

# PRODUCT INFORMATION



## BIO-1211 (trifluoroacetate salt)

Item No. 35007

**Formal Name:** N-[2-[4-[[[(2-methylphenyl)amino]carbonyl]amino]phenyl]acetyl]-L-leucyl-L- $\alpha$ -aspartyl-L-valyl-L-proline, trifluoroacetate salt

**MF:** C<sub>36</sub>H<sub>48</sub>N<sub>6</sub>O<sub>9</sub> • XCF<sub>3</sub>COOH

**FW:** 708.8

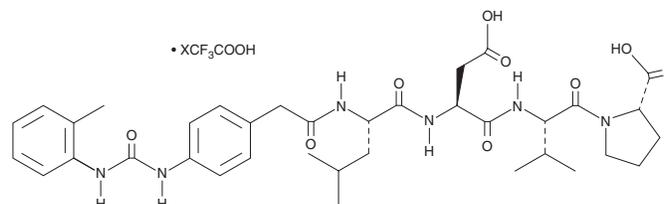
**Purity:**  $\geq$ 98%

**UV/Vis.:**  $\lambda_{\max}$ : 258 nm

**Supplied as:** A solid

**Storage:** -20°C

**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

BIO-1211 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the BIO-1211 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. BIO-1211 (trifluoroacetate salt) is soluble in DMSO.

### Description

BIO-1211 is a peptide inhibitor of  $\alpha$ 4 $\beta$ 1 integrin (IC<sub>50</sub> = 0.004  $\mu$ M).<sup>1</sup> It is selective for  $\alpha$ 4 $\beta$ 1 integrin over  $\alpha$ 4 $\beta$ 7 (IC<sub>50</sub> = 2  $\mu$ M), as well as  $\alpha$ 1 $\beta$ 1,  $\alpha$ 5 $\beta$ 1,  $\alpha$ 6 $\beta$ 1,  $\alpha$ L $\beta$ 2, and  $\alpha$ IIb $\beta$ 3 (IC<sub>50</sub>s = >100  $\mu$ M for all). BIO-1211 binds to  $\alpha$ 4 $\beta$ 1 (IC<sub>50</sub> = 1 nM) and inhibits Jurkat cell adhesion (IC<sub>50</sub> = 4 nM). Inhaled administration of BIO-1211 (3 mg/animal) prevents carbachol-induced airway hyperresponsiveness in an ovine model of allergy-induced asthma.

### Reference

1. Lin, K.c., Ateeq, H.S., Hsiung, S.H., *et al.* Selective, tight-binding inhibitors of integrin  $\alpha$ 4 $\beta$ 1 that inhibit allergic airway responses. *J. Med. Chem.* **42**(5), 920-934 (1999).

#### 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, 10/12/2023

#### 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