

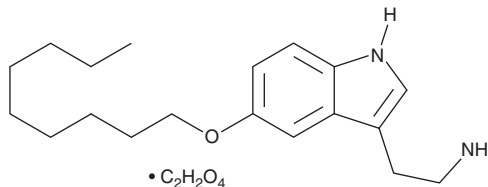
PRODUCT INFORMATION



Nonyloxytryptamine (oxalate)

Item No. 21517

CAS Registry No.: 157798-13-5
Formal Name: 5-(nonyloxy)-1H-indole-3-ethanamine, monoethanedioate
MF: $C_{19}H_{30}N_2O \cdot C_2H_2O_4$
FW: 392.5
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 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

Nonyloxytryptamine (oxalate) is supplied as a crystalline solid. A stock solution may be made by dissolving the nonyloxytryptamine (oxalate) in the solvent of choice. Nonyloxytryptamine (oxalate) is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of nonyloxytryptamine (oxalate) in these solvents is approximately 20 and 30 mg/ml, respectively. Nonyloxytryptamine (oxalate) is also slightly soluble in ethanol.

Nonyloxytryptamine (oxalate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, nonyloxytryptamine (oxalate) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Nonyloxytryptamine (oxalate) has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Nonyloxytryptamine is a potent agonist of the serotonin (5-HT) receptor 5-HT_{1D β} with an EC₅₀ value of 68 nM for adenylate cyclase activity in CHOKM6 cells transfected with the human 5-HT_{1D β} receptor.¹ It is selective for 5-HT_{1D β} over 5-HT_{1A} for its adenylate cyclase activity, with an EC₅₀ greater than 1 μM for 5-HT_{1A}. It binds to 5-HT_{1D β} , 5-HT_{1A}, 5-HT_{1B}, 5-HT_{2A}, and 5-HT_{2C} receptors *in vitro* (K_i s = 1.2-315 nM). Nonyloxytryptamine also inhibits reovirus-mediated cell death in a dose-dependent manner but has no effect on reovirus attachment or internalization.²

References

1. Glennon, R.A., Hong, S.-S., Dukat, M., *et al.* 5-(Nonyloxy)tryptamine: A novel high-affinity 5-HT_{1D β} serotonin receptor agonist. *J. Med. Chem.* **37**(18), 2828-2830 (1994).
2. Mainou, B.A., Ashbrook, A.W., Smith, E.C., *et al.* Serotonin receptor agonist 5-nonyloxytryptamine alters the kinetics of reovirus cell entry. *J. Virol.* **89**(17), 8701-8712 (2015).

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, 11/14/2022

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