

# PRODUCT INFORMATION



## BIO-5192

Item No. 44105

CAS Registry No.: 327613-57-0

Formal Name: (2S)-2-[[[(2S)-1-[(3,5-dichlorophenyl)sulfonyl]-2-pyrrolidiny]carbonyl]amino]-4-[[[(2S)-4-methyl-2-[methyl[2-[4-[[[(2-methyl phenyl)amino]carbonyl]amino]phenyl]acetyl]amino]-1-oxopentyl] amino]-butanoic acid

Synonym: AMD 15057

MF: C<sub>38</sub>H<sub>46</sub>Cl<sub>2</sub>N<sub>6</sub>O<sub>8</sub>S

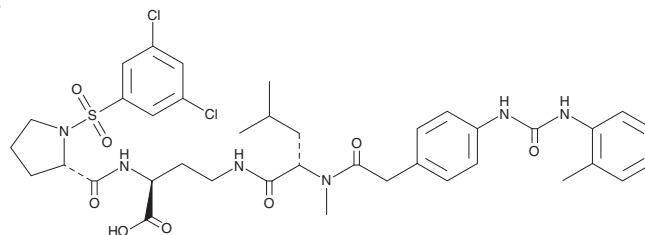
FW: 817.8

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

BIO-5192 is supplied as a solid. A stock solution may be made by dissolving the BIO-5192 in the solvent of choice, which should be purged with an inert gas. BIO-5192 is slightly soluble (0.1-1 mg/ml) in DMSO and a 1:1 solution of acetonitrile:water.

## Description

BIO-5192 is an inhibitor of integrin  $\alpha_4\beta_1$ , also known as very late antigen-4 (VLA-4; IC<sub>50</sub> = 1.8 nM).<sup>1</sup> It is selective for integrin  $\alpha_4\beta_1$  over the integrins  $\alpha_4\beta_7$ ,  $\alpha_9\beta_1$ , and  $\alpha_2\beta_1$  (IC<sub>50</sub>s = >500, 138, and 1,053 nM, respectively). BIO-5192 (1  $\mu$ g/ml) inhibits the adhesion of  $\alpha_4\beta_1$ -expressing A20 cells to fibronectin-coated plates.<sup>2</sup> It improves limb function in a rat model of experimental autoimmune encephalomyelitis (EAE) induced by myelin basic protein (MBP) peptide when administered at a dose of 30 mg/kg twice per day.<sup>1</sup>

## References

1. Leone, D.R., Giza, K., Gill, A., *et al.* An assessment of the mechanistic differences between two integrin  $\alpha_4\beta_1$  inhibitors, the monoclonal antibody TA-2 and the small molecule BIO5192, in rat experimental autoimmune encephalomyelitis. *J. Pharmacol. Exp. Ther.* **305**(3), 1150-1162 (2003).
2. Ramirez, P., Rettig, M.P., Uy, G.L., *et al.* BIO5192, a small molecule inhibitor of VLA-4, mobilizes hematopoietic stem and progenitor cells. *Blood* **114**(7), 1340-1343 (2009).

### 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, 10/22/2025

## 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