

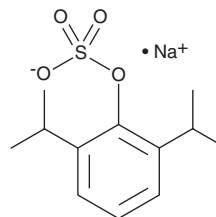
PRODUCT INFORMATION



Propofol Sulfate (sodium salt)

Item No. 9002416

Formal Name: 1-(hydrogen sulfate) 2,6-bis(1-methylethyl)-phenol, monosodium salt
MF: C₁₂H₁₇O₄S • Na
FW: 280.3
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

Propofol is a short-acting anesthetic agent that acts as a potentiator of the GABA_A receptor (EC₅₀ = 10 μM).¹⁻⁴ It also has direct activation effects at the GABA_A receptor, alters receptor desensitization and deactivation, and blocks channel activity.² Propofol shows little or no specificity among the diverse GABA_A receptor subunits.² Intravenous injection of propofol is usually accompanied with pain due to activation of the transient receptor potential ankyrin 1 channel.⁵

References

1. Sooksawate, T. and Simmonds, M.A. Influence of membrane cholesterol on modulation of the GABA_A receptor by neuroactive steroids and other potentiators. *Brit. J. Pharmacol.* **134**, 1303-1311 (2001).
2. Krasowski, M.D., Hong, X., Hopfinger, A.J., et al. 4D-QSAR analysis of a set of propofol analogues: Mapping binding sites for an anesthetic phenol on the GABA_A receptor. *J. Med. Chem.* **45(15)**, 3210-3221 (2002).
3. Lingamaneni, R. and Hemmings, H.C., Jr. Differential interaction of anaesthetics and antiepileptic drugs with neuronal Na⁺ channels, Ca²⁺ channels, and GABA_A receptors. *Br. J. Anaesth.* **90(2)**, 199-211 (2003).
4. Li, M., Zhao, X., Zhang, L., et al. Effects and safety of magnesium sulfate on propofol-induced injection pain, a meta-analysis of randomized controlled trials. *Int. J. Clin. Exp. Med.* **8(5)**, 6813-6821 (2015).
5. Baraldi, P.G., Preti, D., Materazzi, S., et al. Transient receptor potential ankyrin 1 (TRPA1) channel as emerging target for novel analgesics and anti-inflammatory agents. *J. Med. Chem.* **53(14)**, 5085-5107 (2010).

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, 07/25/2023

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