



Product Data Sheet

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

Catalog Number: AS-61206 (1 mg) **Lot Number:** See label on vial

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

Molecular Weight: 1681.0

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized off-white powder

Peptide Reconstitution: Cys(Npys)-(D-Arg)9 peptide is freely soluble in H₂O.

Storage: Cys(Npys)-(D-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 Product:

Name	Cat #	Size
Cys(Npys)-(Arg)9 (C(Npys)RRRRRRRRR-NH2)	AS-61205	1 mg

For Research Use Only