



Product Data Sheet

Product Name: Cyclo(-RGDfK)

Catalog Number: AS-61111 (1 mg) **Lot Number:** See label on vial

Sequence: Cyclo(-Arg-Gly-Asp-D-Phe-Lys) (3-letter code)
Cyclo(-RGDfK) (1-letter code)

Molecular Weight: 603.7

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Cyclo(-RGDfK) peptide is freely soluble in H₂O.

Storage: Cyclo(-RGDfK) 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: In one study where this peptide was labeled with ¹²⁵I, it was found to bind specifically and with high affinity to αvβ3 receptors on neovascular blood vessel sections of different major human cancers. The integrin alpha(IIb)beta(3)-specific cyclic hexapeptide contains an Arg-Gly-Asp (RGD) sequence. Ref: Van Hagen, PM. et al. *Int. J. Cancer* **90**, 186 (2000); Hu, B. et al. *Biochem.* **39**, 2284 (2000).

Published Citation:

Borgman,, MP. et al. *J. Con. Rel.* **132**, 193 (2008).

For Research Use Only