

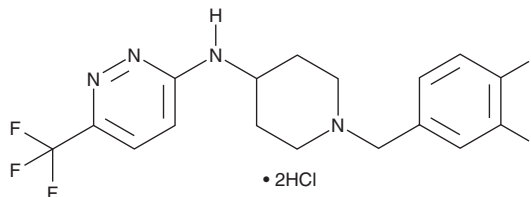
# PRODUCT INFORMATION



## JNJ-37822681 (hydrochloride)

Item No. 37428

**CAS Registry No.:** 2108806-02-4  
**Formal Name:** N-[1-[(3,4-difluorophenyl)methyl]-4-piperidinyl]-6-(trifluoromethyl)-3-pyridazinamine, dihydrochloride  
**MF:**  $C_{17}H_{17}F_5N_4 \cdot 2HCl$   
**FW:** 445.3  
**Purity:**  $\geq 98\%$   
**Supplied as:** A solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

JNJ-37822681 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the JNJ-37822681 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. JNJ-37822681 (hydrochloride) is soluble in the organic solvent DMSO. JNJ-37822681 (hydrochloride) is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

JNJ-37822681 is a dopamine  $D_{2L}$  receptor antagonist ( $K_i = 0.158 \mu M$ ).<sup>1</sup> It is selective for dopamine  $D_{2L}$  receptors over dopamine  $D_1$  and  $D_3$ ,  $\alpha_1$ -adrenergic, and histamine  $H_1$  receptors, and the serotonin (5-HT) receptor subtypes 5-HT<sub>2A</sub> and 5-HT<sub>2C</sub> in radioligand binding assays ( $K_i$ s =  $>3$ , 1.159,  $>5$ , 4.931, 2.896, and  $>4 \mu M$ , respectively). JNJ-37822681 inhibits apomorphine-induced stereotypy and D-amphetamine-induced hyperlocomotion in rat models of psychosis ( $ED_{50}$ s = 0.19 and 1 mg/kg, respectively).

### Reference

1. Langlois, X., Megens, A., Lavreysen, H., *et al.* Pharmacology of JNJ-37822681, a specific and fast-dissociating  $D_2$  antagonist for the treatment of schizophrenia. *J. Pharmacol. Exp. Ther.* **342**(1), 91-105 (2012).

#### 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, 10/25/2022

#### 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