



# Cleaning Log

Maxprep™ Instrument (Cat.# AS9100, AS9101, AS9200, AS9201)

Clean the instrument as described in the Maxprep™ Instrument Operating Manual TM#509 following any spill or at regular intervals as appropriate for instrument usage. Record activities below.

Serial Number: \_\_\_\_\_

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

User Name: \_\_\_\_\_

Please check off steps as they are completed: Comments: \_\_\_\_\_

- Power off the Instrument
- Disconnect from the Power Source
- Remove Deck Trays and any Consumables
- Remove all Carriers with Front Handles and clean (one at a time)
- Clean Deck (as each Carrier is removed)
- Empty and Clean Tip Waste Bin
- Clean Carriers and Tip Holders
- Clean Pipettor Tips
- Clean Other Interior Surfaces
- Clean Exterior Panels and Door
- Reconnect to Power Source

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

User Name: \_\_\_\_\_

Please check off steps as they are completed: Comments: \_\_\_\_\_

- Power off the Instrument
- Disconnect from the Power Source
- Remove Deck Trays and any Consumables
- Remove all Carriers with Front Handles and clean (one at a time)
- Clean Deck (as each Carrier is removed)
- Empty and Clean Tip Waste Bin
- Clean Carriers and Tip Holders
- Clean Pipettor Tips
- Clean Other Interior Surfaces
- Clean Exterior Panels and Door
- Reconnect to Power Source

Download this form at: [www.promega.com/IQOQForms](http://www.promega.com/IQOQForms)