

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 »](#)

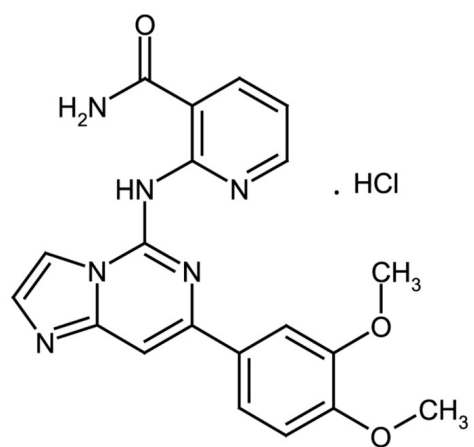
Ordering Information

[Order Online »](#)

ALX-270-479-M001	1mg
ALX-270-479-M005	5mg
ALX-270-479-M025	25mg

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).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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