

PRODUCT INFORMATION



RBI 257 (maleate)

Item No. 17223

CAS Registry No.: 911378-38-6
Formal Name: N-[1-[(4-iodophenyl)methyl]-4-piperidiny]-N-methyl-3-(1-methylethoxy)-2-pyridinamine (2Z)-2-butenedioate

MF: C₂₁H₂₈IN₃O • C₄H₄O₄

FW: 581.4

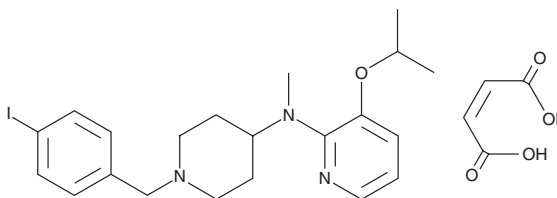
Purity: ≥95%

UV/Vis.: λ_{max}: 240, 295 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

RBI 257 (maleate) is supplied as a crystalline solid. A stock solution may be made by dissolving the RBI 257 (maleate) in the solvent of choice, which should be purged with an inert gas. RBI 257 (maleate) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of RBI 257 (maleate) in these solvents is approximately 30 mg/ml.

RBI 257 (maleate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, RBI 257 (maleate) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. RBI 257 (maleate) 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

RBI 257 is a potent dopamine D₄ receptor antagonist (K_i = 0.3 nM) that is selective over D_{2L}, D₃, D₂-like, and D₁ receptors by 1,721-, 439-, 1,336-, and 8,576-fold, respectively.¹

Reference

1. Kula, N.S., Baldessarini, R.J., Keabian, J.W., *et al.* RBI-257: A highly potent dopamine D₄ receptor-selective ligand. *Eur. J. Pharmacol.* **331**, 333-336 (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, 11/10/2022

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