

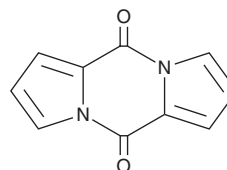
PRODUCT INFORMATION



Pyrocoll

Item No. 26464

CAS Registry No.: 484-73-1
Formal Name: 5H,10H-dipyrrolo[1,2-a:1',2'-d]pyrazine-5,10-dione
MF: $C_{10}H_6N_2O_2$
FW: 186.2
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years
Item Origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

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

Description

Pyrocoll is a bacterial metabolite originally isolated from *Streptomyces*.¹ It inhibits the growth of *A. aureus*, *A. globiformis*, *A. oxydans*, *A. pascens*, and *R. erythropolis* bacteria (MICs = 10, 1, 10, 3, and 10 $\mu\text{g/ml}$, respectively) and HMO2, HepG2, and MCF-7 cancer cells (GI_{50} s = 0.28, 0.42, and 2.2 $\mu\text{g/ml}$, respectively) *in vitro*. Pyrocoll is also active against *P. falciparum* and *T. rhodesiense* (IC_{50} s = 1.19 and 1.97 $\mu\text{g/ml}$, respectively).

Reference

- Dietera, A., Hamm, A., Fiedler, H.P., *et al.* Pyrocoll, an antibiotic, antiparasitic and antitumor compound produced by a novel alkaliphilic *Streptomyces* strain. *J. Antibiot. (Tokyo)* **56(7)**, 639-646 (2003).

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, 12/21/2022

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