

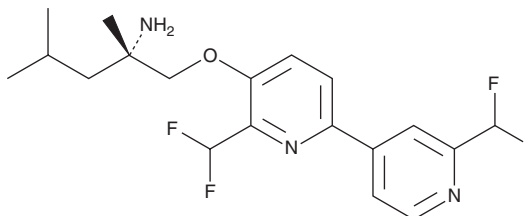
# PRODUCT INFORMATION



**BMS 986176**

Item No. 43668

**CAS Registry No.:** 1815613-42-3  
**Formal Name:** (2S)-1-[[2',6-bis(difluoromethyl)-[2,4'-bipyridin]-5-yl]oxy]-2,4-dimethyl-2-pentanamine  
**Synonym:** LX9211  
**MF:** C<sub>19</sub>H<sub>23</sub>F<sub>4</sub>N<sub>3</sub>O  
**FW:** 385.4  
**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

BMS 986176 is supplied as a solid. A stock solution may be made by dissolving the BMS 986176 in the solvent of choice, which should be purged with an inert gas. BMS 986176 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

## Description

BMS 986176 is an inhibitor of AP2-associated protein kinase 1 (AAK1; IC<sub>50</sub> = 2 nM).<sup>1</sup> It increases the latency to paw withdrawal in a rat model of diabetic neuropathic pain induced by streptozotocin (Item No. 13104) when administered at doses of 0.3 or 3 mg/kg.

## Reference

1. Luo, G., Chen, L., Kostich, W.A., *et al.* Discovery of (S)-1-((2',6-bis(difluoromethyl)-[2,4'-bipyridin]-5-yl)oxy)-2,4-dimethylpentan-2-amine (BMS-986176/LX-9211): A highly selective, CNS penetrable, and orally active adaptor protein-2 associated kinase 1 inhibitor in clinical trials for the treatment of neuropathic pain. *J. Med. Chem.* **65**(6), 4457-4480 (2022).

### 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, 06/26/2025

## 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](http://WWW.CAYMANCHEM.COM)