

PRODUCT INFORMATION

Diaminopropionyl Tripeptide-33 (acetate)

Item No. 45627

Formal Name: (2,3-diaminopropanoyl)-L-alanyl-L-prolyl-L-histidine, acetate

Synonyms: DAP Tripeptide-33,
H-Ala-Pro-His-OH

Peptide Sequence: H-Dap-APH-OH (where Dap = 2,3-diaminopropionic acid)

MF: $C_{17}H_{27}N_7O_5 \cdot XC_2H_4O_2$

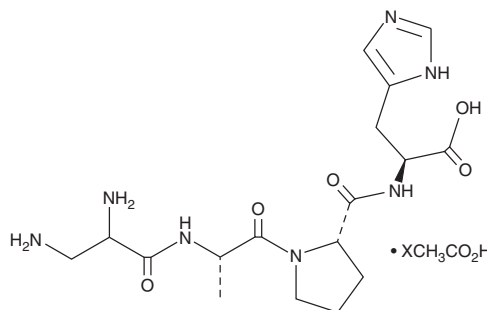
FW: 409.4

Purity: $\geq 95\%$

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

Diaminopropionyl tripeptide-33 (acetate) is supplied as a solid. A stock solution may be made by dissolving the diaminopropionyl tripeptide-33 (acetate) in the solvent of choice, which should be purged with an inert gas. Diaminopropionyl tripeptide-33 (acetate) slightly soluble (0.1-1 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 diaminopropionyl tripeptide-33 (acetate) can be prepared by directly dissolving the solid in aqueous buffers. Diaminopropionyl tripeptide-33 (acetate) is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Diaminopropionyl tripeptide-33 is a peptide containing alanine, proline, and histidine and modified with a 2,3-diaminopropionyl (DAP) group.

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, 06/04/2026

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