

PRODUCT INFORMATION



ACTH (1-39) (trifluoroacetate salt)

Item No. 24257

Synonym:	Adrenocorticotropin Hormone (1-39)	H—Ser—Tyr—Ser—Met—Glu—His—Phe—Arg—Trp—Gly—
MF:	C ₂₀₇ H ₃₀₈ N ₅₆ O ₅₈ S • XCF ₃ COOH	Lys—Pro—Val—Gly—Lys—Lys—Arg—Arg—Pro—Val—
FW:	4,541.1	Lys—Val—Tyr—Pro—Asn—Gly—Ala—Glu—Asp—Glu—
Purity:	≥95%	Ser—Ala—Glu—Ala—Phe—Pro—Leu—Glu—Phe—OH
Supplied as:	A lyophilized powder	• 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

ACTH (1-39) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the ACTH (1-39) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. ACTH (1-39) (trifluoroacetate salt) is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

Description

ACTH (1-39) is a potent agonist of melanocortin receptor 2 (MC2R) with an EC₅₀ value of 57 pM in HeLa cells expressing the mouse receptor.¹ It induces accumulation of cAMP and release of corticosterone from rat adrenal cells in a concentration-dependent manner.² ACTH (1-39) also protects neurons from death induced by staurosporine (Item No. 81590), glutamate, NMDA (Item No. 14581), AMPA (Item No. 14571), kainate, quinolinic acid (Item No. 14941), and reactive oxygen species when used at a concentration of 400 nM.³ Formulations containing ACTH (1-39) have been used in the treatment of systemic lupus erythematosus.

References

1. Kapas, S., Cammas, F.M., Hinson, J.P., *et al.* Agonist and receptor binding properties of adrenocorticotropin peptides using the cloned mouse adrenocorticotropin receptor expressed in a stably transfected HeLa cell line. *Endocrinology* **137(8)**, 3291-3294 (1996).
2. Bristow, A.F., Gleed, C., Fauchère, J.L., *et al.* Effects of ACTH (corticotropin) analogues on steroidogenesis and cyclic AMP in rat adrenocortical cells. Evidence for two different steroidogenically responsive receptors. *Biochem J.* **186(2)**, 599-603 (1980).
3. Lisak, R.P., Nedelkoska, L., Bealmear, B., *et al.* Melanocortin receptor agonist ACTH 1-39 protects rat forebrain neurons from apoptotic, excitotoxic and inflammation-related damage. *Exp. Neurol.* **273**, 161-167 (2015).

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, 10/25/2022

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