

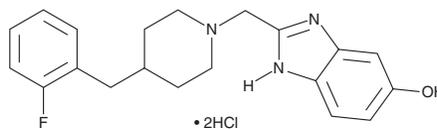
# PRODUCT INFORMATION



## TCN 237 (hydrochloride)

Item No. 42266

**CAS Registry No.:** 1784252-90-9  
**Formal Name:** 2-[[4-[(2-fluorophenyl)methyl]-1-piperidinyl]methyl]-1H-benzimidazol-6-ol, dihydrochloride  
**MF:** C<sub>20</sub>H<sub>22</sub>FN<sub>3</sub>O • 2HCl  
**FW:** 412.3  
**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

TCN 237 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the TCN 237 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. TCN 237 (hydrochloride) is soluble (≥10 mg/ml) in DMSO.

### Description

TCN 237 is an antagonist of NR2B subunit-containing NMDA receptors ( $K_i = 0.85 \text{ nM}$ ).<sup>1</sup> It selectively inhibits glutamate and glycine-induced currents in whole-cell patch-clamp assays using L-M(TK-) cells expressing NR2B subunit-containing NMDA receptors over L-M(TK-) cells expressing NR2A- or NR2D subunit-containing NMDA receptors and HEK293 cells expressing NR2C subunit-containing NMDA receptors. TCN 237 (3 mg/kg) inhibits carrageenan-induced mechanical hyperalgesia in rats.

### Reference

1. McCauley, J.A., Theberge, C.R., Romano, J.J., *et al.* NR2B-selective N-methyl-D-aspartate antagonists: Synthesis and evaluation of 5-substituted benzimidazoles. *J. Med. Chem.* **47**, 2089-2096 (2004).

#### 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, 12/12/2024

#### 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