

# PRODUCT INFORMATION

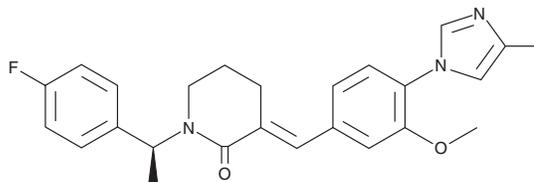


**E2012**

Item No. 29710

**CAS Registry No.:** 870843-42-8  
**Formal Name:** (3E)-1-[(1S)-1-(4-fluorophenyl)ethyl]-3-[[3-methoxy-4-(4-methyl-1H-imidazol-1-yl)phenyl]methylene]-2-piperidinone

**MF:** C<sub>25</sub>H<sub>26</sub>FN<sub>3</sub>O<sub>2</sub>  
**FW:** 419.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 286 nm  
**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

E2012 is supplied as a solid. A stock solution may be made by dissolving the E2012 in the solvent of choice, which should be purged with an inert gas. E2012 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of E2012 in these solvents is approximately 10 mg/ml.

E2012 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, E2012 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. E2012 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

E2012 is a modulator of  $\gamma$ -secretase.<sup>1</sup> It selectively inhibits APP intracellular signaling domain (AICD) cleavage to amyloid- $\beta$  (A $\beta$ ) over Notch cleavage to its signaling effector Notch intracellular domain (NICD), processes both mediated by  $\gamma$ -secretase, when used at a concentration of 1  $\mu$ M in luciferase assays.<sup>2</sup> E2012 also inhibits the activity of the cholesterol synthesis enzyme 3 $\beta$ -hydroxysterol  $\Delta$ <sup>24</sup>-reductase (DHCR24) in primary rat hepatocytes and HepG2 cells (IC<sub>50</sub>s = 11 and 15 nM, respectively).<sup>3</sup> It reduces A $\beta$  (1-42) (A $\beta$ 42) production in primary rat embryonic cerebral cortex neurons with an IC<sub>50</sub> value of 220 nM.<sup>4</sup> E2012 (100 mg/kg) also decreases A $\beta$ 42 levels in guinea pig brain and cerebral spinal fluid (CSF).<sup>5</sup>

## References

1. Hahn, S., Brüning, T., Ness, J., *et al.* Presenilin-1 but not amyloid precursor protein mutations present in mouse models of Alzheimer's disease attenuate the response of cultured cells to  $\gamma$ -secretase modulators regardless of their potency and structure. *J. Neurochem.* **116**(3), 385-395 (2011).
2. Dimitrov, M., Alattia, J.-R., Lemmin, T., *et al.* Alzheimer's disease mutations in APP but not  $\gamma$ -secretase modulators affect epsilon-cleavage-dependent AICD production. *Nat. Commun.* **4**:2246 (2013).
3. Nakano-Ito, K., Fujikawa, Y., Hihara, T., *et al.* E2012-induced cataract and its predictive biomarkers. *Toxicol. Sci.* **137**(1), 249-258 (2014).
4. Kimura, T., Kawano, K., Doi, E., *et al.* Cinnamide compound. *Eisai Co., Ltd.* **WO2005/115990A1** (2005).
5. Lu, Y., Riddell, D.R., Hajos-Korcsok, E., *et al.* Cerebrospinal fluid amyloid- $\beta$  (A $\beta$ ) as an effect biomarker for brain A $\beta$  lowering verified by quantitative preclinical analyses. *J. Pharmacol. Exp. Ther.* **342**(2), 366-375 (2012).

### 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