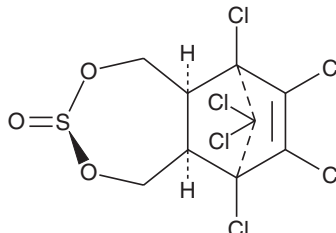


PRODUCT INFORMATION



Endosulfan I Item No. 24253

CAS Registry No.: 959-98-8
Formal Name: (3 α ,5 $\alpha\beta$,6 α ,9 α ,9 $\alpha\beta$)-6,7,8,9,10,10-hexachloro-1,5,5a,6,9,9a-hexahydro-6,9-methano-2,4,3-benzodioxathiepin 3-oxide
Synonyms: α -Endosulfan, Endosulfan A
MF: C₉H₆Cl₆O₃S
FW: 406.9
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.

Laboratory Procedures

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

Description

Endosulfan I is an organochlorine insecticide and a stereoisomer of endosulfan II (Item No. 24254).¹ It is active against a variety of insects, including bollworms (LD₅₀ = 0.63 mg/g).^{1,2} Endosulfan I binds to GABA receptors in rat brain membranes with an IC₅₀ value of 30 nM and is toxic to rats (LD₅₀ = 18 mg/kg).³

References

1. 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).
2. 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).
3. 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).

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