

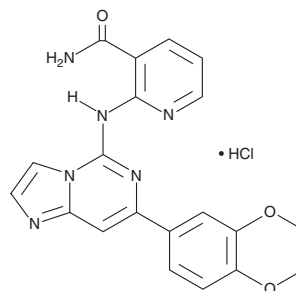
# PRODUCT INFORMATION



## BAY 61-3606 (hydrochloride)

Item No. 11423

**CAS Registry No.:** 1615197-10-8  
**Formal Name:** 2-[[7-(3,4-dimethoxyphenyl)imidazo[1,2-c]pyrimidin-5-yl]amino]-3-pyridinecarboxamide, monohydrochloride  
**MF:** C<sub>20</sub>H<sub>18</sub>N<sub>6</sub>O<sub>3</sub> • HCl  
**FW:** 426.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 331 nm  
**Supplied as:** A crystalline 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

BAY 61-3606 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the BAY 61-3606 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. BAY 61-3606 (hydrochloride) is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

### Description

BAY 61-3606 is an inhibitor of spleen tyrosine kinase (Syk;  $K_i = 7.5$  nM).<sup>1</sup> It is selective for Syk over Lyn, Fyn, Src, Itk, and BTK ( $K_i$ s = >4.7 μM for all). BAY 61-3606 inhibits anti-IgE-induced histamine release in isolated human mast cells sensitized with IgE ( $IC_{50} = 5.1$  nM), as well as reduces anti-IgM-induced proliferation of isolated mouse splenic B cells ( $IC_{50} = 58$  nM). It sensitizes MCF-7 human breast cancer cells to apoptosis induced by TNF-related apoptosis-inducing ligand (TRAIL) when used at concentrations ranging from 0.31 to 2.5 μM.<sup>2</sup> BAY 61-3606 inhibits passive cutaneous anaphylaxis in rats ( $ED_{50} = 8$  mg/kg).<sup>1</sup> It reduces tumor growth in an MCF-7 mouse xenograft model when administered at a dose of 50 mg/kg alone or in combination with TRAIL.<sup>2</sup>

### References

1. Yamamoto, N., Takeshita, K., Shichijo, M., *et al.* The orally available spleen tyrosine kinase inhibitor 2-[7-(3,4-dimethoxyphenyl)-imidazo[1,2-c]pyrimidin-5-ylamino]nicotinamide dihydrochloride (BAY 61-3606) blocks antigen-induced airway inflammation in rodents. *J. Pharmacol. Exp. Ther.* **306**(3), 1174-1181 (2003).
2. Kim, S.Y., Park, S.E., Shim, S.M., *et al.* Bay 61-3606 sensitizes TRAIL-induced apoptosis by downregulating Mcl-1 in breast cancer cells. *PLoS One* **10**(12), (2016).

#### 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, 11/28/2022

#### 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