

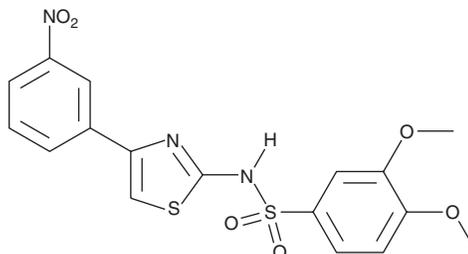
PRODUCT INFORMATION



Ro 61-8048

Item No. 15199

CAS Registry No.: 199666-03-0
Formal Name: 3,4-dimethoxy-N-[4-(3-nitrophenyl)-2-thiazolyl]-benzenesulfonamide
MF: C₁₇H₁₅N₃O₆S₂
FW: 421.4
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 249, 278 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

Ro 61-8048 is supplied as a crystalline solid. A stock solution may be made by dissolving the Ro 61-8048 in the solvent of choice, which should be purged with an inert gas. Ro 61-8048 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of Ro 61-8048 in these solvents is approximately 0.5, 14, and 25 mg/ml, respectively.

Ro 61-8048 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Ro 61-8048 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Ro 61-8048 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Ro 61-8048 is a potent inhibitor of kynurenine 3-hydroxylase (IC₅₀ = 37 nM), preventing the conversion of kynurenine (Item No. 11305) to 3-hydroxy kynurenine.¹ Ro 61-8048 inhibition of kynurenine 3-hydroxylase reduces ischemia-mediated neuronal damage in tissues.^{1,2} Ro 61-8048 can be used *in vivo* by direct injection.³

References

1. Röver, S., Cesura, A.M., Huguenin, P., *et al.* Synthesis and biochemical evaluation of N-(4-phenylthiazol-2-yl)benzenesulfonamides as high-affinity inhibitors of kynurenine 3-hydroxylase. *J. Med. Chem.* **40**(26), 4378-4385 (1997).
2. Carpenedo, R., Meli, E., Peruginelli, F., *et al.* Kynurenine 3-mono-oxygenase inhibitors attenuate post-ischemic neuronal death in organotypic hippocampal slice cultures. *J. Neurochem.* **82**(6), 1465-1471 (2002).
3. Hamann, M., Sander, S.E., and Richter, A. Effects of the kynurenine 3-hydroxylase inhibitor Ro 61-8048 after intrastriatal injections on the severity of dystonia in the dtsz mutant. *Eur. J. Pharmacol.* **586**(1-3), 156-159 (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, 12/20/2022

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