

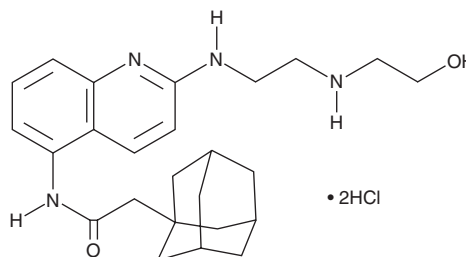
# PRODUCT INFORMATION



## AZ 10606120 (hydrochloride)

Item No. 35820

**CAS Registry No.:** 607378-18-7  
**Formal Name:** N-[2-[[2-[(2-hydroxyethyl)amino]ethyl]amino]-5-quinoliny]-tricyclo[3.3.1.1<sup>3,7</sup>]decane-1-acetamide, dihydrochloride  
**MF:** C<sub>25</sub>H<sub>34</sub>N<sub>4</sub>O<sub>2</sub> • 2HCl  
**FW:** 495.5  
**Purity:** ≥95%  
**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 10606120 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the AZ 10606120 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. AZ 10606120 (hydrochloride) is soluble in DMSO. AZ 10606120 (hydrochloride) is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

AZ 10606120 is an antagonist of the purinergic P2X<sub>7</sub> receptor (IC<sub>50</sub> = 10 nM).<sup>1</sup> It reduces cell migration and proliferation of PANC-1 pancreatic ductal adenocarcinoma cells when used at a concentration of 10 μM.<sup>2</sup> AZ 10606120 (0.01 μg/kg) reduces cecal slurry-induced impairment of endothelial-dependent relaxation but increases spleen bacterial burden in a mouse model of polymicrobial sepsis.<sup>3</sup>

### References

1. Guile, S.D., Alcaraz, L., Birkinshaw, T.N., *et al.* Antagonists of the P2X<sub>7</sub> receptor. From lead identification to drug development. *J. Med. Chem.* **52**(10), 3123-3141 (2009).
2. Giannuzzo, A., Pedersen, S.F., and Novak, I. The P2X<sub>7</sub> receptor regulates cell survival, migration and invasion of pancreatic ductal adenocarcinoma cells. *Mol. Cancer* **14**, 203 (2015).
3. Meegan, J.E., Komalavilas, P., Cheung-Flynn, J., *et al.* Blocking P2X<sub>7</sub> receptor with AZ 10606120 exacerbates vascular hyperpermeability and inflammation in murine polymicrobial sepsis. *Physiol. Rep.* **10**(11), e15290 (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, 12/15/2022

#### 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