

**ORDER FORM
EMISSION FILTER
FLUOR-S MULTI-IMAGER**



Omega series filters (see website www.omegafilters.com for more information)

Mechanical Specifications

Laminated construction edges sealed.
Enclosure: Black anodized aluminum ring (921-0136).
Diameter: 25.4 +0/-0.25mm
Thickness: 4.4 ± 0.1mm without ring
4.8 ± 0.1mm with ring
Clear aperture: 21.4+0.1/-0mm
TWFD: ≤ Lambda / 4 per 25mm

For dual imaging applications filters must be ordered in matched sets

Max. difference in thickness between filters (without ring) in a set: 0.1mm

Check this box for each dual imaging filter ordered.

Optical Specifications

Wavelength Type (Check one)
(>450 nm) BP (Bandpass) WB (Wideband)
 DF (Discriminating) CF (Cut-on) EF (Edge)

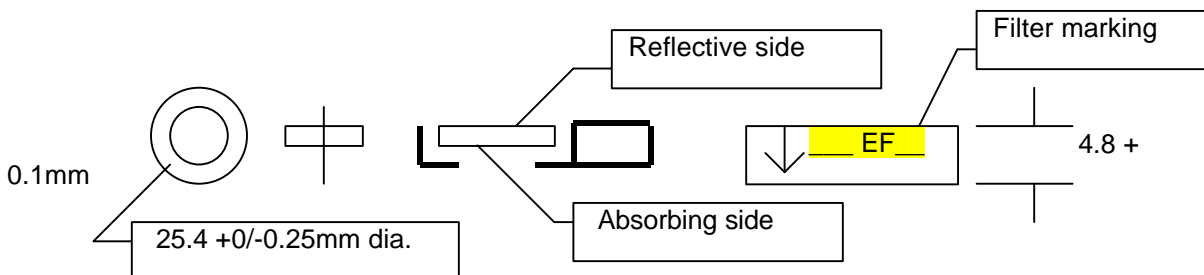
50% Transmission Point: ± 10nm

Transmission Band: Minimum to nm @ > 80% Avg.

Blocking: Minimum O.D. > 5.0 from cuton to UV

Edge slope: Per Omega published data for EF series.

Each filter whose construction incorporates absorbing glass shall be oriented with the absorbing glass facing the rim of the aluminum ring. This orientation shall be clearly marked on the edge of the holder with an arrow. The head shall point toward the absorbing glass. The filter shall be used in an orientation with light going in the direction of the arrow. Thus, the light shall first strike the side furthest away from the absorbing (and possibly fluorescing) glass and autofluorescing shall be minimized.



FAX order form to:
Omega Optical 3 Grove St. P.O. Box 573 Brattleboro, VT 05301
Phone: 802/254-2690 **FAX 802/254-3937**
Bio-Rad order form 410-0167 Rev. A