

Product Data Sheet

Product Name: ACTH (1-24), human

Catalog Number: AS-20613 (1 mg) Lot Number: See label on vial

AS-20614 (5 mg)

Sequence: H-Ser-Tyr-Ser-Met-Glu-His-Phe-Arg-Trp-Gly-Lys-Pro-Val-Gly-Lys-Lys-

Arg-Arg-Pro-Val-Lys-Val-tyr-pro-OH (3-letter code) SYSMEHFRWGKPVGKKRRPVKVYP (1-letter code)

Molecular Weight: 2934.5

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Using H₂O, reconstitute by adding 100 μl to 1 mg ACTH peptide. This

peptide is soluble in 1%NH₄OH and DMSO

Storage: ACTH 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.

References: Baumann, G. and JP. Felber, *J. Clin. Endocrinol. Metab.* **42**, 160 (1976); Burt, DS. and DR. Stanworth, *Biochim. Biophy. Acta* **762**, 458 (1983).

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

We used two other peptides as control peptides in the fusion assay, the oxidized chain B of insulin from bovine pancreas (Sigma, St. Louis, MO) and human adrenocorticotropic hormone (ACTH) residues 1–24 (AnaSpec, San Jose, CA), dissolved in diH_2O without further modification. In the case of oxidized chain B of insulin and ACTH, we added PBS to the diluted solutions to match the osmolarity and ionic strength of the solutions of WT-B5 peptide at a given concentration of peptide-Estes, D. J. et al. *Biophys J.* **91**, 233 (2006).

Published Citations:

Estes, D. J. et al. *Biophys J.* **91**, 233 (2006). O'Donoghue, AJ. et al. *J. Biol. Chem.* **283**, 29186 (2008).

For Research Use Only