

Adjusting the Horizontal Offset for Hamamatsu C4880 Cameras

Document ID
T10155

Created
29-Sep-1998

Abstract

Troubleshooting: Hamamatsu C4880 cameras.

Instructions

Some Hamamatsu C4880 cameras may exhibit a horizontal shift, which appears as a black stripe down the left or right side of the image. This shift will vary in width depending on the binning. To correct this shift, you can adjust some parameters in the c4880xxx.cam file in the mm\vinput\mutech\camera directory.

If you look at the very bottom of the camfile, there is a section named [HOffset] with different values. These are defined for each specific binning parameter. This model of the camera requires different offsets for each binning value. The values listed below are for the standard 4880-10. You'll need to adjust these by trial and error. These will get refreshed each time you change binning - so there is no need to open/close MetaMorph.

To adjust these:

1. Open MetaMorph
2. Open c4880xxx.cam file with Notepad

To adjust 1x1 binning offset

1. Edit 1=X value and Save file in Notepad
2. Define Acquisition Settings - change binning from 1x1 to 2x2 and back to 1x1: this will read the new HOffset parameter.
3. Acquire image and then repeat 3-4 until edge is diminished.

To adjust 2x2 binning offset

1. Edit 2=X value and Save file in Notepad
2. Define Acquisition Settings - change binning from 2x2 to 1x1 and back to 2x2: this will read the new HOffset parameter.
3. Acquire image and then repeat 3-4 until edge is diminished.

Continue until all of the binning ranges that the camera supports have been adjusted. The default values for the C4880-10 are:

[HOffset]

1=68

2=57

4=52

8=50

16=0

Keywords: hamamatsu hardware

Issue Type: Hardware