

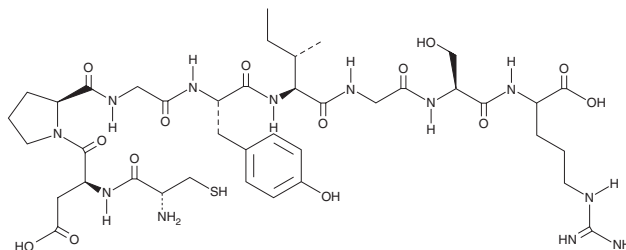
# PRODUCT INFORMATION



## Laminin (925-933)

Item No. 31141

**CAS Registry No.:** 110590-60-8  
**Formal Name:** L-cysteinyl-L- $\alpha$ -aspartyl-L-prolylglycyl-L-tyrosyl-L-isoleucylglycyl-L-seryl-L-arginine  
**Synonyms:** CDPGYIGSR, Cys-Asp-Pro-Gly-Tyr-Ile-Gly-Ser-Arg  
**MF:** C<sub>40</sub>H<sub>62</sub>N<sub>12</sub>O<sub>14</sub>S  
**FW:** 967.1  
**Purity:**  $\geq$ 98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

Laminin (925-933) is supplied as a crystalline solid. A stock solution may be made by dissolving the laminin (925-933) in the solvent of choice, which should be purged with an inert gas. Laminin (925-933) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of laminin (925-933) in these solvents is approximately 30 mg/ml.

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 laminin (925-933) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of laminin (925-933) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Laminin (925-933) is a peptide fragment corresponding to the laminin  $\beta$ 1 chain.<sup>1</sup> It binds to the laminin receptor when used at a concentration of 1 mg/ml. Laminin (925-933) stimulates the attachment of HT-1080 and CHO cells to culture plates when used at concentrations of 100 and 300  $\mu$ g/ml. It induces chemotaxis of, as well as inhibits chemotaxis induced by full-length laminin, but not fibronectin, in B16/F10 murine melanoma cells. Laminin (925-933) conjugated to an agarose hydrogel backbone enhances neurite outgrowth in isolated chick embryo dorsal root ganglia and PC12 cells.<sup>2</sup>

### References

1. Graf, J., Iwamoto, Y., Sasaki, M., *et al.* Identification of an amino acid sequence in laminin mediating cell attachment, chemotaxis, and receptor binding. *Cell* **48(6)**, 989-996 (1987).
2. Bellamkonda, R., Ranieri, J.P., and Aebischer, P. Laminin oligopeptide derivatized agarose gels allow three-dimensional neurite extension in vitro. *J. Neurosci. Res.* **41(4)**, 501-509 (1995).

#### 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, 12/13/2022

#### 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