

PRODUCT INFORMATION



β -Carboline Alkaloids Mixture

Item No. 39405

Supplied as: A solution in methanol

Storage: -20°C

Stability: ≥ 1 year

Special conditions: Vortex or sonicate before use

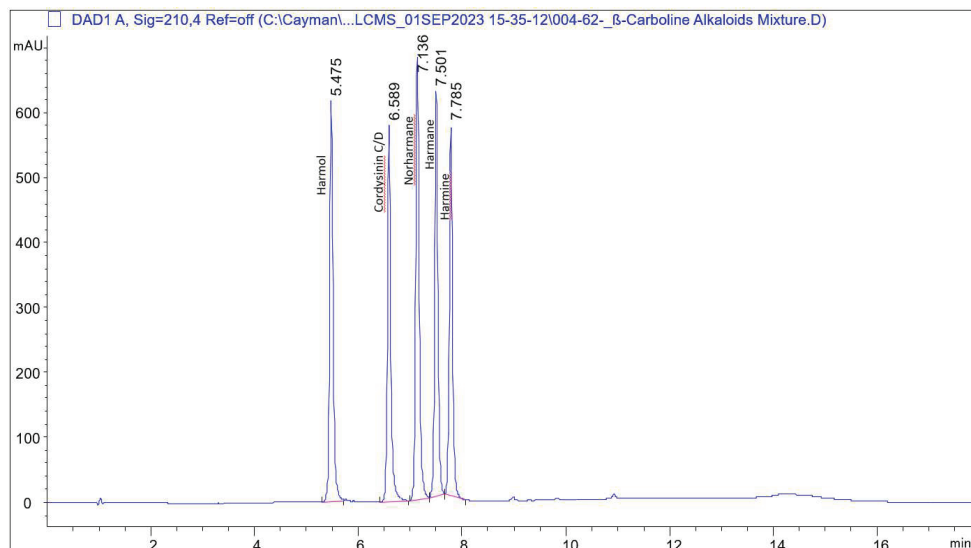
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

β -Carboline Alkaloids Mixture (Item No. 39405) is a mixture of analytical reference standards containing β -carboline alkaloids. The mixture contains harmane, norharmane, harmol, harmine, and cordysinin C/D. The mixture is supplied in an amber ampule in which the headspace has been purged with argon. It is designed for use in method development, system suitability, and as a reference standard for these substances during routine analytical checks using HPLC-UV and LC-MS instrumentation. This product is intended for research and forensic applications.

This mixture contains five β -carboline alkaloids. Included are:

Item No. 39405	β -Carboline Alkaloids Mixture	
Item Name	CAS Number	Concentration (mg/ml)
Harmane	486-84-0	1.0
Norharmane	244-63-3	1.0
Harmol	487-03-6	1.0
Harmine	442-51-3	1.0
Cordysinin C/D	148683-93-6	1.0



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, 09/20/2023

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