

ORDER FORM  
EMISSION FILTER  
Ultramark Microplate Reader



Omega series filters (see website [www.omegafilters.com](http://www.omegafilters.com) for more information)

Mechanical Specifications

Laminated construction edges sealed.  
Enclosure: ABS plastic ring.  
Diameter: 11.9 +0/-0.25mm without ring  
14.0 +0/-0.25mm with ring  
Thickness: 4.9 + 0.1mm without ring  
4.9 ± 0.1mm with ring  
Clear aperture: 7.0+0.1/-0mm  
TWFD:  $\leq \text{Lambda} / 10$

Optical Specifications

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

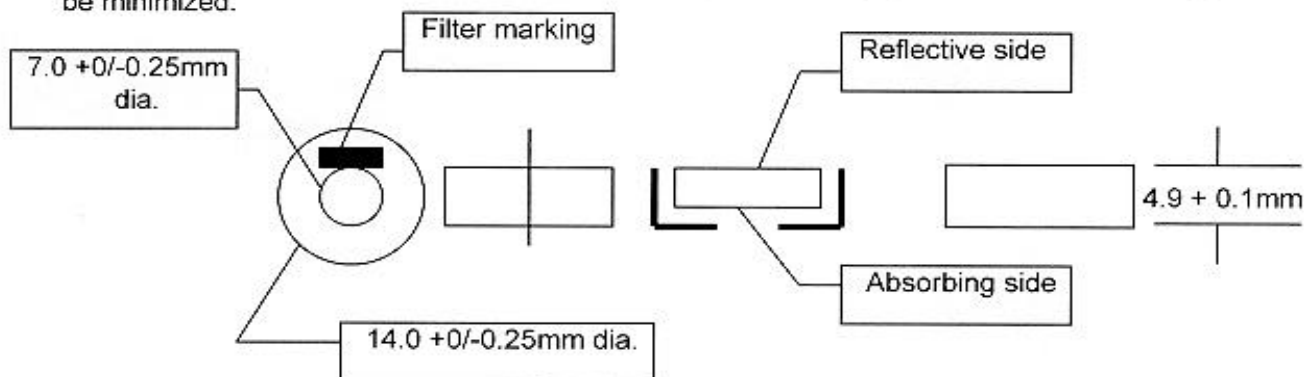
50% Transmission Point: \_\_\_\_\_ ± 10nm

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

Blocking: Minimum O.D. > 5.0 from cut-on 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 410-0170 Rev. A