

PRODUCT INFORMATION



Alsactide (trifluoroacetate salt)

Item No. 41462

Formal Name: 1-β-alanine-17-[N-(4-aminobutyl)-L-lysineamide]-α¹⁻¹⁷-corticotropin, trifluoroacetate salt

Synonyms: [β-Ala¹,Lys¹⁷]ACTH-(1-17)-4-amino-n-butylamide, Hoechst 433

Peptide Sequence: XYSMEHFRWVGKPVGKKK-NH-(CH₂)₄-NH₂ (where X = β-alanine)

MF: C₉₉H₁₅₅N₂₉O₂₁S • XCF₃COOH

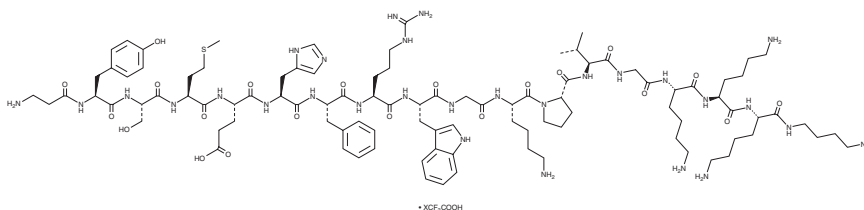
FW: 2,119.6

Purity: ≥95%

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

Alsactide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the alsactide (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Alsactide (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in ethanol and soluble (≥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 alsactide (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. Alsactide (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

Alsactide is a peptide derivative of adrenocorticotrophic hormone (ACTH; Item No. 24257), a peptide hormone found in the brain that is involved in the biological stress response.¹ It increases mean arterial blood pressure (MAP) following bleeding-induced hypotension in a rat model of hemorrhagic shock 15-30 minutes after administration at a dose of 160 µg/kg.

Reference

1. Bertolini, A., Guarini, S., Rompianesi, E., *et al.* α-MSH and other ACTH fragments improve cardiovascular function and survival in experimental hemorrhagic shock. *Eur. J. Pharmacol.* **130**(1-2), 19-26 (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, 09/19/2024

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