



Western Blotting

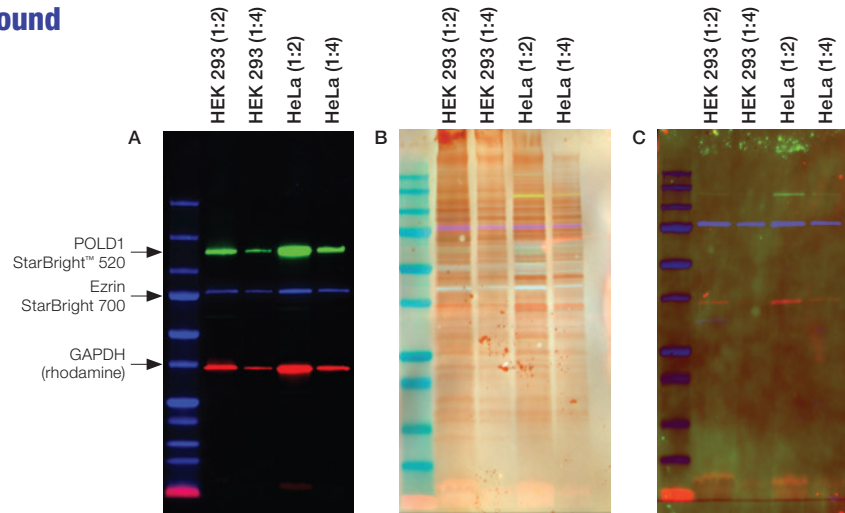
EveryBlot Blocking Buffer

Rapid blocking without compromise
for western blot and ELISA

See Your Target, Not Your Background

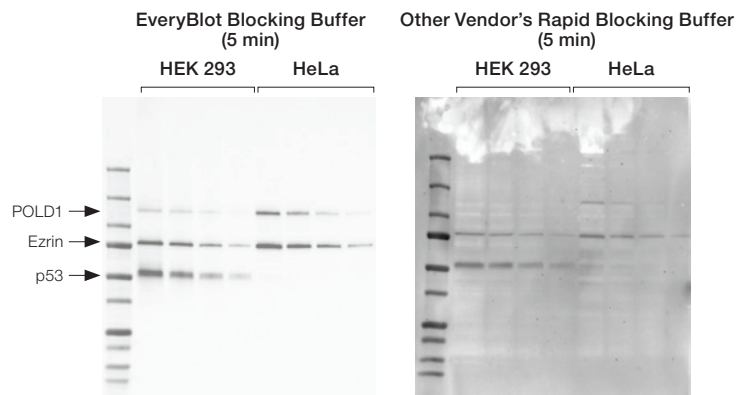
Fluorescence

Better sensitivity and lower background than other vendor's rapid fluorescence-optimized blocker or 3% nonfat dry milk. HEK 293 and HeLa cell lysates at 2-fold dilutions were separated and transferred to a nitrocellulose membrane. Blots were either blocked with EveryBlot Blocking Buffer for 5 min (A), other vendor's rapid fluorescence-optimized blocker for 10 min (B), or 3% nonfat dry milk in Tris buffered saline (TBS) for 30 min (C). After blocking, blots were processed and imaged identically and are displayed with identical contrast settings.



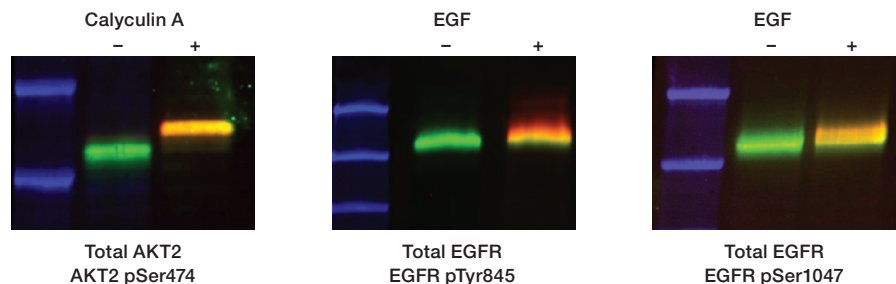
Chemiluminescence

Better sensitivity and lower background with chemiluminescence detection. HEK 293 and HeLa cell lysates were loaded at 2-fold dilutions and transferred to a PVDF membrane. After blocking with either EveryBlot Blocking Buffer (left) or other vendor's rapid blocking buffer (right), blots were processed and imaged identically and are displayed with identical contrast settings for three targets.



Phosphoprotein Detection

Nothing to interfere with phosphoprotein detection. Detection of total (green) and phospho-specific (red) targets after blocking with EveryBlot Blocking Buffer for 5 min. Left, Jurkat cells unstimulated (-) or stimulated (+) with calyculin A. Center and right, A431 cells unstimulated (-) or stimulated (+) with EGF. Other vendor's rapid blocking buffer is not compatible with phosphoprotein detection.

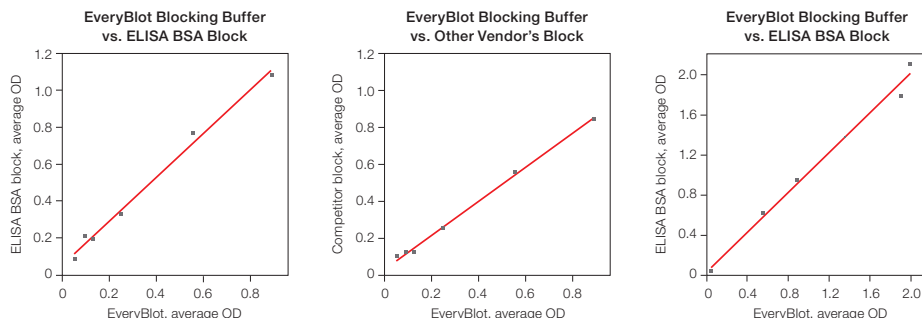


Visit bio-rad.com/EveryBlot for more information.

BIO-RAD

Indirect and Sandwich ELISA Assays

ELISA blocking in seconds. Indirect and sandwich ELISA blocking using EveryBlot Blocking Buffer for <1 min showed comparable results with good correlation coefficient when compared to 1 hr blocking with standard ELISA BSA blocking buffer or other vendor's ELISA blocking buffer. Left and center, indirect ELISA; right, sandwich ELISA.

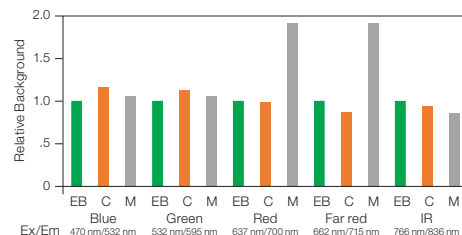


Advanced Formulation for the Best Results

Low Background across the Entire Spectrum

EveryBlot Blocking Buffer resulted in lower fluorescence background compared to other vendor's fluorescence-optimized blocking buffer and nonfat dry milk. Low-fluorescence PVDF membranes without sample or antibodies were blocked in either EveryBlot Blocking Buffer (■),

other vendor's fluorescence-optimized blocking buffer (■), or 3% nonfat dry milk (■). After washing, membranes were imaged side by side on the ChemiDoc™ MP Imaging System. Ex/Em, excitation/emission; IR, infrared.



Better Blocking in Less Time

Speed without compromise. Amphiphilic compounds increase blocking efficiency while maintaining sensitivity and reducing background.

Typical Blockers



1 hr

EveryBlot Blocking Buffer



5 min
(western blot)

<1 min
(ELISA)

Ordering Information

Catalog # Description

EveryBlot Blocking Buffer

12010020 EveryBlot Blocking Buffer, 500 ml

StarBright Blue Secondary Antibodies

12004158 StarBright Blue 700 Goat Anti-Mouse IgG, 400 µl
 12004161 StarBright Blue 700 Goat Anti-Rabbit IgG, 400 µl
 12005866 StarBright Blue 520 Goat Anti-Mouse IgG, 400 µl
 12005869 StarBright Blue 520 Goat Anti-Rabbit IgG, 400 µl

hFAB Rhodamine Housekeeping Protein Fluorescent Primary Antibodies

12004163 hFAB Rhodamine Anti-Actin Primary Antibody, 200 µl
 12004165 hFAB Rhodamine Anti-Tubulin Primary Antibody, 200 µl
 12004167 hFAB Rhodamine Anti-GAPDH Primary Antibody, 200 µl

Clarity™ and Clarity Max Western ECL Blotting Substrates

1705061 Clarity Western ECL Substrate, 500 ml
 1705062 Clarity Max Western ECL Substrate, 100 ml

Catalog # Description

Trans-Blot™ Turbo Transfer System

1704150 Trans-Blot Turbo Transfer System
 1704156 Trans-Blot Turbo Mini 0.2 µm PVDF Transfer Packs, pkg of 10
 1704157 Trans-Blot Turbo Midi 0.2 µm PVDF Transfer Packs, pkg of 10
 1704158 Trans-Blot Turbo Mini 0.2 µm Nitrocellulose Transfer Packs, pkg of 10
 1704159 Trans-Blot Turbo Midi 0.2 µm Nitrocellulose Transfer Packs, pkg of 10
 1704270 Trans-Blot Turbo RTA Mini 0.2 µm Nitrocellulose Transfer Kit, for 40 blots
 1704271 Trans-Blot Turbo RTA Midi 0.2 µm Nitrocellulose Transfer Kit, for 40 blots
 1704272 Trans-Blot Turbo RTA Mini 0.2 µm PVDF Transfer Kit, for 40 blots
 1704273 Trans-Blot Turbo RTA Midi 0.2 µm PVDF Transfer Kit, for 40 blots
 1704274 Trans-Blot Turbo RTA Mini 0.2 µm LF PVDF Transfer Kit, for 40 blots
 1704275 Trans-Blot Turbo RTA Midi 0.2 µm LF PVDF Transfer Kit, for 40 blots

Visit bio-rad.com/EveryBlot for more information.

BIO-RAD, CHEMIDOC, CLARITY, STARBRIGHT, and TRANS-BLOT are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2024 Bio-Rad Laboratories, Inc. Clarity Max Western ECL Substrate is manufactured for Bio-Rad by Cyanagen Srl and is the subject of U.S. and ex-U.S. patents, including US7855287, EP1950207, US9040252, AU2011202658, CA2742025, US8129136, EP1962095, and CN102313732.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23
France 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050
Italy 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 080 007 7373 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670
The Netherlands 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23
Russian Federation 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23
Switzerland 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

