



Product Data Sheet

Product Name: Parathyroid Hormone (1-34)-Lys(Biotin), human

Catalog Number: AS-23647 (1.0 mg) **Lot Number:** See label on vial

Sequence: H-Ser-Val-Ser-Glu-Ile-Gln-Leu-Met-His-Asn-Leu-Gly-Lys-His-Leu-Asn-Ser-Met-Glu-Arg-Val-Glu-Trp-Leu-Arg-Lys-Lys-Leu-Gln-Asp-Val-His-Asn-Phe-Lys(Biotin)-OH (3-letter code)
SVSEIQLMHNLGKHLNSMERVEWLRKKLQDVHNF(Biotin)
(1-letter code)

Molecular Weight: 4472.2

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Parathyroid peptide is freely soluble in water.

Storage: Parathyroid 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
Parathyroid Hormone (1-34), human SVSEIQLMHNLGKHLNSMERVEWLRKKLQDVHNF	AS-20708	1 mg

For Research Use Only