

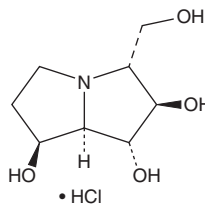
PRODUCT INFORMATION



Australine (hydrochloride)

Item No. 30051

CAS Registry No.: 186766-07-4
Formal Name: (1R,2R,3R,7S,7aR)-hexahydro-3-(hydroxymethyl)-1H-pyrrolizine-1,2,7-triol, monohydrochloride
Synonyms: 7a-*epi*-Alexine, (+)-Australine
MF: C₈H₁₅NO₄ • HCl
FW: 225.7
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years
Item Origin: Synthetic



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

Laboratory Procedures

Australine is supplied as a crystalline solid. A stock solution may be made by dissolving the australine in water. The solubility of australine in water is approximately 20 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Australine is a pyrrolizidine alkaloid originally isolated from *C. australe* that has enzyme inhibitory activities.^{1,2,3} It is an inhibitor of glucoamylase (IC₅₀ = 5.8 μM) that also inhibits glucosidase I, sucrase, maltase, and *A. niger* α-glucosidase (IC₅₀s = 20, 28, 35, and 28 μM, respectively).^{2,3} Australine is selective for these enzymes over glucosidase II, α- and β-mannosidase, and α- and β-galactosidase up to 500 μM, β-glucosidase, with only 5% inhibition at 66 μM, as well as isomaltase and trehalase (IC₅₀ = 97 and 160 μM, respectively). Australine (500 μg/ml) inhibits glycoprotein processing of viral glycoproteins in influenza virus-infected MDCK cells and induces the accumulation of glycoproteins.²

References

1. Molyneux, R.J., Benson, M., Wong, R.Y., *et al.* Australine, a novel pyrrolizidine alkaloid glucosidase inhibitor from *Castanospermum australe*. *J. Nat. Prod.* **51(6)**, 1198-1206 (1988).
2. Tropea, J.E., Molyneux, R.J., Kaushal, G.P., *et al.* Australine, a pyrrolizidine alkaloid that inhibits amyloglucosidase and glycoprotein processing. *Biochemistry* **28(5)**, 2027-2034 (1989).
3. Kato, A., Kano, E., Adachi, I., *et al.* Australine and related alkaloids: Easy structural confirmation by ¹³C NMR spectral data and biological activities. *Tetrahedron Asymmetry* **14(3)**, 325-331 (2003).

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, 04/15/2020

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