

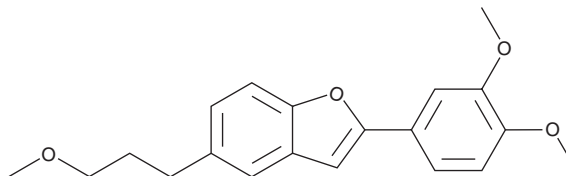
# PRODUCT INFORMATION



## DWK-1339

Item No. 27642

**CAS Registry No.:** 1018946-38-7  
**Formal Name:** 2-(3,4-dimethoxyphenyl)-5-(3-methoxypropyl)-benzofuran  
**MF:** C<sub>20</sub>H<sub>22</sub>O<sub>4</sub>  
**FW:** 326.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 214, 315, 329 nm  
**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

DWK-1339 is supplied as a crystalline solid. A stock solution may be made by dissolving the DWK-1339 in the solvent of choice, which should be purged with an inert gas. DWK-1339 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of DWK-1339 in ethanol is approximately 2 mg/ml and approximately 30 mg/ml in DMSO and DMF.

### Description

DWK-1339 is an inhibitor of amyloid- $\beta$  (A $\beta$ ) aggregation.<sup>1</sup> It inhibits aggregation of monomeric A $\beta$  (1-42) (A $\beta$ 42) and induces disaggregation of A $\beta$ 42 fibrils *in vitro* when used at concentrations ranging from 3.1 to 50  $\mu$ M. DWK-1339 (10  $\mu$ M) reduces A $\beta$ 42-induced toxicity in HT-22 cells. *In vivo*, DWK-1339 (10 mg/kg) increases step-through latency in a passive avoidance test and increases spontaneous alteration in the Y-maze compared with vehicle control mice in a mouse model of A $\beta$ 42-induced acute Alzheimer's disease. It also decreases brain levels of A $\beta$ 42 and increases spontaneous alteration in the Y-maze in the APP/PS1 transgenic mouse model of Alzheimer's disease.

### Reference

1. Ha, H., Kang, D.W., Kim, H.-M., *et al.* Discovery of an orally bioavailable benzofuran analogue that serves as a  $\beta$ -amyloid aggregation inhibitor for the potential treatment of Alzheimer's disease. *J. Med. Chem.* **61**(1), 396-402 (2018).

**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, 12/22/2022

### 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