

P505-15

Highly selective Syk inhibitor

P505-15 is a highly selective and potent inhibitor of spleen tyrosine kinase (Syk). It potently inhibited B cell antigen receptor-mediated B cell signaling/activation and basophil degranulation. It displayed anti-inflammatory activity in rodent models of rheumatoid arthritis. P505-15 displayed efficacy *in vitro* and *in vivo* in non-Hodgkin lymphoma and chronic lymphocytic leukemia models. It protected against toxin-induced hepatic fibrosis, associated hepatocellular injury, and hepatocarcinogenesis. P505-15 was also found to inhibit PINK1 in HeLa cells and dopaminergic neurons.

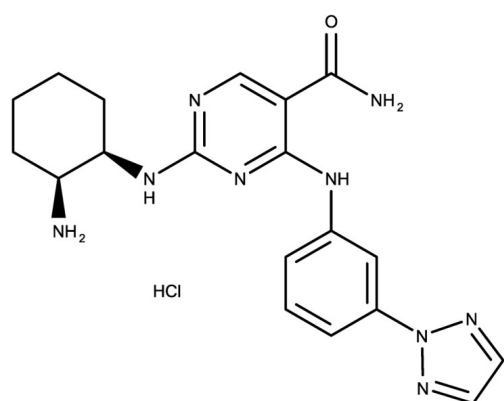
Ordering Information

[Order Online »](#)

ENZ-CHM454-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



HCl

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 2-[[[(1R,2S)-2-Aminocyclohexyl]amino]-4-[3-(triazol-2-yl)anilino]pyrimidine-5-carboxamide hydrochloride

Appearance White solid.

CAS 1370261-97-4

Couple Type Inhibitor

Formula $C_{19}H_{23}N_9O \cdot HCl$

Identity Determined by NMR.

MW 429.91

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (25 mg/ml) or water (>25 mg/ml).



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-
fr@enzolifesciences.com](mailto:info-fr@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)