

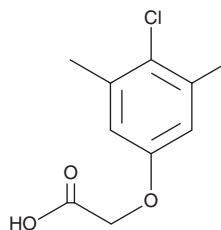
PRODUCT INFORMATION



(4-Chloro-3,5-dimethylphenoxy)acetic Acid

Item No. 16877

CAS Registry No.: 19545-95-0
Formal Name: 2-(4-chloro-3,5-dimethylphenoxy)-acetic acid
MF: C₁₀H₁₁ClO₃
FW: 214.6
Purity: ≥98%
UV/Vis.: λ_{max}: 230, 280 nm
Supplied as: A crystalline 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-Chloro-3,5-dimethylphenoxy)acetic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the (4-chloro-3,5-dimethylphenoxy)acetic acid in the solvent of choice, which should be purged with an inert gas. (4-Chloro-3,5-dimethylphenoxy)acetic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of (4-chloro-3,5-dimethylphenoxy)acetic acid in these solvents is approximately 30 mg/ml.

(4-Chloro-3,5-dimethylphenoxy)acetic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (4-chloro-3,5-dimethylphenoxy)acetic acid should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. (4-Chloro-3,5-dimethylphenoxy)acetic acid has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(4-Chloro-3,5-dimethylphenoxy)acetic acid is a synthetic auxin, a class of plant growth regulators, that has been used to promote hypocotyl elongation during seedling emergence.¹

Reference

1. Savaldi-Goldstein, S., Baiga, T.J., Pojer, F., *et al.* New auxin analogs with growth-promoting effects in intact plants reveal a chemical strategy to improve hormone delivery. *Proc. Natl. Acad. Sci. USA* **105(39)**, 15190-15195 (2008).

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, 11/02/2022

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