

Calcium Chloride Solution

Instruction Manual

Product	Size	Catalog Number
Calcium Chloride Solution (0.5 M)	2 ml	C-34006

Product Description

Some types of primary human cells, especially Normal Human Epidermal Keratinocytes (NHEK), respond sensitively to the calcium concentration in the culture medium. Calcium Chloride solution allows the adjustment of the optimal amount of calcium required by the cells. Our Calcium Chloride (CaCl₂) Solution is prepared as a solution from highly pure tissue culture tested Calcium Chloride and cell culture-grade water. It is a clear colorless and sterile solution. The concentration of CaCl₂ is 0.5 M.

Storage and Stability

Store at -20°C in the dark immediately after arrival. If stored properly, the product is stable until the expiry date stated on the label.

Quality Control

All lots of PromoCell Calcium Chloride Solution are subjected to comprehensive quality control tests. Each lot is routinely tested for the use with human primary cells, 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.

If you require special media modifications, we offer a custom media service starting at 10 bottles per order.
Contact us at info@promocell.com to find out more.

PromoCell GmbH
Sickingenstr. 63/65
69126 Heidelberg
Germany

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

info@promocell.com
www.promocell.com

© PromoCell GmbH