



Product Data Sheet

Product Name: Biotin - Bradykinin

Catalog Number: AS-23974 (1mg) **Lot Number:** See label on vial

Sequence: Biotin-Arg-Pro-Pro-Gly-Phe-Ser-Pro-Phe-Arg-OH (3-letter code)
Biotin-RPPGFSPFR (1-letter code)

Molecular Weight: 1286.5

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Biotin-Bradykinin is freely soluble in water.

Storage: Biotin-Bradykinin 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.

For Research Use Only