

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 ⁺⁺ / Mg ⁺⁺	500 ml	C-40232

Description

Buffers and Salt Solutions are frequently used in tissue culture, cell isolation procedures and molecular biology. For convenience, PromoCell offers 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. Dul-

becco's PBS should be stored at 4–8°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.

Concentration [g/l]	HEPES-BSS C-40000 C-40010 C-40020
---------------------	--

■ Inorganic Salts

CaCl ₂ anhydrous	-
KCl	0.22
MgSO ₄ x 7H ₂ O	-
NaCl	7.7
NaHCO ₃	-
Na ₂ HPO ₄	0.14
NaH ₂ PO ₄ x H ₂ O	-

■ Other Components

D-Glucose	1.8
HEPES	7.15
Phenol Red	0.001

Concentration [g/l]	Dulbecco's PBS C-40230	Dulbecco's PBS C-40232 w/o Ca ⁺⁺ /Mg ⁺⁺
---------------------	---------------------------	---

■ Inorganic Salts

CaCl ₂ x 2H ₂ O	0.132	-
KCl	0.2	0.2
KH ₂ PO ₄	0.2	0.2
MgCl ₂ x 6H ₂ O	0.1	-
NaCl	8	8
Na ₂ HPO ₄	1.15	1.15

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