

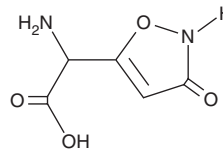
PRODUCT INFORMATION



Ibotenic Acid

Item No. 14584

CAS Registry No.: 2552-55-8
Formal Name: α-amino-2,3-dihydro-3-oxo-5-isoxazoleacetic acid
Synonym: (±)-Ibotenic Acid, IBO, NSC 204850
MF: C₅H₆N₂O₄
FW: 158.1
Purity: ≥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

Ibotenic acid is supplied as a solid. A stock solution may be made by dissolving the ibotenic acid in water. The solubility of ibotenic acid in water is approximately 10 mM. We do not recommend storing the aqueous solution for more than one day.

Description

Ibotenic acid is a neuroexcitatory amino acid originally isolated from *Amanita* species that functions as a NMDA and metabotropic glutamate receptor agonist.¹ As a neurotoxin, ibotenic acid is often used to induce brain lesions in animals that model cognitive dysfunctions resulting from neurodegenerative diseases, traumatic brain injury, and stroke.²

References

1. Curtis, D.R., Lodge, D., and McLennan, H. The excitation and depression of spinal neurones by ibotenic acid. *J. Physiol.* **291**, 19-28 (1979).
2. Tayebati, S.K. Animal models of cognitive dysfunction. *Mech. Ageing Dev.* **127(2)**, 100-108 (2005).

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, 02/22/2023

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