



Product Data Sheet

Peptide Name:	GRGESP	
Catalog Number:	AS-22950 (5 mg)	Lot Number: See label on vial
Sequence:	H-Gly-Arg-Gly-Glu-Ser-Pro-OH (3-letter code) GRGESP (1-letter code)	
Molecular Weight:	601.6	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: GRGESP peptide is freely soluble in water.

Storage: GRGESP 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 inactive control for the integrin-binding peptide, GRGDTP.

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 prepare T3₅₋₇ CTD C974S/N1049K with different calcium loadings, protein samples were diluted 1:200 from a 9.8 mg/ml stock into TBS containing either 2, 0.5 or 0.1 mM CaCl₂, or 1 mM EDTA, each in the absence or presence of 1 mM dithiothreitol. Blocking and washing buffers contained the same additions. Cells were added for the 1 h assay in standard serum-free DMEM. GRGDSP or GRGESP synthetic peptides (>98% purity; AnaSpec, San Jose, CA) were added to final concentrations of 1 mM-[Kvansakul, M. et al. *Embo J* **23**, 1223 \(2004\).](#)

Control cells were resuspended without blocking antibody or with nonblocking mouse IgG antibody. Additionally, small peptides containing the RGD sequence (1 mM GRGDSP, Anaspec, San Jose, CA) or control RGE sequence (GRGESP, Anaspec) were used to investigate whether cell attachment was dependent on this common integrin-binding peptide sequence contained in matrix proteins such as fibronectin-[Gilchrist, C. et al. *J Orthopaedic Res* **25**, 829 \(2007\).](#)

Published Citations:

- Rajotte, D. and E. Ruoslahti *J Biol Chem* **274**, 11593 (1999).
- Nadanaka, S. et al. *J Biol Chem* **276**, 33657 (2001).
- Kvansakul, M. et al. *Embo J* **23**, 1223 (2004).
- Huang, J. and F. Grinnell *Mol Biol Cell* **16**, 5070 (2005).
- Gilchrist, C. et al. *J Orthopaedic Res* **25**, 829 (2007).
- Sawai, H. et al. *JSR* **144**, 117 (2008).
- Bunch, TA. *JBC* **285**, 1841 (2010).

Related Products:

Name	Cat #	Size
Cyclo-[GRGESP]	AS-64447	1 mg
Cyclo-[GRGESP]		

For Research Use Only