

# PRODUCT INFORMATION



**PF-3084014**

Item No. 27061

**CAS Registry No.:** 1290543-63-3  
**Formal Name:** (2S)-2-[[[(2S)-6,8-difluoro-1,2,3,4-tetrahydro-2-naphthalenyl]amino]-N-[1-[2-[(2,2-dimethylpropyl)amino]-1,1-dimethylethyl]-1H-imidazol-4-yl]-pentanamide

**MF:** C<sub>27</sub>H<sub>41</sub>F<sub>2</sub>N<sub>5</sub>O

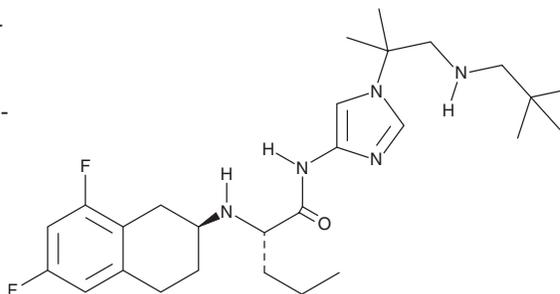
**FW:** 489.6

**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

PF-3084014 is supplied as a solid. A stock solution may be made by dissolving the PF-3084014 in the solvent of choice, which should be purged with an inert gas. PF-3084014 is soluble in the organic solvent DMSO.

## Description

PF-3084014 is a  $\gamma$ -secretase inhibitor (IC<sub>50</sub>s = 6.2 and 1.2 nM in an enzyme assay and in whole cells, respectively).<sup>1</sup> It inhibits  $\gamma$ -secretase activity without inducing Notch-related toxicity (EC<sub>50</sub>s = 1,300-3,000 nM for B cells in fetal thymic organ cultures). It reduces the level of amyloid- $\beta$  (1-40) (A $\beta$ 40) fragments in mouse brain three hours following administration of 1 to 18 mg/kg doses and reduces A $\beta$ 42 fragment levels when administered at doses of 9 or 18 mg/kg. It also reduces A $\beta$  fragments in guinea pig brain when administered at doses ranging from 0.3 to 3 mg/kg per day for five days.

## Reference

1. Lanz, T.A., Wood, K.M., Richter, K.E.G., *et al.* Pharmacodynamics and pharmacokinetics of the  $\gamma$ -secretase inhibitor PF-3084014. *J. Pharmacol. Exp. Ther.* **334**(1), 269-277 (2010).

### 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