

PRODUCT INFORMATION



Cosyntropin (acetate)

Item No. 23912

CAS Registry No.: 60189-34-6

Formal Name: α^{1-24} -corticotropin, acetate

Synonyms: ACTH (1-24), Adrenocorticotrophic Hormone 1-24, Tetracosactide

MF: $C_{136}H_{210}N_{40}O_{31}S \cdot XC_2H_4O_2$

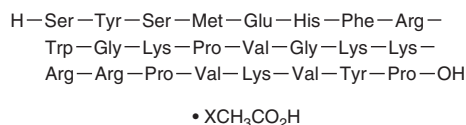
FW: 2,993.5

Purity: $\geq 95\%$

Supplied as: A crystalline solid

Storage: $-20^\circ C$

Stability: ≥ 4 years



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

Laboratory Procedures

Cosyntropin (acetate) is supplied as a crystalline solid. Aqueous solutions of cosyntropin (acetate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of cosyntropin (acetate) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Cosyntropin is a synthetic peptide analog of adrenocorticotrophic hormone (ACTH) residues 1-24.¹ It increases the plasma concentration of hydrocortisone (Item No. 20739) in mares more rapidly and consistently than natural ACTH 1-24. Formulations containing cosyntropin have been used to diagnose adrenal deficiencies in humans, dogs, and horses.

Reference

1. Elier, H., Goble, D., and Oliver, J. Adrenal gland function in the horse: Effects of cosyntropin (synthetic) and corticotropin (natural) stimulation. *Am. J. Vet. Res.* **40(5)**, 724-726 (1979).

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, 10/24/2022

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