

PRE-INSTALLATION GUIDE

# CloneSelect Imager



Welcome to the family of Molecular Devices CloneSelect® Imager users. This document provides you with the information required to prepare your facility for the installation of your new CloneSelect Imager.

This document will cover the following topics:

- Certifications
- Environment requirements
- Space and table requirements
- Power requirements
- Facility receiving requirements
- Utility, water, other service, and safety precautions and hazard identification sections
- Service and support

## Certifications

CloneSelect Imagers meet all current requirements for CE approval and is ISO9001:2008 certified.

## Environment requirements

As with any precision optical instrument, take care to maintain a low-dust, low-vibration environment.

Below are the environmental requirements:

### Indoor use only

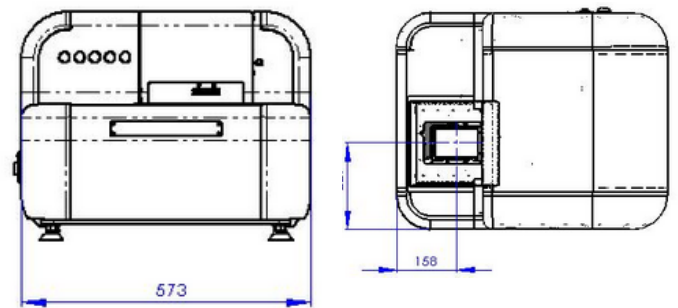
Temperature	10°C to 40°C
Humidity	20% to 80% non-condensing
Altitude	Up to 2,000M
Transient overvoltage	Installation category (Overvoltage category) II
Rated pollution	Pollution degree 2

## Space and table requirements

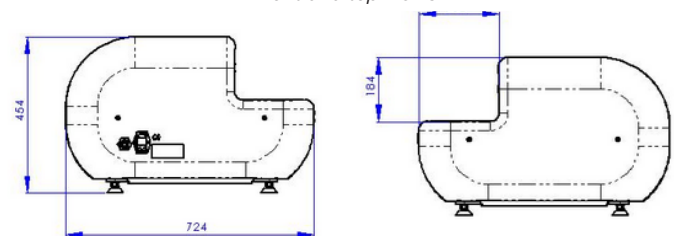
The CloneSelect Imager is a benchtop unit. The support surface needs to be capable of supporting more than 61 kg (131 lbs). The instrument comes with an external computer, monitor, keyboard, and mouse, which requires additional space next to the instrument.

### Dimensions and weight

Instrument Dimensions (height x width x depth)	45.4 cm (17.9 in.) x 57.4 cm (22.6 in.) x 72.4 cm (28.5 in.)
Instrument weight	45 kg (99 lbs)
Computer Dimensions (height x width x depth)	41.79 cm (16.45 in.) x 17.65 cm (6.94 in.) x 51.83 cm (20.4 in.)
Computer weight	15.9 kg (32 lbs)
Peripheral Dimensions (height x width x depth)	50 cm (19.7 in.) x 60 cm (23.6 in.) x 40 cm (15.75 in.)



Front and top views



Side views

## Power requirements

At least three power sockets are required on the wall (instrument, computer, monitor). The CloneSelect Imager's power is auto-switching.

Voltage: 100 – 240V AC 50-60 Hz single phase

Power: 100 VA

Max Amp: 2.5 A

Connections: IEC input

Fuses: F1 - F3A

Fluctuations must be within  $\pm 10\%$  of the nominal voltage.

## Facility receiving requirements

**Only Molecular Devices personnel are authorized to open the crate(s).**

Equipment arrives in one crate including the computer system.

Dimensions (cm/in.)	L	W	H	Weight (kg/lbs)
Instrument crate	157.5/62	94/37	86.4/34	155/342

All boxes should be stored indoors and protected from environmental extremes until the equipment is removed.

Crate storage	Specification
Ambient storage conditions	Temperature: -20°C to +60°C (-4°F to 140°F) Relative humidity: 15% to 75% relative humidity (non-condensing)

## Utility, other service, safety precautions, and hazard identification sections

### Transport and storage

CloneSelect Imager must be stored and transported in temperatures within the range -25°C to +55°C.

The instrument should not be moved after installation. If relocation is necessary, it should be lifted by two people and placed on a rolling cart capable of supporting the weight. Call Molecular Devices Technical Support for more information.

### Lifting points

CloneSelect Imager is intended for benchtop operation. Care should be taken when moving the instrument. It is recommended that two people lift it.

### Electrical safety

CloneSelect Imager must be connected to a properly grounded power outlet to protect users from the risk of electric shock. The main chassis of the instrument is grounded together with all associated electrical components. Do not remove any of the fixed covers, as there are no user serviceable parts inside. All internal work should be referred to Molecular Devices approved service personnel.

## Support and service

**Only a CloneSelect Imager system-trained Molecular Devices engineer is authorized to install or uninstall the instrument.**

For support or service, please contact us at:

Web: [www.moleculardevices.com/support](http://www.moleculardevices.com/support)  
[www.moldev.com/support](http://www.moldev.com/support)

### Contact Us

Phone: +1.800.635.5577

Web: [www.moleculardevices.com](http://www.moleculardevices.com)

Email: [info@moldev.com](mailto:info@moldev.com)

Check our website for a current listing of worldwide distributors.

### Regional Offices

USA and Canada +1.800.635.5577

United Kingdom +44.118.944.8000

Europe\* 00800.665.32860

China (Beijing) +86.10.6410.8669

China (Shanghai) +86.21.3372.1088

Hong Kong +852.3971.3530

Japan +81.3.6362.9109

South Korea +82.2.3471.9531

\*Austria, Belgium, Denmark, Finland, France, Germany, Ireland, Netherlands, Spain, Sweden and Switzerland