

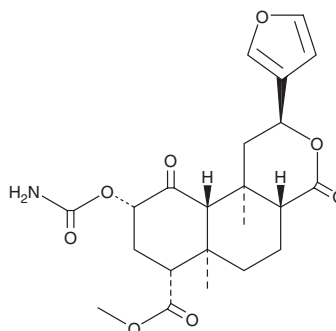
PRODUCT INFORMATION



Salvinorin A Carbamate

Item No. 22245

CAS Registry No.: 858345-57-0
Formal Name: (2S,4aR,6aR,7R,9S,10aS,10bR)-9-[(aminocarbonyloxy)-2-(3-furanyl)dodecahydro-6a,10b-dimethyl-4,10-dioxo-2H-naphtho[2,1-c]pyran-7-carboxylic acid, methyl ester
Synonyms: Divinorin A Carbamate, Sal A Carbamate
MF: C₂₂H₂₇NO₈
FW: 433.5
Purity: ≥90%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Salvinorin A carbamate is supplied as a crystalline solid. A stock solution may be made by dissolving the salvinorin A carbamate in the solvent of choice. Salvinorin A carbamate is soluble in organic solvents such as acetonitrile, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of salvinorin A carbamate in acetonitrile and DMF is approximately 1 mg/ml and approximately 2 mg/ml in DMSO.

Salvinorin A carbamate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, salvinorin A carbamate should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Salvinorin A carbamate has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Salvinorin A carbamate is a potent κ -opioid receptor (KOR) full agonist that is almost as potent as salvinorin A (Item No. 11487) with EC₅₀ values of 6.2 and 4.5 nM, respectively, for activation of the human KOR to enhance binding of [³⁵S]GTPγS.¹ The addition of a carbamate group to salvinorin A increases biological stability by decreasing deacetylation.

Reference

1. Béguin, C., Richards, M.R., Wang, Y., *et al.* Synthesis and *in vitro* pharmacological evaluation of salvinorin A analogues modified at C(2). *Bioorg. Med. Chem. Lett.* **15(11)**, 2761-2765 (2005).

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, 10/23/2025

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