

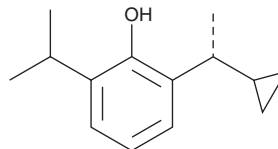
PRODUCT INFORMATION



Cipecfol

Item No. 40951

CAS Registry No.: 1637741-58-2
Formal Name: 2-[(1R)-1-cyclopropylethyl]-6-(1-methylethyl)-phenol
Synonyms: Ciprofol, HSK3486
MF: C₁₄H₂₀O
FW: 204.3
Purity: ≥98%
Supplied as: A neat liquid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Cipecfol is supplied as a neat liquid. A stock solution may be made by dissolving the cipecfol in the solvent of choice, which should be purged with an inert gas. Cipecfol is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

Cipecfol is a general anesthetic.¹ It binds to the GABA_A receptor when used at a concentration of 10 μM. Cipecfol inhibits decreases in superoxide dismutase (Sod) levels, increases in NADPH oxidase (Nox) levels, and apoptosis induced by the β₁- and β₂-adrenergic receptor agonist isoproterenol (Item No. 15592) in H9c2 rat cardiomyocytes and reduces infarct size in a mouse model of isoproterenol-induced myocardial infarction.² It induces loss of righting reflex (LORR) in mice (ED₅₀ = 1.5 mg/kg).¹

References

1. Qin, L., Ren, L., Wan, S., *et al.* Design, synthesis, and evaluation of novel 2,6-disubstituted phenol derivatives as general anesthetics. *J. Med. Chem.* **60**(9), 3606-3617 (2017).
2. Yang, Y., Xia, Z., Xu, C., *et al.* Ciprofol attenuates the isoproterenol-induced oxidative damage, inflammatory response and cardiomyocyte apoptosis. *Front. Pharmacol.* **13**, 1037151 (2022).

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, 05/11/2026

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