



Product Data Sheet

Product Name:	Cyclo (-GRGDSP)	
Catalog Number:	AS-61110 (1 mg) AS-61110-5 (5 mg)	Lot Number: See label on vial
Sequence:	Cyclo(-Gly-Arg-Gly-Asp-Ser-Pro) (3-letter code) GRGDSP, N to C cyclized (1-letter code)	
Molecular Weight:	569.6	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Cyclo (-GRGDSP) peptide is freely soluble in water.

Storage: Cyclo (-GRGDSP) peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at -20°C or lower. Reconstituted peptide can be aliquoted and stored at -20°C or lower.

Description: This cyclic peptide is a potent vasodilator. It is more powerful than the linear GRGDSP (Cat# 22945, 22946) in changing the vascular tone of arterioles isolated from rat cremaster muscle. Ref: Mogford, JE. et al. *Circulation Res.* **79**, 821 (1996).

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

To block integrin binding in the microvessel constructs, 10 mM cyclo-GRGDSP peptide (AnaSpec, San Jose, CA) was added to the construct before polymerization. Vessel constructs were then polymerized and imaged on *day 0* and *day 1* and compared with untreated control constructs. [Kirkpatrick, ND. et al. *Am. J. Physiol. Heart Circ. Physiol.* **292**, H3198 \(2007\).](#)

The Petri dish was again rinsed three times with TBS and filled with 5 ml TBS plus 1% BSA and the indicated cations. In some experiments, 10 $\mu\text{g}/\text{ml}$ HFN7.1, 1 mg/ml cyclo(GRGDSP) (AnaSpec, Inc.), or 10 $\mu\text{g}/\text{ml}$ TS2/16 or 12G10 (Fig. 6 D) were added to the buffer-[Kong, F. et al. *J. Cell Biol.* **185**, 1275 \(2009\).](#)

Published Citations:

- Nadanaka, S. et al. *J Biol Chem* **276**, 33657 (2001).
- Kirkpatrick, ND. et al. *Am. J. Physiol. Heart Circ. Physiol.* **292**, H3198 (2007).
- Kong, F. et al. *J. Cell Biol.* **185**, 1275 (2009).

Related Products:

Name	Cat #	Size
GRGDSP	AS-22945	1 mg
GRGDSP	AS-22946	5 mg
Prototype of RGD-containing peptide, FITC-labeled FITC-LC-GRGDSP	AS-60619	1 mg
GRGDSPK	AS-61109	1 mg
GRGDSPK		

For Research Use Only