

# PRODUCT INFORMATION



## Calcineurin Autoinhibitory Peptide (acetate)

Item No. 44133

**Formal Name:** L-isoleucyl-L-threonyl-L-seryl-L-phenylalanyl-L- $\alpha$ -glutamyl-L- $\alpha$ -glutamyl-L-alanyl-L-lysylglycyl-L-leucyl-L- $\alpha$ -aspartyl-L-arginyl-L-isoleucyl-L-asparaginyl-L- $\alpha$ -glutamyl-L-arginyl-L-methionyl-L-prolyl-L-prolyl-L-arginyl-L-arginyl-L- $\alpha$ -aspartyl-L-alanyl-L-methionyl-L-proline, acetate

**Peptide Sequence:** Ac-ITSFEEAKGLDRINERMPPRRDAMP-OH

**MF:**  $C_{124}H_{205}N_{39}O_{39}S_2 \cdot XC_2H_4O_2$

**FW:** 2,930.4

**Purity:**  $\geq 95\%$

**Supplied as:** A solid

**Storage:**  $-20^\circ C$

**Stability:**  $\geq 4$  years

H-Ile-Thr-Ser-Phe-Glu-Glu-Ala-Lys-Gly-Leu-  
Asp-Arg-Ile-Asn-Glu-Arg-Met-Pro-Pro-Arg-  
Arg-Asp-Ala-Met-Pro-OH  
 $\cdot XCH_3CO_2H$

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

### Laboratory Procedures

Calcineurin autoinhibitory peptide (acetate) is supplied as a solid. A stock solution may be made by dissolving the calcineurin autoinhibitory peptide (acetate) in the solvent of choice, which should be purged with an inert gas. Calcineurin autoinhibitory peptide (acetate) is soluble ( $\geq 10$  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 calcineurin autoinhibitory peptide (acetate) can be prepared by directly dissolving the solid in aqueous buffers. Calcineurin autoinhibitory peptide (acetate) is soluble ( $\geq 10$  mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

Calcineurin autoinhibitory peptide is a peptide fragment corresponding to residues 457 to 482 of calcineurin A that inhibits the phosphatase activity of calcineurin A in a cell-free assay ( $IC_{50} = 10 \mu M$ ).<sup>1</sup>

### Reference

1. Hashimoto, Y., Perrino, B.A., and Soderling, T.R. Identification of an autoinhibitory domain in calcineurin. *J. Biol. Chem.* **265**(4), 1924-1927 (1990).

#### 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, 03/12/2026

#### 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