



Product Data Sheet

Product Name: Oxytocin

Catalog Number: AS-24275 (5 mg) Lot Number: See label on vial
AS-24276 (25 mg)

Sequence: H-Cys-Tyr-Ile-Gln-Asn-Cys-Pro-Leu-Gly-NH₂ (Disulfide bridge: 1-6)
(3-letter code)
CYIQNCPLG-NH₂ (Disulfide bridge: 1-6) (1-letter code)

Molecular Weight: 1007.2

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Oxytocin peptide is freely soluble in water.

Storage: Oxytocin 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.

Reference: Ludwig, M. *Am. J. Physio.* **268**, E537 (1995).

Related Products:

Name	Cat #	Size
Biotin-Oxytocin	AS-23994	1 mg
Biotin-CYIQNCPLG-NH ₂ (Disulfide bridge:1-6)		

For Research Use Only