

PRODUCT INFORMATION



ACTH (1-4) (trifluoroacetate salt)

Item No. 43818

Formal Name:	L-seryl-L-tyrosyl-L-seryl-L-methionine	
Synonyms:	Adrenocorticotrophic hormone (1-4), Corticotropin (1-4), Ser-Tyr-Ser-Met-OH	
Peptide Sequence:	SYSM-OH	H-Ser-Tyr-Ser-Met-OH
MF:	$C_{20}H_{30}N_4O_8S \cdot XCF_3COOH$	$\bullet XCF_3COOH$
FW:	486.5	
Purity:	≥98%	
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

ACTH (1-4) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the ACTH (1-4) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. ACTH (1-4) (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in acetonitrile. It is also slightly soluble (0.1-1 mg/ml) in water. We do not recommend storing the aqueous solution for more than one day.

Description

ACTH (1-4) is an N-terminal peptide fragment of adrenocorticotrophic hormone (ACTH; Item No. 24257).¹ It potentiates natriuresis induced by arginine vasopressin (argipressin; Item No. 24154) in cats when administered at a dose of 20 nmol/kg. Intracerebroventricular administration of ACTH (1-4) (3 µg/animal) potentiates excessive grooming induced by α-melanocyte-stimulating hormone (α-MSH; Item No. 29923) in rats.²

References

1. Cort, J.H., Cort, J., Nováková, J., *et al.* Interaction of vasopressins and linear N-terminal ACTH fragments in the induction of natriuresis. *Eur. J. Clin. Invest.* **4(5)**, 293-298 (1974).
2. De Graan, P.N.E., Spruijt, B.M., Eberle, A.N., *et al.* ACTH₁₋₄ potentiates α-MSH-induced melanophore dispersion and excessive grooming. *Peptides* **7(1)**, 1-4 (1986).

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, 09/05/2025

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