



EveryBlot Blocking Buffer

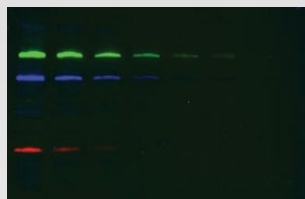
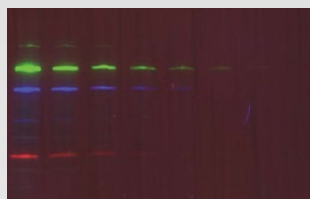
5 minute blocking for ALL western blots...without compromises

See your target, not your background

Fluorescence

Other vendor's new fluorescence-optimized blocker, 1 hour

EveryBlot Blocking Buffer, 5 minutes



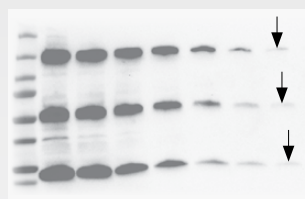
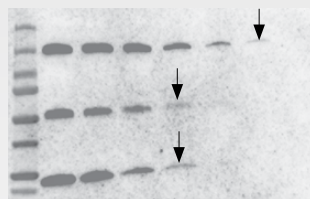
← RNA Helicase A (DyLight 800)
 ← ALIX (StarBright 700)
 ← Tubulin (hFAB Rhodamine)

Better sensitivity and lower background than other vendor's new fluorescence-optimized blocker. HEK 293 cell proteins were separated and transferred to low fluorescence PVDF. Left, blot blocked in the other vendor's new fluorescence-optimized blocker for 1 hour. Right, blot blocked in EveryBlot Blocking Buffer for 5 minutes. After blocking, both blots were processed and imaged identically and are displayed with identical contrast settings.

Chemiluminescence

Milk, 30 minutes

EveryBlot Buffer, 5 minutes



← RNA Helicase A
 ← p53
 ← PRDX 3

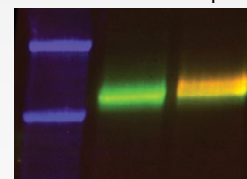
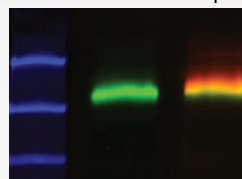
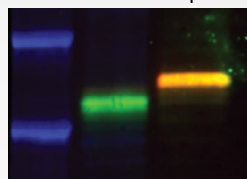
Better sensitivity and lower background with chemiluminescence detection. Three targets were detected in HEK 293 cells. After blocking, both blots were processed and imaged identically and are displayed with identical contrast settings.

Phosphoprotein Detection

Calyculin A

EGF

EGF



Total AKT2
 AKT2 pSer474

Total EGFR
 EGFR pTyr845

Total EGFR
 EGFR pSer1047

Nothing to interfere with phosphoprotein detection. Detection of total (green) and phospho-specific (red) targets. Left, Jurkat cells unstimulated (-) or stimulated (+) with calyculin A. Center and right, A431 cells unstimulated (-) or stimulated (+) with EGF.

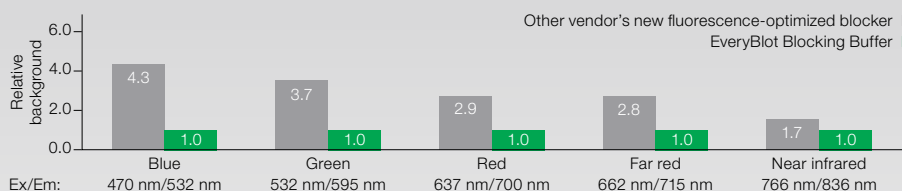


Visit bio-rad.com/EveryBlot for more information.



Advanced formulation for the best results

Low background across the entire spectrum



Lower fluorescence background than other vendor's new fluorescence-optimized blocker. Low fluorescence PVDF membranes without sample or antibodies were blocked in either the other vendor's new fluorescence-optimized blocking buffer or EveryBlot Blocking Buffer. After washing, membranes were imaged side by side on the ChemiDoc MP Imaging System.

Better blocking in less time

Typical Blockers



1 hour

EveryBlot Blocking Buffer



5 minutes

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

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
 12004159 StarBright Blue 700 Goat Anti-Mouse IgG, 80 µl
 12004161 StarBright Blue 700 Goat Anti-Rabbit IgG, 400 µl
 12004162 StarBright Blue 700 Goat Anti-Rabbit IgG, 80 µl
 12005866 StarBright Blue 520 Goat Anti-Mouse IgG, 400 µl
 12005867 StarBright Blue 520 Goat Anti-Mouse IgG, 80 µl
 12005869 StarBright Blue 520 Goat Anti-Rabbit IgG, 400 µl
 12005870 StarBright Blue 520 Goat Anti-Rabbit IgG, 80 µl

hFAB Rhodamine Housekeeping Protein Fluorescent Primary Antibodies

12004164 hFAB Rhodamine Anti-Actin Primary Antibody, 40 µl
 12004163 hFAB Rhodamine Anti-Actin Primary Antibody, 200 µl
 12004166 hFAB Rhodamine Anti-Tubulin Primary Antibody, 40 µl
 12004165 hFAB Rhodamine Anti-Tubulin Primary Antibody, 200 µl
 12004168 hFAB Rhodamine Anti-GAPDH Primary Antibody, 40 µl
 12004167 hFAB Rhodamine Anti-GAPDH Primary Antibody, 200 µl

Catalog # Description

Clarity and Clarity Max Western ECL Blotting Substrates

1705060 Clarity Western ECL Substrate, 200 ml
 1705061 Clarity Western ECL Substrate, 500 ml
 1705062 Clarity Max Western ECL Substrate, 100 ml

Trans-Blot Turbo Transfer System

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

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

BIO-RAD and TRANS-BLOT are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. Clarity Max Western ECL Substrate is manufactured by Cyanagen Srl and is the subject of patent application numbers US7855287, EP1950207, US9040252, AU2011202658, CA2742025, US8129136, and EP1962095, together with other equivalent granted patents and patent applications in other countries like CN102313732. All trademarks used herein are the property of their respective owner.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00
Germany 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600
Japan 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30
Poland 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580
Sweden 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300
United Kingdom 44 01923 47 1301

