

How to setup a programmable electronic high power crimper

Author

Thermo Fisher Scientific

Keywords

Electronic crimper, vial caps

Introduction

This protocol explains the setup and testing of the Thermo Scientific™ Programmable Electronic High Power Crimper (60180-ECRHB, 60180-ECRH11KIT, and the 60180-ECRH20KIT). The crimper can be used with different size jaw sets (i.e., ECMJ-11, ECDJ-11, ECMJ-20, etc.) depending on the vial to be crimped.

Important notes

- The ECRH series of crimpers allows the use of different size jaws for work in a mixed vial environment
- The unit must be programmed and tested for each different type of vial used

Materials required

- ECRH
- Appropriate size jaws: crimping jaws, flip off jaws or decappers
- Test vial and caps
- Charger

Protocol

1. Assemble the stand according to the manufacturer's instructions provided with the unit.
2. Insert the correct jaw set into the unit from the bottom using a counter clockwise turn to lock the jaws into position. The button on the jaws should click into position and the jaws should be firmly locked into place.
3. Power up the unit by plugging in the cord from the power supply to the top of the crimper. When inserted properly it should click into place.
4. The configured jaw type last used on the crimper will appear in the display on the top of the unit.
5. Use the minus (-) button to scroll through the available types and choose the correct size of jaws. Use the plus (+) button to scroll through and select the correct jaw type (i.e., 20C for a 20 mm decapper) jaw set.
6. Click the 'crimp' button on the front of the unit to save the settings.
7. Place a crimp top on your vial and lift under the unit. Place the seal and cap on the vial and rest the tool on top of the cap. Squeeze the go-button lightly to engage the motor. This button must be held down until the crimp is complete. Hold the 'crimp' button until the crimping process completes and releases.
8. Check that the crimp is done. Crimp at least 5 vials to make sure the settings are correct. The top should be able to be slightly turned with a high force. If the septum has been deformed or shifted then over-crimping has been performed. If the top spins freely then the crimper needs to be adjusted to a higher setting.
9. To adjust the crimp setting, use the plus or minus button to increase or decrease the distance the crimp head moves. Retest with a new vial to determine if the correct crimping power has been applied.



Related products

| Description | Part Number |
|---|-------------|
| Accessory Base for Programmable High Powered Electronic Crimper | ECRH-STAND |
| Crimper Jaw Set, 8 mm, standard seals | ECMJ-8 |
| Crimper Jaw Set, 11 mm, standard seals | ECMJ-11 |
| Crimper Jaw Set, 13 mm, standard seals | ECMJ-13 |
| Crimper Jaw Set, 20 mm, standard seals | ECMJ-20 |
| Decrimper Jaw Set, 11 mm | ECDJ-11 |
| Decrimper Jaw Set, 13 mm | ECDJ-13 |
| Decrimper Jaw Set, 20 mm | ECDJ-20 |
| Flip-Off Crimper Jaw Set, 13 mm (for circular rim flip-off caps*) | ECMJ-13FO |
| Flip-Off Crimper Jaw Set, 20 mm (for circular rim flip-off caps*) | ECMJ-20FO |
| Flip-Off Crimper Jaw Set, no overhang, 20 mm | ECMJ-20FONO |

*Flip-Off caps need to be used with Flip-Off Jaw sets

Current versions of product instructions are available at separatedbyexperience.com/chromexpert

Learn more at

separatedbyexperience.com/chromexpert

ThermoFisher
SCIENTIFIC