

PRODUCT INFORMATION



Goralatide (acetate)

Item No. 36207

Formal Name: N-acetyl-L-seryl-D- α -aspartyl-L-lysyl-L-proline, acetate

Synonyms: N-Acetyl-Ser-Asp-Lys-Pro, AcSDKP

MF: $C_{20}H_{33}N_5O_9 \cdot XC_2H_4O_2$

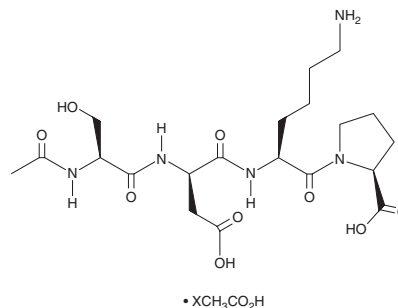
FW: 487.5

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

Goralatide (acetate) is supplied as a solid. A stock solution may be made by dissolving the goralatide (acetate) in the solvent of choice. Goralatide (acetate) is soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Goralatide is a tetrapeptide regulator of hematopoiesis.¹ It decreases the proportion of mouse high proliferative potential colony forming cells (HPP-CFCs) in S phase when cultured using regenerating femoral marrow conditioned medium at a concentration of 1 ng/ml. Goralatide (4 $\mu\text{g/kg}$ twice per day) reduces decreases in the number of HPP-CFCs, burst-forming unit erythroid (BFU-E) cells, and colony forming unit granulocyte-macrophages (CFU-GMs) induced by 5-fluorouracil (5-FU; Item No. 14416) in mice.² It also reduces mortality induced by doxorubicin (Item No. 15007) in mice.³

References

1. Robinson, S., Lenfant, M., Wdzieczak-Bakala, J., *et al.* The molecular specificity of action of the tetrapeptide acetyl-N-Ser-Asp-Lys-Pro (AcSDKP) in the control of hematopoietic stem cell proliferation. *Stem Cells* **11**(5), 422-427 (1993).
2. Aidoudi, S., Guigon, M., Lebeurier, I., *et al.* *In vivo* effect of platelet factor 4 (PF4) and tetrapeptide AcSDKP on haemopoiesis of mice treated with 5-fluorouracil. *Br. J. Haematol.* **94**(3), 443-448 (1996).
3. Massé, A., Ramirez, L.H., Bindoula, G., *et al.* The tetrapeptide acetyl-N-Ser-Asp-Lys-Pro (Goralatide) protects from doxorubicin-induced toxicity: Improvement in mice survival and protection of bone marrow stem cells and progenitors. *Blood* **91**(2), 441-449 (1998).

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, 12/19/2022

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