

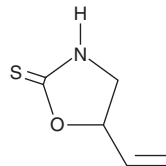
PRODUCT INFORMATION



(±)-Goitrin

Item No. 26382

CAS Registry No.: 13190-34-6
Formal Name: 5-ethenyl-2-oxazolidinethione
Synonyms: DL-Goitrin, (R,S)-Goitrin
MF: C₅H₇NOS
FW: 129.2
Purity: ≥98%
UV/Vis.: λ_{max}: 246 nm
Supplied as: A crystalline 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

(±)-Goitrin is supplied as a crystalline solid. A stock solution may be made by dissolving the (±)-goitrin in the solvent of choice, which should be purged with an inert gas. (±)-Goitrin is slightly soluble in ethanol and DMSO.

Description

(±)-Goitrin is a goitrogen and antithyroid compound that has been found in plants of the *Brassica* family.^{1,2} Dietary administration of (±)-goitrin (0.06-0.18% w/w) increases thyroid weight and decreases thyroid radioactive iodine uptake in rats fed a low-iodine diet.¹ It also increases thyroid weight in chicks in a dose-dependent manner.² (±)-Goitrin inhibits norepinephrine formation by bovine dopamine-β-hydroxylase *in vitro* when used at a concentration of 0.2 mM.³

References

1. Faiman, C., Ryan, R.J., and Eichel, H.J. Effect of goitrin analogues and related compounds on the rat thyroid gland. *Endocrinology* **81(1)**, 88-92 (1967).
2. Akiba, Y. and Matsumoto, J. Antithyroid activity of goitrin in chicks. *Poult. Sci.* **55(2)**, 716-719 (1976).
3. Zenker, N., Hubbard, L.S., and Wright, J. Inhibition of dopamine β-hydroxylase by goitrin, a natural antithyroid compound. *J. Nat. Prod.* **51(5)**, 862-865 (1988).

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, 12/06/2022

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