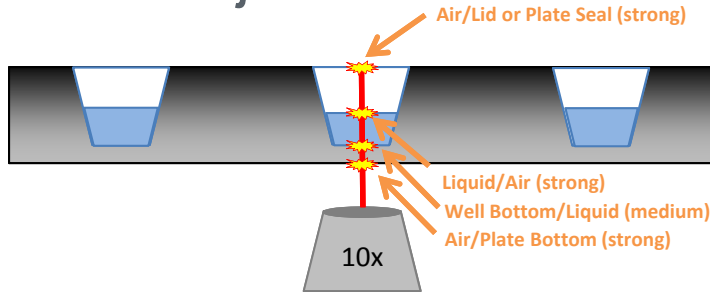


Laser Autofocus on Plates and Slides with Air and Oil Objectives

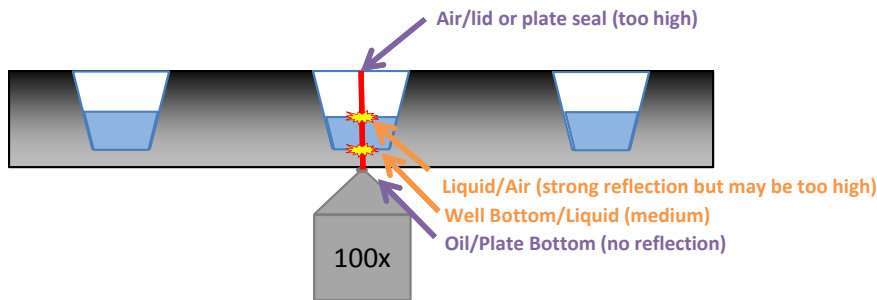
The laser autofocus mechanism detects a reflection of the laser spot where there is a change in refractive index (e.g. glass-air transition)

Typical Multi-well Plate

- Air objective

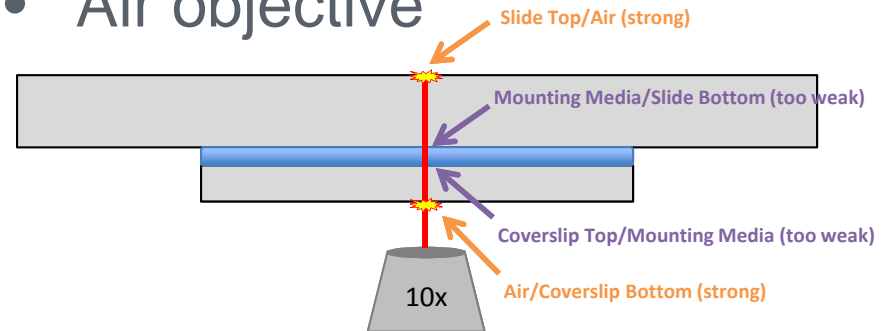


- Oil objective

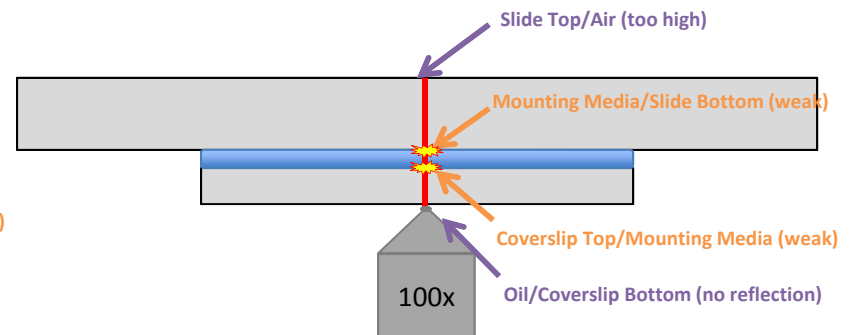


Typical Slide / Coverslip

- Air objective



- Oil objective



Diagrams are not to scale

