

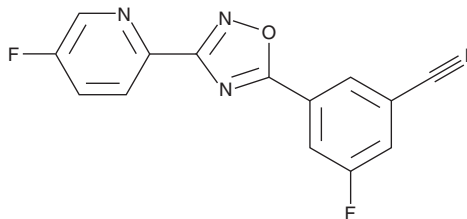
# PRODUCT INFORMATION



**AZD 9272**

Item No. 29717

CAS Registry No.: 327056-26-8  
Formal Name: 3-fluoro-5-[3-(5-fluoro-2-pyridinyl)-1,2,4-oxadiazol-5-yl]-benzonitrile  
MF:  $C_{14}H_6F_2N_4O$   
FW: 284.2  
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

## Laboratory Procedures

AZD 9272 is supplied as a solid. A stock solution may be made by dissolving the AZD 9272 in the solvent of choice, which should be purged with an inert gas. AZD 9272 is soluble in the organic solvent DMSO at a concentration of approximately 50 mM.

## Description

AZD 9272 is a brain-penetrant negative allosteric modulator of metabotropic glutamate receptor 5 (mGluR5;  $IC_{50} = 7.6$  nM in a FLIPR calcium mobilization assay).<sup>1</sup> It is selective for mGluR5 over mGluR1-4 and 6-8 in a FLIPR calcium mobilization assay ( $IC_{50}s = >30$   $\mu\text{M}$ ) and a panel of 134 receptors, ion channels, nuclear hormone receptors, transporters, and enzymes at 10  $\mu\text{M}$  in radioligand binding assays.

## Reference

1. Raboisson, P., Breitholtz-Emanuelsson, A., Dahllöf, H., *et al.* Discovery and characterization of AZD9272 and AZD6538—Two novel mGluR5 negative allosteric modulators selected for clinical development. *Bioorg. Med. Chem. Lett.* **22**(22), 6974-6979 (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, 10/07/2025

## 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