

# Buffers and salt solutions

## Instruction Manual

Product	Size	Catalog Number
HEPES Buffered Saline Solution	30 ml	C-40000
	125 ml	C-40010
	250 ml	C-40020
Dulbecco's PBS	500 ml	C-40230
Dulbecco's PBS, w/o Ca <sup>++</sup> / Mg <sup>++</sup>	500 ml	C-40232

### Product Description

Buffers and Salt Solutions are frequently used in tissue culture, cell isolation procedures and molecular biology. For convenience, we offer ready-to-use Buffers and Salt Solutions. They are prepared from raw materials of the highest quality using tissue culture grade water.

### Storage and Stability

Store HEPES Buffered Saline Solution at -20°C immediately after arrival. Dulbecco's PBS

should be stored at 4 to 8°C and should not be frozen. Please note: C-40230 is light-sensitive and should be stored in the dark. If stored properly, the product is stable until the expiry date stated on the label.

### Quality Control

All lots of PromoCell buffers and salt solutions are subjected to comprehensive quality control tests. Each lot is routinely tested for absence of cytotoxicity, and physical parameters such as osmolality and pH level. Approved

in-house lots are used as a reference.

In addition, all lots have been tested for the absence of microbial contaminants (fungi, bacteria).

### 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.

## HEPES-BSS

Concentration [g/l]	HEPES-BSS C-40000   C-40010   C-40020
CaCl <sub>2</sub> anhydrous	—
KCl	0.22
MgSO <sub>4</sub> x 7H <sub>2</sub> O	—
NaCl	7.7
NaHCO <sub>3</sub>	—
Na <sub>2</sub> HPO <sub>4</sub>	0.14
NaH <sub>2</sub> PO <sub>4</sub> x H <sub>2</sub> O	—
D-Glucose	1.8
HEPES	7.15
Phenol Red	0.001

## Dulbecco's PBS

Concentration [g/l]	Dulbecco's PBS C-40230	Dulbecco's PBS C-40232 w/o Ca <sup>++</sup> /Mg <sup>++</sup>
CaCl <sub>2</sub> x 2H <sub>2</sub> O	0.132	—
KCl	0.2	0.2
KH <sub>2</sub> PO <sub>4</sub>	0.2	0.2
MgCl <sub>2</sub> x 6H <sub>2</sub> O	0.1	—
NaCl	8	8
Na <sub>2</sub> HPO <sub>4</sub>	1.15	1.15

**PromoCell GmbH**  
Sickingerstr. 63/65  
69126 Heidelberg  
Germany

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