# 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.

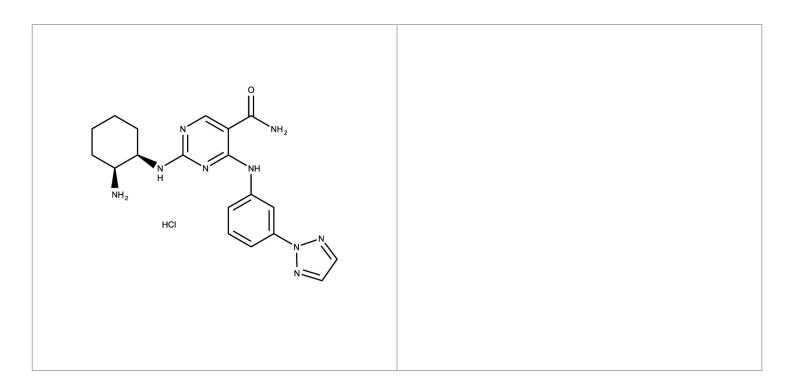
### **Ordering Information**

**Order Online** »

**ENZ-CHM454-0025** 25mg

Manuals, SDS & CofA

View Online »



### **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  $C_{19}H_{23}N_9O$  . HCI

**Identity** Determined by NMR.

MW 429.91

Purity ≥98% (HPLC)

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

Last modified: January 24, 2025

