

## FluoroQuest™ Fluorescence Quantum Yield Determination Kit

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 20	Keep in freezer Avoid exposure to light	Fluorescence Spectrophotometer Fluorescent microplate readers

### Introduction

When a fluorophore absorbs a photon of light, an energetically excited state is formed. The fate of this species is varied, depending upon the exact nature of the fluorophore and its surroundings, but the end result is deactivation (loss of energy) and return to the ground state. The main deactivation processes which occur are fluorescence (loss of energy by emission of a photon), internal conversion and vibrational relaxation (non-radiative loss of energy as heat to the surroundings), and intersystem crossing to the triplet manifold and subsequent non-radiative deactivation. The fluorescence quantum yield is the ratio of photons absorbed to photons emitted through fluorescence. In other words the quantum yield gives the probability of the excited state being deactivated by fluorescence rather than by another, non-radiative mechanism. The kit provides all the essential components for fluorescence quantum yield determination for biological conjugates. It is optimized for determining the fluorescence quantum yields of fluorescent conjugates of proteins, peptides, nucleotides and nucleic acids.

### Kit Components

Components	Amount	Fluorescence Quantum Yield in Water ( $\Phi_s$ value)	Conjugates Labeled with the Following Dyes or the Dyes of Similar Wavelength (for the Reference Standard Selection)
Reference A	1 ml	0.98	FAM, 6-TET, 6-HEX, 6-JOE, FITC, Cy®2, Alexa Fluor® 488 and 514, iFluor™ 488 and 514, DyLight™ 488 or other dyes that have emission of $500 \pm 50$ nm
Reference B	1 ml	0.20	Cy3®, Alexa Fluor® 514, 532, 546 and 555, iFluor™ 514, 532 and 555, DyLight™ 555, TRITC or other dyes that have emission of $550 \pm 50$ nm
Reference C	1 ml	0.44	Texas Red®, Texas Red®-X, Alexa Fluor® 594, iFluor™ 594, California Red™, DyLight™ 594 or other dyes that have emission of $600 \pm 50$ nm
Reference D	1 ml	0.24	Cy5®, Cy5.5®, Cy7®, Alexa Fluor® 633, 647, 700 and 750, iFluor™ 633, 647, 700 and 750, DyLight™ 650, 680 and 755 or other dyes that have emission of $650 \pm 50$ nm

### General Experimental Considerations:

The most reliable method for recording fluorescence quantum yield is the comparative method of Williams *et al* (Ref 1). It involves the use of well characterized standard samples with known fluorescence quantum yield values. Essentially, solutions of the standard and test samples with identical absorbance at the same excitation wavelength can be assumed to be absorbing the same number of photons. Hence, a simple ratio of the integrated fluorescence intensities of the two solutions (recorded under identical conditions) will yield the ratio of the quantum yield values. Since the fluorescence quantum yield for the standard sample ( $\Phi_s$ ) is known, it is trivial to calculate the fluorescence quantum yield for the sample ( $\Phi_x$ ).

The standard reference should be chosen to have certain absorbance at the excitation wavelength of choice for the test sample, and, if possible, emit in a similar region to the test sample. Standard 10 mm path length fluorescence cuvettes are sufficient for running the fluorescence measurements. For small amount samples, either a microplate reader or nanodrop device can be used. The absorbance in the 10 mm cuvette should be  $< 0.2$ . Above this level, non-linear effects may be observed due to inner filter effects, and the resulting quantum yield values may be inaccurate.

This kit is only designed for a one-point quick estimation of the fluorescence quantum yield of a biological conjugate. For more accurate deamination of fluorescence quantum yield please refer to the references cited since there are many factors that have a significant effect on fluorescence quantum yields.

**Procedure:**

- 1) Make 1 mM stock solutions of the selected standard and test samples using proper solvent (such as water or DMSO).
- 2) Record the absorbance spectra of the selected standard and test samples (0.1 to 5  $\mu$ M) using the diluted solutions prepared from their stock solutions (from Step #1) with an aqueous buffer (such as PBS or TRIS).
- 3) Select the excitation wavelength based on the absorption maxima (from Step#2). We recommend to use this following formula to select the desired excitation wavelength:

$$\text{Desired excitation (nm)} = \text{the shorter absorption maximum} - 50 \text{ nm}$$

(\*Selected from either the reference standard or test sample. For example, if your test sample has absorption maximum at 500 nm, and the selected reference standard has the absorption maximum at 490 nm, you should select 440 nm (490nm -50 nm) as excitation wavelength)

- 4) Adjust the concentrations of the selected reference standard and test samples to have the same absorbance (between 0.2 to 0.8) at the selected excitation wavelength.  
*Note: It might be easier to choose the isosbestic point of the reference standard and test sample absorption spectra.*
- 5) Dilute both the selected reference standard and test sample by 10 times (must be diluted by the **same** proportion) using the same buffer or pure water.
- 6) Record the fluorescence spectra of the same diluted solutions (from Step #5) in the 10 mm fluorescence cuvette using a fluorescence spectrophotometer or a microplate reader that can generate corrected fluorescence spectrum.  
*Note: that all the fluorescence spectra must be recorded with the same instrument settings. Changing instrument settings between samples will invalidate the quantum yield measurement.*
- 7) Calculate the integrated fluorescence intensity (the area of the fluorescence spectrum) from the fully corrected fluorescence spectrum (see your instrument instruction).
- 8) Calculate the fluorescence quantum yield using the following equation:

$$\Phi_x = \Phi_s \frac{A_x}{A_s}$$

$\Phi_x$  and  $\Phi_s$  are the fluorescence quantum yields of the selected reference standard and test sample respectively.  $A_x$  and  $A_s$  are the integrated fluorescence intensity (the curve area) of selected reference standard and test sample respectively.

**References:**

- 1) A. T. R. Williams, S. A. Winfield and J. N. Miller, Relative fluorescence quantum yields using a computer controlled luminescence spectrometer, *Analyst*, 1983, **108**, 1067.
- 2) S. Dhimi, A. J. de Mello, G. Rumbles, S. M. Bishop, D. Phillips and A. Beeby, Phthalocyanine fluorescence at high concentration: dimers or reabsorption effect? *Photochem. Photobiol.*, 1995, 61, 341.
- 3) J. R. Lakowicz, *Principles of Fluorescence Spectroscopy*, Kluwer Academic/Plenum Press, New York, 1999, Second Edition.
- 4) J. C. Scaiano (Ed.), *Handbook of Organic Photochemistry*, CRC Press, 1989.