



Product Data Sheet

Product Name: Cys(Npys)-(Arg)9

Catalog Number: AS-61205 (1 mg) **Lot Number:** See label on vial

Sequence: H-Cys(Npys)-Arg-Arg-Arg-Arg-Arg-Arg-Arg-Arg-NH2 (3-letter code)
C(Npys)RRRRRRRRR-NH2 (1-letter code)

Molecular Weight: 1681.1

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Using H₂O, reconstitute by adding 100 µl to 1 mg Cys(Npys)-(Arg)9 peptide.

Storage: Cys(Npys)-(Arg)9 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
Cys(Npys)-(D-Arg)9 (C(Npys)rrrrrrrr-NH2)	AS-61206	1 mg
(Arg)9 (RRRRRRRRR)	AS-61204	1 mg
(Arg)9, FAM-labeled, Abs/Em=492/518 nm (FAM-RRRRRRRRR)	AS-61207	1 mg
(Arg)9, TAMRA-labeled, Abs/Em=544/572 nm (TAMRA-RRRRRRRRR)	AS-61208	1 mg

For Research Use Only