

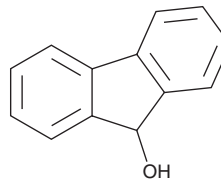
PRODUCT INFORMATION



Fluorenol

Item No. 19799

CAS Registry No.: 1689-64-1
Formal Name: 9H-fluoren-9-ol
Synonyms: 9-Fluorenol Alcohol,
9-Hydroxyfluorene, NSC 5320
MF: C₁₃H₁₀O
FW: 182.2
Purity: ≥98%



UV/Vis.: λ_{max}: 209, 221, 227, 235, 271 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly

Description

Fluorenol (Item No. 19799) is an analytical reference standard that is classified as a nootropic. It is a dopamine reuptake inhibitor (IC₅₀ = 9 μM) and a major metabolite of a compound developed as a wakefulness-promoting agent.¹ At 100 mg/kg, fluorenol is reported to promote wakefulness in mice with 39% increased effectiveness over the structurally related compound, modafinil (Item No. 15417).¹ This product is intended for forensic and research applications.

Reference

1. Dunn, D., Hostetler, G., Igbal, M., *et al.* Wake promoting agents: Search for next generation modafinil, lessons learned: part III. *Bioorg. Med. Chem. Lett.* **22(11)**, 3751-3753 (2016).

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, 07/29/2016

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