

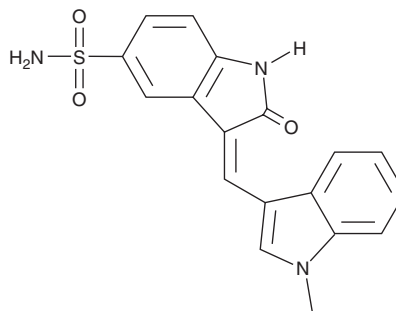
# PRODUCT INFORMATION



## Syk Inhibitor

Item No. 14829

**CAS Registry No.:** 622387-85-3  
**Formal Name:** 2,3-dihydro-3-[(1-methyl-1H-indol-3-yl)methylene]-2-oxo-1H-indole-5-sulfonamide  
**Synonyms:** OXSI 2, Spleen Tyrosine Kinase  
**MF:** C<sub>18</sub>H<sub>15</sub>N<sub>3</sub>O<sub>3</sub>S  
**FW:** 353.4  
**Purity:** ≥95% (mixture of isomers)  
**UV/Vis.:** λ<sub>max</sub>: 223, 279, 428 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Syk Inhibitor is supplied as a crystalline solid. A stock solution may be made by dissolving the Syk inhibitor in the solvent of choice, which should be purged with an inert gas. Syk Inhibitor is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

### Description

Syk is a non-receptor tyrosine kinase that upon phosphorylation binds to immunoreceptor tyrosine-based activation motifs (ITAMs) of Fcγ chains and mediates downstream signaling related to platelet function and inflammation.<sup>1</sup> Syk inhibitor is an oxindole compound that potently blocks Syk activity with an IC<sub>50</sub> value of 14 nM and inhibits FcεRI-mediated rat RBL-2H3 basophil cell degranulation (EC<sub>50</sub> = 313 nM).<sup>2</sup> At 2 μM, this compound completely abolishes convulxin-induced platelet aggregation and shape change.<sup>3</sup>

### References

1. Singh, R., Masuda, E.S., and Payan, D.G. Discovery and development of spleen tyrosine kinase (SYK) inhibitors. *J. Med. Chem.* **55**(8), 3614-3643 (2012).
2. Lai, J.Y.Q., Cox, P.J., Patel, R., *et al.* Potent small molecule inhibitors of spleen tyrosine kinase (Syk). *Bioorg. Med. Chem. Lett.* **13**(18), 3111-3114 (2003).
3. Bhavaraju, K., Kim, S., Daniel, J.L., *et al.* Evaluation of [3-(1-methyl-1H-indol-3-yl-methylene)-2-oxo-2,3-dihydro-1H-indole-5-sulfonamide] (OXSI-2), as a Syk selective inhibitor in platelets. *Eur. J. Pharmacol.* **580**(3), 285-290 (2008).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/15/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM