

# PRODUCT INFORMATION



## Salvinorin B

Item No. 23582

**CAS Registry No.:** 92545-30-7  
**Formal Name:** (2S,4aR,6aR,7R,9S,10aS,10bR)-2-(3-furanyl) dodecahydro-9-hydroxy-6a,10b-dimethyl-4,10-dioxo-2H-naphtho[2,1-c]pyran-7-carboxylic acid, methyl ester

**Synonyms:** Divinorin B, Sal B

**MF:** C<sub>21</sub>H<sub>26</sub>O<sub>7</sub>

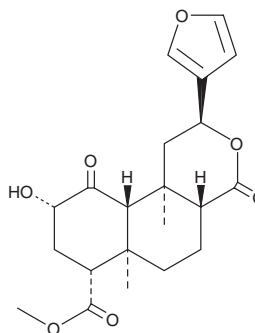
**FW:** 390.4

**Purity:** ≥95%

**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 B (Item No. 23582) is an analytical reference standard categorized as a diterpene.<sup>1</sup> Salvinorin B is a presumptive metabolite of salvinorin A (Item Nos. 38208 | 11487).<sup>2,3</sup> It has been found in *S. divinorum*, a plant with hallucinogenic properties.<sup>3,4</sup> This product is intended for research and forensic applications.

### References

1. Vortherms, T.A. and Roth, B.L. Salvinorin A: From natural product to human therapeutics. *Mol. Interv.* **6(5)**, 257-265 (2006).
2. Schmidt, M.D., Schmidt, M.S., Butelman, E.R., et al. Pharmacokinetics of the plant-derived  $\kappa$ -opioid hallucinogen salvinorin A in nonhuman primates. *Synapse* **58(3)**, 208-210 (2005).
3. Schmidt, M.S., Prisinzano, T.E., Tidgewell, K., et al. Determination of Salvinorin A in body fluids by high performance liquid chromatography-atmospheric pressure chemical ionization. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **818(2)**, 221-225 (2005).
4. 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).

#### 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, 06/20/2023

#### 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