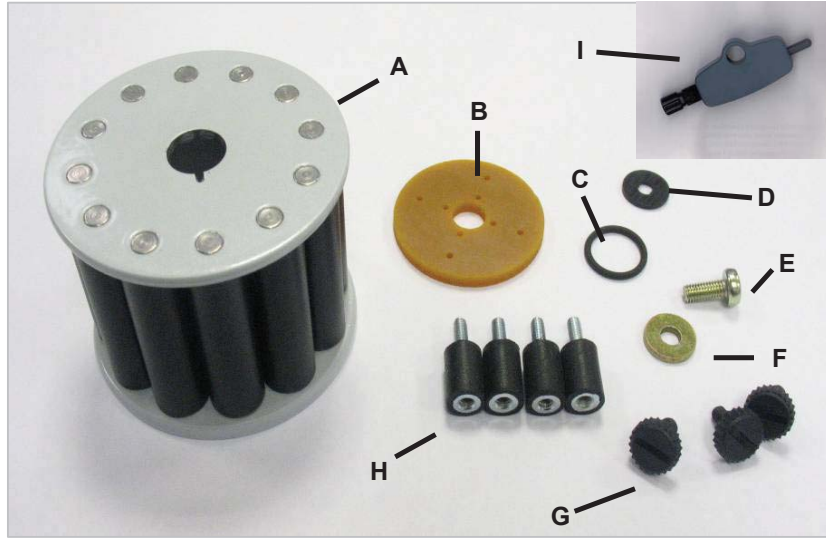


**3.10 iCAP 6000 & 7200 Pump Service Spares Kit**

3 channel pump kit	8423 155 50241
4 channel pump kit	8423 155 50251
Set of 4 rubber motor mounts	8423 155 50671

**Contents of the kits:**



- A** pump head (3 or 4 channel)
- B** rubber motor coupling
- C** o-ring for pump head
- D** nylon washer
- E** metal screw
- F** metal washer
- G** plastic screws x 3
- H** rubber motor mounts x 4
- I** pump tubing clamps x 2

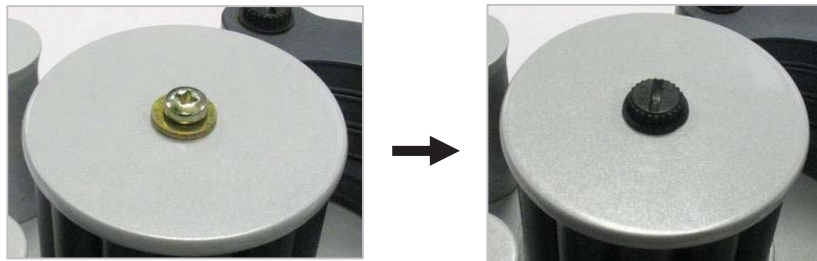
**How to change the pump head**



**VERY IMPORTANT:**

**Ensure that the cut-out on the pump head (1) is aligned with the key on the drive shaft (2).**

A rubber o-ring (C) should sit on the pump head mount before fitting the pump head.



Use the metal screw and washer (E & F) to pull the pump head into position.

**Warning:** Using the plastic screw at this stage may damage the plastic thread.

Once the pump head is in place remove the metal screw and washer and fit the plastic screw (G) with the nylon washer (D).

This page (1/1) is part numbered 4301 460 00141