

Product Data Sheet

Product Name: Cyclo (-GRGDSP)

Catalog Number: AS-61110 (1 mg) Lot Number: See label on vial

AS-61110-5 (5 mg)

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 μ g/ml HFN7.1, 1 mg/ml cyclo(GRGDSP) (AnaSpec, Inc.), or 10 μ g/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 GRGDSPK	AS-61109	1 mg

For Research Use Only