

Clone Select Imager Training Guide

Clone Select Imager Hardware Overview



Date Revised 03/22/2017 Version C

Index

- Index
- [Chapter Purpose](#)
- [Clone Select Imager Specifications](#)
- [Indicator Lights](#)
- [Plate Carrier](#)
- [Computer Workstation](#)
- [Instrument Power & Ethernet Connections](#)
- [Instrument Power Up Procedure](#)
- [Support Resources](#)



Chapter Purpose

The purpose of this chapter is to orient the user to features of the CloneSelect Imager hardware and to the power up procedure.

This guide does not include detailed descriptions around sample imaging or analysis. Please refer to corresponding chapters for details on these topics.



CloneSelect Imager Specifications

- Accommodates plate formats from 384-well to single well
- Windows 7 and XP compatible software
- Illumination: Xenon flash
- Imaging: 16-bit cooled CCD camera
- Standard pixel resolution: 3.6 μ m
- 4x objective lens
- Integrated barcode reader
- Automation compatible



CloneSelect Imager Hardware Orientation: Indicator Lights



Indicator Lights

The indicator lights show the status of the instrument.



On



Ready



Imaging



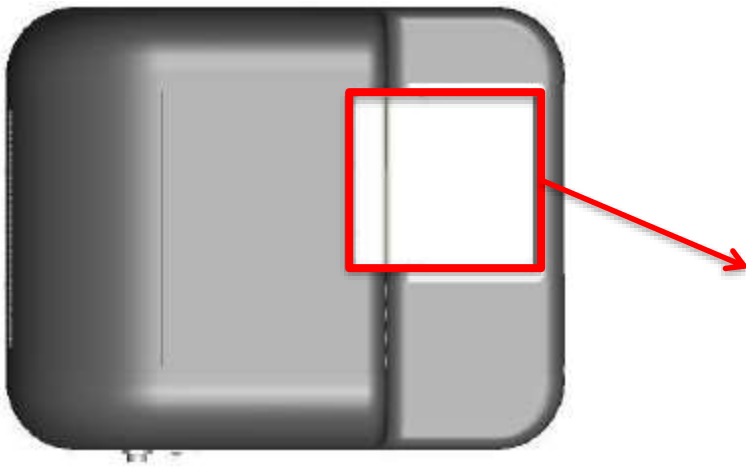
Vacuum On



Warning



CloneSelect Imager Hardware Orientation: Plate Carrier



- The **plate carrier** consists of a glass plate surrounded by a vacuum bed seal.
- The **microplate** is placed on the plate carrier so that the **skirt is over the vacuum bed seal**.
- When the plate carrier retracts into the instrument for imaging, two **pushers** gently push the microplate into the **back right corner** of the plate carrier so that the plate is always in the correct position for imaging.
- Before imaging, a **vacuum** is applied under the plate so that it is held down flat on the glass plate so that **all the wells are in the same focal plane**. Therefore there is no need to autofocus on every well – accelerating imaging!

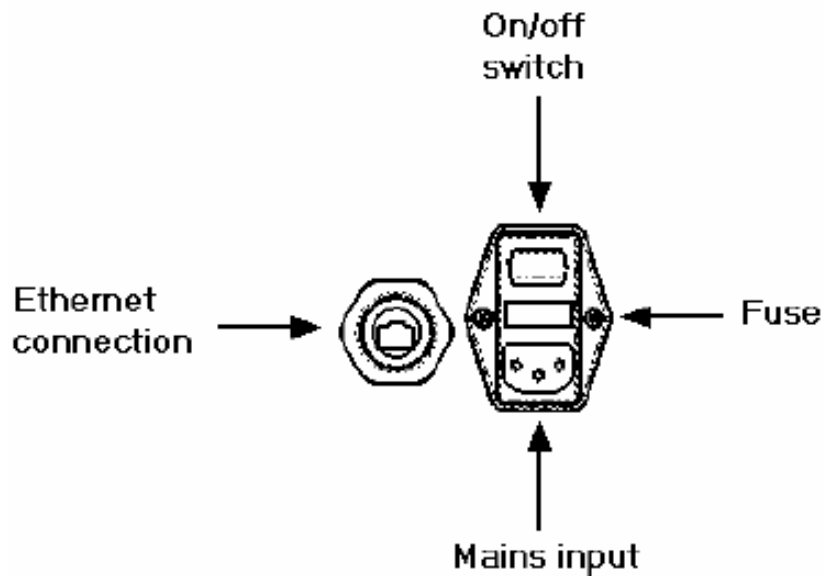
CloneSelect Imager Hardware Orientation: Computer



- CloneSelect Imager is supplied with a **Dell Tower 5810 Workstation with special hardware and software** to support the imaging function.
- **Do not** attempt to use any other workstation to operate CloneSelect Imager.
- The workstation is supplied with **Microsoft Windows 7™** and will require **security configuration** if it is to be connected to a **network**.
 - If the workstation , ensure that **no changes** are made to the configuration of the **private network connection** to the CloneSelect Imager since that will stop it operating.



CloneSelect Imager Hardware Orientation: Instrument Power & Ethernet Connection

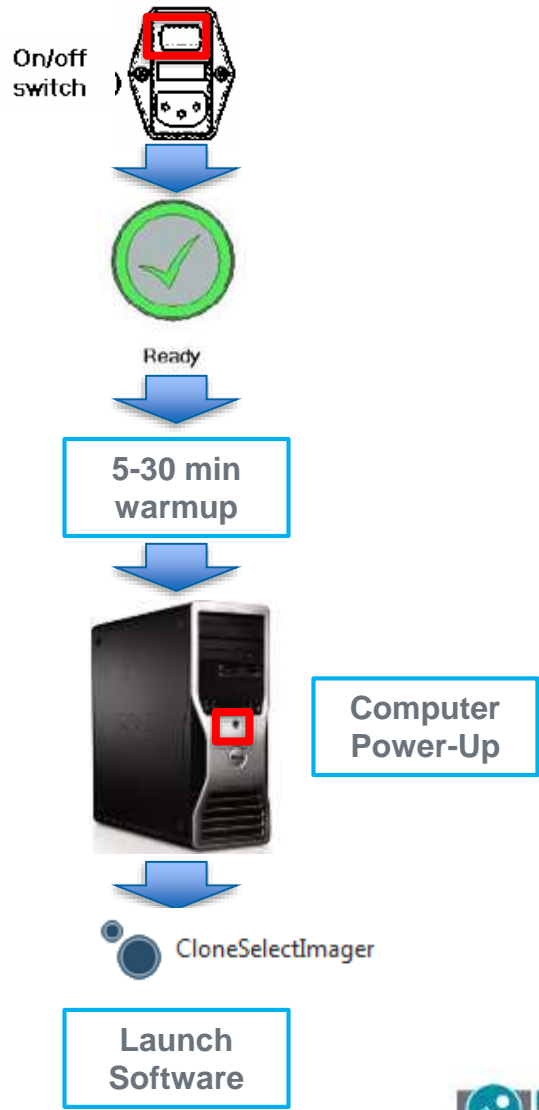


- The **power & ethernet connections** to CloneSelect Imager are located on the **connections panel** on the **left hand side** of the instrument.
- Ensure that the **mains cables** to the **CloneSelect Imager**, the **workstation** and to the **monitor** are connected **before making the network connections**.
- Note that the connection to CloneSelect Imager is a **private Gigabit Ethernet link** and the connection **must** be made to the workstation **add-on card Gigabit Ethernet port.****

****NOTE: This will have been performed on install by your Molecular Devices Field Service Engineer.**

CloneSelect Imager Hardware Orientation: Instrument Power-Up Procedure

- 1) Switch **on** CloneSelect Imager.
- 2) Wait for the **Ready light** to illuminate.
- 3) Leave CloneSelect Imager in the **'Ready'** state for at least **5 minutes** prior to use.
 - ✓ In cases where the unit is **cold** to the touch, wait for **30 minutes** prior to use.
- 4) Switch **on** the **computer workstation** and launch the **CloneSelect Imager software** by clicking on the **desktop shortcut**.
- 5) The instrument is now ready for use.



Support Resources

- Go to the HELP menu within CSI Software
- Support and Knowledge Base: <http://mdc.custhelp.com/>
- Request Support: <http://mdc.custhelp.com/app/ask> or via email support@moldev.com
- Technical Support can also be reached by telephone:
 - 1 (800) 635-5577
 - Select options for Tech Support → Biotherapeutics Products → Clone Select Imager





MOLECULAR DEVICES

ADVANCING PROTEIN AND CELL BIOLOGY