub>18</sub>H<sub>15</sub>OP  
**FW:** 278.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 224 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

TPPO is supplied as a crystalline solid. A stock solution may be made by dissolving the TPPO in an organic solvent purged with an inert gas. TPPO is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of TPPO in ethanol is approximately 20 mg/ml and approximately 3 mg/ml in DMSO and DMF.

TPPO is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TPPO should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. TPPO has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

TPPO is a coordinating solvent used to activate crystallization of chemical compounds. It has been used in flame retardant applications, as an epoxy cure catalyst, and more recently, to produce nanostructures.

#### 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, 06/17/2020

#### 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