

PRODUCT INFORMATION



Salvinorin A

Item No. 11487

CAS Registry No.: 83729-01-5
Formal Name: (2S,4aR,6aR,7R,9S,10aS,10bR)-9-(acetyloxy)-2-(3-furanyl)dodecahydro-6a,10b-dimethyl-4,10-dioxo-2H-naphtho[2,1-c]pyran-7-carboxylic acid, methyl ester

Synonyms: Divinorin A, Sal A

MF: C₂₃H₂₈O₈

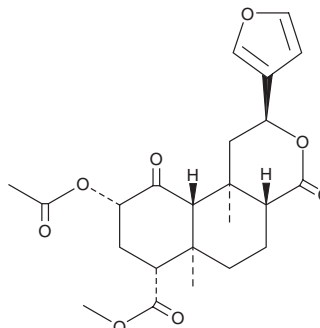
FW: 432.5

Purity: ≥95%

Supplied as: A neat solid

Storage: -20°C

Stability: ≥5 years



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

Description

Salvinorin A (Item No. 11487) is an analytical reference material categorized as an opioid and a diterpene.¹ Salvinorin A has been found in *S. divinorum*, a plant with hallucinogenic properties.¹⁻⁴ This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

References

1. Roth, B.L., Baner, K., Westkaemper, R., *et al.* Salvinorin A: A potent naturally occurring nonnitrogenous κ opioid selective agonist. *Proc. Natl. Acad. Sci. USA* **99**(18), 11934-11939 (2002).
2. Harding, W.W., Tidgewell, K., Schmidt, M., *et al.* Salvinicins A and B, new neoclerodane diterpenes from *Salvia divinorum*. *Org. Lett.* **7**(14), 3017-3020 (2005).
3. Tsujikawa, K., Kuwayama, K., Miyaguchi, H., *et al.* Determination of salvinorin A and salvinorin B in *Salvia divinorum*-related products circulated in Japan. *Forensic Sci. Int.* **180**(2-3), 105-109 (2008).
4. Munro, T.A., Duncan, K.K., Xu, W., *et al.* Standard protecting groups create potent and selective κ opioids: Salvinorin B alkoxymethyl ethers. *Bioorg. Med. Chem.* **16**(3), 1279-1286 (2008).

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, 04/19/2023

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