

# PRODUCT INFORMATION



**SUN11602**

Item No. 36572

**CAS Registry No.:** 704869-38-5  
**Formal Name:** 4-[[4-[[2-[[4-amino-2,3,5,6-tetramethylphenyl)amino]acetyl]methylamino]-1-piperidinyl]methyl]-benzamide

**MF:** C<sub>26</sub>H<sub>37</sub>N<sub>5</sub>O<sub>2</sub>

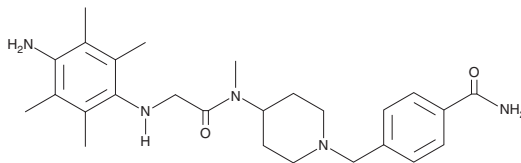
**FW:** 451.6

**Purity:** ≥98%

**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

SUN11602 is supplied as a crystalline solid. A stock solution may be made by dissolving the SUN11602 in the solvent of choice, which should be purged with an inert gas. SUN11602 is soluble in DMSO.

## Description

SUN11602 is a mimetic of FGF2, also known as basic FGF (bFGF).<sup>1</sup> It protects against glutamate-induced toxicity in primary rat cortical neurons when used at concentrations of 0.1, 0.3, and 1 μM, an effect that can be reversed by the FGFR1 inhibitor PD 166866 (Item No. 22464). SUN11602 (1 μM) stimulates neurite outgrowth in primary rat hippocampal neurons.<sup>2</sup> It reduces cognitive deficits induced by injection of aggregated amyloid-β (1-40) (Aβ40) and ibotenic acid (Item No. 14584) into the hippocampus in a rat model of Alzheimer's disease when administered at a dose of 10 mg/kg.

## References

1. Murayama, N., Kadoshima, T., Takemoto, N., *et al.* SUN11602, a novel aniline compound, mimics the neuroprotective mechanisms of basic fibroblast growth factor. *ACS Chem. Neurosci.* **4**(2), 266-276 (2012).
2. Ogino, R., Murayama, N., Noshita, T., *et al.* SUN11602 has basic fibroblast growth factor-like activity and attenuates neuronal damage and cognitive deficits in a rat model of Alzheimer's disease induced by amyloid β and excitatory amino acids. *Brain Res.* **1585**, 159-166 (2014).

### 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, 11/21/2022

## 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