

# PRODUCT INFORMATION



**AZ 13705339**

Item No. 42087

**CAS Registry No.:** 2016806-57-6  
**Formal Name:** 2-[[[2-[[3-(ethylsulfonyl)-4-(4-methyl-1-piperazinyl)phenyl]amino]-5-fluoro-4-pyrimidinyl][5-(hydroxymethyl)-2-methylphenyl]amino]methyl]-benzonitrile

**MF:** C<sub>33</sub>H<sub>36</sub>FN<sub>7</sub>O<sub>3</sub>S

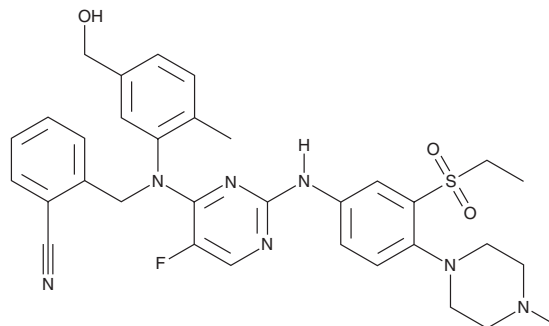
**FW:** 629.8

**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

AZ 13705339 is supplied as a solid. A stock solution may be made by dissolving the AZ 13705339 in the solvent of choice, which should be purged with an inert gas. AZ 13705339 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

## Description

AZ 13705339 is an inhibitor of p21-activated kinase 1 (PAK1; IC<sub>50</sub> = 0.33 nM).<sup>1</sup> It is selective for PAK1 over a panel of 117 kinases at 100 nM. AZ 13705339 binds to PAK1 and PAK2 (K<sub>d</sub> = 0.28 and 0.32 nM, respectively). It inhibits *S. enterica* infection in HAP1 chronic myeloid leukemia cells when used at a concentration of 10 μM.<sup>2</sup> AZ 13705339 inhibits cell death induced by an antibody targeting sialic acid binding Ig-like lectin 8 (Siglec-8) in IL-5-stimulated primary human eosinophils (IC<sub>50</sub> = 59 nM).<sup>3</sup>

## References

1. McCoull, W., Hennessy, E.J., Blades, K., *et al.* Optimization of highly kinase selective bis-anilino pyrimidine PAK1 inhibitors. *ACS Med. Chem. Lett.* **7**(12), 1118-1123 (2016).
2. Davidson, A., Tyler, J., Hume, P., *et al.* A kinase-independent function of PAK is crucial for pathogen-mediated actin remodelling. *PLoS Pathog.* **17**(8), e1009902 (2021).
3. Carroll, D.J., Cao, Y., Bochner, B.S., *et al.* Siglec-8 signals through a non-canonical pathway to cause human eosinophil death *in vitro*. *Front. Immunol.* **12**, 737988 (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, 12/04/2024

## 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