

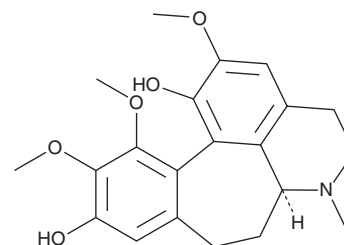
PRODUCT INFORMATION



Floramultine

Item No. 31419

CAS Registry No.: 1354-66-1
Formal Name: 4,5,6,6aS,7,8-hexahydro-2,11,12-trimethoxy-6-methylbenzo[6,7]cyclohept[1,2,3-ij]isoquinoline-1,10-diol
Synonyms: (S)-Floramultine, (S)-Bechuanine, Merenderine
MF: C₂₁H₂₅NO₅
FW: 371.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Tripladenia cunninghamii*



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

Laboratory Procedures

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

Description

Floramultine is an isoquinoline alkaloid that has been found in *C. speciosum*.¹ It inhibits butyrylcholinesterase (BChE) and acetylcholinesterase (AChE; K_s = 23.44 and 151.3 μM, respectively).

Reference

1. Basova, N.E., Rozengart, E.V., and Suvorov, A.A. Cholinergic activity of isoquinoline alkaloids from the showy autumn crocus (*Colchicum speciosum* Stev.). *Dokl. Biochem. Biophys.* **406**, 27-31 (2006).

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, 11/09/2022

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