

PRODUCT INFORMATION



N-Acetyl ACTH (1-17) (human) (trifluoroacetate salt)

Item No. 27404

Formal Name:	N-acetyl- α^{1-17} -corticotropin, trifluoroacetate salt
Synonyms:	Ac-SYSMEHFRWGKPVGKKR, N-Acetyl Adrenocorticotrophic Hormone (1-17) $\text{Ac-Ser-Tyr-Ser-Met-Glu-His-Phe-Arg-Trp-Gly-Lys-Pro-Val-Gly-Lys-Lys-Arg-OH}$
MF:	$\text{C}_{97}\text{H}_{147}\text{N}_{29}\text{O}_{24}\text{S} \cdot \text{XCF}_3\text{COOH}$
FW:	2,135.5
Purity:	$\geq 95\%$
Supplied as:	A solid
Storage:	-20°C
Stability:	≥ 4 years

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

Laboratory Procedures

N-Acetyl ACTH (1-17) (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the N-acetyl ACTH (1-17) (human) (trifluoroacetate salt) in water. The solubility of N-acetyl ACTH (1-17) (human) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

N-Acetyl ACTH (1-17) is an N-terminal peptide fragment of adrenocorticotrophic hormone (ACTH), a peptide hormone found in the brain that is involved in the biological stress response.^{1,2} It has been found in pituitary extracts from *Cpe^{fat/fat}* mice that lack the carboxypeptidase E active form, which is required for processing prohormones.³

References

1. Fenger, M. α -Melanocyte-stimulating-hormone precursors in the pig pituitary. *Biochem. J.* **235**(3), 715-722 (1986).
2. Strand, F.L., Lee, S.J., Zuccarelli, L.A., *et al.* Non-corticotrophic ACTH peptides modulate nerve development and regeneration. *Rev. Neurosci.* **4**(4), 321-363 (1993).
3. Che, F.-Y. and Fricker, L.D. Quantitative peptidomics of mouse pituitary: Comparison of different stable isotopic tags. *J. Mass. Spectrom.* **40**(2), 238-249 (2005).

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, 12/22/2022

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