
UNPACKING INSTRUCTIONS

The VersaDoc components come packaged in smaller boxes that are then packed in a large reusable container. Please follow the steps below to unpack the system:

1. Remove top of the container.
2. Remove the packing material and smaller boxes from within the container.
3. Remove the separator sheet and top foam cap.
4. Remove the outer sleeve.
5. Remove the plastic bag.
6. Lift the system out of the bottom the container and place it on the floor.
7. Carefully open all the boxes and take an inventory of what has been received.

The following is a list of items that will be found in the package(s):

1. VersaDoc System.
2. Camera Assembly.
3. Camera Electronics Unit [CEU] (for VersaDoc 5000 systems) with power cord.
4. Camera Power Supply(for VersaDoc 3000 systems) with power cord.
5. UV EPI lamp modules (left and right).
6. White Light EPI lamp modules (left and right).
7. Chemi Sample Tray.
8. Fluorescence Reference Plate
9. UV Shield.
10. Nikon 50mm Lens.
11. Tamron Zoom Lens.
12. Accessories Box:
 - a) Filters:
 - Large (660nm)
 - Filter # 1 (520nm)
 - Filter # 2 (530nm)
 - Filter # 3 (610nm)
 - Filter # 4 (clear)
 - 52mm to 58mm Adapter Ring
 - b) Cables:
 - 8000247 (AIA Cable)
 - 9210780 (DC Power Cable – To Camera)
 - Serial Cable (DB9 to DB9)
 - USB Cable (type A to type B)
 - Mac Serial Cable (USB to Serial)
 - Power Cable (US, Canada)
 - c) Miscellaneous:
 - Sample Holders (4)
 - Lens and Filter Cleaning Kit

- Focusing Target

13. Documentation:

- Warranty Card
- Users Manual
- Declaration of Conformity
- Hardware and Software Acceptance Sheet
- CD (Camera Specification File)

We suggest the customer store and reuse the packaging if there is a need to move or ship the instrument again.

Refer to the Users Manual for assembly instructions.