

*e*Blot™

Protein Transfer System

For fast semi-dry electroblotting of proteins from mini polyacrylamide gels to nitrocellulose or PVDF membranes

Version No. 08152014

User Manual



GenScript USA Inc.

860 Centennial Ave.

Piscataway, NJ 08854

Tel: 732-885-9188, 732-885-9688

Fax: 732-210-0262, 732-885-5878

Email: product@genscript.com



Table of Contents

Product Contents	1
Product Specifications	3
eBlot™ Protein Transfer Device	5
Accessory Products	7
Introduction	8
Quick Reference Guide	11
Protocols	12
Troubleshooting	21
Examples of Results	22
Technical Support	23
Warranty	24
Appendix	25

Product Contents

Type of products

This manual is supplied with the eBlot™ Protein Transfer Device (L03010).

eBlot™ Protein Transfer Device Contents

The contents of the eBlot™ Protein Transfer Device are listed below

Component	Quantity
eBlot™ Protein Transfer Device	1 each
eBlot™ Graphite Electrode (Installed inside the device)	1 each
Sponge Cushion (Installed inside the device)	2 each
Regional specific power cord	1 each
Forceps	1 each
Shovel	1 each
Shallow Tray	2 each

Upon Receiving the Instrument

Examine the unit carefully for any damage incurred during transit. File any damage claims with the carrier. The warranty does not cover in-transit damage.

eBlot™ Protein Transfer Pads

The following eBlot™ Protein Transfer Pads are available from GenScript:

Product	Cat. No.
eBlot™ Protein Transfer Pads (Basic, without membrane, 20-pak)	L03011
eBlot™ Protein Transfer Pads (Nitrocellulose, 20-pak)	L03013
eBlot™ Protein Transfer Pads (PVDF, 20-pak)	L03014

The eBlot™ Protein Transfer Pads come with the following components.

Component	Basic	Nitrocellulose	PVDF
eBlot™ Protein Transfer Pad	20	20	20
Nitrocellulose membranes (20 sheets)	-	1	-
PVDF membranes (20 sheets)	-	-	1
eBlot™ Equilibration Buffer (125 ml)	2	2	2
Gel Window (66 mmx56 mm)	1	1	1
Gel Window (76 mmx64 mm)	1	1	1
Gel Window (88 mmx78 mm)	1	1	1
Absorbent Filter Paper	1	1	1
Sponge Cushion	2	2	2

Product Contents, continued

Components for eBlot™ Protein Transfer Pad are as follows:

Component	Quantity
1x eBlot™ Cathode Pad	1
1x eBlot™ Anode Pad	1

Store the eBlot™ Protein Transfer Pads at room temperature. For best results, use the eBlot™ Protein Transfer Pads before the expiration date printed on the package.

eBlot™ Graphite Electrode

The eBlot™ Graphite Electrode (L03012) is available separately from GenScript:

Product	Cat. No.
eBlot™ Graphite Electrode	L03012

For best results, when a graphite electrode has been used for 100 times of protein transfer, replace it with a new one.

eBlot™ Equilibration Buffer

The eBlot™ Equilibration Buffer (M01078, 125 ml) is available separately from GenScript.

Product	Cat. No.
eBlot™ Equilibration Buffer	M01078

⚠ Important

DO NOT start a run without properly assembled Transfer Stacks in place!

Product Specifications

Intended Use For research use only. Not intended for human or animal diagnostic or therapeutic uses.

eBlot™ Protein Transfer Device Specifications

Weight:	1.6 kg
Dimensions:	325 mm (l) x 195 mm (w) x 70 mm (h)
Electrical Parameters:	100-120 V, 220-240 V, 50/60 Hz, 3.2 A
Built-in Features:	Digital Display, Alarm, Light LED
Compatibility:	Suitable for fast electroblotting of proteins from mini polyacrylamide gels to PVDF or nitrocellulose membranes
Materials:	Acrylonitrile Butadiene Styrene, Polycarbonate, Aluminum, Titanium, Plasticized silicone.
Operating Temperature:	15 - 40 °C
Forceps:	Stainless steel
Shovel:	Polycarbonate
Shallow Tray:	Polycarbonate

Avoid acetone, dimethyl sulfoxide, and acetic acid. These reagents can erode or damage the device.

eBlot™ Protein Transfer Pads Specifications

The eBlot™ Protein Transfer Pads are used with the eBlot™ Protein Transfer Device.

The specifications of the eBlot™ Protein Transfer Pads are listed below:

eBlot™ Protein Transfer Pad

eBlot™ Anode Pad:	90 mm (l) x 80 mm (w) x 2.5 mm (thickness)
eBlot™ Cathode Pad:	90 mm (l) x 80 mm (w) x 2.5 mm (thickness)
Materials:	Blotting filter paper presoaked with proprietary anode or cathode buffer

Membrane

Nitrocellulose:	90 mm (l) x 80 mm (w)
PVDF:	90 mm (l) x 80 mm (w)

eBlot™ Equilibration Buffer

Size:	125 ml x 2
Formulation:	Proprietary

Product Specifications, continued

Gel Window

Inter frame size:	66 mm x 56 mm
	76 mm x 64 mm
	88 mm x 78 mm
Materials:	Polycarbonate

Absorbent Filter Paper

Size:	80 mm x 70 mm
Materials:	Vegetable fiber

eBlot™ Graphite Electrode Specifications

The eBlot™ Electrode is used as the replaceable anode electrode of eBlot™ Protein Transfer Device and available separately from GenScript. The specifications for eBlot™ Graphite Electrode are listed below:

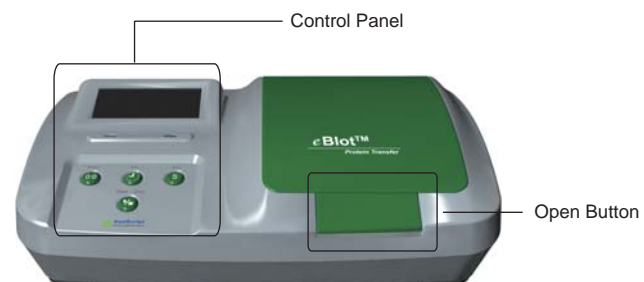
eBlot™ Graphite Electrode

Dimensions:	105 mm (l) x 95 mm (w) x 10 mm (h)
Weight:	140 g
Materials:	Powdered carbon, Clay, and Stainless steel

eBlot™ Protein Transfer Device

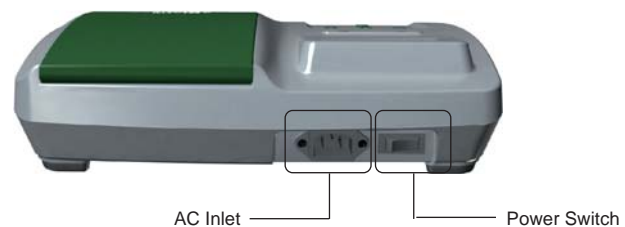
Front View of eBlot™ Device

The front-top view showing various parts of the eBlot™ Protein Transfer Device is shown below.



Rear View of eBlot™ Device

The rear view showing various parts of the eBlot™ Protein Transfer Device is shown below.



eBlot™ Protein Transfer Device, continued

Control Panel of eBlot™ Device

The control panel of the eBlot™ Protein Transfer Device is described below.

The **Digital Display** shows two rows of multi-digits that specify the transfer conditions as follows:

The upper three digits after text <PN> indicate how many times of transfer the graphite electrode has been used for.

The lower four digits specify the running time of protein transfer in minute and second.

The two **Status Lights** show the working mode of the eBlot™ Protein Transfer Device. When the right status light is on, the device is switched on and working at transferring mode; when both left and right status lights are on, the device is working at numbering mode.

The **Reset** button is used to clear parameters.

The **Min.** button is used to shift between transferring and numbering mode, and to set running time. Each short press will increase one minute. Each long press (2 seconds) will toggle working mode from transferring to numbering or the opposite.

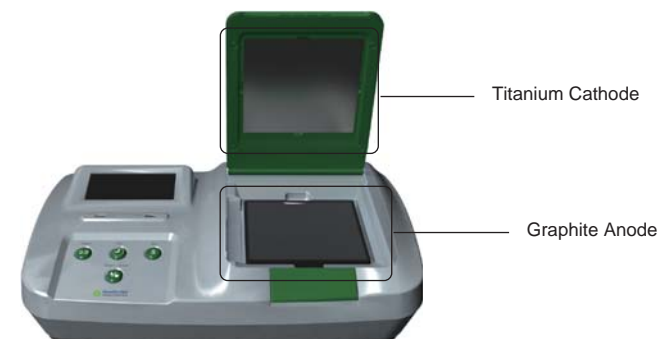
The **Sec.** button is also used to set running time, each press will increase 5 seconds.

The **Start/Stop** button is used to activate/stop the transfer program.



Top View of Open eBlot™ Device

The top view of open eBlot™ Protein Transfer Device identifying various parts.



Accessory Products

Precast Gels and Premade Buffers

The precast Express™ PAGE Gels as well as premade buffers and buffer powders are available from GenScript. For details, contact Technical Support or visit www.genscript.com.

ONE-HOUR Western™ Detection Kits

The ONE-HOUR Western™ Detection Kits used for Western blotting analysis are available from GenScript. Ordering information is provided below. For more information, visit www.genscript.com or call Technical Support.

Product	Quantity	Cat. No.
ONE-HOUR Western™ Basic Kit (Rabbit)	1 Kit (5 Assays)	L00204
ONE-HOUR Western™ Basic Kit (Mouse)	1 Kit (5 Assays)	L00205
ONE-HOUR Western™ Basic Kit (Goat)	1 Kit (5 Assays)	L00399
ONE-HOUR Western™ Standard Kit (Rabbit)	1 Kit (5 Assays)	L00204C
ONE-HOUR Western™ Standard Kit (Mouse)	1 Kit (5 Assays)	L00205C
ONE-HOUR Western™ Standard Kit (Goat)	1 Kit (5 Assays)	L00228
ONE-HOUR Western™ Standard Kit with TMB (Rabbit)	1 Kit (5 Assays)	L00204T
ONE-HOUR Western™ Standard Kit with TMB (Mouse)	1 Kit (5 Assays)	L00205T
ONE-HOUR Western™ Standard Kit with TMB (Goat)	1 Kit (5 Assays)	L00228T
ONE-HOUR Western™ Advanced Kit (Rabbit)	1 Kit (5 Assays)	L00241
ONE-HOUR Western™ Advanced Kit (Mouse)	1 Kit (5 Assays)	L00242
ONE-HOUR Western™ Advanced Kit (Goat)	1 Kit (5 Assays)	L00243
ONE-HOUR IP-Western Kit (Rabbit)	1 Kit (5 Assays)	L00231
ONE-HOUR IP-Western Kit (Mouse)	1 Kit (5 Assays)	L00232
ONE-HOUR IP-Western Kit (Goat)	1 kit (5 Assays)	L00233
ONE-HOUR Western™ Fluorescent Kit	1 Kit (10 Assays)	L00397
ONE-HOUR Western™ Multiplex Fluorescent Kit	1 Kit (10 Assays)	L00398

Introduction

System Overview

Semi-dry Western blotting is a common technique applied in protein research. Conventional semi-dry blotting is a cumbersome process, requiring time-consuming reagent preparation and setup, followed by an electrophoretic transfer that could take one hour or more. GenScript's eBlot™ Protein Transfer System accelerates the semi-dry blotting process without sacrificing performance. The eBlot™ Protein Transfer System, consisting of the eBlot™ Protein Transfer Device and eBlot™ Protein Transfer Pads, enables researchers to quickly, reliably perform electrophoretic transfer of proteins from various types of mini polyacrylamide gels to membranes in 7 to 10 minutes without the need to prepare additional buffers. The proteins transferred using the eBlot™ Protein Transfer System exhibit high detection sensitivity as to proteins transferred using other existing blotting methods.

System Components

The eBlot™ Protein Transfer System consists of:

eBlot™ Protein Transfer Device

The eBlot™ Protein Transfer Device is a self-contained electroblotting unit with a power supply built-in that allows for fast transfer of proteins.

eBlot™ Protein Transfer Pads

As the consumable part of eBlot™ Protein Transfer System, the eBlot™ Protein Transfer Pads are the mixed assortment of eBlot™ Protein Transfer Pad, nitrocellulose or PVDF membranes, Equilibration Buffer and Gel Window. Each pack of eBlot™ Protein Transfer Pad contains an 1×eBlot™ Cathode Pad and an 1×eBlot™ Anode Pad presoaked with proprietary cathode buffer and anode buffer respectively, allowing for rapid, convenient and reliable protein blotting without the need to prepare additional buffers.

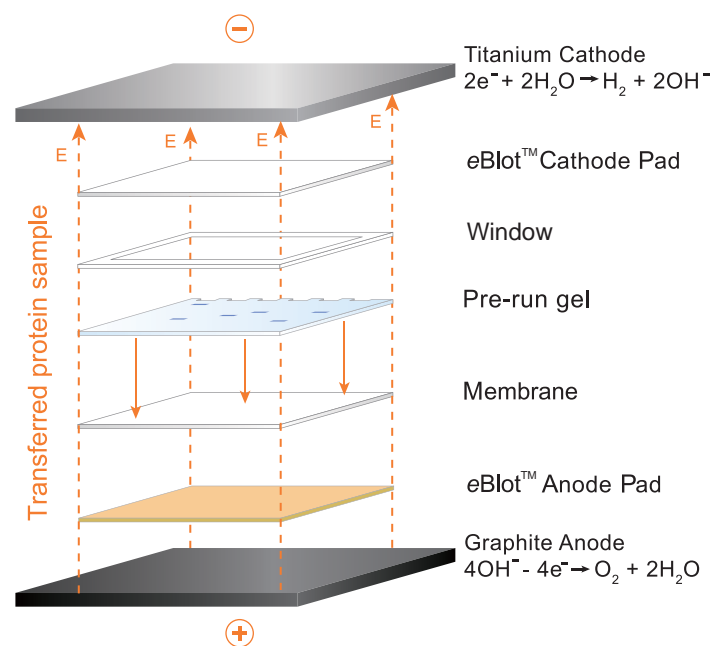
Introduction, continued

System Mechanism

The eBlot™ Protein Transfer System is based on the proprietary fast semi-dry blotting technique developed by GenScript.

To use the eBlot™ Protein Transfer System for rapid protein blotting, assemble the eBlot™ Protein Transfer Pad with your pre-run gel and the membrane on the eBlot™ Protein Transfer Device. The eBlot™ Cathode Pad and eBlot™ Anode Pad act as ion reservoirs that contain the appropriate anode and cathode buffers. The design of the eBlot™ Protein Transfer Device reduces the distance between electrodes and the integrated power supply enables the system to generate a certain definite voltage allowing for rapid and directional movement of negatively charged protein molecules from the gel matrix onto the membrane within 7-10 minutes. Meanwhile, the membrane presoaked with eBlot™ Equilibration Buffer shows good combination ability with proteins that prevents the flow through of low molecular weight proteins from the membrane.

Schematic mechanism of eBlot™ Protein Transfer System showing the flow of current:



Introduction, continued

System Features

Important features of the eBlot™ Protein Transfer System are listed below:

- Unique semi-dry electroblotting technique created for fast, reliable protein transfer within 7-10 minutes.
- Self-contained electroblotting unit with an integrated power supply for easy and convenient procedures.
- Consumable transfer pads offering convenience without the need for additional buffers.
- Compatible for use with various types of mini polyacrylamide gels.
- Proprietary formula without methanol .
- High transfer efficiency as compared to other existing blotting methods.

Quick Reference Guide

Introduction

A quick reference guide for operating the eBlot™ Protein Transfer Device is provided below.

Mode	Action	Sound	Light	Display
eBlot™ Device plugged in	Connect eBlot™ Device to an electrical outlet and power switch is on	–	Steady right light	Default running time (00:00)
eBlot™ Device and transfer stack assembled	Assemble transfer stack on the device and close lid	–	Steady right light	Default running time (00:00)
Time selection	Press Min. and Sec. button to select desired running time	–	Steady right light	User specified running time (00:00)
Run	Press Start/Stop button	–	Flashing right light	Counting down time
End of run	Automatic	Continuous beeping for 2 minutes	Steady right light	Default running time (00:00)
Checking the number of uses of the graphite anode	Press and hold Min. button for 2 seconds	–	Steady left and right lights	Times of the graphite anode has been used for transfer
Replacement of worn graphite anode	Switch off the device and replace the worn graphite anode with a new one	–	–	–

Protocols

Recommendations

To obtain the best results, follow these recommendations:

1. Wear gloves at all times during the entire blotting procedures to prevent contamination of pads, gels and membranes.
2. Do not touch the gel or membrane with bare or gloved hands. This may contaminate the gel or membrane and interfere with further analysis. If needed, always use forceps to adjust the membrane or gel.
3. Avoid using expired eBlot™ Protein Transfer Pads and eBlot™ Equilibration Buffer. Always use the pads and buffer before the specified expiration date printed on the package.
4. Remove any trapped air bubbles between the gel and membrane during the assembly of the transfer stack using the small shovel supplied with the device.

Installing the eBlot™ Device

1. Check the Power Cord supplied with the unit to ensure that the cord is compatible with the local socket format.
2. Place the eBlot™ Protein Transfer Device on a levelled laboratory bench. Keep the area around the device clear to ensure proper ventilation of the unit.
3. For your safety: Position the device properly such that the Power Switch and the AC inlet located on the rear of the unit are easily accessible.
4. Ensure the Power Switch is in the **Off** position.
5. Open the closed lid of the eBlot™ Protein Transfer Device by pressing the **Open** button. Place one or two pieces of Sponge Cushion in the anode tank depending on gel thickness. For 0.75 and 1.0mm gel, use one or two pieces of Sponge Cushion; for 1.5mm gel, use one or zero piece of Sponge Cushion.

Note: After 20 times of protein transfer, replace the used Sponge Cushions with new ones. A pair of new Sponge Cushions are included in each box of eBlot™ Protein Transfer Pads.

6. Insert the eBlot™ Graphite Electrode into the anode tank as described in Section "Replacing the eBlot™ Graphite Electrode", then close the lid of the device.

Protocols, continued

Installing the eBlot™ Device , continued

- Pull out the Waste Tray from the right side of the device. Place a new Absorbent Filter Paper inside the tray and then push the tray in.

Note: Change absorbent paper with a new one when opening a new box of pads.



- Attach the power cord to the AC inlet and then to the electrical outlet. Use only properly grounded AC outlets and power cords.
- When the electrophoresis of your samples is almost complete, press the Power Switch (located on the rear of the device) to turn **ON** the eBlot™ Protein Transfer Device. The right Status Light is on indicating you are using transfer mode. The lower four digits of the Digital Display show the default running time (00:00).



You are ready to use the eBlot™ Protein Transfer Device for blotting application.

Assembling the Transfer Stack

- Open the closed lid by pressing the **Open** button.



Protocols, continued

Assembling the Transfer Stack, continued

- Remove one package labeled as eBlot™ Protein Transfer Pad from the eBlot™ Protein Transfer Pads box and tear the laminated sealing of the package. Remove the two small packages respectively labeled 1xeBlot™ Cathode Pad and 1xeBlot™ Anode Pad.



- Tear the sealing of the 1xeBlot™ Anode Pad package. Remove the eBlot™ Anode Pad from the package and place it on the anode plate of the eBlot™ Protein Transfer Device.

Note: The pad must be placed in the middle of the graphite electrode position, can not contact on both sides of the metal frame.



- Pour 10 ml eBlot™ Equilibration Buffer into the shallow tray supplied with the eBlot™ Device.
- Tear the sealing of the Nitrocellulose or PVDF Membranes package. Remove one sheet of membrane and soak it in the eBlot™ Equilibration Buffer for 1 minute.

Note: If a PVDF membrane is used, the membrane must be pre-wetted with methanol before equilibrating in eBlot™ Equilibration Buffer. For users of eBlot™ Protein Transfer Pads (Basic, 20-pak) (L03011), the pre-cut membrane should be prepared by themselves.

Protocols, continued

Assembling the Transfer Stack, continued

- Place the equilibrated membrane on the eBlot™ Anode Pad. Gently remove air bubbles between the membrane and the anode pad using the small shovel supplied with the device.



- Carefully remove the pre-run gel containing your protein samples from the gel cassette and briefly rinse the gel with distilled water in another shallow tray to remove any small gel pieces attached to the gel and facilitate easy positioning of the gel on the membrane.
- Place the gel on the membrane. Gently remove air bubbles between the gel and the membrane.



Protocols, continued

Assembling the Transfer Stack, continued

- Select appropriate Gel Window according to actual size of the gel (see the table below). Place Gel Window on the gel. Ensure that the Gel Window fully covers the margin of the membrane.

Pre-run gel	Gel Window
Gel with size > 80 × 60 mm	Gel Window 88 × 78 mm
Gel with size 66 × 56 mm to 80 × 60 mm	Gel Window 76 × 64 mm
Gel with size < 66 × 56 mm	Gel Window 66 × 56 mm

Note: Gel Window is used as the spacer between anode pad and cathode pad to prevent short circuit.



- Tear the sealing of the 1×eBlot™ Cathode Pad package. Remove the eBlot™ Cathode Pad from the package and place it on the gel.



Note: During assembling of the transfer stack, make sure to remove all the air bubbles trapped between the transfer pads, pre-run gel and the membrane, which may prevent the transfer of proteins and cause empty spots on the transferred membrane.

Protocols, continued

Assembling the Transfer Stack, continued

11. Press the **Open** button, and then push back and close the lid of eBlot™ Protein Transfer Device.

Performing Blotting

Perform protein blotting as described below within 15 minutes of assembling the transfer stack.

1. Press the **Min.** and **Sec.** buttons to set appropriate running time based on the protein size (see table below). If an undesired running time is set by mistake, press **Reset** button to clear the wrong time, and then press the **Min.** and **Sec.** buttons to choose the desired running time.



Protein size (kDa)	Recommended start running time (min)
< 80	7
80-160	8-9
> 160	10

Note: Based on the initial results, the transfer time may need to be optimized to make best transfer results by pressing the Time button in 5-second increment. For proteins greater than 200 kDa, we recommend to start with 11 minutes.

2. Press the **Start/Stop** button to start the transfer. The running time begins to count down and right Status Light keeps flashing during the whole transfer program.



Protocols, continued

Performing Blotting, continued

3. At the end of the transfer, current automatically shuts off and the eBlot™ Protein Transfer Device signals the end of transfer with repeated beeping sound. The right status light stops flashing and the lower five digits show text (00:00).
4. Press any button on the control panel to stop the beeping.
5. Proceed to disassemble the stack and clean the device.

Disassembling and Cleaning the eBlot™ Device

To obtain good transfer and detection results, disassemble the transfer stack right away after ending the blotting procedure.

1. Open the closed lid by pressing the **Open** button.
2. Carefully separate the transferred membrane from the transfer stack and proceed with further protein detection procedures.

Note: If you are using PVDF membrane, place the membrane immediately into the blocking solution (or water) as PVDF membrane dries quickly. If the PVDF membrane is dried, rewet the membrane with methanol and then rinse it with distilled water to wet it completely before use.

3. Discard the gel and the used eBlot™ Protein Transfer Pad.

Note: Do not re-use the eBlot™ Protein Transfer Pad after blotting. Discard after each use.

4. Clean the titanium cathode plate, graphite anode plate and its surrounding area with a dry cloth or paper tissue.
5. Replace the Absorbent Filter Paper in the Waste Tray with a new one when it has soaked up the waste from 20 times of transfer. Pull out the Waste Tray from the right side of the device to perform the replacement.

Note: After about 40 runs, take out the graphite electrode, soak in distilled water for 30 minutes, dry the surface with a clean paper towel, then leave air dry overnight with graphite side up.

At this point, the eBlot™ Protein Transfer Device is ready for another run.

If you are not using the device, turn off the Power Switch located on the back of the device.

For any other repairs and service, contact Technical Support. Do not perform any repairs or service on the eBlot™ Protein Transfer Device by yourself to avoid any possible damages to the device.

Protocols, continued

Replacing the Graphite Electrode

During the transfer process, the Graphite Electrode will absorb ions from anode pad as well as lose carbon composition of the anode buffer. For best blotting results, after having been used for **100 times** of transfer, the worn Graphite Electrode should be replaced by a new one.

1. If the eBlot™ Protein Transfer Device works at transferring mode, press and hold **Min.** button for 2 seconds to toggle to numbering mode. If the upper three digits show "100" or a number greater than "100", perform the replacing protocol as describe below.



2. Switch **Off** the eBlot™ Protein Transfer Device. Open the lid of the device and take the worn Graphite Electrode out of the device.



3. Tear the sealing of a new Graphite Electrode package and take the new Graphite Electrode out of the package. Place the new Graphite Electrode into the anode tank and close the lid of the device.

Protocols, continued

Replacing the Graphite Electrode, continued

4. Switch **On** the eBlot™ Device. Press and hold **Min.** button for 2 seconds to toggle to numbering mode. When the upper three digits are flashing, press **Reset** button to zero the transferring times.



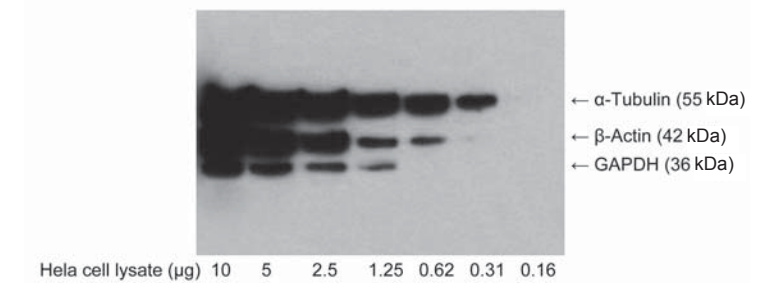
After successfully installing Graphite Electrode into the eBlot™ Device, you are ready to use the eBlot™ Device for another blotting application.

Troubleshooting

Problem	Cause	Solution
The right Status Light doesn't flash during blotting process.	Incomplete electric circuit due to improper assembly of the transfer pads.	Ensure the transfer stack is assembled correctly: use the eBlot™ Anode Pad first followed by the membrane, the pre-run gel, Gel Window and eBlot™ Cathode Pad.
The left and right Status Lights flash simultaneously.	Excessive current is flowing through the Device.	Check the transfer stack and ensure Gel Window covered correctly on the gel.
Inefficient transfer	1. Salt built-up on plate electrodes	1. Clean the titanium cathode plate, graphite anode plate with a wet cloth or paper tissue followed by a dry one to remove any insoluble salts.
	2. Membrane insufficiently equilibrated in eBlot™ Equilibration buffer	2. Equilibrate membrane in eBlot™ Equilibration buffer before transfer.
	3. Incorrect transfer conditions or insufficient transfer time	3. Use a gel of lower concentration to separate high molecular weight proteins. Increase the transfer time in 5-second increments.
	4. PVDF membrane was not prewet with methanol	4. Pre-wet PVDF membrane with methanol before transfer.
	5. Confusion of the eBlot™ Anode Pad and Cathode Pad	5. Ensure the transfer stack is assembled correctly: <i>Bottom-eBlot™</i> Anode Pad (yellow), <i>Top-eBlot™</i> Cathode Pad (white).
Empty spots on the membrane	Air bubbles trapped between gel and membrane prevent the transfer of proteins.	When assembling transfer stack, use the small shovel supplied with the device to remove any air bubbles between the gel and the membrane.

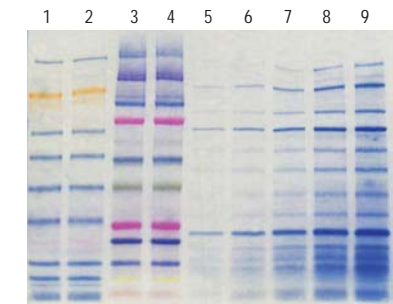
Examples of Results

Western blot result after using the eBlot™ for protein transfer



An Express™ PAGE Gels 8-16% (GenScript MG816W12) was blotted using the eBlot™. 10.0, 5.0, 2.5, 1.25, 0.62, 0.31 and 0.16 μg of HeLa cell lysate were loaded respectively. Proteins on the Nitrocellulose membrane were detected using THE™ alpha Tubulin Antibody (GenScript, A01410), THE™ beta Actin Antibody (GenScript, A00702) and GAPDH Antibody (GenScript, A01632). The secondary antibody was Anti-MOUSE IgG (H & L) (GOAT) Antibody Peroxidase Conjugated (Rockland, 610-1302). The signal was developed with LumiSensor™ HRP Substrate Kit (GenScript, L00221V500).

Result Using Nitrocellulose



Good transfer of protein standard bands onto the nitrocellulose membrane. Protein Standards were separated on a 4-20% GenScript Express™ Plus PAGE Gel (Bis-Tris). After electrophoresis, the gel was blotted using the eBlot™ protein transfer system for 11 min as described in this manual. Lane 1 and 2, Invitrogen Pre-Stain protein marker (5 μL, MW: 4-250 kDa), Lane 3 and 4, Bio-rad Pre-Stain protein marker (10 μL, MW: 10-250 kDa), and Lane 5-9, NEB All Blue Pre-Stain protein marker (0.3, 0.6, 1.3, 2.5, and 5 μL, MW: 10-230 kDa).

Technical Support

Web Resources

Visit the GenScript Web site at www.genscript.com for:

1. Technical resources, including manuals, vector maps and sequences, application notes, MSDSs, FAQs, formulations, citations, handbooks, *etc.*
2. Complete technical support contact information
3. Access to the GenScript Online Catalog
4. Additional product information and special offers

Contact Us

For more information or technical assistance, call, write, fax, or email.

GenScript USA Inc.

860 Centennial Ave.

Piscataway, NJ 08854

Tel: 732-885-9188, 732-885-9688

Fax: 732-210-0262, 732-885-5878

Email: product@genscript.com

Warranty

eBlot™ Protein Transfer Device

GenScript warrants that eBlot™ Protein Transfer Device will be free from defects in material and workmanship for a period of one year from date of purchase. If any defects occur in the product during this warranty period, GenScript will, at its option, repair, replace, or refund the purchase price of this product at no charge to you. The following defects, however, are specifically excluded:

1. Defects caused by improper operation.
2. Repair or modification done by anyone other than GenScript or an authorized agent.
3. Use of fittings or other spare parts supplied by anyone other than GenScript.
4. Damage caused by accident or misuse.
5. Damage caused by disaster.
6. Corrosion due to use of improper solvent or sample.

For any inquiry or request for repair service, contact GenScript after confirming the model and serial number of your instrument. For your protection, items being returned must be insured against possible damage or loss. This warranty shall be limited to the replacement of defective products. It is expressly agreed that this warranty will be in lieu of all warranties of fitness and in lieu of the warranty of merchantability.

Appendix

CECS Center Europe Certification Service

Certificate No. **CECS/0020110523A**

VERIFICATION OF EMC COMPLIANCE

EU COUNCIL DIRECTIVE 2004/108/EC

Applicant : Nanjing Genscript Co.,LTD
Address : No.78 Shuangbai Road,Xuanwu District,Nanjing China.
Manufacturer : Nanjing Genscript Co.,LTD
Address : No.78 Shuangbai Road,Xuanwu District,Nanjing China.
Sample Name : Multifunction Gel Processor
Model : GS-01, GS-02
Test Report No. : BZT082110520/EMC
Codes/Standards Applied : EN 55014-1:2006; EN 55014-2:1997+A1: 2001, EN 61000-3-2:2006,; EN 61000-3-3:1995+A1:2001+A2:2005
Remarks : The sample meets the requirements of the above standards.
Date Of Issuance : Jul 4,2011

Conclusion : We Confirm That The Technical Construction File And Manufacturing, Inspection And Testing Processes For Above Mentioned Sample Comply With The Essential Safety Requirements of EU COUNCIL DIRECTIVE OF 2004/108/EC Applied Codes And Standards.

President of CECS : Robert Owen

 Signature

CE

Center Europe Certification Service Co., Ltd
 Abbey Orchard St, Westminster, Greater London SW1P
 info@cecsuk.com

Appendix, continued

CECS Center Europe Certification Service

Certificate No. **CECS/0020110523B**

VERIFICATION OF LVD COMPLIANCE

EU COUNCIL DIRECTIVE 2006/95/EC

Applicant : Nanjing Genscript Co.,LTD
Address : No.78 Shuangbai Road,Xuanwu District,Nanjing China.
Manufacturer : Nanjing Genscript Co.,LTD
Address : No.78 Shuangbai Road,Xuanwu District,Nanjing China.
Sample Name : Multifunction Gel Processor
Model : GS-01, GS-02
Test Report No. : BZT082110520 /LVD
Codes/Standards Applied : EN 60335-1:2001+A1:2004+A2:2006
Remarks : The sample meets the requirements of the above standards.
Date Of Issuance : Jul 4,2011

Conclusion : We confirm that the technical construction file and manufacturing, inspection and testing processes for above mentioned sample comply with the essential safety requirements of EU COUNCIL LOW VOLTAGE DIRECTIVE OF 2006/95/EC applied codes and standards.

President of CECS : Robert Owen

 Signature

CE

Center Europe Certification Service Co., Ltd
 Abbey Orchard St, Westminster, Greater London SW1P
 info@cecsuk.com

Appendix, continued

VERIFICATION OF CONFORMITY
According to FCC Part 15B

Certificate No.: SEM11058591

Responsible Party's Name : NanJing Genscript Co., Ltd.
Address : No.78 Shuangbai Road, Xuanwu District, Nanjing, China

Manufacturer : NanJing Genscript Co., Ltd.
Address : No.78 Shuangbai Road, Xuanwu District, Nanjing, China

Description of Product : Multifunction Gel Processor
Model No. : GS-01, GS-02

Trade Name : eStain, eBlot

Report No. : STR11058051E-3

Compliance With Part 15B of FCC Rules.



This device complies with Part 15 of the FCC Rules. Operation is subject to the following two conditions: (1) This device may not cause harmful interference, and (2) This device must accept any interference received, including interference that may cause undesired operation.

Responsible Party:

Responsible Signature: _____

Name / Title: _____

Date: _____

Tested By:



SEM.Test Compliance Service Co., Ltd.
3/F, Jinbao Commerce Building, Xin'an Fanshen Road,
Bao'an District, Shenzhen, P.R.C.

Issued By: _____

Name / Title: Jandy So / PQ Manager

Date of Issue: Jul 12, 2011



The Certification of Verification shows that the tested sample technically compliances with the FCC Part 15. The certification applies to the tested sample above mentioned only and should not implied an assessment of the whole.

Tel.: +86-755-33663308 Fax.: +86-755-33663309 E-mail: sem@semtest.com.cn Website: www.semtest.com.cn