

PRODUCT INFORMATION

ACTH (3-24) (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt)
Item No. 41464

Synonyms:	Adrenocorticotrophic Hormone (3-24), Corticotropin (3-24)	H—Ser—Met—Glu—His—Phe—Arg—Trp—Gly—Lys—Pro— Val—Gly—Lys—Lys—Arg—Arg—Pro—Val—Lys—Val— Tyr—Pro—OH
Peptide Sequence:	SMEHFRWGKPVGKKRRPVKVYP-OH	
MF:	C ₁₂₄ H ₁₉₆ N ₃₈ O ₂₇ S • XCF ₃ COOH	
FW:	2,683.2	
Purity:	≥95%	
Supplied as:	A solid	• XCF ₃ COOH
Storage:	-20°C	
Stability:	≥4 years	

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

Laboratory Procedures

Adrenocorticotrophic hormone (ACTH) (3-24) (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the ACTH (3-24) (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. ACTH (3-24) (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) is sparingly soluble (1-10 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 (3-24) (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. ACTH (3-24) (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

ACTH (3-24) is a potential impurity found in commercial preparations of tetracosactide, also known as cosyntropin, a synthetic form of ACTH (1-24).¹

Reference

1. Taichrib, A., Scriba, G.K., and Neusüss, C. Identification and characterization of impurities of tetracosactide by capillary electrophoresis and liquid chromatography coupled to time-of-flight mass spectrometry. *Anal. Bioanal. Chem.* **401**(4), 1365-1375 (2011).

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, 10/04/2024

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