

Changing the 16-Pin Bayonet Cartridge

Introduction

The 16-pin bayonet cartridge (reorder number 5065-4413), see [Figure 1](#), contains 16 electrodes. These are configured to fit in to the wells of an analysis chip.

The electrodes make contact with the liquid in the wells when the lid of the Agilent 2100 Bioanalyzer instrument is closed.

The cartridge, which includes the pin set, can be removed if the electrodes become contaminated or damaged.

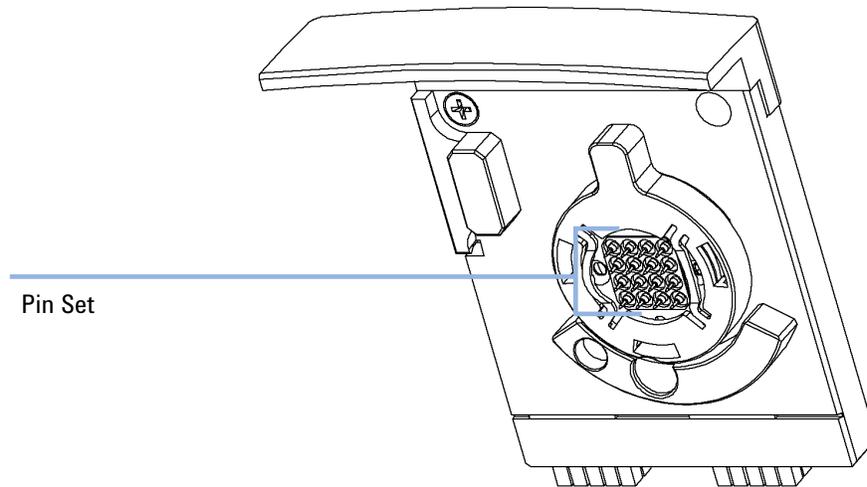


Figure 1 16-Pin Cartridge and Pin Set of Cartridge



Removing the 16-Pin Cartridge

CAUTION

Touching the electrodes

Possible damage to the electrodes and high voltage power supplies

→ Do not touch the electrodes while the cartridge is in the 2100 Bioanalyzer instrument.

The line switch is located at the rear of the 2100 Bioanalyzer instrument.

- 1 Turn off line power to the 2100 Bioanalyzer instrument.
- 2 Open the lid.
- 3 Pull the metal lever on the inside left of the lid to the vertical position as shown in [Figure 2](#).

When the lever is in the vertical position, the cartridge is released from the lid by about 10 mm.

- 4 Gently pull the cartridge out of the lid as shown [Figure 2](#).

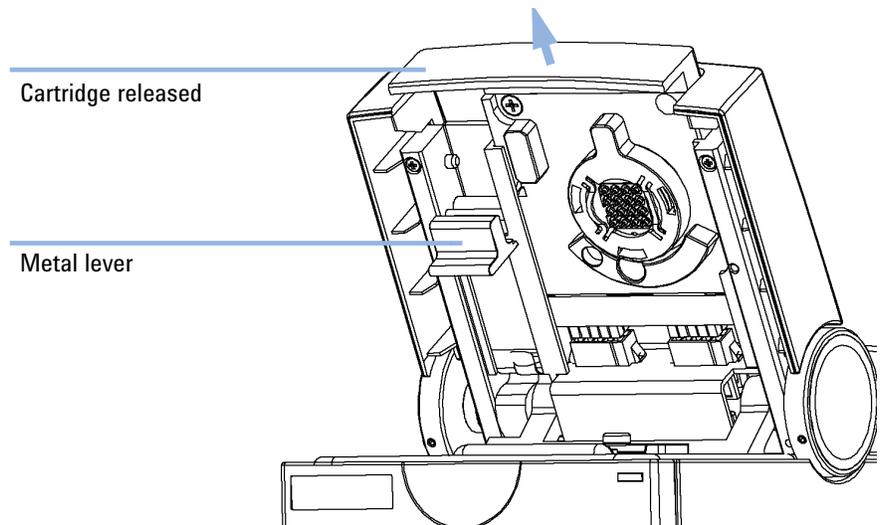


Figure 2 Removing the 16-Pin Cartridge

Removing the Pin Set of the 16-Pin Cartridge

- 1 Remove the 16-pin cartridge as described above.
- 2 Open the bayonet socket of the pin set by turning the plastic lever to the left as described in [Figure 3](#).

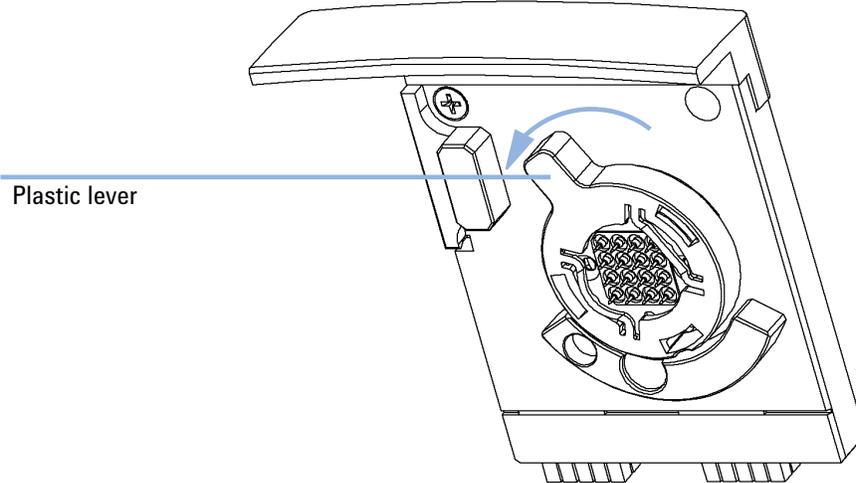


Figure 3 Opening the Bayonet Socket of the Pin Set

Removing the Pin Set of the 16-Pin Cartridge

- 3 Remove the cover of the bayonet socket by gently pulling the plastic lever as shown in [Figure 4](#). The pin set may stick to the electrode base. Remove it by carefully pulling it off. See [Figure 4](#).

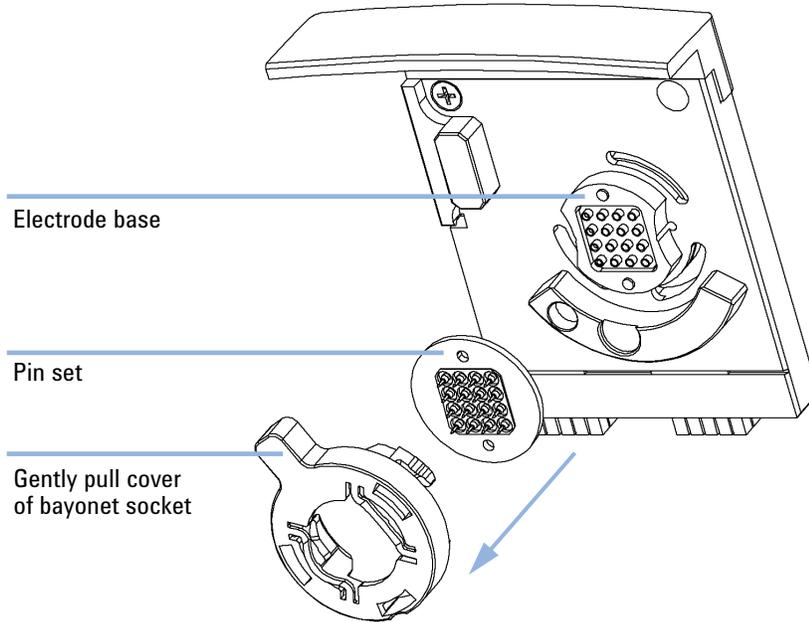


Figure 4 Releasing the Pin Set

NOTE

For hints on how to clean the pin set, refer to [“Cleaning the Pin Set of the 16-pin Cartridge”](#) on page 8.

Inserting the Pin Set of the 16-Pin Cartridge

CAUTION

Liquids (even small amounts) on the pin set

Possible damage to the high voltage power supply

→ Make sure that the pin set is completely dry before placing it back into the electrode base.

- 1 Put the pin set on the cartridge base and the bayonet cover on the pin set (see [Figure 5](#)).

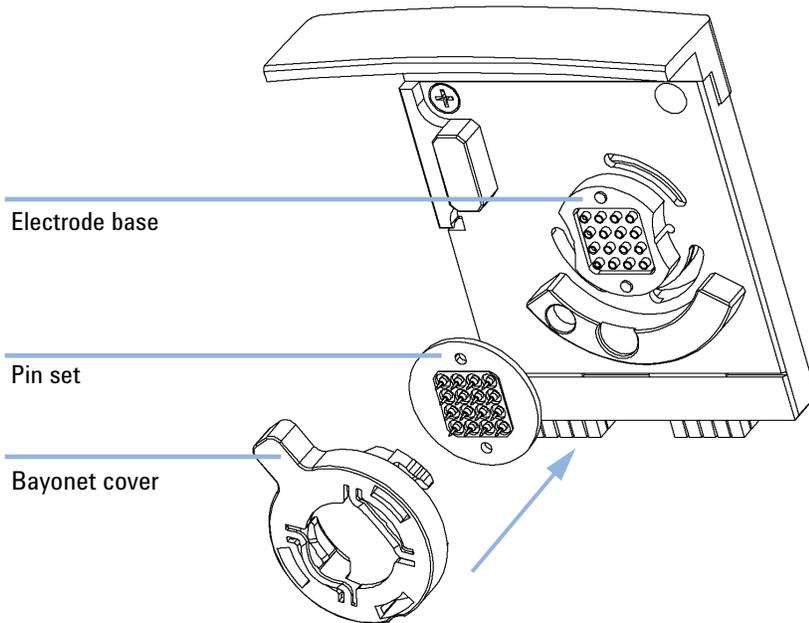


Figure 5 Inserting the Pin Set

Inserting the Pin Set of the 16-Pin Cartridge

- 2 Lock the pin set to the electrode base by pushing the plastic lever of the bayonet cover to the right as shown in [Figure 6](#).

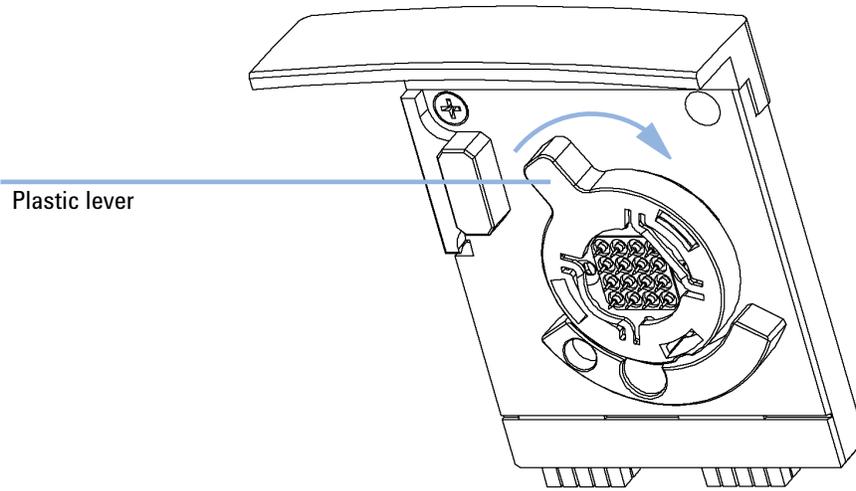


Figure 6 Closing the Socket of the Pin Set

Inserting the 16-Pin Cartridge

CAUTION

Liquids (even small amounts) on the pin set

Possible damage to the high voltage power supply

→ Make sure the pin set is be completely dry before putting in the 16-pin cartridge.

- 1 Slide the 16-pin cartridge into the 2100 Bioanalyzer lid as shown in [Figure 7](#).
- 2 Move the metal lever in the flat (closed) position.
- 3 Push the metal front of the 16-pin cartridge to ensure a tight connection to the 2100 Bioanalyzer.

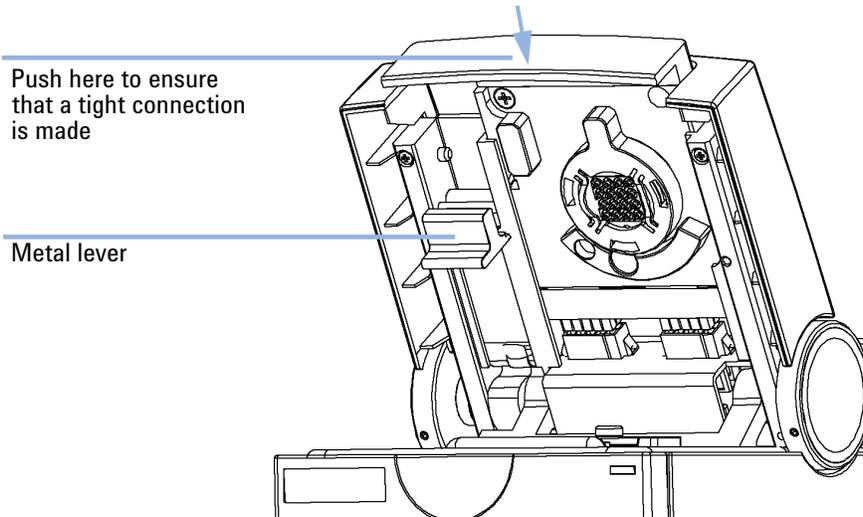


Figure 7 Inserting the 16-Pin Cartridge

Cleaning the Pin Set of the 16-pin Cartridge

After removing the pin set from the 16-pin cartridge you can clean it by using

- de-ionized water, or
- isopropanol, or
- RNase Zap.

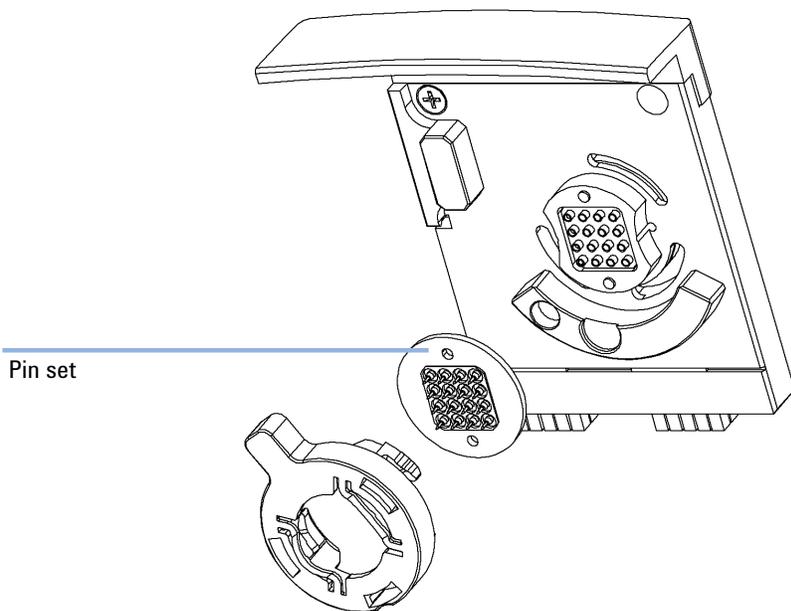


Figure 8 Pin Set

On a regular quarterly basis, or after contamination, gently clean the pin set with a lint-free surgical cotton swab damped in de-ionized water. In case of highly contaminated or dirty pins you may autoclave the pin set.

NOTE

The pins of the pin set should not be bent or misaligned. Both will lead to poor quality results or pre-terminated assay runs.

NOTE

For autoclaving the pin set, follow your standard procedures for plastic material.

CAUTION

- 1 Sonicate the pin set for 10 min.
- 2 Gently clean the pin set with a soft tooth brush.

Liquids (even small amounts) on the pin set
Possible damage to the high voltage power supply

→ Make sure the pin set is completely dry before putting in the 16-pin cartridge.

- 3 Insert the 16-pin cartridge, see [“Inserting the 16-Pin Cartridge”](#) on page 7.

