

Product Name:	GRGDSP	
Catalog Number:	AS-22945 (1 mg) AS-22946 (5 mg)	Lot Number: See label on vial
Sequence:	H-Gly-Arg-Gly-Asp-Ser-Pro-OH (3-letter code) GRGDSP (1-letter code)	
Molecular Weight:	588.6	
% Peak Area by HPLC	: ≥ 95	

Appearance: Lyophilized white powder

Peptide Reconstitution: Using H₂O, reconstitute by adding 50-60 µl to 1 mg GRGDSP peptide.

Storage: 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: An inhibitor of cell attachment to salmosin and vitronectin. Ref: Kang, I-C. et al. *Cancer Res.* **59**, 3754 (1999).

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.

GRGDSP or GRGESP synthetic peptides (>98% purity; AnaSpec, San Jose, CA) were added to final concentrations of 1 mM, based on previous titration experiments (Adams and Lawler, 1993b, 1994). In some experiments, cells were fixed and processed for phalloidin staining of F-actin as described-Kvansakul, M. et al. *The Embo Journal.* **23**, 1223 (2004).

Additionally, small peptides containing RGD sequence (1 mM GRGDSP, Anaspec, San Jose, CA) or control RGE sequence (GRGESP, Anaspec) were used to investigate whether cells attachment was dependent on this common integrin-binding peptide sequence contained in matrix proteins such as fibronectin-<u>Gilchrist, CL. et al. J. Orthopaedic Res. **25**, 829 (2007).</u>

Published Citations:

Chan, BP. et al. *Biotech. Prog.* **20**, 566 (2004). Jiang, H. and F. Grinnell *Mol. Biol. Cell* **16**, 5070 (2005). Kvansakul, M. et al. *The Embo Journal.* **23**, 1223 (2004). Lee, J-W. et al. *Biomacromolecules* **7**, 1695 (2006). Fukushima, Y. et al. *J. Neurosurgery* **107**, 578 (2007). Gilchrist, CL. et al. *J. Orthopaedic Res.* **25**, 829 (2007). Lee, H. et al. *Lab Chip* **7**, 331 (2007). Sawai, H. et al. *J. Surg. Res.* **144**, 117 (2008).

Related Products:

Name Cyclo(-GRGDSP) (GRGDSP, N to C cyclized)	Cat # AS-61110 AS-61110-5	Size 1 mg 5 mg
Prototype of RGD-containing peptide, FITC-labeled (FITC-LC-GRGDSP)	AS-60619	1 mg
GRGESP (GRGESP)	AS-22950	5 mg

For Research Use Only

© AnaSpec, Inc. 34801 Campus Drive, Fremont, CA 94555 Tel: (800) 452-5530 | (510) 791-9560 | service@anaspec.com | www.anaspec.com