

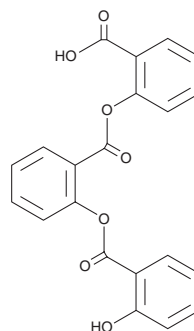
# PRODUCT INFORMATION



## Tri-Salicylic Acid

Item No. 37348

**CAS Registry No.:** 85531-17-5  
**Formal Name:** 2-[(2-hydroxybenzoyl)oxy]-benzoic acid,  
2-carboxyphenyl ester  
**MF:** C<sub>21</sub>H<sub>14</sub>O<sub>7</sub>  
**FW:** 378.3  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Tri-salicylic acid is supplied as a solid. A stock solution may be made by dissolving the tri-salicylic acid in the solvent of choice, which should be purged with an inert gas. Tri-salicylic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of tri-salicylic acid in ethanol, DMSO, and DMF is approximately 10 mg/ml in DMSO and approximately 25 mg/ml in ethanol and DMF.

### Description

Tri-salicylic acid is a degradation product of aspirin.<sup>1</sup> It is formed under conditions of thermal stress.

### Reference

1. Reepmeyer, J.C. Thermal decomposition of aspirin: formation of linear oligomeric salicylate esters. *J. Pharm. Sci.* **72(3)**, 322-323 (1983).

#### 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, 10/04/2023

#### 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