

Recommended Media for Endothelial Cells



ENDOTHELIAL CELL TYPE	RECOMMENDED ENDOTHELIAL CELL MEDIA
HUAEC (C-12202)	Endothelial Cell Growth Medium (C-22010, C-22110) or
HPAEC (C-12241)	
HUVEC (C-12200, C-12203, C-12205)	Endothelial Cell Growth Medium 2 (C-22011, C-22111)
HSaVEC (C-12231)	
HUVEC (C-12206, C-12208)	Endothelial Cell Growth Medium 2 (C-22011, C-22111)
HAoEC (C-12271)	Endothelial Cell Growth Medium MV (C-22020, C-22120) or
HCAEC (C-12221)	
HCMEC (C-12285)	Endothelial Cell Growth Medium MV2 (C-22022, C-22121)
HUtMEC (C-12295)	
HPMEC (C-12281)	
HDMEC (C-12210, C-12212, C-12215)	
HDBEC (C-12211, C-12225)	Endothelial Cell Growth Medium MV (C-22020, C-22120)
HDLEC (C-12216, C-12217)	Endothelial Cell Growth Medium MV2 (C-22022, C-22121)

Recommended Media for Endothelial Cells

Frequently Asked Questions:

What's the difference between Endothelial Cell Growth Medium and Endothelial Cell Growth Medium 2?

The two media differ in their supplementation: Endothelial Cell Growth Medium contains Endothelial Cell Growth Supplement = ECGS (bovine hypothalamus extract). In Endothelial Cell Growth Medium 2, ECGS is replaced by VEGF, IGF and ascorbic acid and the concentrations of the other supplements were adapted.

When should I use Endothelial Cell Growth Medium, when should I use Endothelial Cell Growth Medium 2?

Endothelial Cell Growth Medium can be used for all endothelial cells that were isolated in Endothelial Cell Growth Medium and for most standard applications. If you prefer a more defined media composition, Endothelial Cell Growth Medium 2 is the medium of choice (2% FCS, no ECGS). If you plan to do VEGF stimulation, then Endothelial Cell Growth Medium is the recommended medium because the cells have not been in contact with stimulating concentrations of VEGF before.

What's the difference between Endothelial Cell Growth Medium MV and Endothelial Cell Growth Medium MV2?

The two media differ in their supplementation: Endothelial Cell Growth Medium MV contains Endothelial Cell Growth Supplement = ECGS (bovine hypothalamus extract). In Endothelial Cell Growth Medium MV2, ECGS is replaced by VEGF, IGF, basic FGF, EGF and ascorbic acid. A few other adaptations were made.

When should I use Endothelial Cell Growth Medium MV, when should I use Endothelial Cell Growth Medium MV2?

Endothelial Cell Growth Medium MV can be used for all (microvascular) endothelial cells that were isolated in Endothelial Cell Growth Medium MV and for most standard applications. If you prefer a more defined media composition, Endothelial Cell Growth Medium MV2 is the medium of choice (5% FCS, no ECGS). If you plan to do VEGF stimulation, then Endothelial Cell Growth Medium MV is the recommended medium because the cells have not been in contact with stimulating concentrations of VEGF before.

PromoCell GmbH

Sickingenstr. 63/65
69126 Heidelberg
Germany

Email: info@promocell.com
www.promocell.com

USA/Canada

Phone: 1 – 866 – 251 – 2860 (toll free)
Fax: 1 – 866 – 827 – 9219 (toll free)

Deutschland

Telefon: 0800 – 776 66 23 (gebührenfrei)
Fax: 0800 – 100 83 06 (gebührenfrei)

France

Téléphone: 0800 – 90 93 32 (ligne verte)
Téléfax: 0800 – 90 27 36 (ligne verte)

United Kingdom

Phone: 0800 – 96 03 33 (toll free)
Fax: 0800 – 169 85 54 (toll free)

Other Countries

Phone: +49 6221 – 649 34 0
Fax: +49 6221 – 649 34 40

© PromoCell GmbH