

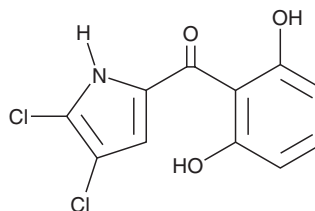
PRODUCT INFORMATION



Pyoluteorin

Item No. 19457

CAS Registry No.: 25683-07-2
Formal Name: (4,5-dichloro-1H-pyrrol-2-yl)
(2,6-dihydroxyphenyl)-methanone
Synonym: NSC 143092
MF: C₁₁H₇Cl₂NO₃
FW: 272.1
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

Pyoluteorin is supplied as a solid. A stock solution may be made by dissolving the pyoluteorin in the solvent of choice, which should be purged with inert gas. Pyoluteorin is soluble in the organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Pyoluteorin is a bichlorinated pyrrole produced by several species of *Pseudomonas* that demonstrates antibiotic, antifungal, and herbicidal activity. Pyoluteorin production is influenced by positive autoregulation and has been shown to function as a signal molecule negatively regulating additional secondary metabolites produced by *Pseudomonas*.¹

Reference

1. Brodhagen, M., Henkels M. D., and Loper, J. E. Positive autoregulation and signaling properties of pyoluteorin, an antibiotic produced by the biological control organism *Pseudomonas fluorescens* Pf-5. *Appl. Environ. Microbiol.* **70(3)**, 1758-1766 (2004).

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, 12/19/2022

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