

# PRODUCT INFORMATION

## ACTH (11-24) (human, mouse, rat, porcine, bovine, ovine) (acetate) Item No. 41559

**Formal Name:**  $\alpha^{11-24}$ -corticotropin, acetate  
**Synonyms:**  $\alpha^{11-24}$ -ACTH, Adrenocorticotropin  
 Hormone (11-24), Corticotropin (11-24)  
**Peptide Sequence:** KPVGKKRRPVKVYP-OH H—Lys—Pro—Val—Gly—Lys—Lys—Arg—Arg—Pro—Val—  
**MF:**  $C_{77}H_{134}N_{24}O_{16} \cdot XC_2H_4O_2$  Lys—Val—Tyr—Pro—OH  
**FW:** 1,652.0  $\cdot XCH_3CO_2H$   
**Purity:**  $\geq 98\%$   
**Supplied as:** A solid  
**Storage:**  $-20^\circ C$   
**Stability:**  $\geq 4$  years

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

ACTH (11-24) (human, mouse, rat, porcine, bovine, ovine) (acetate) is supplied as a solid. A stock solution may be made by dissolving the ACTH (11-24) (human, mouse, rat, porcine, bovine, ovine) (acetate) in the solvent of choice, which should be purged with an inert gas. ACTH (11-24) (human, mouse, rat, porcine, bovine, ovine) (acetate) is slightly soluble (0.1-1 mg/ml) in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of ACTH (11-24) (human, mouse, rat, porcine, bovine, ovine) (acetate) can be prepared by directly dissolving the solid in aqueous buffers. ACTH (11-24) (human, mouse, rat, porcine, bovine, ovine) (acetate) is soluble ( $\geq 10$  mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

Adrenocorticotropin hormone (ACTH) (11-24) is a peptide fragment of ACTH (Item No. 24257), a peptide hormone found in the brain that is involved in the biological stress response, and an antagonist of melanocortin receptor 2 (MC2R), also known as the ACTH receptor ( $IC_{50} = \sim 1$  nM in HeLa cells expressing the mouse receptor).<sup>1,2</sup> It induces corticosterone and aldosterone production in isolated rat zona glomerulosa and zona fasciculata cells, respectively, with minimum effective concentrations of 4 and 3  $\mu M$ , respectively.<sup>3</sup> Unlike ACTH (4-11) (Item No. 27430), ACTH (11-24) does not induce differentiation of primary mouse melanocytes.<sup>4</sup> *In vivo*, ACTH (11-24) (1.69  $\mu g$ /animal, i.c.v.) inhibits ACTH-induced stretching, yawning, and penile erections in rats.<sup>5</sup>

### References

1. Strand, F.L., Lee, S.J., Zuccarelli, L.A., et al. *Rev. Neurosci.* **4**(4), 321-363 (1993).
2. Kapas, S., Cammas, F.M., Hinson, J.P., et al. *Endocrinology* **137**(8), 3291-3294 (1996).
3. Szalay, K.S., De Wied, D., and Stark, E. *J. Steroid Biochem.* **32**(2), 259-262 (1989).
4. Hirobe, T. and Hiroyuki, A. *J. Exp. Zool.* **286**(6), 632-640 (2000).
5. Genedani, S. and Bernardi, M. *Neuropeptides* **26**(4), 241-244 (1994).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 09/11/2024

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM