



The IVIS® will take a snapshot of the bioluminescent or fluorescent source and record the counts. Sensitivity settings used for the snapshot image will be adjusted in order to obtain the desired user definable **Target Count Minimum**. You can define separate targets for bioluminescence, epi-fluorescence and trans-fluorescence images. It is advantageous to raise the Minimum Target Count at or around 30,000 counts as to bring out low intensity signals (e.g., finding small metastases in a mouse with a large primary tumor, imaging multiple mice for high throughput). Raising the target above 30,000 counts, though, will increase the chance of oversaturation.

Sensitivity settings will be adjusted beginning with your **First Preference** until the target count level has been attained. If the target level cannot be closely approximated by adjusting the first preference, the software also uses the second and then third preferences to attempt to reach the target level during image acquisition using the user defined settings on the control panel as a starting point.

Default settings are, Exposure time of 0.5 – 60 sec. (max. of 600 sec.) for First Preference and Binning of 1-8 (Small binning = 4, Medium = 8, Large = 16) for the Second Preference. Third Preference defaults to f/Stop ranging from 1 to 8.

