

# EndoGo™ XF Culture Media

A Defined, Xeno-Free Culture  
Medium for the Expansion  
of Endothelial Cells



## Benefits

- 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

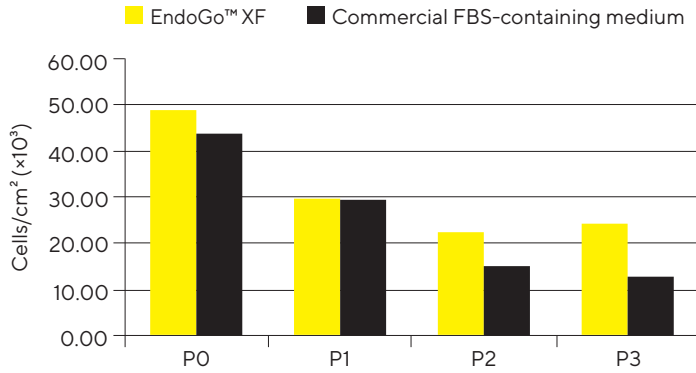
## Product Information

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.

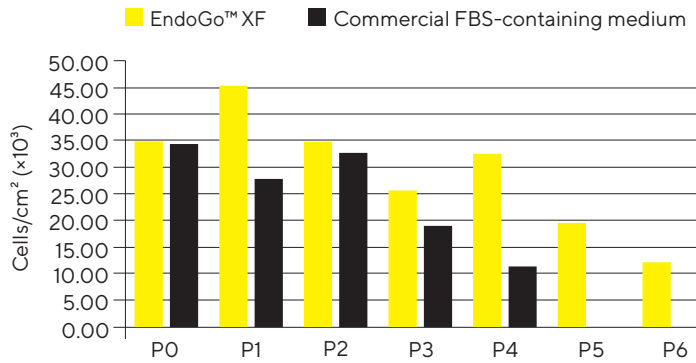
## Cell Proliferation

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

### A. HDMEC proliferation



### B. HPMEC proliferation



### C.

	P0	P1	P2	P3	PDL
EndoGo™ XF	2.07	2.56	2.15	2.26	9.04
Commercial FBS-containing medium	1.42	2.07	1.59	1.35	6.43

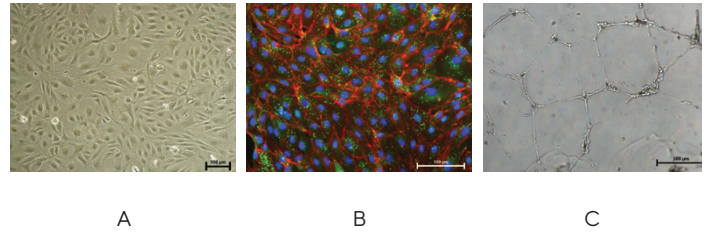
**Figure 1:** Cell counts and population doubling level (PDL) of HDMEC (A and C), HPMEC (B) expanded for several passages in EndoGo™ XF in comparison to commercial FBS-containing medium. Viable cells were counted using ChemoMetec Viability and Cell Count Assay.

\*HPMEC did not survive P5 in the FBS-containing medium.

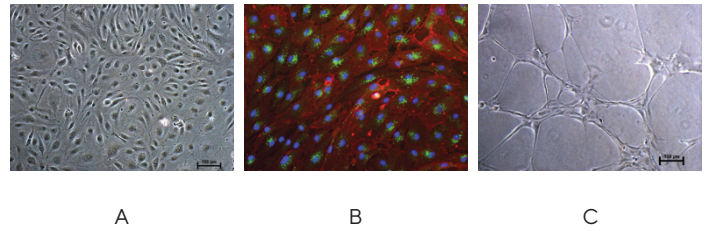
## Cell Morphology and Characteristics

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
NutriFreez® D10 Cryopreservation Medium without Phenol Red	05-714-1B	100 mL
	05-714-1E	50 mL
	05-714-1D	10 mL
NutriFreez® D5 Salt-based Cryopreservatiom Solution	05-715-1A	500 mL
	05-715-1B	100 mL
	05-715-1D	10 mL


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).

**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

 For further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)