

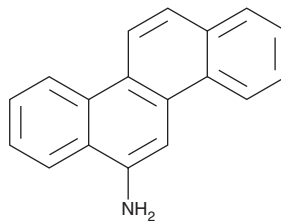
# PRODUCT INFORMATION



## 6-Chrysenamine

Item No. 35780

CAS Registry No.: 2642-98-0  
Synonyms: 6-Aminochrysene, NSC 80186  
MF:  $C_{18}H_{13}N$   
FW: 243.3  
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

6-Chrysenamine is supplied as a solid. A stock solution may be made by dissolving the 6-chrysenamine in the solvent of choice, which should be purged with an inert gas. 6-Chrysenamine is slightly soluble in chloroform and methanol.

### Description

6-Chrysenamine is a carcinogen and an anticancer agent.<sup>1,2</sup> Neonatal administration of 6-chrysenamine (600  $\mu\text{g}/\text{animal}$ ; s.c.) induces the formation of hepatomas and pulmonary adenocarcinomas in adult mice.<sup>1</sup> 6-Chrysenamine (3-110 mg/kg) inhibits tumor growth in an EO771 murine mammary carcinoma model.<sup>2</sup>

### References

1. Roe, F.J., Carter, R.L., and Adamthwaite, S. Induction of liver and lung tumours in mice by 6-aminochrysene administered during the 1st 3 days of life. *Nature* **221(5185)**, 1063-1064 (1969).
2. Gelzer, J. and Loustalot, P. Chrysenex<sup>®</sup> in experimental advanced mammary cancer. *Eur. J. Cancer* **3(1)**, 79-80 (1967).

#### 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, 12/13/2022

#### 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