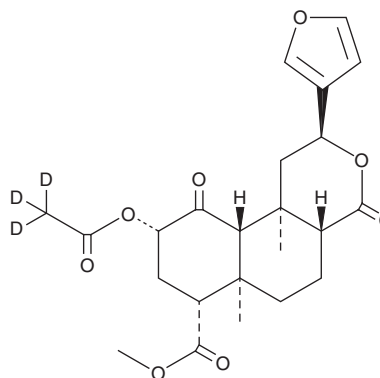


PRODUCT INFORMATION



Salvinorin A-d₃ Item No. 36410

CAS Registry No.: 791114-98-2
Formal Name: (2S,4aR,6aR,7R,9S,10aS,10bR)-9-(acetyl-d₃-oxy)-2-(3-furanyl)dodecahydro-6a,10b-dimethyl-4,10-dioxo-2H-naphtho[2,1-c]pyran-7-carboxylic acid, methyl ester
Synonyms: Divinorin A-d₃, Sal A-d₃
MF: C₂₃H₂₅D₃O₈
FW: 435.5
Chemical Purity: ≥98% (Salvinorin A)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
Supplied as: A crystalline 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-d₃ (Item No. 36410) is intended for use as an internal standard for the quantification of salvinorin A (Item No. 11487) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Salvinorin A is categorized as an opioid and a diterpene.¹ It has been found in *S. divinorum*, a plant with hallucinogenic properties.¹⁻⁴ This product is intended for research and forensic applications.

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, 07/07/2023

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