

Product Data Sheet

Product Name: Cyclo(-RGDyK)

Catalog Number: AS-61183-5 (5 mg) Lot Number: See label on vial

Sequence: Cyclo(-Arg-Gly-Asp-D-Tyr-Lys) (3-letter code)

Cyclo(-RGDyK) (1-letter code)

Molecular Weight: 620.7

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Using H₂O, reconstitute by adding 100 μl to 1 mg Cyclo(-RGDyK)

peptide.

Storage: Cyclo(-RGDyK) 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.

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.

Cyclo(RGDyK) peptide was from AnaSpec (CA, USA). In order to introduce a thiol group into cyclo(RGDyK) peptide (potent integrin $\alpha_V\beta_3$ antagonist), 1 mg of cyclo(RGDyK) was dissolved in 500 µL of dry DMSO, to which was added 1 mg of S-acetylthioglycolic acid N-hydroxysuccinimide ester (SATA), and stirred over night at room temperature Then 1 mL of 10% hydroxylamine was added and stirred for 3 h to deprotect a thiol group and to yield the thiolated RGD peptide cyclo(RGDy(ϵ -acetylthiol)K), denoted as RGD-SH. Zako, T. et al. *Biochem. Biophys. Res. Comm.* **381**, 54 (2009).

Published Citation:

Zako, T. et al. Biochem. Biophys. Res. Comm. 381, 54 (2009).

Related Products:

| Name | Cat # | Size |
|--|------------------|--------|
| Cyclo [-RGDyK(HiLyte Fluor™ 750)], Abs/Em = 754/778 | 3 nm AS-62333-01 | 0.1 mg |
| Cyclo (-RGDfK) (Cyclo(-RGDfK)) | AS-61111 | 1 mg |
| Cyclo (-RADfK-), RGD negative control (Cyclo(-RADfK-)) | AS-62351 | 1 mg |

For Research Use Only