

S3e™ Individual Filter Holder

Catalog #	Description
144-1007	S3e Individual Filter Holder , pkg of 1, empty filter holder with lens retaining ring

For research purposes only.

Intended Use

The S3e Individual Filter Holder is used with the S3™ or S3e Cell Sorter to customize filter combinations for experimental fluorophore panels. The filter holders are designed to be easily accessible on the instrument for the quick exchange of band pass filters or for adding neutral density filters to block light. They are compatible with commercially available filters and installation is quick and easy. Unmounted 1 inch or 25 mm filters are recommended.

Contents

- Filter holder
- Lens retaining ring

Tools Required

- Spanner wrench (included in S3e Accessory Kit, catalog #145-1065)



Warning and Precautions

The filters are very delicate and require the utmost care in handling. These filters are made of glass and can cause injury, especially if they are cracked or shattered during the installation procedure.

If the filters are incorrectly installed, light transmission will not be directed to the correct channel. **In most cases, the labeled filter side is installed facing the incoming light.**

Note: Please refer to the filter vendor instructions for filter orientation.

Directions for Filter Installation

1. Set the spanner wrench into the S3e Individual Filter Holder and ensure that the notches align with the retaining ring inside the holder.
2. Unscrew counterclockwise until the retaining ring disengages from the filter holder.



S3e Individual Filter Holder and lens retaining ring.

3. Carefully handling the filter only by its edges, install the filter by placing it in the holder with the arrow indicator or lettering facing the source of emitted light.
4. Secure the retaining ring back in place by using the spanner wrench and rotating clockwise. (It may help to set the ring in place lying flat in the opening before using the spanner wrench.)
5. Rotate the spanner wrench until the retaining ring fits gently against the filter without overtightening.
6. Verify that everything is tight and does not rattle.

Visit [bio-rad.com/S3eFilterHolder](https://www.bio-rad.com/S3eFilterHolder) for more information.