

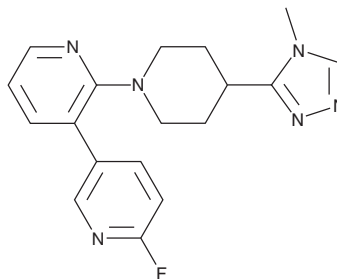
# PRODUCT INFORMATION



## SEN-177

Item No. 40449

CAS Registry No.: 2117405-13-5  
Formal Name: 6'-fluoro-2-[4-(4-methyl-4H-1,2,4-triazol-3-yl)-1-piperidinyl]-3,3'-bipyridine  
MF:  $C_{18}H_{19}FN_6$   
FW: 338.4  
Purity:  $\geq 98\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

SEN-177 is supplied as a solid. A stock solution may be made by dissolving the SEN-177 in the solvent of choice, which should be purged with an inert gas. SEN-177 is soluble in methanol, acetonitrile, and DMSO.

### Description

SEN-177 is a glutamyl cyclase inhibitor ( $IC_{50} = 53 \text{ nM}$ ).<sup>1</sup> It reduces polyglutamine aggregation in HeLa cells expressing EGFP-HTT-Q74, an *in vitro* model of Huntington's disease when used at concentrations of 12.5 and 25  $\mu\text{M}$ .<sup>1</sup> SEN-177 (50  $\mu\text{M}$ ) reduces polyglutamine aggregation in primary mouse neurons expressing EGFP-Q80.

### Reference

1. Jimenez-Sanchez, M., Lam, W., Hannus, M., et al. siRNA screen identifies QPCT as a druggable target for Huntington's disease. *Nat. Chem. Biol.* **11**(5), 347-354 (2015).

#### 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, 03/06/2024

#### 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