



## Electroporation cuvettes

Technical Data Sheet

References: CE-0001-50  
CE-0002-50  
CE-0004-50

### General Specifications:

<b>Material</b>	Polished Aluminium
<b>Sterility</b>	Gamma irradiated sterile
<b>Gap tolerance</b>	+/-2%
<b>Max electric field</b>	25kV/cm
<b>Packaging</b>	50, individually sealed

References	CE-0001-50	CE-0002-50	CE-0004-50
<i>Cap color</i>	Black	Yellow	Blue
<i>Gap (mm)</i>	1mm	2mm	4mm
<i>Min Volume (µl)</i>	10	20	80
<i>Max Volume (µl)</i>	90	400	800

### Shipping conditions

Shipping at ambient temperature

### Storage conditions

Storage at Room temperature.

### For any further information required please contact our Customer Help Desk:

*For Europe:*

E-mail: [info@eurogentec.com](mailto:info@eurogentec.com)

Tel: +32 4 372 76 65

*For USA:*

E-mail: [info.usa@eurogentec.com](mailto:info.usa@eurogentec.com)

Tel: +1 858 793 26 61



#### Eurogentec Headquarters

LIEGE SCIENCE PARK • 4102 SERAING • Belgium • Tel: +32 4 372 76 65  
Fax: +32 4 372 75 00 • [info@eurogentec.com](mailto:info@eurogentec.com) • [www.eurogentec.com](http://www.eurogentec.com)

#### Eurogentec North America

3347 Industrial Ct • Suite A • San Diego, CA 92121 • USA • Tel: +1 858 793 26 61  
858 793 26 66 • [info.usa@eurogentec.com](mailto:info.usa@eurogentec.com) • [www.eurogentec.com](http://www.eurogentec.com)