

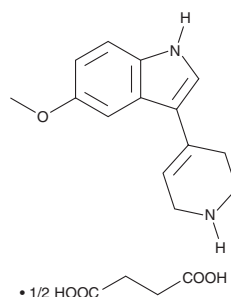
PRODUCT INFORMATION



RU-24969 (succinate)

Item No. 33142

CAS Registry No.: 66611-27-6
Formal Name: 5-methoxy-3-(1,2,3,6-tetrahydro-4-pyridinyl)-1H-indole, butanedioic acid (2:1)
MF: C₁₄H₁₆N₂O • 1/2C₄H₆O₄
FW: 287.3
Purity: ≥95%
UV/Vis.: λ_{max}: 223, 264 nm
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

RU-24969 (succinate) is supplied as a crystalline solid. A stock solution may be made by dissolving the RU-24969 (succinate) in the solvent of choice, which should be purged with an inert gas. RU-24969 (succinate) is slightly soluble in ethanol, DMSO, and dimethyl formamide.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of RU-24969 (succinate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of RU-24969 (succinate) in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

RU-24969 is an agonist of the serotonin (5-HT) receptor subtypes 5-HT_{1A} and 5-HT_{1B}.¹ It binds to 5-HT_{1A} and 5-HT_{1B} receptors (IC₅₀s = 9.5 and 6.6 nM, respectively) and is selective for these receptors over 5-HT₂ receptors in cell-free assays (IC₅₀ = 5,120 nM). RU-24969 inhibits tritium efflux induced by the 5-HT antagonist methiothepin (Item No. 23138) in tritium-preloaded rat frontal cortex slices (pA₂ = 6.27). *In vivo*, RU-24969 (3 mg/kg) potentiates cocaine-induced increases in nucleus accumbens dopamine levels, as well as the reinforcing effects of cocaine self-administration, in rats.²

References

1. Middlemiss, D.N. The putative 5-HT₁ receptor agonist, RU 24969, inhibits the efflux of 5-hydroxytryptamine from rat frontal cortex slices by stimulation of the 5-HT autoreceptor. *J. Pharm. Pharmacol.* **37(6)**, 434-437 (1985).
2. Parsons, L.H., Koob, G.F., and Weiss, F. RU 24969, a 5-HT_{1B/1A} receptor agonist, potentiates cocaine-induced increases in nucleus accumbens dopamine. *Synapse* **32(2)**, 132-135 (1999).

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, 10/25/2022

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