

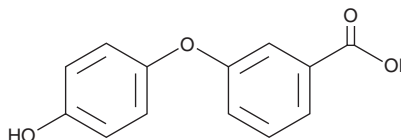
PRODUCT INFORMATION



(4'-Hydroxy)phenoxybenzoic Acid

Item No. 29891

CAS Registry No.: 35065-12-4
Formal Name: 3-(4-hydroxyphenoxy)-benzoic acid
MF: C₁₃H₁₀O₄
FW: 230.2
Purity: ≥95%
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

(4'-Hydroxy)phenoxybenzoic acid is supplied as a solid. A stock solution may be made by dissolving the (4'-hydroxy)phenoxybenzoic acid in the solvent of choice, which should be purged with an inert gas. (4'-Hydroxy)phenoxybenzoic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (4'-hydroxy)phenoxybenzoic acid in these solvents is approximately 5, 10, and 15 mg/ml, respectively.

(4'-Hydroxy)phenoxybenzoic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (4'-hydroxy)phenoxybenzoic acid should first be dissolved in DMF and then diluted with the aqueous buffer of choice. (4'-Hydroxy)phenoxybenzoic acid has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(4'-Hydroxy)phenoxybenzoic acid is a pyrethroid insecticide metabolite.^{1,2}

References

1. Akhtar, M.H., Mahadevan, S., and Russell, F. Cleavage of 3-phenoxybenzoic acid by chicken microsomal preparations. *J. Environ. Sci. Health B.* **28(5)**, 527-543 (1993).
2. Maloney, S.E., Maule, A., and Smith, A.R. Microbial transformation of the pyrethroid insecticides: Permethrin, deltamethrin, fastac, fenvalerate, and fluvalinate. *Appl. Environ. Microbiol.* **54(11)**, 2874-2876 (1988).

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, 12/08/2022

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