

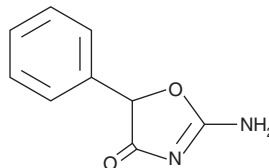
PRODUCT INFORMATION



Pemoline

Item No. 14208

CAS Registry No.: 2152-34-3
Formal Name: 2-amino-5-phenyl-4(5H)-oxazolone
Synonyms: NSC 25159, NSC 169499
MF: $C_9H_8N_2O_2$
FW: 176.2
Purity: $\geq 98\%$
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.

Description

Pemoline (Item No. 14208) is an analytical reference standard categorized as a stimulant.^{1,2} Pemoline induces hyperactivity and repetitive self-biting in mice and rats.¹ Pemoline prevents memory formation in the passive avoidance test in rats.² Pemoline is regulated as a Schedule IV compound in the United States. This product is intended for research and forensic applications.

References

1. Genovese, E., Napoli, P.A., and Bolego-Zonta, N. Selfaggressiveness: A new type of behavioral change induced by pemoline. *Life Sci.* **8(9)**, 513-515 (1969).
2. Gurowitz, E.M., Lubar, J.F., Ain, B.R., *et al.* Disruption of passive avoidance learning by magnesium pemoline. *Psychon. Sci.* **8(1)**, 19-20 (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, 01/13/2026

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