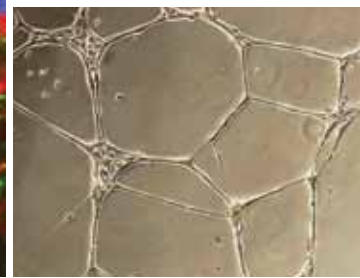
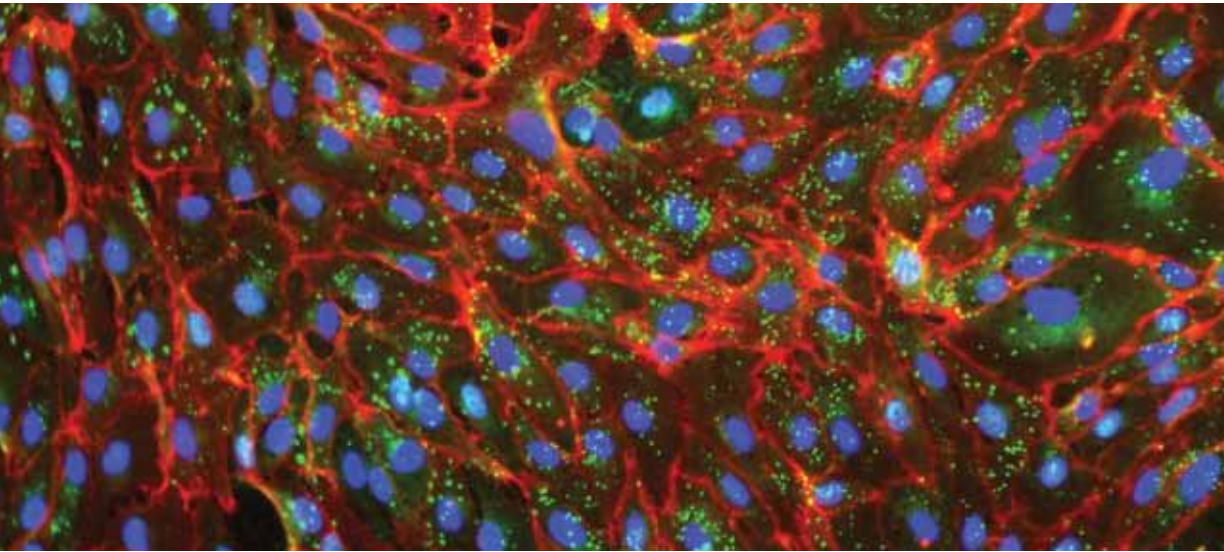




**BI**  
Biological Industries  
*Culture of Excellence*

# EndoGo™ XF Culture Media

A defined, xeno-free culture medium for the expansion of endothelial cells



## Expanding cell culture excellence and advancing clinical applications

- **Defined, xeno-free, serum-free medium**
- **Supports long-term expansion of large and small vessels endothelial cells**
- **Maintains high proliferation potential, typical morphology and EC marker expression**

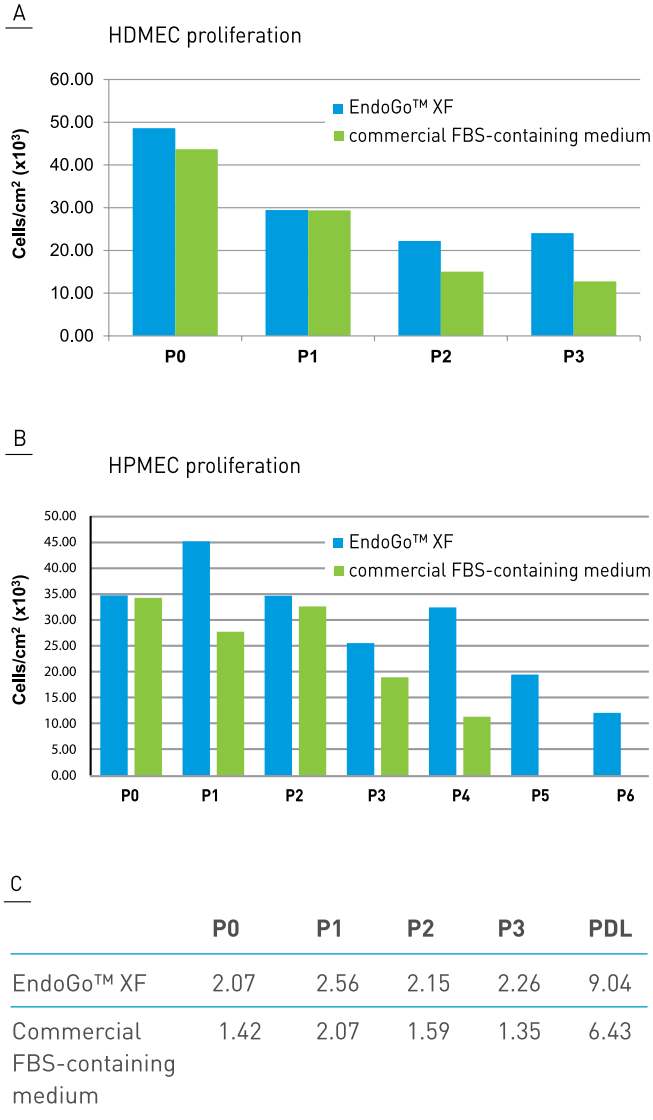
EndoGo™ XF is a novel XF culture medium specially designed for long-term expansion of large and small vessels EC from various sources. The medium provides optimally balanced nutritional environment that selectively promotes proliferation of normal human EC, while maintaining typical cobblestone-like cell morphology, phenotypic surface marker profile, and angiogenic differentiation potential. EndoGo™ XF supports microvascular EC (MVEC) from blood and lymph vessel as well as EC derived from: dermal, cardiac, lung, bladder and adipose tissues. In addition, EndoGo™ XF supports EC from arterial or venous (e.g. HUVEC).

No adaptation is required for the transition from bovine serum-containing medium to EndoGo™ XF.

Before use, supplementation with 2-5% of human AB serum (off the clot) or human platelet lysate (hPL) is required. It does not contain any non-human origin ingredients e.g., Bovine Brain Extract (BBE). It is recommended for use with Human Fibronectin Solution for optimal attachment, spreading and proliferation of cells. For optimal cell passage and long term culture of cells, it is recommended to use Recombinant Trypsin Solution with EDTA. For inhibition it is recommended to use Soybean Trypsin Inhibitor (SBTI).

## Cell Proliferation

Superior cell number and PDL of human microvascular endothelial cells in EndoGo™ XF.



**Figure 1:** Cell counts and population doubling level (PDL) of HDMEC (**A and C**), HPMEC (**B**) expanded for several passages in EndoGo™ XF in comparison to commercial FBS-containing medium. Viable cells were counted using ChemoMetec Viability and Cell Count Assay. \*HPMEC did not survive P5 in the FBS-containing medium.

## How to Order

**Biological Industries** | T. 972-4-996-0595 F. 972-4-996-8896 | [info@bioind.com](mailto:info@bioind.com)

**Biological Industries USA** | T. 860.316.2702 F. 860.269.0596 | [orders@bioindusa.com](mailto:orders@bioindusa.com)

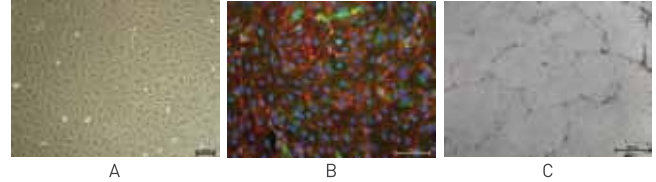
©2018 Biological Industries. All rights reserved. The trademarks mentioned herein are the property of Biological Industries and/or its affiliates or their respective owners.

*Culture of Excellence*

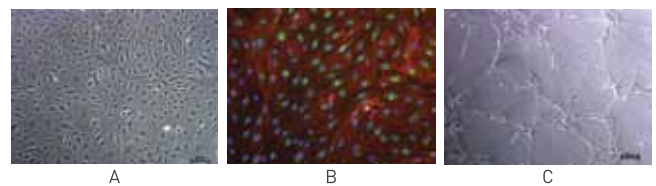
## Cell Morphology and Characteristics

EndoGo™ XF promotes proliferation of both micro and macro EC from a variety of sources while maintaining classical EC morphology, classical profile of EC markers, similar gene expression profile pattern, and similar angiogenic features of EC.

Macro vascular EC HUVEC



Micro vascular EC HDMEC



**Figure 2:** Microvascular EC (upper panel) and macrovascular EC (lower panel) maintain classic cobblestone like morphology after expansion for several sequential passages with equal seeding (5000 cells/cm²) in EndoGo™ XF +2% OTC human AB serum on hFN pre-coated dishes (A), expanded cells preserved endothelial cell features (EC markers expression) (B) and angiogenic potential to form capillary-like tubes (C).

## Ordering Information

Product	Cat. #	Qty
EndoGo™ XF	05-400-1A	500 mL
EndoGo™ XF Supplement Mix	05-410-1-25	2.5 mL
Human Fibronectin Solution	05-750-1H	5 mL
	05-750-1F	1 mL
Soybean Trypsin Inhibition	03-048-1C	20 mL
Recombinant Trypsin EDTA Solution	03-079-1B	100 mL
NutriFreez™ D10 Cryopreservation Medium	05-713-1A	500 mL
	05-713-1B	100 mL
	05-713-1C	20 mL
	05-713-1D	10 mL
	05-713-1E	50 mL



[www.bioind.com](http://www.bioind.com)