

# PRODUCT INFORMATION



## NTNCB (hydrochloride)

Item No. 43216

**CAS Registry No.:** 191931-56-3  
**Formal Name:** 2-nitro-N-[[trans-4-[[[(1,2,3,4-tetrahydro-2-naphthalenyl)methyl]amino]methyl]cyclohexyl]methyl]-benzenesulfonamide, monohydrochloride

**MF:** C<sub>25</sub>H<sub>33</sub>N<sub>3</sub>O<sub>4</sub>S • HCl

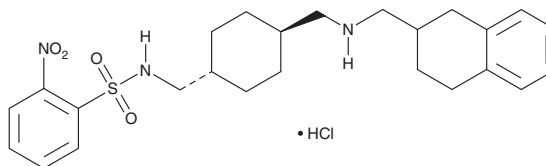
**FW:** 508.1

**Purity:** ≥98%

**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

NTNCB (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the NTNCB (hydrochloride) in the solvent of choice, which should be purged with an inert gas. NTNCB (hydrochloride) is sparingly soluble (1-10 mg/ml) in DMSO.

### Description

NTNCB is an antagonist of neuropeptide Y (NPY) receptor Y<sub>5</sub> (K<sub>i</sub> = 8 nM).<sup>1</sup> It is selective for neuropeptide Y<sub>5</sub> over Y<sub>1</sub> receptors (K<sub>i</sub> = >10,000 nM). NTNCB (10 and 20 μM) suppresses colony formation by PC3 and DU145 prostate cancer cells.<sup>2</sup> *In vivo*, NTNCB (1, 3, and 10 mg/kg per day) reduces tumor volume in a PC3 mouse xenograft model.

### References

1. Islam, I., Dhanoa, D., and Finn, J., *et al.* Discovery of potent and selective small molecule NPY Y<sub>5</sub> receptor antagonists. *Bioorg. Med. Chem. Lett.* **12(13)**, 1767-1769 (2002).
2. Tian, S., Liao, X. and Cao, W., *et al.* GSFM: A genome-scale functional module transformation to represent drug efficacy for *in silico* drug discovery. *Acta Pharm. Sin. B.* **15(1)**, 133-150 (2025).

#### 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, 04/14/2025

#### 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