

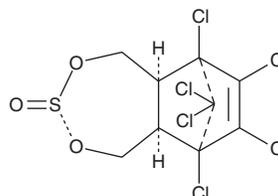
PRODUCT INFORMATION



Endosulfan II

Item No. 24254

CAS Registry No.: 33213-65-9
Formal Name: (3 α ,5 α ,6 β ,9 β ,9 α)-6,7,8,9,10,10-hexachloro-1,5,5a,6,9,9a-hexahydro-6,9-methano-2,4,3-benzodioxathiepin 3-oxide
Synonyms: β -Endosulfan, Endosulfan B
MF: C₉H₆Cl₆O₃S
FW: 406.9
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

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

Description

Endosulfan II is an organochlorine insecticide and a stereoisomer of endosulfan I.¹ It is active against a variety of insects including bollworms and tobacco budworms (LD₅₀s = 4.14 and 4.95 mg/g, respectively).^{2,3} Endosulfan II binds to GABA receptors in rat brain membranes with an IC₅₀ value of 60 nM and is less toxic to rats than endosulfan I (LD₅₀s = 240 and 18 mg/kg, respectively).

References

1. Abalis, I.M., Eldefrawi, M.E., and Eldefrawi, A.T. High-affinity stereospecific binding of cyclodiene insecticides and γ -hexachlorocyclohexane to γ -aminobutyric acid receptors of rat brain. *Pest. Biochem. Phys.* **24(1)**, 95-102 (1985).
2. Wolfenbarger, D.A. and Guerra, A.A. Toxicity of endosulfan and its isomers to the bollworm and tobacco budworm. *J. Econ. Entomol* **65(4)**, 1122-1123 (1972).
3. Busvine, J.R. The insecticidal potency of γ -BHC and the chlorinated cyclodiene compounds and the significance of resistance to them. *Bull. Entom. Res.* **55(2)**, 271-288 (1964).

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, 09/19/2023

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