

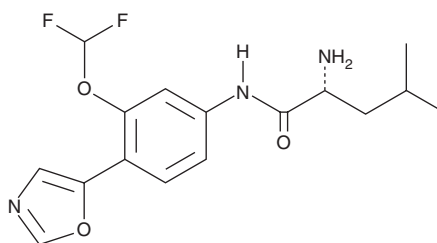
# PRODUCT INFORMATION



## BMS 911172

Item No. 45308

**CAS Registry No.:** 1644248-18-9  
**Formal Name:** 2R-amino-N-[3-(difluoromethoxy)-4-(5-oxazolyl)phenyl]-4-methylpentanamide  
**MF:** C<sub>16</sub>H<sub>19</sub>F<sub>2</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 339.3  
**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 911172 is supplied as a solid. A stock solution may be made by dissolving the BMS 911172 in the solvent of choice, which should be purged with an inert gas. BMS 911172 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

### Description

BMS 911172 is an AP2-associated kinase 1 (AAK1) inhibitor that reduces the phosphorylation of AP-2 complex subunit mu2 (AP2M2) in HEK293 cells overexpressing AAK1 and AP2M2 (IC<sub>50</sub> = 0.051 μM).<sup>1</sup> It is selective for AAK1 over a panel of 219 kinases, including BMP2-inducible kinase (BIKE), cyclin G-associated kinase (GAK), and serine/threonine kinase 16 (STK16; IC<sub>50</sub>s = 0.120, 1.06, and 15.8 μM, respectively). BMS 911172 also inhibits AP2M2 phosphorylation in mouse brain lysates (ED<sub>50</sub> = 6 mg/kg). *In vivo*, BMS 911172 (30 mg/kg) reduces formalin-induced paw flinching and biting in a rat model of persistent pain. It also reduces thermal and mechanical hyperalgesia, as well as cold and mechanical allodynia, in a rat model of chronic constriction injury when administered at a dose of 60 mg/kg.

### Reference

1. Hartz, R.A., Ahuja, V.T., Nara, S.J., *et al.* Discovery, structure-activity relationships, and *in vivo* evaluation of novel aryl amides as brain penetrant adaptor protein 2-associated kinase 1 (AAK1) inhibitors for the treatment of neuropathic pain. *J. Med. Chem.* **64**(15), (2021).

**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, 03/12/2026

### 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