



## Product Data Sheet

**Product Name:** Corticotropin Releasing Factor, CRF, human, rat  
**Catalog Number:** AS-24254 (1 mg) **Lot Number:** See label on vial

**Sequence:** H-Ser-Glu-Glu-Pro-Pro-Ile-Ser-Leu-Asp-Leu-Thr-Phe-His-Leu-Leu-Arg-Glu-Val-Leu-Glu-Met-Ala-Arg-Ala-Glu-Gln-Leu-Ala-Gln-Gln-Ala-His-Ser-Asn-Arg-Lys-Leu-Met-Glu-Ile-Ile-NH<sub>2</sub> (3-letter code)  
SEPPISLDLTFHLLREVLEMARAEQLAQQAHSNRKLMEII-NH<sub>2</sub> (1-letter code)

**Molecular Weight:** 4757.5

**% 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:** The peptide, corticotropin releasing factor (CRF) was first isolated in the mammalian brain that regulates the hypothalamic-pituitary-adrenocortical axis. It also plays a key role in modulating the endocrine, autonomic, and behavioral responses to stress and cardiovascular, gastrointestinal, and immune activities. Ref: Lewis, K. et al. *PNAS* **98**, 7570 (2001); Chen, R. et al. *PNAS* **90**, 8967 (1993).

*For Research Use Only*