

PRODUCT INFORMATION



Benzoic Acid

Item No. 43179

CAS Registry No.: 65-85-0

Synonyms: Benzenecarboxylic Acid,
Carboxybenzene, NSC 149

MF: $C_7H_6O_2$

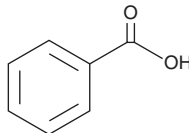
FW: 122.1

Purity: $\geq 99\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

Benzoic acid is supplied as a solid. A stock solution may be made by dissolving the benzoic acid in the solvent of choice, which should be purged with an inert gas. Benzoic acid is soluble (≥ 10 mg/ml) in ethanol and DMSO.

Description

Benzoic acid is an aromatic carboxylic acid and a synthetic intermediate.¹ It has been used in industrial phenol production. Formulations containing benzoic acid have been used as preservatives and flavoring agents in food, hygiene, cosmetic, and pharmaceutical products.

Reference

1. Del Olmo, A., Calzada, J., and Nuñez, M. Benzoic acid and its derivatives as naturally occurring compounds in foods and as additives: Uses, exposure, and controversy. *Crit. Rev. Food Sci. Nutr.* **57(14)**, 3084-3103 (2017).

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, 03/03/2025

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