



## Product Data Sheet

---

**Product Name:** Orexin A, bovine, human, mouse, rat  
**Catalog Number:** AS-24470 (1 mg) **Lot Number:** See label on vial

**Sequence:** Pyr-Pro-Leu-Pro-Asp-Cys-Cys-Arg-Gln-Lys-Thr-Cys-Ser-Cys-Arg-Leu- Tyr-Glu-Leu-Leu-His-Gly-Ala-Gly-Asn-His-Ala-Ala-Gly-Ile-Leu-Thr-Leu- NH<sub>2</sub> (Disulfide bridge: 6-12 and 7-14) (3-letter code)  
Pyr-PLPDCCRQKTCSCRLYELLHGAGNHAAGILTL-NH<sub>2</sub> (Disulfide bridge: 6-12 and 7-14) (1-letter code)

**Molecular Weight:** 3561.2

**% Peak Area by HPLC:** ≥ 95

**Appearance:** Lyophilized white powder

**Peptide Reconstitution:** Use distilled or higher quality water. Add water directly to the lyophilized peptide powder to obtain a final concentration of approximately 0.5 mg/mL to 1mg/mL or less. Gently vortex to mix. For peptides that have poor solubility in the suggested solvent, brief sonication may increase solubility in some cases.

**Storage:** Peptide is shipped at ambient temperature. Upon receipt, store lyophilized powder at –20°C or lower. Reconstituted peptide should be aliquoted into several freezer vials and stored at –20°C or lower. Do not freeze thaw.

**Description:** This hypothalamic neuropeptide has been shown to regulate feeding behavior. Ref: Sakurai, T. et al. *Cell* **92**, 573 (1998).

***For Research Use Only***