

ProStain™

simple, sensitive fluorescent protein quantification

Protein quantification is an essential step in many cell biology procedures, such as Western blot and ELISA. Active Motif's ProStain™ Protein Quantification Kit simplifies protein quantification by providing highly sensitive detection reagents in a convenient and easy-to-use format. The kit's detection reagent is also resistant to many commonly found contaminants such as detergent and salt (Figure 1), which can limit the usefulness of other protein quantification systems. In addition, the large Stokes shift (Figure 2), fast reaction times and limited free dye quantum yield make ProStain a significant improvement over other commonly used detection systems such as the Bradford or BCA assay. The ProStain Kit is ideal for use with many of Active Motif's Sample Preparation products, including the popular Nuclear Extract Kit and Mitochondrial Fractionation Kit.

The ProStain™ advantage

- Large Stokes shift for reduced background
- Fast and simple – results in less than 1 hour
- Robust – limited effect from contaminants like detergent and salt
- Highly sensitive
- Wide dynamic range

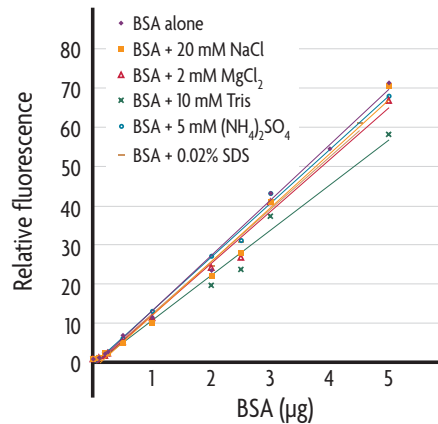


Figure 1: Linear, reproducible quantification in the presence of contaminants.

Increasing amounts of BSA protein were quantified using the ProStain Protein Quantification Kit in the presence of a variety of contaminants.

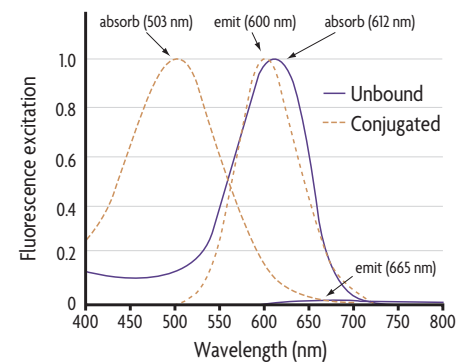


Figure 2: Absorption/emission spectra of free/bound dye.

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

Product	Format	Catalog No.
ProStain™ Protein Quantification Kit	1000 rxns	15001

CONTENTS

Each ProStain Protein Quantification Kit supplies sufficient reagents for 1000 fluorescent protein quantification reactions.