

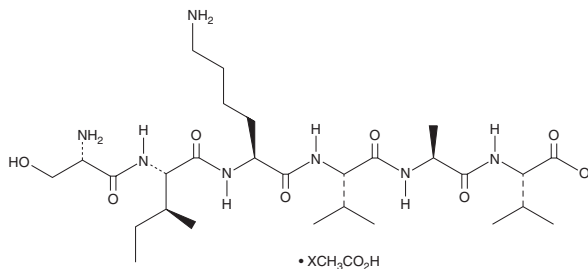
# PRODUCT INFORMATION



## SIKVAV (acetate)

Item No. 35159

**Formal Name:** L-seryl-L-isoleucyl-L-lysyl-L-valyl-L-alanyl-L-valine, acetate  
**Synonyms:** Hexapeptide-10, Ser-Ile-Lys-Val-Ala-Val  
**MF:** C<sub>28</sub>H<sub>53</sub>N<sub>7</sub>O<sub>8</sub> • XC<sub>2</sub>H<sub>4</sub>O<sub>2</sub>  
**FW:** 615.8  
**Purity:** ≥98%  
**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

SIKVAV (acetate) is supplied as a solid. Aqueous solutions of SIKVAV (acetate) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of SIKVAV (acetate) in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

SIKVAV is a laminin  $\alpha$ 1-derived peptide.<sup>1</sup> It binds to  $\alpha$ 3,  $\alpha$ 6, and  $\beta$ 1 integrins and induces adhesion of CAC2 adenoid cystic carcinoma cells when used at a concentration of 1  $\mu$ g/ml. SIKVAV (100  $\mu$ g/ml) stimulates the proliferation and migration of primary mouse fibroblasts, as well as induces the secretion of EGF, VEGF, TGF- $\beta$ , and FGF in the same cells.<sup>2</sup> Topical administration of a chitosan hydrogel containing SIKVAV promotes wound healing and angiogenesis in mice.

### References

- Freitas, V.M., Vilas-Boas, V.F., Pimenta, D.C., *et al.* SIKVAV, a laminin  $\alpha$ 1-derived peptide, interacts with integrins and increases protease activity of a human salivary gland adenoid cystic carcinoma cell line through the ERK 1/2 signaling pathway. *Am. J. Pathol.* **171(1)**, 124-138 (2007).
- Chen, X., Zhang, M., Chen, S., *et al.* Peptide-modified chitosan hydrogels accelerate skin wound healing by promoting fibroblast proliferation, migration, and secretion. *Cell Transplant.* **26(8)**, 1331-1340 (2017).

#### 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, 10/26/2022

#### 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