

# PRODUCT INFORMATION

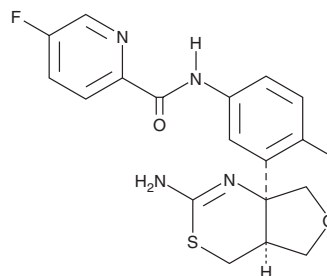


**LY2886721**

Item No. 21599

**CAS Registry No.:** 1262036-50-9  
**Formal Name:** N-[3-[(4aS,7aS)-2-amino-4a,5-dihydro-4H-furo[3,4-d][1,3]thiazin-7a(7H)-yl]-4-fluorophenyl]-5-fluoro-2-pyridinecarboxamide

**Synonym:**  $\beta$ -Secretase  
**MF:** C<sub>18</sub>H<sub>16</sub>F<sub>2</sub>N<sub>4</sub>O<sub>2</sub>S  
**FW:** 390.4  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 279 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

## Laboratory Procedures

LY2886721 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY2886721 in the solvent of choice, which should be purged with an inert gas. LY2886721 is soluble in organic solvents such as ethanol and DMSO. It is also soluble in water. The solubility of LY2886721 in ethanol and water is approximately 1 mg/ml and approximately 9 mg/ml in DMSO. We do not recommend storing the aqueous solution for more than one day.

## Description

LY2886721 selectively inhibits the  $\beta$ -site amyloid protein cleaving enzyme (BACE) ( $IC_{50}$  = 20 nM for recombinant hBACE1).<sup>1</sup> BACE performs the first cleavage step in the A $\beta$  cascade; thus, its inhibition by LY2886721 decreases production of all forms of the A $\beta$  peptide.<sup>2</sup> A Phase II clinical study for the use of LY2886721 in the treatment of Alzheimer's disease was terminated early due to abnormal liver biochemistry values in several patients. The toxic effects did not appear to be due to BACE inhibition.

## References

1. May, P.C., Boggs, L.N., Brier, R., *et al.* Preclinical characterization of LY2886721: A BACE1 inhibitor in clinical development for early Alzheimer's disease. *J. Alzheim. Assoc.* **8(4)**, 95 (2003).
2. Vassar, R., Kovacs, D.M., Yan, R., *et al.* The beta-secretase enzyme BACE in health and Alzheimer's disease: regulation, cell biology, function, and therapeutic potential. *J. Neurosci.* **29(41)**, 12787-12794 (2009).

### 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, 12/15/2022

## 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