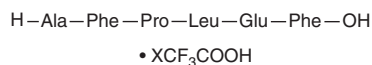


# PRODUCT INFORMATION

## ACTH (34-39) (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt)

Item No. 43820

<b>Formal Name:</b>	L-alanyl-L-phenylalanyl-L-prolyl-L-leucyl-L- $\alpha$ -glutamyl-L-phenylalanine
<b>Synonyms:</b>	Adrenocorticotrophic Hormone (34-39), Corticotropin (34-39), H-Ala-Phe-Pro-Leu-Glu-Phe-OH
<b>Peptide Sequence:</b>	H-AFPLEF-OH
<b>MF:</b>	$C_{37}H_{50}N_6O_9 \cdot XCF_3COOH$
<b>FW:</b>	722.8
<b>Purity:</b>	$\geq 98\%$
<b>Supplied as:</b>	A solid
<b>Storage:</b>	-20°C
<b>Stability:</b>	$\geq 4$ years



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

### Laboratory Procedures

ACTH (34-39) (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the ACTH (34-39) (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. ACTH (34-39) (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in DMSO and a 1:1 solution acetonitrile:water.

### Description

Adrenocorticotrophic hormone (ACTH) (34-39) is a peptide fragment of ACTH, a peptide hormone produced by the anterior pituitary gland that is involved in the biological stress response.<sup>1</sup>

### Reference

1. Lightman, S.L., Birnie, M.T., and Conway-Campbell, B.L. Dynamics of ACTH and cortisol secretion and implications for disease. *Endocr. Rev.* **41**(3), 470-490 (2020).

#### 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, 09/19/2025

#### 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](http://WWW.CAYMANCHEM.COM)