

ProStain™ Protein Quantification Kit

(version D)

Catalog No. 15001

Active Motif North America

1914 Palomar Oaks Way, Suite 150

Carlsbad, California 92008, USA

Toll free: 877 222 9543

Telephone: 760 431 1263

Fax: 760 431 1351

Active Motif Europe

104 Avenue Franklin Roosevelt

B-1330 Rixensart, Belgium

UK Free Phone: 0800 169 31 47

France Free Phone: 0800 90 99 79

Germany Free Phone: 0800 181 99 10

Telephone: +32 (0)2 653 0001

Fax: +32 (0)2 653 0050

Active Motif Japan

Azuma Bldg, 7th Floor

2-21 Ageba-Cho, Shinjuku-Ku

Tokyo, 162-0824, Japan

Telephone: +81 3 5225 3638

Fax: +81 3 5261 8733

Copyright 2008 Active Motif, Inc.

Information in this manual is subject to change without notice and does not constitute a commitment on the part of Active Motif, Inc. It is supplied on an “as is” basis without any warranty of any kind, either explicit or implied. Information may be changed or updated in this manual at any time.

This documentation may not be copied, transferred, reproduced, disclosed, or duplicated, in whole or in part, without the prior written consent of Active Motif, Inc. This documentation is proprietary information and protected by the copyright laws of the United States and international treaties.

The manufacturer of this documentation is Active Motif, Inc.

© 2008 Active Motif, Inc., 1914 Palomar Oaks Way, Suite 150; Carlsbad, CA 92008. All rights reserved.

All trademarks, trade names, service marks or logos referenced herein belong to their respective companies.

The purification column used in this kit is covered under US patent 005618418A.

TABLE OF CONTENTS	Page
Introduction	1
Kit Components and Storage	2
Additional Materials Required	2
Preparation of Reagents	2
ProStain™ Protein Quantification Protocol	3
Appendix	
Section A: Troubleshooting Guide	5
Section B: Related Products	6
Technical Services	8

Introduction

Active Motif's ProStain™ Protein Quantification Kit simplifies protein quantification by providing highly sensitive detection reagents in a convenient, easy-to-use format. The kit's detection reagent is also resistant to many effects, which can limit the usefulness of other protein quantification systems, such as pH and many commonly found contaminants; for example, detergents and salts. In addition, the large Stokes shift, fast reaction time and limited free dye quantum yield make this kit a significant improvement over other photometric or fluorescent-based detection systems.

The kit provides lyophilized dye reagent, dilution buffer and BSA for preparation of standards. Simply resuspend the lyophilized dye in methanol to create the concentrated stock solution, dilute the stock solution, load 100 µl into the wells of a microplate, add 100 µl of sample, mix, then read the fluorescence. The assay is performed at room temperature, and the signal is stable for up to 2 hours. Common contaminants, such as pH, salts, solvents and some detergents are well tolerated in this assay, but buffers containing high amounts of free amines will affect sensitivity.

Advantages

- Large Stokes shift for reduced background
- Fast and simple procedure
- Robust – limited effect from contaminating substances
- Increased quantum yield for improved sensitivity and wide dynamic range

product	format	catalog no.
ProStain™ Protein Quantification Kit	1000 assays	15001

* Sufficient components are provided for performing 1000 assays using fluorescent-based detection. This assay can also be easily adapted for use in smaller or larger formats such as 384-well plates or cuvettes.

ProStain™ is for research use only. Not for use in diagnostic procedures.

Kit Components and Storage

The ProStain Kit is for research use only. Not for use in diagnostic procedures. Kit components arrive at room temperature. We recommend storing each component at the temperatures recommended in the table below:

Reagents	Quantity	Storage / Stability
Dye Reagent AM1	0.5 mg	4°C for 6 months
Dilution Buffer	100 ml	4°C for 6 months
BSA	1 mg	4°C for 6 months

Additional materials required

- Multi-channel pipettor & pipettor reservoirs
- Fluorescent detector
- Distilled water
- Black microtiter plates or cuvettes
- Methanol (for reconstituting the Dye Reagent)

Preparation of Reagents

Dye Reagent Stock Solution

The Dye Reagent is supplied lyophilized. Prepare the Dye Reagent Stock Solution by resuspending the lyophilized Dye Reagent in 12.5 ml methanol in the provided amber bottle. This stock solution can be stored at 4°C for 6 months.

Dye Reagent Working Solution

Prepare the Dye Reagent Working Solution by diluting the Dye Reagent Stock Solution 1:10 with distilled water. This solution should be prepared fresh on the day the assay is performed.

Reagent	10 rxns	50 rxns	100 rxns
Dye Reagent Stock Solution	110 µl	550 µl	1.1 ml
Distilled H ₂ O	990 µl	4.95 ml	9.9 ml

Dilution Buffer

This is supplied ready to use.

Stock BSA Solution

The BSA is supplied lyophilized. Prepare the Stock BSA Solution by resuspending the lyophilized BSA in 1 ml of dH₂O in the provided tube to make a 1 mg/ml solution. This stock solution can be stored at -20°C for 6 months.

Protein Quantification Kit Protocol

PLEASE READ THE ENTIRE PROTOCOL BEFORE STARTING!

1. Remove kit contents from 4°C and bring all components to room temperature before use.
2. Set up a BSA Standard Curve in duplicate using the following concentrations: 10.0, 5.0, 2.5, 1.25, 0.63, 0.32, 0.15 and 0.0 µg/ml. (See table below for suggested layout.)
3. Add 198 µl of Dilution Buffer to wells A1 and A2.
4. Add 100 µl of Dilution Buffer to wells B1 through H1 and B2 through H2
5. Pipette 2 µl stock BSA solution (1 mg/ml) into wells A1 and A2.
6. Mix wells A1 and A2 by pipetting.
7. Transfer 100 µl from well A1 to B1 and A2 to B2.
8. Mix wells B1 and B2 by pipetting.
9. Transfer 100 µl from well B1 to C1 and B2 to C2.
10. Continue this procedure to wells G1 and G2. After mixing, discard 100 µl of solution from wells G1 and G2.
11. Wells H1 and H2 are blanks and should contain only 100 µl of Dilution Buffer.

	1	2	3	4	5	6	7	8
A	10.0 µg/ml	10.0 µg/ml	–	–	–	–	–	–
B	5.0 µg/ml	5.0 µg/ml	–	–	–	–	–	–
C	2.5 µg/ml	2.5 µg/ml	–	–	–	–	–	–
D	1.25 µg/ml	1.25 µg/ml	–	–	–	–	–	–
E	0.63 µg/ml	0.63 µg/ml	–	–	–	–	–	–
F	0.32 µg/ml	0.32 µg/ml	–	–	–	–	–	–
G	0.15 µg/ml	0.15 µg/ml	–	–	–	–	–	–
H	Blank	Blank	–	–	–	–	–	–

12. **Sample wells:** For protein determination of unknown samples, prepare a series of dilutions with the Dilution Buffer, for example: 1:50, 1:100 and 1:200. Pipette 100 μ l into each well. Duplicates of each sample are recommended.
13. Add 100 μ l/well of Dye Reagent Working Solution (see Preparation of Reagents section on page 2 for directions on preparing the Dye Reagent Working Solution) and mix by pipetting up and down.
14. Incubate for 30 minutes at room temperature (20-25°C) without agitation.

Note: The signal intensity is stable for up to 1.5 hours after the 30-minute incubation; read sample before 1.5 hours have elapsed.
15. Measure the fluorescence (excitation: 488 nm, emission: 635 nm).

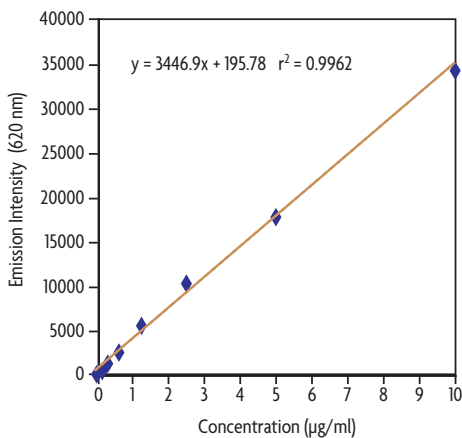
Note: When measuring fluorescence, ensure that the gain settings are set to optimal (~ 140 gain), the number of flashes are set to 3 with no lag time, and the integration time is ~ 40 μ s.
16. Use the standard curve to calculate the protein concentrations of the unknown samples. For the BSA standard curve, fit with the function $y = Ax + B$. The r^2 should be over 0.95.

Calculation of protein concentration using the BSA standard curve

Average the duplicate readings for each standard and sample and subtract the value obtained from the zero standard. Plot the fluorescence for the standards against the quantity of the standards and draw the best fit curve. To quantify the amount of protein in the samples, find the fluorescence value for the samples on the y-axis and extend a horizontal line to the standard curve. At the intersection point extend a vertical line to the x-axis and read the corresponding standard value. Note: If the samples have been diluted, the value read from the standard curve must be multiplied by the dilution factor.

Example standard curve.

The standard curve at right is provided for demonstration only. You should make a standard curve should be made every time an experiment is performed.



Appendix

Section A: Troubleshooting Guide

I. Buffer Compatibility and Contaminating Substances

A number of common contaminants have been tested with ProStain, and most are well tolerated; however, samples containing high concentrations of free amines are not recommended.

Contaminant	Final Concentration in the assay	Concentration in 100 μ l of sample	Result
Sodium chloride	20 mM	20 mM	OK
Sodium chloride	2 mM	4 mM	OK
Tris	10 mM	20 mM	OK
Ammonium sulfate	5 mM	10 mM	OK
Tween	0.001%	0.002%	Not recommended
Triton	0.001%	0.002%	Not recommended
SDS	0.04%	0.08%	OK

* BSA standards were assayed in the presence or absence of contaminants at the indicated final concentrations. Equivalent concentrations in 100 μ l sample volumes are also listed. This is not a complete list of contaminants. To determine your buffer compatibility, prepare two BSA curves – one in the same buffer as your samples and one with the supplied Dilution Buffer to determine if there is any interference.

II. Protein-to-Protein Variation

ProStain determines the protein concentration of a sample relative to a BSA standard curve. If you are quantitating recombinant or purified protein rather than a protein extract, accuracy may be improved by using the same protein at a known concentration to make the standard curve, if available.

III. Excitation and Emission Filters

The excitation and emission maxima for Dye Reagent bound to protein are 503 and 602 nm, respectively. We recommend using filters with the following ranges: 485-525 nm Excitation and 575-650 nm Emission.

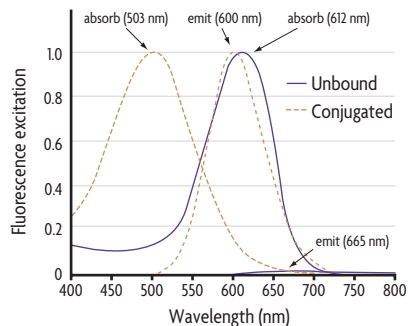


Figure 1: Absorption/emission spectra of free/bound dye.

Normalized absorption and emission spectra of free (solid lines) and conjugated dye (dotted lines) in phosphate buffer of pH 7.2.

Section B. Related Products

Sample Preparation	Format	Catalog No.
Nuclear Extract Kit	100 reactions	40010
	400 reactions	40410
Mitochondrial Fractionation Kit	100 reactions	40015
GAPDH Whole-cell Normalization Kit	1 x 96-well plate	48007
	5 x 96 well-plates	48507

Fluorescent Detection	Format	Catalog No.
CE Dye 503	1 kit	15101
CE Dye 540	1 kit	15102
Albumin Blue Fluorescent Assay Kit	1 kit	15002
ToxCount™ Cell Viability Assay	20 x 96 rxns	18010
LavaCell™ Membrane Stain	200 µg	15004

Fluorescent Dyes	Format	Catalog No.
Chromeo™ 488 Biotin	1 mg	15512
Chromeo™ 488 Carboxylic Acid	1 mg	15510
Chromeo™ 488 NHS-Ester	1 mg	15511
Chromeo™ 488 Streptavidin	1 mg	15513
Chromeo™ 494 Biotin	1 mg	15112
Chromeo™ 494 Carboxylic Acid	1 mg	15110
Chromeo™ 494 NHS-Ester	1 mg	15111
Chromeo™ 494 Streptavidin	1 mg	15113
Chromeo™ 546 Biotin	1 mg	15212
Chromeo™ 546 Carboxylic Acid	1 mg	15210
Chromeo™ 546 NHS-Ester	1 mg	15211
Chromeo™ 546 Streptavidin	1 mg	15213
Chromeo™ 642 Biotin	1 mg	15312
Chromeo™ 642 Carboxylic Acid	1 mg	15310
Chromeo™ 642 NHS-Ester	1 mg	15311
Chromeo™ 642 Streptavidin	1 mg	15313

Fluorescent Secondary Antibodies	Format	Catalog No.
Chromeo™ 488 Goat anti-Mouse IgG	1 mg	15031
Chromeo™ 488 Goat anti-Rabbit IgG	1 mg	15041
Chromeo™ 494 Goat anti-Rabbit IgG	1 mg	15042
Chromeo™ 546 Goat anti-Mouse IgG	1 mg	15033
Chromeo™ 546 Goat anti-Rabbit IgG	1 mg	15043
Chromeo™ 642 Goat anti-Mouse IgG	1 mg	15034
Chromeo™ 642 Goat anti-Rabbit IgG	1 mg	15044
ATTO 594 Goat anti-Mouse IgG	250 µl	15037
ATTO 594 Goat anti-Rabbit IgG	250 µl	15047
ATTO 647N (STED) Goat anti-Mouse IgG	250 µl	15038
ATTO 647N (STED) Goat anti-Rabbit IgG	250 µl	15048
ATTO 655 (STED) Goat anti-Mouse IgG	250 µl	15039
ATTO 655 (STED) Goat anti-Rabbit IgG	250 µl	15049

Fluorescent Secondary Antibodies	Format	Catalog No.
Chromo™ 488 Antibody Labeling Kit	1 kit	15090
Chromo™ 494 Antibody Labeling Kit	1 kit	15091
Chromo™ 546 Antibody Labeling Kit	1 kit	15092
Chromo™ 642 Antibody Labeling Kit	1 kit	15093

Immunofluorescence	Format	Catalog No.
MAX Stain™ Universal Immunofluorescence System	1 kit	15250
MAXpack™ Immunostaining Media Kit	1 kit	15251
MAXblock™ Blocking Medium	150 ml	15252
MAXbind™ Staining Medium	250 ml	15253
MAXwash™ Washing Medium	1000 ml	15254
MAXfluor™ TRIO Mounting Media Kit	1 kit	15255
MAXfluor™ Mounting Medium	2 ml	15256
MAXfluor™ DAPI Mounting Medium	2 ml	15257
MAXfluor™ PI Mounting Medium	2 ml	15258

Chromatin Immunoprecipitation	Format	Catalog No.
ChIP-IT™ Express	25 rxns	53008
ChIP-IT™ Express Enzymatic	25 rxns	53009
ChIP-IT™ Express HT	96 rxns	53018
ChIP-IT™ Protein G Magnetic Beads	25 rxns	53014
Re-ChIP-IT™	25 rxns	53016
ChIP-IT™	25 rxns	53001
ChIP-IT™ Shearing Kit	10 rxns	53002
ChIP-IT™ Enzymatic	25 rxns	53006
Enzymatic Shearing Kit	10 rxns	53005
Salmon Sperm DNA/Protein G agarose	25 rxns	53003
ChIP-IT™ Control Kit – Human	5 rxns	53010
ChIP-IT™ Control Kit – Mouse	5 rxns	53011
ChIP-IT™ Control Kit – Rat	5 rxns	53012
Ready-to-ChIP HeLa Chromatin	10 rxns	53015

Co-Immunoprecipitation	Format	Catalog No.
Universal Magnetic Co-IP Kit	25 rxns	54002
Nuclear Complex Co-IP Kit	50 rxns	54001

Histone Purification	Format	Catalog No.
Histone Purification Kit	10 rxns	40025
Histone Purification Mini Kit	10 rxns	40026

Chromatin Assembly	Format	Catalog No.
Chromatin Assembly Kit	10 rxns	53500

DNA Methylation	Format	Catalog No.
MethylDetector™	50 rxns	55001
MethylCollector™	25 rxns	55002
Fully Methylated Jurkat DNA	10 µg	55003

Technical Services

If you need assistance at any time, please call Active Motif Technical Service at one of the numbers listed below.

Active Motif North America

1914 Palomar Oaks Way, Suite 150
Carlsbad, CA 92008
USA

Toll Free: 877 222 9543
Telephone: 760 431 1263
Fax: 760 431 1351
E-mail: tech_service@activemotif.com

Active Motif Europe

104 Avenue Franklin Roosevelt
B-1330 Rixensart, Belgium
UK Free Phone: 0800 169 31 47
France Free Phone: 0800 90 99 79
Germany Free Phone: 0800 181 99 10
Telephone: +32 (0)2 653 0001
Fax: +32 (0)2 653 0050
E-mail: eurotech@activemotif.com

Active Motif Japan

Azuma Bldg, 7th Floor
2-21 Ageba-Cho, Shinjuku-Ku
Tokyo, 162-0824, Japan
Telephone: +81 3 5225 3638
Fax: +81 3 5261 8733
E-mail: japantech@activemotif.com

Visit Active Motif on the worldwide web at <http://www.activemotif.com>

At this site:

- Read about who we are, where we are, and what we do
- Review data supporting our products and the latest updates
- Enter your name into our mailing list to receive our catalog, *MotifVations* newsletter and notification of our upcoming products
- Share your ideas and results with us
- View our job opportunities

Don't forget to bookmark our site for easy reference!