



# StrataCooler Cryo Preservation Module

## Instruction Manual

**Catalog #400005**

Revision B.0

**For Research Use Only. Not for use in diagnostic procedures.**

IN# 70081-00



**Agilent Technologies**

## LIMITED PRODUCT WARRANTY

This warranty limits our liability to replacement of this product. No other warranties of any kind, express or implied, including without limitation, implied warranties of merchantability or fitness for a particular purpose, are provided by Agilent. Agilent shall have no liability for any direct, indirect, consequential, or incidental damages arising out of the use, the results of use, or the inability to use this product.

## ORDERING INFORMATION AND TECHNICAL SERVICES

### Email

techservices@agilent.com

### World Wide Web

www.genomics.agilent.com

### Telephone

Location	Telephone
United States and Canada	800 227 9770
Austria	01 25125 6800
Benelux	02 404 92 22
Denmark	45 70 13 00 30
Finland	010 802 220
France	0810 446 446
Germany	0800 603 1000
Italy	800 012575
Netherlands	020 547 2600
Spain	901 11 68 90
Sweden	08 506 4 8960
Switzerland	0848 8035 60
UK/Ireland	0845 712 5292
All Other Countries	Please visit <a href="http://www.agilent.com/genomics/contactus">www.agilent.com/genomics/contactus</a>

# StrataCooler Cryo Preservation Module

## SPECIFICATIONS

Preservation module	Catalog #	Specifications		
		Capacity	Dimensions	Weight
StrataCooler Cryo preservation module	400005	32 standard cryovials	12 cm W × 21 cm D × 8 cm H	1 kg

## APPLICATION

---

Most cell lines of higher eukaryotes can be preserved indefinitely by storage at  $-185^{\circ}\text{C}$  in liquid nitrogen. This temperature ensures that the survival rate of the cells will not significantly change over years of cryopreservation.<sup>1</sup> Cells are prepared for freezing by suspension in special, tissue culture freeze media. These freeze media have a high protein content provided by fetal calf serum (FCS) and dimethylsulfoxide (DMSO) to prevent water crystal formation.

Prior to permanent storage in liquid nitrogen, the temperature of the cells must be lowered to approximately  $-80^{\circ}\text{C}$ . The proportion of cells that survive this process is determined by the freezing method used. Most cell types show  $>50\%$  survival when cooling rates are kept at  $0.2\text{--}1^{\circ}\text{C}/\text{minute}$ . The StrataCooler Cryo preservation module has been designed to freeze mammalian and insect cells at a controlled rate of  $0.4\text{--}0.6^{\circ}\text{C}/\text{minute}$  to achieve an  $80\text{--}90\%$  survival rate.

## PROTOCOL

---

**Note** *Before use, the StrataCooler Cryo preservation module must be cooled to 4°C by overnight refrigeration.*

1. Harvest the cells by spinning in a centrifuge. Resuspend the cells in ice-cold freeze medium (see *Medium*). Transfer the cell suspension to labeled cryovials.

**Note** *Most cell lines freeze well at 10<sup>6</sup> cells/ml. Freeze in 1-ml aliquots.*

2. Transfer the cryovials to the prechilled (4°C) StrataCooler Cryo preservation module.
3. Place the StrataCooler Cryo preservation module in a –80°C freezer.
4. After overnight incubation at –80°C, transfer the vials to liquid nitrogen for long-term storage.

## MEDIUM

---

### Freeze Medium

95% FCS

5% DMSO

Combine, filter sterilize and store at –20°C

## REFERENCE

---

1. Simione, F. P., and Brown, E. M., eds. (1991) "ATCC Preservation Methods," Second Ed. American Type Culture Collection, Rockville, Maryland.

## MSDS INFORMATION

---

Material Safety Data Sheets (MSDSs) are provided online at <http://www.genomics.agilent.com>. MSDS documents are not included with product shipments.