

Dual LED Blue/White Light Transilluminator

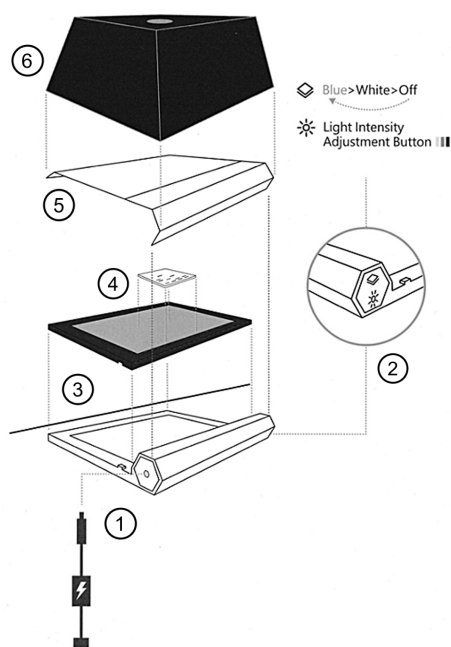
Catalog Number LB0100

Doc. Part No. 2162228 Pub. No. MAN0018667 Rev. B.0

Contents

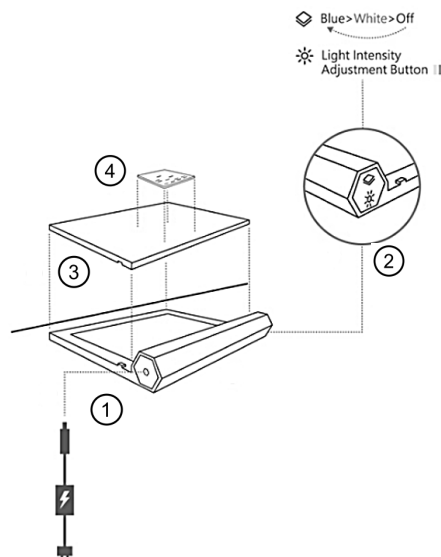
- Dual LED blue/white transilluminator base
- Amber filter cover (filters to ~580 nm)
- Power cord and adapter
- Imaging hood
- White uniform plate
- Blue uniform plate

Visualize protein and DNA gels using Blue Light mode



- ① Connect the power cord to the power socket.
- ② Press the ON/OFF switch to the Blue Light mode.
- ③ Place the blue uniform plate in position.
- ④ Place the protein- or DNA-containing gel on the glass.
- ⑤ Position the amber filter (580 nm) to magnetically connect with the transilluminator's base.
- ⑥ *(Optional)* Place the imaging hood on top of the amber filter for observing stained protein gels, membranes, or nucleic acid gels previously stained with fluorescent dyes excited in the range of 430-490 nm. For example, gels stained with SYPRO™ red stain, SYPRO™ Ruby, SYPRO™ Orange, SYPRO™ Tangerine, SYBR™ Green, and SYBR™ SAFE can be observed using the blue filter.
- ⑦ *(Optional)* Press the light intensity adjustment button for optimal viewing.
- ⑧ *(Optional)* Place a camera or cellular device on top of the imaging hood opening to capture images.
- ⑨ Once complete, turn the transilluminator off.

Visualize protein gels using White Light mode



- ① Connect the power cord to the power socket.
- ② Press the ON/OFF switch to the White Light mode.
- ③ Place the white uniform plate in position.
- ④ Place the protein gel, membrane, or X-ray film on the glass.
- ⑤ Capture images as needed.
- ⑥ Once complete, turn the transilluminator off.

Specifications

Parameter	Specification
LED source	Blue and white LED modules
LED life	> 30,000 hours
Emission maxima	470 nm (blue LED)
Input voltage	100-240 VAC
Input current	2.0 A
Automatic power-off	5 minutes
Recommended storage temperature	15-30°C
Gel viewing dimensions	12 × 18 cm
Unit dimensions (L × W × H)	22 × 18.5 × 3 cm

Troubleshooting

Observation	Possible cause	Recommended action
Low signal.	Suboptimal labeling of proteins or nucleic acids.	Optimize the labeling protocol.
	Too much ambient light.	Use the imaging hood cover with the viewing hole card on top.
No signal.	The fluorescent reagent used to label proteins or DNA was not excitable using blue light illumination.	Use fluorescent stains that can be excited at 470nm.
Cell phone image is pixelated.	The glass uniform plate was not in place.	Use the provided blue glass uniform plate while using a 470 nm excitable staining protocol or a white glass uniform plate when taking pictures of a non-fluorescent gel or membrane.

Clean the transilluminator

Use a clean, wet cotton cloth to wipe the surfaces of the transilluminator and blue/white uniform plates.

Safety information

The Dual LED Blue/White Light Transilluminator is an electrical device. The included power cord has a universal transformer compatible with 110 V to 220 V. Only the power cord supplied with the Dual LED Blue/White Light Transilluminator should be used to power the device.

- Never touch the power cord or outlet with wet hands.
- Do not use this device in damp areas or while standing on damp floors.
- Do not use the device if a crack is apparent in illumination surface.
- Do not attempt to open the transilluminator.
- The device has an automatic shut-off after 5 minutes.
- Always disconnect the transilluminator from the electrical outlet before cleaning the device.

- The transilluminator does not produce UV light. However, the intense blue light emitted by the transilluminator may promote macular degeneration upon prolonged exposure, especially in individuals prone to such problems (e.g., older individuals or those with fair complexion and blue eyes or nutritional or endocrine defects). Use the transilluminator amber filter unit to protect your eyes.

Note: The amber filter unit is not designed as protection for UV emission, and will NOT protect your eyes when viewing gels on UV transilluminators.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Thermo Fisher Scientific | 3747 N. Meridian Road | Rockford, Illinois 61101 USA
For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.