

Cleaning Procedure:

This general-purpose optics cleaning kit can be used to clean the following items in a Bio-Rad Fluor-S and GEL DOC, Insta Doc systems:

1. The Sample Platen, Filters on the UV Lamps/transilluminator
2. Emission filters the IR filter on the lens and Lenses and Lamps

This kit consists of:

1. A general purpose Industrial Optical Cleaner for Plastic Optics
2. Lens Cleaning paper.

PROCEDURE:

1. To clean Quartz platen, Transilluminator filter, UV filter on lamps:

Squirt some cleaner on the surface and then wipe it dry with a folded lens tissue.

2. To clean filters, lenses and lamps:

WARNING: These are coated optics and can be scratched easily. All precautions must be exercised in cleaning these optics. Clean both surfaces of these optics in the same way.

- This procedure requires use of powderless gloves.
- Fold the lens tissue four times using a pair of hemostats or forceps.
- Put some cleaner on the tissue. Shake off the extra cleaner.
- Hold the optical surface to be cleaned at an angle so that its surface can be easily seen under room lights.
- Gently clean the surface using the wet soft cushion made from the folded tissue and the cleaner and wiping in strokes in the same direction every time.
- Allow the glass surface to dry. Make sure that no streaks are visible on the surface of the filter.

After cleaning reinstall the optics and close the cap on the optic cleaner and store both the Cleaner and the Lens Tissue in a clean and dry location.

NOTE: A 50/50 mixture of Reagent/HPLC grade Methanol and Acetone can be used in place of the Plastic Optic Cleaner. The only item you can not clean with this Acetone Methanol mixture is the PLASTIC filter on the Scanning White light bulb.