

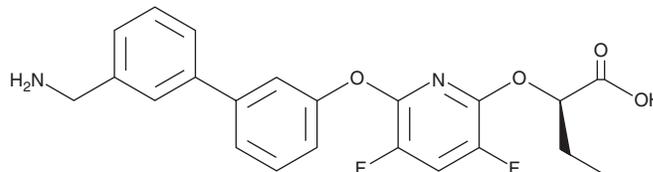
# PRODUCT INFORMATION



**ZK 824859**

Item No. 37789

**CAS Registry No.:** 2254001-81-3  
**Formal Name:** 2-[[6-[[3'-(aminomethyl)[1,1'-biphenyl]-3-yl]oxy]-3,5-difluoro-2-pyridinyl]oxy]-butanoic acid  
**MF:** C<sub>22</sub>H<sub>20</sub>F<sub>2</sub>N<sub>2</sub>O<sub>4</sub>  
**FW:** 414.4  
**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

ZK 824859 is supplied as a solid. A stock solution may be made by dissolving the ZK 824859 in the solvent of choice, which should be purged with an inert gas. ZK 824859 is soluble in methanol and DMSO.

## Description

ZK 824859 is an inhibitor of urokinase-type plasminogen activator (uPA; IC<sub>50</sub>s = 79 and 410 nM for the human and mouse proteins, respectively).<sup>1</sup> It is selective for human uPA over human tissue-type plasminogen activator (tPA) and plasmin (IC<sub>50</sub>s = 1,580 and 1,330 nM, respectively) but not selective for mouse uPA over mouse tPA and plasmin (IC<sub>50</sub>s = 910 and 1,600 nM, respectively). ZK 824859 (25 or 50 mg/kg per day) reduces the incidence of limb weakness and paralysis in mice in a model of experimental autoimmune encephalomyelitis (EAE).

## Reference

1. Islam, I., Yuan, S., West, C.W., et al. Discovery of selective urokinase plasminogen activator (uPA) inhibitors as a potential treatment for multiple sclerosis. *Bioorg. Med. Chem. Lett.* **28**(20), 3372-3375 (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, 03/06/2024

## 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