

# BAY 61-3606 . hydrochloride

Syk inhibitor

Cell permeable, potent, ATP-competitive and reversible inhibitor of spleen tyrosine kinase (Syk) ( $IC_{50}=10nM$ ). Highly selective if compared to other kinases; shows no inhibitory effect against Btk, Fyn, Itk, Lyn, and Src up to concentrations of  $4.7\mu M$ . Anti-inflammatory and orally active. Exhibits good *in vivo* efficacy in the treatment of various allergy and asthma conditions in rat models.

Citations: 5

[View Online »](#)

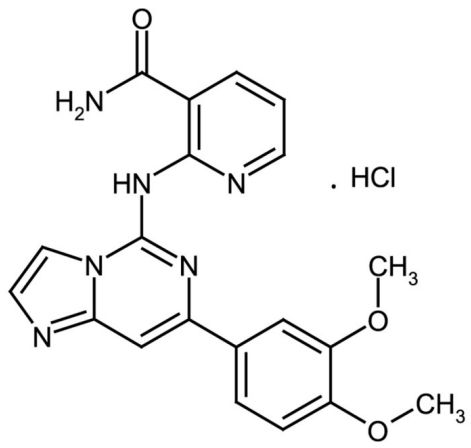
Ordering Information

[Order Online »](#)

|                  |     |
|------------------|-----|
| ALX-270-479-M001 | 1mg |
|------------------|-----|

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

|                   |  |
|-------------------|--|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. |
| Handling          | Protect from light. Hygroscopic.                                 |
| Long Term Storage | -20°C  |
| Shipping          | Ambient Temperature  |

**Regulatory Status** RUO - Research Use Only

## Product Details

|                  |  |
|------------------|--|
| Alternative Name | 2-[[7-(3,4-Dimethoxyphenyl)imidazo[1,2-c]pyrimidin-5-yl]amino]pyridine-3-carboxamide . HCl |
| Appearance       | Yellow solid.  |
| CAS              | 648903-57-5  |
| Couple Target    | Syk  |
| Couple Type      | Inhibitor  |
| Formula          | $C_{20}H_{18}N_6O_3 \cdot HCl$   |
| Identity         | Determined by MS.  |
| MW               | 390.4 . 36.5   |
| Purity           | ≥98% (HPLC(UV))  |
| Solubility       | Soluble in water (30mg/ml), DMSO (2mg/ml) or methanol (1mg/ml).                            |

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)