

Vitronectin ACF

(Human Recombinant)

Manufactured Under
Animal-Free Conditions



Description

Vitronectin is a secreted glycoprotein that can support cell adhesion through binding to various integrins and proteoglycans. Recombinant vitronectin ACF can function as a chemically-defined matrix component for the attachment of human embryonic stem cell and IPS cells in a feeder-free culture system. Recombinant Vitronectin ACF is a 459 amino acid, single-chain, monomeric protein, which migrates at an apparent molecular weight of 75 kDa by SDS-PAGE under reducing conditions. The calculated molecular weight of Vitronectin ACF is 52.2 kDa.

Storage & Preparation

Vitronectin ACF is a lyophilized protein and should be stored at -20 to -80°C up to expiration date.

Preparation of Vitronectin ACF, 0.5mg/ml

Reconstitution procedure should be performed on ice.

1. Spin down the vial before reconstitution.
2. To the original 0.2mg vial add 0.4ml of sterile tissue culture water. **Do not vortex.**
3. Incubate on ice for a few minutes.
4. Gently mix by pipetting up and down.
5. Keep on ice for immediate use (up to 1 week at 2-8°C)

For long-term storage (3 months) aliquot and freeze at -80°C. Additional freeze-thaw cycles are not recommended.

Specifications

Sterile filtered and lyophilized.

Authenticity:

Verified by N-terminal and Mass Spectrometry analyses.

Purity:

≥ 95% by SDS-PAGE gel and HPLC analyses.

Endotoxin:

Endotoxin level is < 0.1 EU/μg of protein.

Protein Content:

Verified by UV Spectroscopy and | or SDS-PAGE gel.

Biological Activity:

Promotes attachment of hPSC in serum-free conditions.

Germany


Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

Israel

Biological Industries Israel Beit Haemek Ltd.
2511500 Kibbutz Beit Haemek
Phone: 972 4 9960595

 For further contacts, visit
www.sartorius.com