

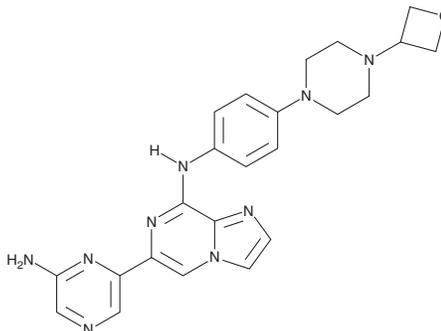
# PRODUCT INFORMATION



## Lanraplenib

Item No. 32518

**CAS Registry No.:** 1800046-95-0  
**Formal Name:** 6-(6-amino-2-pyrazinyl)-N-[4-[4-(3-oxetanyl)-1-piperazinyl]phenyl]-imidazo[1,2-a]amine  
**Synonyms:** GS-9876, GS-SYK  
**MF:** C<sub>23</sub>H<sub>25</sub>N<sub>9</sub>O  
**FW:** 443.5  
**Purity:** ≥98%  
**Supplied as:** A 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

Lanraplenib is supplied as a solid. A stock solution may be made by dissolving the lanraplenib in the solvent of choice, which should be purged with an inert gas. Lanraplenib is soluble in the organic solvent DMSO (sonicated) at a concentration of approximately 40 mg/ml.

### Description

Lanraplenib is an inhibitor of spleen tyrosine kinase (Syk; IC<sub>50</sub> = 6.2 nM).<sup>1</sup> It is selective for Syk over a panel of 395 kinases but does inhibit JAK2 (IC<sub>50</sub> = 120 nM). Lanraplenib reduces anti-IgM- and anti-CD40-induced proliferation of isolated human B cells (EC<sub>50</sub> = 108 nM). It also inhibits immune complex-induced production of TNF-α, IL-1β, and IL-6 in human macrophages (EC<sub>50</sub>s = 180, 90, and 700 nM, respectively). It reduces proteinuria and increases survival in the NZBWF1 mouse model of systemic lupus erythematosus (SLE).

### Reference

1. Blomgren, P., Chandrasekhar, J., Di Paolo, J.A., *et al.* Discovery of lanraplenib (GS-9876): A once-daily spleen tyrosine kinase inhibitor for autoimmune diseases. *ACS Med. Chem. Lett.* **11(4)**, 506-513 (2020).

#### 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, 12/16/2022

#### 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