

Instruction Manual

Product	Size	Catalog Number
BPE-15 (7.5 mg/ml Protein)	2 ml	C-30020
	10 x 2 ml	C-30040
BPE-26 (13 mg/ml Protein)	2 ml	C-30021
	10 x 2 ml	C-30041
ECGS (3 mg/ml Protein)	2 ml	C-30160
	10 x 2 ml	C-30180
ECGS / H (3 mg/ml Protein + Heparin)	2 ml	C-30120
	10 x 2 ml	C-30140

Description

PromoCell Growth Supplements are produced from bovine brain tissue exclusively originating from New Zealand and thus classified as BSE-free. The Growth Supplements are aqueous protein extracts from mixed-sex bovine pituitary glands (BPE) or hypothalami (ECGS and ECGS/H). They are especially useful for the cultivation of different types of primary human cells under serum-reduced or serum-free conditions.

Bovine Pituitary Extract (BPE) -15 and -26 are standardized to a protein concentration of 7.5 mg/ml or 13 mg/ml, respectively. For Normal Human Epidermal Keratinocytes (NHEK) the optimal concentration of BPE-15 is 0.4% (v/v) in the medium supplemented with the appropriate growth factors and hormones. Other epithelial and non-epithelial cells need the higher concentrated BPE-26.

Endothelial Cell Growth Supplement (ECGS) is standardized to a protein content of

3 mg/ml. ECGS/H is additionally supplemented with 22.5 mg/ml Heparin, which enhances the growth promoting effect on endothelial cells. For Human Umbilical Vein Endothelial Cells (HUVEC) the optimal concentration of ECGS/H is 0.4% (v/v) in the medium supplemented with the appropriate growth factors and hormones.

Storage and Stability

Store Growth Supplements at -20°C immediately after arrival. After thawing, store at 4°C and use within one week. Avoid repeated freeze-thaw cycles. If stored properly, the product is stable until the expiry date stated on the label.

Note: Some precipitation might be observed after addition of PromoCell Growth Supplements to the medium. This does not affect the biological activity.

Quality Control

All lots of PromoCell Growth Supplements are subjected to comprehensive quality control tests. Each lot is routinely tested for its growth promoting activity using primary human cells, the absence of cytotoxicity, and physical parameters such as protein content. Approved in-house lots are used as a reference.

In addition, all lots of PromoCell Growth Supplements have been tested for the absence of microbial contaminants (fungi, bacteria and mycoplasma).

Intended Use

The products are for *in vitro* research use only and not for diagnostic or therapeutic procedures. For safety precautions please see appropriate MSDS.

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