

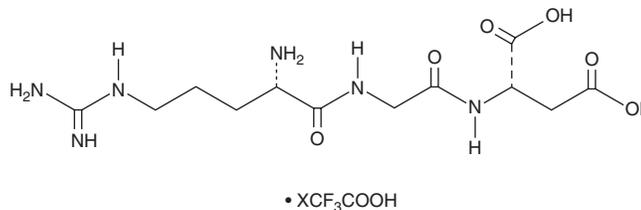
# PRODUCT INFORMATION



## Arg-Gly-Asp (trifluoroacetate salt)

Item No. 31369

**CAS Registry No.:** 120103-89-1  
**Formal Name:** N-(N-L-arginylglycyl)-L-aspartic acid, trifluoroacetate salt  
**Synonym:** RGD  
**MF:** C<sub>12</sub>H<sub>22</sub>N<sub>6</sub>O<sub>6</sub> • XCF<sub>3</sub>COOH  
**FW:** 346.3  
**Purity:** ≥95%  
**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

Arg-Gly-Asp (RGD) (trifluoroacetate salt) is supplied as a crystalline solid. Aqueous solutions of RGD (trifluoroacetate salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of RGD (trifluoroacetate salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

RGD is a synthetic peptide corresponding to the cell adhesion sequence for the integrins  $\alpha$ V $\beta$ 1,  $\alpha$ V $\beta$ 3,  $\alpha$ V $\beta$ 5,  $\alpha$ V $\beta$ 6,  $\alpha$ V $\beta$ 8,  $\alpha$ 8 $\beta$ 1, and  $\alpha$ 5 $\beta$ 1.<sup>1,2</sup> It is a component of radioligands and peptidomimetics that have been used in the study of integrin expression and activity *in vivo*.<sup>2,3</sup>

### References

1. Ruoslahti, E. RGD and other recognition sequences for integrins. *Annu. Rev. Cell Dev. Biol.* **12**, 697-715 (1996).
2. Danhier, F., Le Breton, A., and Pr  at, V. RGD-based strategies to target alpha(v) beta(3) integrin in cancer therapy and diagnosis. *Mol. Pharm.* **9(11)**, 2961-2973 (2012).
3. Schottelius, M., Laufer, B., Kessler, H., *et al.* Ligands for mapping  $\alpha$ <sub>v</sub> $\beta$ <sub>3</sub>-integrin expression *in vivo*. *Acc. Chem. Res.* **42(7)**, 969-980 (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, 12/09/2022

#### 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