
UNPACKING INSTRUCTIONS

The Fluor-S MultiImager components come packaged in smaller boxes which are then packed in a large reusable container. Please follow the steps below to unpack:

1. Remove top and sides of container. Remove securing nails if possible to allow for reuse of shipping container.
2. Remove the packing material and smaller boxes from within the container.
3. Remove the straps holding the brace which straddles the Fluor-S. Carefully remove the brace.
4. 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:

1. Fluor-S main case
2. Protective foam pad (located inside of main case)
3. White plastic platen (located inside of main case)
4. Fluor-S Tower Assembly with camera
5. Filters:
 - Large (660 nm)
 - Filter #1 (520 nm)
 - Filter #2 (530 nm)
 - Filter #3 (610 nm)
 - Filter #4 (clear)
6. Tamron Zoom Lens
7. Cables:
 - 8004952 (8 pin DIN cable)
 - 8004938 (15 pin motor control cable) Note Warning Label
 - 8004953 (9 pin camera power cable)
 - 8004951 (15 pin EPI power cable)
 - 8004950 (68 pin HD camera cable)
 - 9007622 (Power cable)Note: Two EPI cables (8004952) and one ground cable (8004974) are attached to Fluor-S
8. Documentation:
 - Warranty Card
 - Instruction Manual
 - Declaration of Conformity
 - Hardware and Software Acceptance Sheet
 - Diskette (Camera dark pixel map)
9. Two empty boxes

Advise the customer to store and reuse the packaging if there is a need to move or ship the instrument again. A part number for the entire shipping container is included in the spare parts section of the Service Manual.