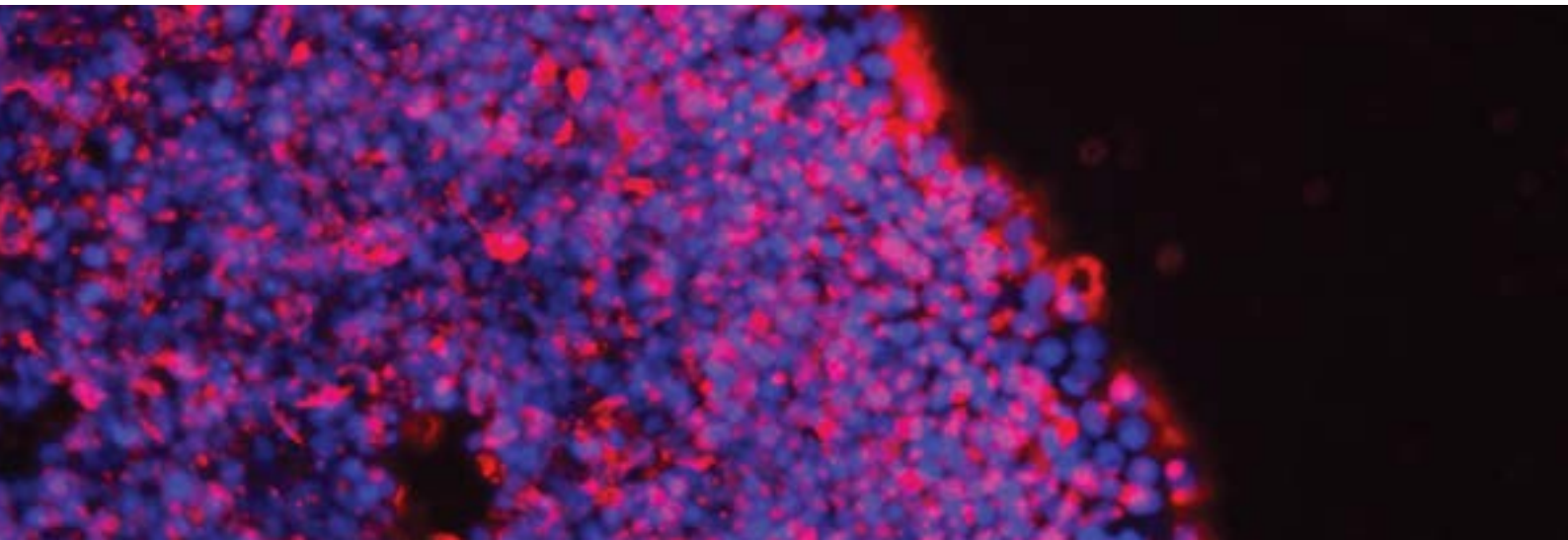




NutriStem[®] V9 XF

A superior, defined, xeno-free, serum-free culture medium for hPSC cells on vitronectin



- Superior proliferation rates in long-term culture on recombinant vitronectin
- Precoating-free protocol
- ROCK inhibitor-free protocol for seeding, passaging and thawing.
- Produced under cGMP
- Maintains high pluripotency at high passages
- Weekend-free feeding
- Cytokine-free basal medium, applicable for EBs formation, reprogramming, and differentiation

NutriStem[®] V9 XF is a defined, xeno-free, serum-free medium designed to support the growth and expansion of human pluripotent stem cells (hPSCs), on recombinant vitronectin matrice and enzyme-free passaging as small aggregates. NutriStem[®] V9 XF medium contains only the essential components required for long-term maintenance of hPSCs. This medium shows superior proliferation rates during long-term cultures, while maintaining the pluripotency of the cells. NutriStem[®] V9 XF medium allows the culture of hPSCs in vitronectin pre-coated culture, as well as the direct addition of vitronectin to the medium (no need for precoating).

NutriStem[®] V9 XF medium shows superior proliferation rates in long-term culture

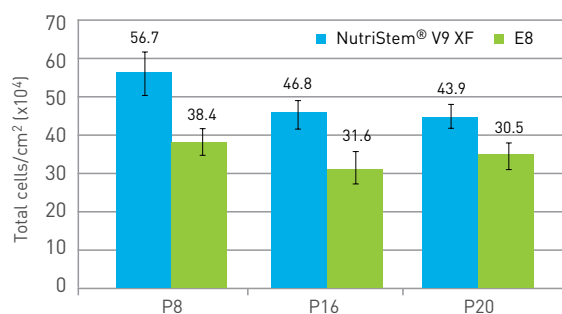


Figure 1: Nucleocounts performed on equal volume of cell suspension during long-term expansion of H1 cultured in NutriStem[®] V9 XF medium and competitor medium using 0.5µg/cm² Vitronectin ACF.

Normal cell morphology and assessment of pluripotency

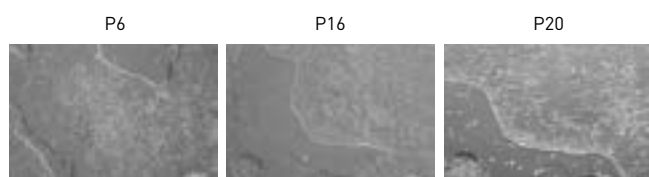


Figure 2: Phase contrast images (x100) of H1 hESC culture maintained in NutriStem® ACF medium, using 0.5µg/cm² Vitronectin ACF. Representative images from culture at P6, P16 and P20.

Undifferentiated colonies maintained typical hPSC colony morphology during long-term culture of H1 hPSC.

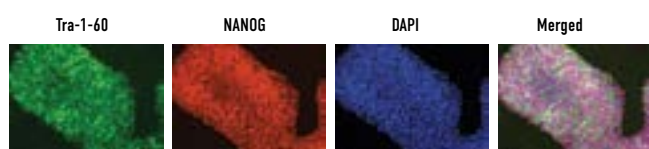


Figure 3: Immunofluorescence analysis of human pluripotent markers of H1 hPSC expanded in NutriStem® V9 XF medium using Vitronectin ACF. Cells from P8 were fixed and stained for the classic pluripotent surface markers: TRA 1-60 (green) and nuclear conjugated markers: NANOG (red) both counterstained with DAPI (blue). Scale bar 200µm.

Cells cultured in NutriStem® V9 XF medium on vitronectin express high levels of pluripotent markers.

Ordering Information

Cat.No.	Product Name	Size
05-105-1A	NutriStem® V9 XF basal medium	500 mL
05-106-1F	NutriStem® V9 XF supplement mix	1 mL
05-754-0002	Vitronectin ACF	200 µg
01-862-1B	EDTA Solution 0.5M	100 mL
05-713-1A	NutriFreez™ D10 Cryopreservation Medium	500 mL
05-713-1B	NutriFreez™ D10 Cryopreservation Medium	100 mL
05-713-1C	NutriFreez™ D10 Cryopreservation Medium	20 mL
05-713-1D	NutriFreez™ D10 Cryopreservation Medium	10 mL
05-713-1E	NutriFreez™ D10 Cryopreservation Medium	50 mL

How to Order

Biological Industries | T. 972-4-996-0595 F. 972-4-996-8896 | info@bioind.com

Biological Industries USA | T. 860.316.2702 F. 860.269.0596 | orders@bioindusa.com

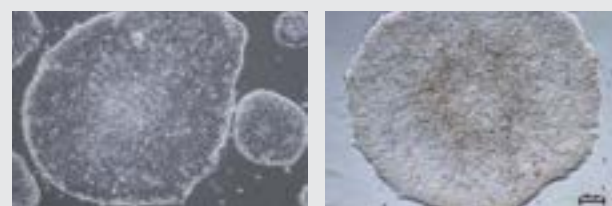
©2019 Biological Industries. All rights reserved. The trademarks mentioned herein are the property of Biological Industries and/or its affiliates or their respective owners.

Culture of Excellence

Precoating-Free Procedure

The use of a recombinant protein matrix, such as vitronectin, markedly improves stable maintenance of hPSCs. However, its use is time-consuming and laborious. Therefore, a user-friendly protocol has been developed to eliminate the precoating procedure.

While seeding, Vitronectin ACF is added directly into NutriStem® V9 XF medium, making pre-coating unnecessary.



A- H1

B- iPSC

Figure 4: H1 cultured in NutriStem® V9 XF medium for 6 sequential passages using a pre-coating free protocol. Vitronectin ACF was added directly to NutriStem® V9 XF medium before cell seeding. Representative images of H1 colony (A) iPSC ACS 1019 (ATCC) (B) colony morphology (x100). The pre-coating-free protocol supports classic colony morphology while maintaining high levels of pluripotent markers expression.

Immunofluorescence staining

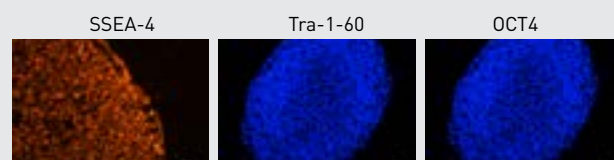


Figure 5: Immunofluorescence analysis of human pluripotent markers of H1 hESC expanded in NutriStem® V9 XF medium using a pre-coating-free protocol. Cells from P8 were fixed and stained for the classic pluripotent surface markers: SSEA4 (red), TRA 1-60 (green) and nuclear conjugated markers: OCT-4 (blue). Scale bar 200µm.

Cells cultured in NutriStem® V9 XF medium using the pre-coating-free protocol express high levels of pluripotent markers.



www.bioind.com