

# PRODUCT INFORMATION



## Naltrindole (hydrochloride)

Item No. 9000705

**CAS Registry No.:** 111469-81-9  
**Formal Name:** (4bS,8R,8aS,14bR)-7-(cyclopropylmethyl)-5,6,7,8,14,14b-hexahydro-4,8-methanobenzofuro[2,3-a]pyrido[4,3-b]carbazole-1,8a(9H)-diol, monohydrochloride

**MF:** C<sub>26</sub>H<sub>26</sub>N<sub>2</sub>O<sub>3</sub> • HCl  
**FW:** 451.0

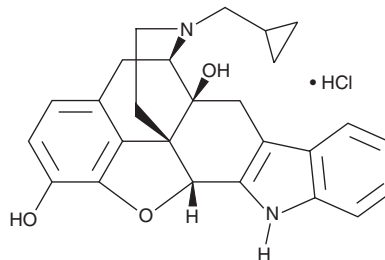
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 210, 284 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

Naltrindole (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the naltrindole (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Naltrindole (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of naltrindole (hydrochloride) in ethanol is approximately 10 mg/ml and approximately 5 mg/ml in DMSO and DMF.

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

### Description

Naltrindole is a potent antagonist of the human  $\delta$ -opioid receptor ( $K_i = 0.02-0.3$  nM), with much lower affinities for  $\kappa$ - and  $\mu$ -opioid receptors ( $K_{iS} = 10-66$  and  $6-64$  nM, respectively).<sup>1,2</sup> Naltrindole is commonly used to investigate the role of the  $\delta$ -opioid receptor in signaling responses to test compounds.<sup>3-5</sup>

### References

1. Meng, F., Wei, Q., Hoversten, M.T., *et al.* Switching agonist/antagonist properties of opiate alkaloids at the  $\delta$  opioid receptor using mutations based on the structure of the orphanin FQ receptor. *J. Biol. Chem.* **275(29)**, 21939-21945 (2000).
2. Raynor, K., Kong, H., Chen, Y., *et al.* Pharmacological characterization of the cloned  $\kappa$ -,  $\delta$ -, and  $\mu$ -opioid receptors. *Mol. Pharm.* **45(2)**, 330-334 (1994).
3. Steinmiller, C.L. and Young, A.M. Pharmacological selectivity of CTAP in a warm water tail-withdrawal antinociception assay in rats. *Psychopharmacology (Berl)* **195(4)**, 497-507 (2008).
4. Shannon, H.E. and Lutz, E.A. Comparison of the peripheral and central effects of the opioid agonists loperamide and morphine in the formalin test in rats. *Neuropharmacology* **42(2)**, 253-261 (2002).
5. Mendes, G.L., Santos, A.R.S., Malheiros, A., *et al.* Assessment of mechanisms involved in antinociception caused by sesquiterpene polygodial. *J. Pharmacol. Exp. Ther.* **292(1)**, 164-172 (2000).

#### 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