

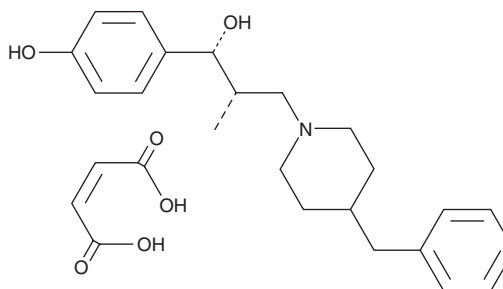
PRODUCT INFORMATION



Ro 25-6981 (maleate)

Item No. 14578

CAS Registry No.: 1312991-76-6
Formal Name: αR-(4-hydroxyphenyl)-βS-methyl-4-(phenylmethyl)-1-piperidinepropanol, 2Z-butenedioate
MF: C₂₂H₂₉NO₂ • C₄H₄O₄
FW: 455.6
Purity: ≥95%
Supplied as: A 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

Ro 25-6981 (maleate) is supplied as a solid. A stock solution may be made by dissolving the Ro 25-6981 (maleate) in the solvent of choice. Ro 25-6981 (maleate) is soluble in the organic solvent DMSO, which should be purged with an inert gas. It is also soluble in water. The solubility of Ro 25-6981 (maleate) in DMSO and water is approximately 100 and 25 mM, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

Ro 25-6981 (maleate) is a potent, selective activity-dependent blocker of NMDA receptors containing the NR2B subunit (IC_{50S} = 9 nM and 52 μM for cloned receptor subunit combinations NR1C/NR2B and NR1C/NR2A, respectively).¹ Ro 25-6981 (maleate) displays neuroprotectant effects *in vitro*, protecting cortical neurons against glutamate toxicity and combined oxygen and glucose deprivation with IC₅₀ values of 0.4 and 0.04 μM, respectively.¹

Reference

1. Fischer, G., Mutel, V., Trube, G., *et al.* Ro 25-6981, a highly potent and selective blocker of N-methyl-D-aspartate receptors containing the NR2B subunit. Characterization *in vitro*. *J. Pharmacol. Exp. Ther.* **283**(3), 1285-1292 (1997).

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, 01/16/2023

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