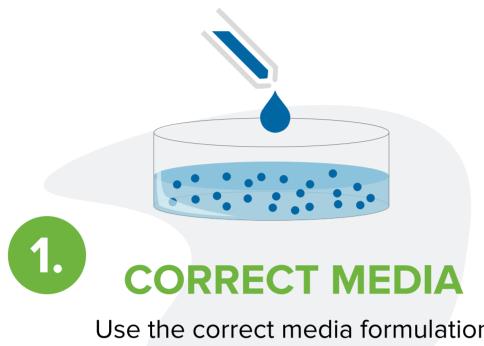
# Tips for running a successful live cell imaging experiment

Matthew Hammer | Cellular Imaging Applications Scientist







Use the correct media formulation when preparing your samples



Control and maintain proper environmental conditions when imaging your samples



# **HCA SOFTWARE**

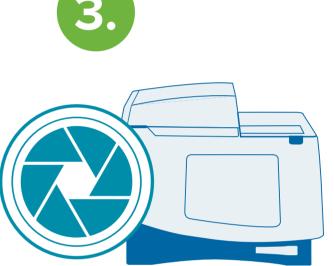
Utilize powerful high-content analysis (HCA) software and machine learning to obtain reproducible and meaningful data



Tips & techniques to maintain

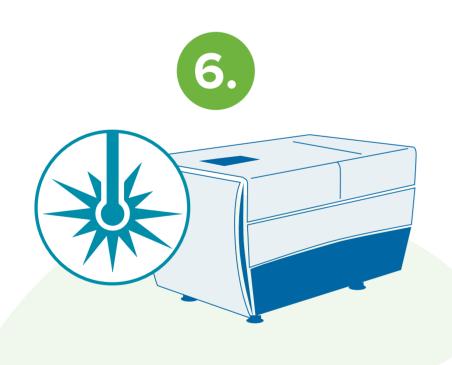
sample integrity and acquire

high resolution cell images



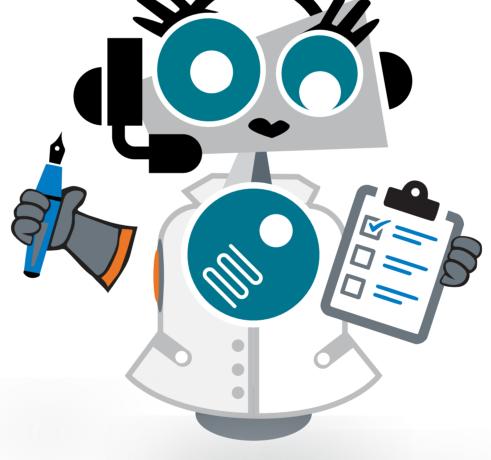
### **AUTOFOCUS**

Use robust autofocus modes during image acquisition



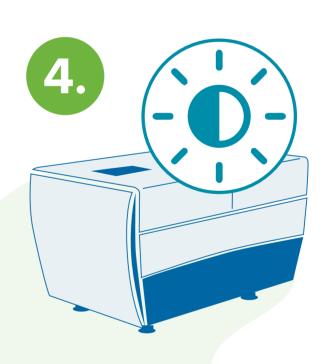
## LASER ILLUMINATION & BINNING

Use confocal imaging systems that allow for the use of high-power laser light sources and binning



#### **DECONVOLUTION**

If appropriate, employ image deconvolution algorithms when using widefield microscopy systems



### **EXPOSURE**

Minimize the power of the illuminating light source as well as your exposure time during image acquisition

Read the full post on our blog

