

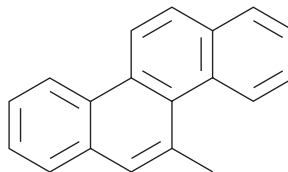
PRODUCT INFORMATION



5-methyl Chrysene

Item No. 18677

CAS Registry No.: 3697-24-3
Formal Name: 5-methyl-chrysene
Synonym: NSC 407620
MF: C₁₉H₁₄
FW: 242.3
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 269, 300, 312, 326 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

5-methyl Chrysene is supplied as a crystalline solid. A stock solution may be made by dissolving the 5-methyl chrysene in the solvent of choice, which should be purged with an inert gas. 5-methyl Chrysene is soluble in organic solvents such as chloroform and methanol.

5-methyl Chrysene is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

5-methyl Chrysene is a carcinogenic polycyclic aromatic hydrocarbon consisting of four fused rings and acts as an aryl hydrocarbon receptor agonist.¹ It is produced by the incomplete combustion of organic matter and found primarily in gasoline exhaust and tobacco smoke.¹ 5-methyl Chrysene can be activated to mutagenic metabolites by cytochrome P450 1A1 and 1B1.¹

Reference

1. Ahmad, S., Kabler, S.I., Rudd, L., *et al.* Cytotoxicity and mutagenicity of 5-methylchrysene and its 1,2-dihydrodiol in V79MZ cells modified to express human CYP1A1 or CYP1B1, in the presence or absence of human GSTP1 coexpression. *Toxicol. Lett.* **183(1-3)**, 99-104 (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, 10/20/2022

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