



Product Data Sheet

Product Name: Cyclo [-RGDyK(HiLyte Fluor™ 750)]

Catalog Number: AS-62333-01 (0.1 mg) **Lot Number:** See label on vial

Sequence: Cyclo[-Arg-Gly-Asp-D-Tyr-Lys(HiLyte Fluor™ 750)] (3-letter code)
Cyclo[-RGDy-K(HiLyte Fluor™ 750)] (1-letter code)

Molecular Weight: 1630.5

Peptide Purity: >95%

Appearance: Lyophilized blue powder

Peptide Reconstitution: Cyclo [-RGDyK(HiLyte Fluor™ 750)] peptide is freely soluble in water.

Storage: Cyclo [-RGDyK(HiLyte Fluor™ 750)] 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.

Related Products:

Name	Cat #	Size
Cyclo(-RGDyK)	AS-61183-5	5 mg

For Research Use Only