

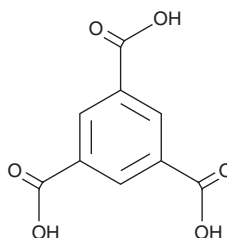
# PRODUCT INFORMATION



## Trimesic Acid

Item No. 19198

**CAS Registry No.:** 554-95-0  
**Formal Name:** 1,3,5-benzenetricarboxylic acid  
**Synonyms:** NSC 3998, TMA  
**MF:** C<sub>9</sub>H<sub>6</sub>O<sub>6</sub>  
**FW:** 210.1  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 212 nm  
**Supplied as:** A crystalline solid  
**Storage:** Room temperature  
**Stability:** ≥4 years



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

### Laboratory Procedures

Trimesic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the trimesic acid in the solvent of choice, which should be purged with an inert gas. Trimesic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of trimesic acid in ethanol is approximately 0.2 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Trimesic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, trimesic acid should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Trimesic acid has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Trimesic acid is a planar benzoic acid that has been used as a non-immunoglobulin protein scaffold for studies of protein-protein interactions.<sup>1</sup> It also has various industrial uses in synthesizing adhesive and coating materials, plastics engineering, and as a pharmaceutical intermediate.

### Reference

1. Lian, W., Upadhyaya, P., Rhodes, C.A., *et al.* Screening bicyclic peptide libraries for protein-protein interaction inhibitors: Discovery of a tumor necrosis factor-α antagonist. *J. Am. Chem. Soc.* **135**(32), 11990-11995 (2013).

#### 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, 11/11/2022

#### 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](http://WWW.CAYMANCHEM.COM)