

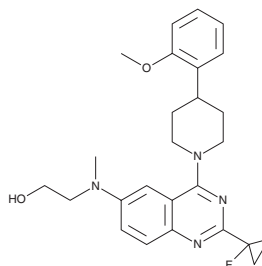
# PRODUCT INFORMATION



## SBI-553

Item No. 36782

**CAS Registry No.:** 1849603-72-0  
**Formal Name:** 2-[[2-(1-fluorocyclopropyl)-4-[4-(2-methoxyphenyl)-1-piperidinyl]-6-quinazoliny]methylamino]-ethanol  
**MF:** C<sub>26</sub>H<sub>31</sub>FN<sub>4</sub>O<sub>2</sub>  
**FW:** 450.6  
**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

SBI-553 is supplied as a solid. A stock solution may be made by dissolving the SBI-553 in the solvent of choice, which should be purged with an inert gas. SBI-553 is soluble in the organic solvent dimethyl formamide at a concentration of approximately 3 mg/ml. SBI-553 is slightly soluble in ethanol and DMSO.

### Description

SBI-553 is a  $\beta$ -arrestin-biased positive allosteric modulator (PAM) of neurotensin receptor 1 (NTSR1).<sup>1</sup> It binds to NSTR1 and increases neurotensin-1 affinity for NSTR1 in HEK293 cells expressing the human receptor when used at concentrations ranging from 0.01 to 10  $\mu$ M. SBI-553 (0.03-30  $\mu$ M) induces  $\beta$ -arrestin recruitment and NTSR1 internalization without stimulating Gq protein activation, 1,4,5-triphosphate (IP<sub>3</sub>) generation, or calcium mobilization in cell-based assays. *In vivo*, SBI-553 (12 mg/kg, i.p.) attenuates cocaine- or methamphetamine-induced hyperlocomotion in wild-type but not  $\beta$ -arrestin-2 knockout mice. It also reduces cocaine self-administration in mice.

### Reference

1. Slosky, L.M., Bai, Y., Toth, K., *et al.*  $\beta$ -Arrestin-biased allosteric modulator of NTSR1 selectively attenuates addictive behaviors. *Cell* **181**(6), 1364-1379.e14 (2018).

#### 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, 09/30/2022

#### 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