



## Product Data Sheet

---

**Product Name:**  $\beta$ -Endorphin, human

**Catalog Number:** AS-24319 (1 mg)                      **Lot Number:** See label on vial

**Sequence:** H-Tyr-Gly-Gly-Phe-Met-Thr-Ser-Glu-Lys-Ser-Gln-Thr-Pro-Leu-Val-Thr-Leu-Phe-Lys-Asn-Ala-Ile-Ile-Lys-Asn-Ala-Tyr-Lys-Lys-Gly-Glu-OH  
(3-letter code)  
YGGFMTSEKSQTPLVTLFKNAIIKNAYKKGE (1-letter code)

**Molecular Weight:** 3465.1

**Peptide Purity:** >95%

**Appearance:** Lyophilized white powder

**Peptide Reconstitution:**  $\beta$ -Endorphin peptide is freely soluble in water.

**Storage:**  $\beta$ -Endorphin peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^{\circ}\text{C}$  or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}\text{C}$  or lower.

*For Research Use Only*