

# PRODUCT INFORMATION

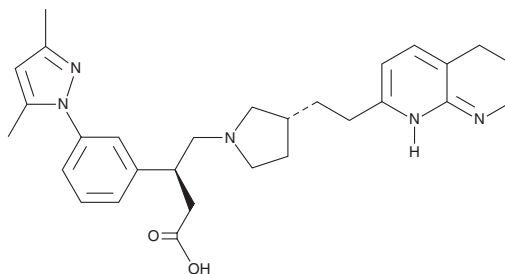


**GSK3008348**

Item No. 43471

**CAS Registry No.:** 1629249-33-7  
**Formal Name:** (βS,3R)-β-[3-(3,5-dimethyl-(1H-pyrazol-1-yl)phenyl)-3-([2-(5,6,7,8-tetrahydro-(1,8-naphthyridin-(2-yl)ethyl]-1-pyrrolidinebutanoic acid

**MF:** C<sub>29</sub>H<sub>37</sub>N<sub>5</sub>O<sub>2</sub>  
**FW:** 487.6  
**Purity:** ≥95%  
**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

GSK3008348 is supplied as a solid. A stock solution may be made by dissolving the GSK3008348 in the solvent of choice, which should be purged with an inert gas. GSK3008348 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

## Description

GSK3008348 is an integrin  $\alpha\text{V}\beta_6$  inhibitor ( $K_i = 0.01 \text{ nM}$ ).<sup>1</sup> It is selective for integrin  $\alpha\text{V}\beta_6$  over integrins  $\alpha\text{V}\beta_1$ ,  $\alpha\text{V}\beta_3$ ,  $\alpha\text{V}\beta_5$ , and  $\alpha\text{V}\beta_8$  ( $\text{IC}_{50}\text{s} = 3.9, 50, 1,000, 126, \text{ and } 16 \text{ nM}$ , respectively, in cell adhesion assays). GSK3008348 induces internalization and degradation of integrin  $\alpha\text{V}\beta_6$  and reduces the phosphorylation of SMAD2 in normal human bronchial epithelial (NHBE) cells.<sup>2</sup> Intranasal administration of GSK3008348 (3 mg/kg) decreases lung collagen levels in a mouse model of lung fibrosis induced by bleomycin (Item No. 13877).

## References

- Procopiou, P.A., Anderson, N.A., Barrett, J., *et al.* Discovery of (S)-3-(3-(3,5-dimethyl-1H-pyrazol-1-yl)phenyl)-4-((R)-3-(2-(5,6,7,8-tetrahydro-1,8-naphthyridin-2-yl)ethyl)pyrrolidin-1-yl)butanoic acid, a nonpeptidic  $\alpha\text{V}\beta_6$  integrin inhibitor for the inhaled treatment of idiopathic pulmonary fibrosis. *J. Med. Chem.* **61**(18), 8417-8443 (2018).
- John, A.E., Graves, R.H., Pun, K.T., *et al.* Translational pharmacology of an inhaled small molecule  $\alpha\text{V}\beta_6$  integrin inhibitor for idiopathic pulmonary fibrosis. *Nat. Commun.* **11**(1), 4659 (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, 07/09/2025

## 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