

# R406\*

Syk inhibitor - InvitroFit™

Catalog code: inh-r406n, inh-r406n-5

<https://www.invivogen.com/r406>

For research use only

Version 23L08-MM

## PRODUCT INFORMATION

Contents R406\* (R406 besylate) is available in two quantities:

- **inh-r406n:** 2 mg R406\* - InvitroFit™
- **inh-r406n-5:** 5 x 2 mg R406\* - InvitroFit™

### Storage and stability

- R406\* is provided lyophilized and shipped at room temperature. Upon receipt, store at -20°C.
- Upon resuspension, store at -20°C. Resuspended product is stable for 6 months when properly stored. Avoid repeated freeze-thaw cycles.

### Quality Control:

- Inhibitory activity has been confirmed using cellular assays.
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

## DESCRIPTION

R406 besylate (also known as Tamatinib besylate) is an R406 salt. R406 is the active metabolite of Fostamatinib. It was initially identified as a potent inhibitor of the spleen tyrosine kinase (Syk). R406 binds to the ATP binding pocket of Syk and inhibits its kinase activity as an ATP-competitive inhibitor<sup>1</sup>. It has also been cited as an inhibitor of the Flt-3 and Ret tyrosine kinases<sup>2</sup>.

Because Syk plays an important role in FcγR-mediated signal transduction and inflammatory propagation, it is considered a good target for the inhibition of various autoimmune conditions, including rheumatoid arthritis and lymphoma. R406 was shown to attenuate autoantibody-induced arthritis in mice<sup>1</sup>, and to induce significant apoptosis in cancers displaying Syk over-expression<sup>3,4</sup>. Interestingly, R406 selectively abrogates Syk-dependent NLRP3 inflammasome activation by *C. albicans* but not by the bacterial toxin nigericin<sup>5</sup>.

1. **Braselmann S. et al., 2006.** R406, an orally available spleen tyrosine kinase inhibitor blocks fc receptor signaling and reduces immune complex-mediated inflammation. *J Pharmacol Exp Ther.* 319(3):998. 2. **Jakus Z et al., 2010.** Genetic deficiency of Syk protects mice from autoantibody-induced arthritis. *Arthritis Rheum.* 62(7):1889. 3. **Chen L. et al., 2008.** SYK-dependent tonic B-cell receptor signaling is a rational treatment target in diffuse large B-cell lymphoma. *Blood.* 111(4):2230. 4. **Zhang J. et al., 2012.** A novel retinoblastoma therapy from genomic and epigenetic analyses. *Nature.* 481(7381):329. 5. **Gross O. et al., 2009.** Syk kinase signalling couples to the Nlrp3 inflammasome for anti-fungal host defence. *Nature* 459, 433. 6. **Mócsai A. et al., 2010.** The SYK tyrosine kinase: a crucial player in diverse biological functions. *Nat Rev Immunol.* 10(6):387.

## CHEMICAL PROPERTIES

CAS number: 841290-81-1

Chemical name: 2H-Pyrido[3,2-b]-1,4-oxazin-3(4H)-one, 6-[[[5-fluoro-2-[(3,4,5-trimethoxyphenyl)amino]-4-pyrimidinyl]amino]-2,2-dimethyl-, benzenesulfonate (1:1)

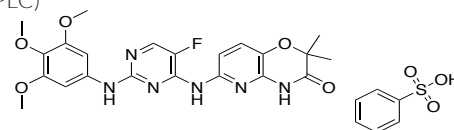
Formula: C<sub>28</sub>H<sub>29</sub>FN<sub>6</sub>O<sub>8</sub>S

Molecular weight: 628.63 g/mol

Solubility: 20 mg/ml in DMSO

Purity: ≥97% (UHPLC)

Structure:



## METHODS

### Preparation of 2 mg/ml stock solution

1. Add 1 ml of DMSO to 2 mg of R406
2. Vortex until completely resuspended.
3. Prepare aliquots and store at -20°C. Once R406 has been resuspended, dilutions can be prepared with aqueous buffers.

## PROTOCOLS (For reference only)

### Cell Culture Assay<sup>1</sup>

Cells: DLBCL (diffuse large B-cell lymphoma) cell lines

Working concentration: 0.3-5 μM

Incubation time: 72 hours

Method: Cell proliferation and apoptosis assays.

### Animal Study<sup>6</sup>

Animal model: Female C57BL/6 mice

Dose: 5 mg/kg/day

Administration: Intragastrically

Solubility: 1% DMSO/30% polyethylene glycol/1% Tween 80

## RELATED PRODUCTS

Product	Description	Cat.Code
HEK-Blue™ hDectin-1a Cells	Reporter cells	hkb-hdect1a
QUANTI-Blue™ Solution	SEAP detection reagent	rep-qbs

## TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)