

## QuantiFluo™ Protein Assay Kit (QFPR-200)

### Quantitative Fluorimetric Determination of Protein/Peptide Concentration

#### DESCRIPTION

The protein or peptide is known as the "building blocks of life" and is one of the most important macromolecules in life science. Protein determination is a very common practice. Simple, direct and automation-ready procedures for measuring protein or peptide concentration are very desirable. BioAssay Systems' QuantiFluo™ protein assay kit is based on an improved o-phthalaldehyde method. This reagent reacts with primary amines in protein or peptide and forms a blue fluorescent product, allowing detection of nanograms of proteins. The fluorescence intensity ( $\lambda_{ex/em} = 360/450\text{nm}$ ) is proportional to the protein concentration in the sample.

#### KEY FEATURES

*Fast and sensitive.* Assay is completed within a few minutes. Linear detection range of 1.3 - 1000  $\mu\text{g/mL}$  BSA.

*Convenient and high-throughput.* Homogeneous "mix-incubate-measure" type assay. No wash and reagent transfer steps are involved. Can be readily automated on HTS liquid handling systems for processing thousands of samples per day.

#### APPLICATIONS

For quantitative determination of protein in various biological samples.

#### KIT CONTENTS

**Reagent:** 20 mL

**Standard:** 1 mL 1mg/mL BSA

*Storage conditions:* This product is shipped at room temperature. For long-term storage, keep kit at  $-20^{\circ}\text{C}$ . Shelf life of 6 months after receipt.

*Precautions:* Reagents are for research use only. Normal precautions for laboratory reagents should be exercised while using the reagents. Please refer to Material Safety Data Sheet for detailed information.

#### ASSAY PROCEDURE FOR 96-WELL PLATE READER

Use black flat-bottom 96-well plates. Prior to assay, bring all reagents to room temperature.

*Sample preparation.* Cultured cells can be assayed directly in a black culture plate. Remove culture media, wash cells 3 times with 1x phosphate-buffered saline (PBS). Alternatively cells (e.g.  $2 \times 10^6$ ) can be lysed by homogenization, followed by centrifugation at 14,000 rpm for 5 min. Use clear supernatant for assay.

Tissue (20 mg) can be homogenized in 200  $\mu\text{L}$  ice-cold water, followed by centrifugation at 14,000 rpm for 5 min.

Samples not measured on the same day can be stored frozen at  $-20^{\circ}\text{C}$ .

*Note: (1). This assay is compatible with most detergents, chelators and buffer components. Primary amine-containing buffers (e.g. Tris, glycine) should be avoided, if possible. For best results, include the same concentration of the sample buffer in the standards and blank. (2). If the sample protein concentration is higher than 1000  $\mu\text{g/mL}$ , dilute sample in water and repeat the assay. Multiply result by the dilution factor.*

1. *Standards.* Dilute standards in  $\text{H}_2\text{O}$  as follows.

No	Standard + $\text{H}_2\text{O}$	Standard ( $\mu\text{g/mL}$ )
1	100 $\mu\text{L}$ + 0 $\mu\text{L}$	1000
2	60 $\mu\text{L}$ + 40 $\mu\text{L}$	600
3	30 $\mu\text{L}$ + 70 $\mu\text{L}$	300
4	0 $\mu\text{L}$ + 100 $\mu\text{L}$	0

Transfer 10  $\mu\text{L}$  standards into separate wells of the plate.

Transfer 10  $\mu\text{L}$  of each sample in separate wells of the plate.

2. *Assay.* Add 90  $\mu\text{L}$  Reagent to all wells. Immediately tap plate to mix. Incubate 5 min at room temperature. Measure fluorescence intensity at 360/450nm on a plate reader. It is best to read all samples at the same time interval after mixing.

*Note: 1. The standard protocol uses a Sample:Reagent ratio of 1:9. Higher sensitivity can be achieved by using higher sample volume (e.g. 1:1 or 5:1).*

*2. For cells cultured in 384-well plates, add 40  $\mu\text{L}$  Reagent to assay wells.*

#### CALCULATION

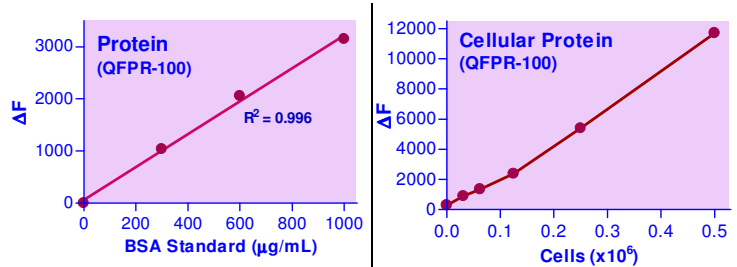
Plot the protein standard curve and determine its Slope. The protein concentration of a Sample is calculated as

$$[\text{Protein}] = \frac{F_{\text{SAMPLE}} - F_{\text{BLANK}}}{\text{Slope}} \quad (\mu\text{g/mL})$$

where  $F_{\text{SAMPLE}}$  and  $F_{\text{BLANK}}$  are the fluorescence intensity values of the Sample and the blank (i.e. #4  $\text{H}_2\text{O}$ ), respectively.

#### MATERIAL REQUIRED BUT NOT PROVIDED

Pipetting devices, centrifuge tubes, black flat bottom 96-well plates and plate reader.



*Left:* Standard curve performed on a 96-well plate reader (Spectramax M2);

*Right:* Human liver HepG2 cells were plated in a black 96-well tissue culture plate.

After overnight culture, cells were washed 3 times with 200  $\mu\text{L}$  PBS. 90  $\mu\text{L}$  Reagent were added to each well. Fluorescence was read after a 5-min incubation.

#### LITERATURE

- Kutchai H, Geddis LM (1977). Determinations of protein in red cell membrane preparations by o-phthalaldehyde fluorescence. *Anal Biochem.* 77:315-9.
- Robrish SA, et al (1978). The use of the o-phthalaldehyde reaction as a sensitive assay for protein and to determine protein in bacterial cells and dental plaque. *Anal Biochem.* 84:196-204.
- Joys TM, Kim H (1979). o-Phthalaldehyde and the fluorogenic detection of peptides. *Anal Biochem.* 94:371-7.

