

final volume of 10 μ L. Electrophoresis was carried out at 90 V for two hours. After separation, proteins were transferred to PVDF membrane (Immobilon-FL, Millipore IPVH00010) overnight at 15V.

The membrane was washed with PBS containing 0.05% Tween-20 (PBST) for 5 minutes. The membrane was then blocked for one hour at room temperature in a 1:10 dilution of BSA blocker (Pierce 18587) as per manufacturers instructions.

The blot was incubated with the primary antibody, Rabbit-anti-Transferrin (Abcam ab1223-1) at 5 μ g/mL for two hours at room temperature.

The blot was washed five times with Pierce wash buffer (1860498) diluted 1:30 as per manufacturers instructions.

The blot was incubated one hour at room temperature with the secondary antibody, Cy5-labeled Goat-anti-Rabbit IgG (GE Healthcare PA45012V) diluted 1:2500 in wash buffer.

The blot was then washed six times in Pierce wash buffer for five minutes each. After a final rinse in PBS, the blot was imaged while still damp, approximately one hour after the PBS rinse. The blot was placed on a low-fluorescent plastic sheet on the sample holder for imaging.

Experiment 2: All samples were loaded in equal volumes (10 μ L). Each lane contained 30 ng of Human IgG, corresponding to approximately 20 ng of heavy chain. Two-fold serial dilutions of Transferrin, and of α 1-fetoprotein (AFP), were loaded on the gel as follows:

Lane	Transferrin (ng)	α 1-Fetoprotein (ng)
1	63	0.004
2	31	0.008
3	16	0.015
4	8	0.03
5	4	0.06
6	2	0.12
7	1	0.24
8	0.49	0.49
9	0.24	1
10	0.12	2
11	0.06	4
12	0.03	8
13	0.015	16
14	0.008	31
15	0.004	63

Separation was carried out at 150 V for one hour,

After separation, proteins were transferred to a PVDF membrane (Immobilon-FL, Millipore IPVH00010). Transfer was conducted at 100 V for one hour with an ice pack inserted in the tank.

The membrane was washed with PBS containing 0.05% Tween-20 (PBST) for 5 minutes. The membrane was then blocked for one hour at room temperature in a 2% solution of ECL Advance blocking agent (GE Healthcare CPK1075) dissolved in PBST.

After blocking, the membrane was washed with PBST for 5 minutes and incubated with a mixture of primary antibodies, containing Rabbit-anti-Transferrin (Abcam ab1223-1) and Mouse-anti-AFP (Meridian Life Sciences H45301M) antibodies, each diluted at 1:5000 in PBST.

The membrane was then subjected to two brief, one 15 minute, and four five-minute washes in PBST.

The membrane was then incubated with a mixture of fluorescently labeled secondary antibodies for one hour at room temperature. The secondary antibodies used were Cy3-Goat-anti-Mouse IgG at 1:2500 dilution, and Cy5-Goat-anti-Rabbit IgG at 1:2500 dilution (both from GE Healthcare). A Cy2-Goat-anti-Human IgG antibody was also added at 1:2500 dilution, to detect Human IgG directly. The Cy2-Goat-anti-Human IgG antibody was prepared from non-labeled antibody (Vector Labs AI-3000) using the Cy2 labeling kit from GE (GE Healthcare PA32000).

After hybridization, the membrane was washed three times with PBST for five minutes each, once with PBS for five minutes, then briefly rinsed with water and dried at room temperature.

The dry membrane was imaged with the FluorChem Q imaging system at 1.7x magnification. The membrane was positioned on the sample holder.

Single channel images of each of the three fluorescent dyes were acquired on the FluorChem Q. The spectral characteristics of the filter sets used are shown in Table 1. The exposure times were adjusted manually for each of the three channels.

Channel	Excitation	Emission Filter	Compatible Dyes
Blue	475/42 nm	537/35 nm	Cy2, Alexa 488, FITC
Green	534/30 nm	614/62 nm	Cy3, Alexa 546
Red	632/22 nm	699/62 nm	Cy5, Alexa 680

Table 1. Spectral characteristics of FluorChem Q filter sets, and compatible fluorescent dyes.

Cell Biosciences products used

- FluorChem Q

Other materials required

- 10-20% linear gradient Ready Gel (Bio-Rad 161-1124)
- Immobilon-FL PVDF membrane (Millipore IPVH00010)
- ECL Advance blocking agent (GE Healthcare CPK1075)
- BSA blocker (Pierce 1858753)

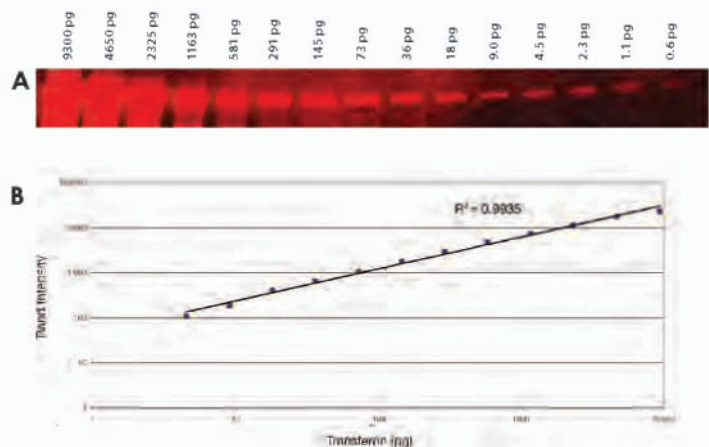


Figure 2. Serial dilutions of Transferrin detected by Western blot and imaged on the FluorChem Q. (a) Image acquired with a 16 second exposure. A lower limit of 1.1 pg Transferrin was detected, and the fluorescence intensity was linear from 4.5 pg to 9300 pg (b), providing a dynamic range over 3.3 orders of magnitude.

- Pierce wash buffer (Pierce 1860498)
- Rabbit-anti-Transferrin antibody (Abcam ab1223-1)
- Mouse-anti-AFP (Meridian Life Sciences H45301M)
- ECL Plex goat- α -mouse IgG-Cy3 (GE Healthcare PA43009)
- ECL Plex goat- α -rabbit IgG-Cy5 (GE Healthcare PA45011)
- Goat-anti-Human IgG (Vector Labs AI-3000)
- Cy2 Ab Labeling Kit (GE Healthcare PA32000)
- ECL Advance blocking agent (GE Healthcare CPK1075)
- Mini PROTEAN Tetra cell (Bio-Rad 165-8004)
- Tris-Glycine running buffer
- Tris-Glycine transfer buffer
- PBS buffer
- PBST buffer

Results and Discussion

Experiment 1: The performance of the FluorChem Q was tested by preparing a Western blot of a dilution series of Transferrin, a protein normally present in Human serum. An image of the Western blot, acquired on the FluorChem Q with a 16 second exposure is shown in Figure 2. A measurable linear dynamic range of over 2000-fold was achieved (Figure 2). Exposing a duplicate blot for 256 seconds gave a lower limit of detection of 0.6 pg (Figure 3)

Experiment 2: The compatibility of the FluorChem Q and ECL Plex fluorescent western blotting system when imaging multicolor Western blots was tested by preparing a Western blot containing serial dilutions of two proteins, Transferrin and AFP, as well as a constant level of a third protein, Human IgG. The three proteins were imaged in one imaging protocol, each in a different fluorescent channel. Though each of the proteins can be imaged independently, the images captured in each of the three fluorescent channels can be merged to allow the three proteins to be visualized simultaneously. Such a merged image depicting each of the three fluorescent channels is shown in Figure 4a. The power of multiplexing is demonstrated by this image, in which Transferrin (73-76 kDa) and AFP (70kDa) are easily distinguished.

The single channel images captured in the red channel (Transferrin, Figure 4b), and green channel (AFP, Figure 4c), demonstrate the large dynamic range and low limit of detection achieved for each protein and dye combination in a true Western blot experiment. The blue channel, IgG, can be used in later analysis as a loading control (Figure 4d).

The linear range and detection limit obtained for a given Western blot experiment will depend on multiple experimental factors including the nature of the protein of interest, the efficiency of transfer of the protein from gel to membrane, the selection of the primary and secondary antibodies, and optimization of the antibody hybridization conditions. The results presented in this report are specific to the materials and conditions described herein.



Figure 3. A lower limit of detection of 0.6 pg Transferrin was achieved with a 256 second exposure time.

Conclusion

The FluorChem Q provides a powerful means to image and quantify multicolor fluorescent Western blots labeled using the ECL Plex system. The sensitivity of the system is reflected in a detection limit of 0.6 pg of Transferrin. With an optimized exposure time, the fluorescence intensity of Transferrin staining was linear over three orders of magnitude, covering a range of protein amounts from 4.5 pg to 9.6 ng, which encompasses the range commonly seen in Western blotting.

The FluorChem Q also performed in multiplex analysis, acquiring images of three proteins, detected with three unique fluorophores on a single blot.

The linearity of the response over a large dynamic range, combined with the ability to detect multiple proteins simultaneously on the same blot, demonstrates the potential for accurate protein quantification, with a minimization of time and sample usage.

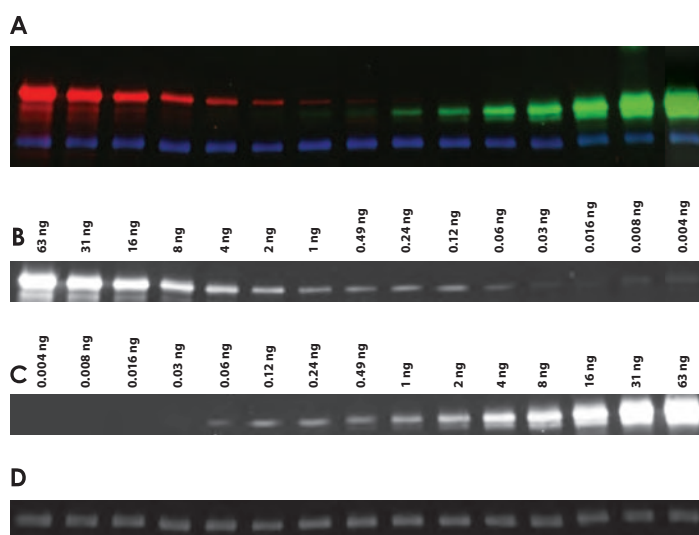


Figure 4. FluorChem Q generated image of ECL Plex labeled western. The Western blot was probed with three different fluorescent secondary antibodies. Panel (a) shows the composite 3-color image, which is 16 bit in all three color channels. Blue: Human IgG. Red: Transferrin. Green: a1-Fetoprotein. Single channel images are shown for (b) the red channel, detecting Transferrin, and (c) the green channel, detecting a1-Fetoprotein and (d) the blue channel detecting IgG.

References

1. Patton, Wayne F. (2000) "A thousand points of light: The application of fluorescence detection technologies to two-dimensional gel electrophoresis and proteomics" *Electrophoresis*. 21: 1123-1144.
2. GE Healthcare Application Note 28-4015-40 AB "Multiplex protein detection using the ECL Plex fluorescent Western blotting system"

Acknowledgement

Cell Biosciences gratefully acknowledges the researchers at Advansta Corporation in Menlo Park, California, for the information contained in this Tech Note.

Copyright © 2008 Cell Biosciences. All rights reserved. The Cell Biosciences logo and the wordmark FluorChem are registered trademarks of the Company. ECL Plex is a trademark of GE Healthcare. All other trademarks, service marks and tradenames appearing in this brochure are the property of their respective owners.

The FluorChem Q Imaging System provides the sensitivity of chemiluminescence and the quantitative power of fluorescence in one easy-to-use instrument. Designed with fast lens technology, the peltier cooled camera captures high-resolution images with a linear dynamic range that outperforms film, and with speeds 10X faster than a laser scanner. Equipped with three integrated excitation sources for multicolor westerns, the FluorChem Q is compatible with Cy dyes, Alexa dyes, Qdots, and chemiluminescent western blotting kits. Additionally, the FluorChem Q provides the flexibility to image fluorescently stained DNA gels, as well as Coomassie and silver stained protein gels.

For more information, visit us at: <http://www.cellbiosciences.com>



Cell  Biosciences