

Fluor-S Multimager Quick Reference Guide

Step 1. Focus lens by aligning the two focusing arrows on the lens.

Note: When using the zoom lens at a zoom setting lower than 28 mm, the lens must be focused manually.

Step 2. Position sample on the platten.

Step 3. Refer to guide below for recommended f-stop, light source, filter and integration time settings.

<u>Sample</u>	<u>Recommended f-stop¹</u>		<u>Light source</u>	<u>Filter</u>	<u>Int. Time²</u>
	<u>Zoom lens</u>	<u>50mm lens</u>			
Ethidium Bromide	f-4	f-2.8	Scan:UV	520 LP	5-30 sec.
SYBR Green	f-4	f-2.8	Scan:UV	530 BP	5-30 sec.
SYPRO Orange	f-4	f-2.8	Scan:UV	520 LP	5-30 sec.
Radiant Red	f-4	f-2.8	Scan:UV	610 LP	5-30 sec.
Fluorescein	f-4	f-2.8	Scan:UV	530 BP	30 sec. - 2 min.
Texas Red ³	f-4	f-2.8	Scan:UV	610 LP	30 sec.- 2 min.
Colorimetric Stains ⁴ e.g. Coomassie or Silver	f-11	f-11	Scan:White light	Clear	1-5 sec.
X-ray film ⁴	f-11	f-11	Scan:White light	Clear	1-5 sec.
Photographs	f-11	f-11	Epi:White light	Clear	1-5 sec.
Chemiluminescence ⁵		f1.8	Chemi:No light	Clear	1-10 min.

Notes:

1. A slight improvement in image sharpness may be noted by closing the f-stop down one or two stops from full open.
2. Increase integration time two fold for each step increase in f-stop.
3. When using Multi-color labels (FITC, Texas Red), the 610 filter requires approximately twice the integration to obtain the same signal as the other filters.
4. For stained gels or x-ray film use the white backing plate.
5. Chemiluminescent samples- The 50 mm lens is recommended. Always remove the 660 filter.