



Agilent 7100 CE

Short Electrode

This technical note provides general and usage information about the CE short electrode (p/n G7100-60033).

General Information

This short electrode is 5 mm shorter than the standard electrode (p/n G7100-60007) and is fitted into the instrument in exactly the same way as the standard electrode.

Instructions for removing and re-installing the electrodes can be found in the *Agilent 7100 Capillary Electrophoresis System User Manual* in the section "*Preparing Vials for Buffer and Sample*", chapter 5 "*Using the Agilent 7100 Capillary Electrophoresis System*".



Using the Short Electrode

- Fill the sample vials and ensure not to exceed 6 mm in the sample vials.
- Fill the run buffer vials to at least 1.8 cm.

Run buffer vials must be filled to at least 1.8 cm in order to ensure good electrical contact between the electrode and the buffer.

NOTE

No electrokinetic injection is possible when using the short electrode with the sample vials (filled to no more than 6 mm). This should not be attempted, since the electrode does not contact the sample under these conditions.



G7100-90118

Part Number:
G7100-90118

Edition: 11/2012
Printed in Germany

© Agilent Technologies, Inc 2012

Agilent Technologies, Inc
Hewlett-Packard-Strasse 8
76337 Waldbronn, Germany