

## **Product Data Sheet**

Product Name: GRGDS, amide

Catalog Number: AS-61108 (5 mg) Lot Number: See label on vial

Sequence: H-Gly-Arg-Gly-Asp-Ser-NH2 (3-letter code)

GRGDS-NH2 (1-letter code)

Molecular Weight: 489.5

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: GRGDS peptide is freely soluble in water.

Storage: GRGDS 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 peptide inhibits the adhesion of human ovarian carcinoma OVCAR-3 cells to fibronectin, but not to laminin. Ref: Kanemoto, T. et al. *Invasion Metastasis* **11**, 84 (1991).

## **Published Citations:**

Vartak, DG. et al. Mol. Pharmaceutics 6, 1856 (2009).

## Related Products:

 Name
 Cat #
 Size

 GRGDS
 AS-61107
 5 mg

**GRGDS** 

GRGDS, LC-biotin labeled AS-62347 1 mg

Biotin-LC-GRGDS

For Research Use Only