

PRODUCT INFORMATION



Benzoylhypaconine

Item No. 34046

CAS Registry No.: 63238-66-4
Formal Name: (1 α ,6 α ,14 α ,15 α ,16 β)-1,6,16-trimethoxy-4-(methoxymethyl)-20-methyl-aconitane-8,13,14,15-tetrol, 14-benzoate

Synonyms: Benzoylhypacoitine, BHA

MF: C₃₁H₄₃NO₉

FW: 573.7

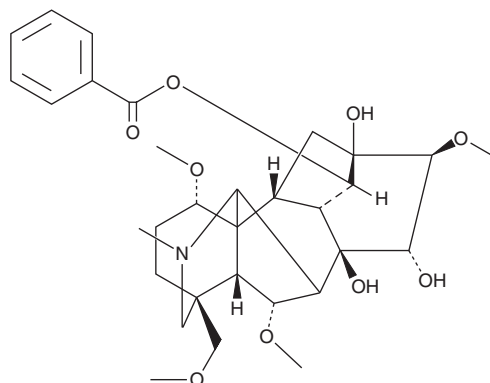
Purity: \geq 98%

UV/Vis.: λ_{max} : 231 nm

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years



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

Laboratory Procedures

Benzoylhypaconine is supplied as a solid. A stock solution may be made by dissolving the benzoylhypaconine in the solvent of choice, which should be purged with an inert gas. Benzoylhypaconine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of benzoylhypaconine in these solvents is approximately 30 mg/ml.

Description

Benzoylhypaconine is a monoester diterpene alkaloid and hydrolyzed product of hypaconitine (Item No. 29656) that has been found in *Aconitum* and has anti-inflammatory activity.¹⁻³ It inhibits COX-2 activity in a cell-free assay (IC₅₀ = 20.5 μ M).² Benzoylhypaconine (200 mg/kg) reduces histamine-induced vascular permeability in rats.³

References

1. Ng, S.W., Ching, C.K., Chan, A.Y.W., *et al.* Simultaneous detection of 22 toxic plant alkaloids (aconitum alkaloids, solanaceous tropane alkaloids, sophora alkaloids, strychnos alkaloids and colchicine) in human urine and herbal samples using liquid chromatography-tandem mass spectrometry. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **942-943**, 63-69 (2013).
2. Zhu, H., Liu, S., Li, X., *et al.* Bioactivity fingerprint analysis of cyclooxygenase-2 ligands from *radix Aconiti* by ultrafiltration-UPLC-MSn. *Anal. Bioanal. Chem.* **405(23)**, 7437-7445 (2013).
3. Hikino, H., Konno, C., Takata, H., *et al.* Antiinflammatory principles of *Aconitum* roots. *J. Pharmacobiodyn.* **3(10)**, 514-525 (1980).

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, 11/17/2022

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