

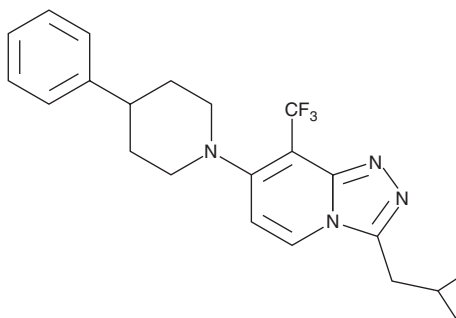
# PRODUCT INFORMATION



**JNJ-42153605**

Item No. 21984

**CAS Registry No.:** 1254977-87-1  
**Formal Name:** 3-(cyclopropylmethyl)-7-(4-phenyl-1-piperidinyl)-8-(trifluoromethyl)-1,2,4-triazolo[4,3-a]pyridine  
**MF:** C<sub>22</sub>H<sub>23</sub>F<sub>3</sub>N<sub>4</sub>  
**FW:** 400.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 260, 336 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

JNJ-42153605 is supplied as a crystalline solid. A stock solution may be made by dissolving the JNJ-42153605 in the solvent of choice, which should be purged with an inert gas. JNJ-42153605 is soluble in the organic solvent dimethyl formamide (DMF) at a concentration of approximately 30 mg/ml.

JNJ-42153605 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JNJ-42153605 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. JNJ-42153605 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

JNJ-42153605 is a positive allosteric modulator of metabotropic glutamate receptor 2 (mGluR2; EC<sub>50</sub> = 17 nM in CHO cells expressing the human receptor).<sup>1</sup> *In vivo*, JNJ-42153605 (3 mg/kg) inhibits mGluR2-mediated rapid eye movement (REM) sleep in rats. It also reverses phencyclidine-induced hyperlocomotion in mice (ED<sub>50</sub> = 5.4 mg/kg).

## Reference

1. Cid, J.M., Tresadern, G., Vega, J.A., *et al.* Discovery of 3-cyclopropylmethyl-7-(4-phenylpiperidin-1-yl)-8-trifluoromethyl[1,2,4]triazolo[4,3-a]pyridine (JNJ-42153605): A positive allosteric modulator of the metabotropic glutamate 2 receptor. *J. Med. Chem.* **55**(20), 8770-8789 (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, 10/04/2022

## 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