

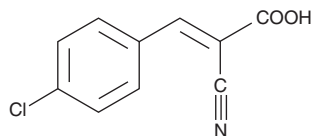
PRODUCT INFORMATION



4-chloro- α -Cyanocinnamic Acid

Item No. 9000981

CAS Registry No.: 20374-46-3
Formal Name: 3-(4-chlorophenyl)-2E-cyano-2-propenoic acid
Synonym: Cl-CCA
MF: C₁₀H₆ClNO₂
FW: 207.6
Purity: \geq 98%
UV/Vis.: λ_{max} : 203, 225, 301 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

4-chloro- α -Cyanocinnamic acid (Cl-CCA) is supplied as a crystalline solid. A stock solution may be made by dissolving the Cl-CCA in the solvent of choice, which should be purged with an inert gas. Cl-CCA is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Cl-CCA in these solvents is approximately 14 mg/ml.

Cl-CCA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Cl-CCA should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Cl-CCA has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Cl-CCA is a phenylpropanoid, related to cinnamic acid, which is synthesized by plants from phenylalanine. Cl-CCA readily co-crystallizes with various biomolecules (e.g., peptides, nucleic acids, etc.).

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, 10/19/2022

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