

Cleaning Log

Serial Number: _____

Maxwell® FSC Instrument (Cat.# AS4600)

Clean the instrument as described in the *Maxwell® FSC Instrument Operating Manual* TM#462 following any spill or at regular intervals as appropriate for instrument usage. Record activities below.

Today's Date (DD/MM/YYYY): _____ **Time:** _____

User Name: _____

Please check off steps as they are completed:

- Disconnect from the Power Source.
- Remove Deck Tray.
- Clean Magnetic Rod Assembly.
- Clean Magnet Bar.
- Clean Plunger Bar.
- Clean Deck.
- Clean Interior.
- Clean Exterior and Door.
- Reinsert Deck Tray.
- Reconnect to Power Source.

Comments: _____

Today's Date (DD/MM/YYYY): _____ **Time:** _____

User Name: _____

Please check off steps as they are completed:

- Disconnect from the Power Source.
- Remove Deck Tray.
- Clean Magnetic Rod Assembly.
- Clean Magnet Bar.
- Clean Plunger Bar.
- Clean Deck.
- Clean Interior.
- Clean Exterior and Door.
- Reinsert Deck Tray.
- Reconnect to Power Source.

Comments: _____
