

## Technical Bulletin #443

# Calculating Oxygen Concentration from Fluorescence Data on the BD™ Oxygen Biosensor System

Mark Timmins, Ph.D. and Tariq Haq  
BD Biosciences Discovery Labware, Bedford, MA 01730

## Introduction

The oxygen-sensitive fluorophore in the BD™ Oxygen Biosensor System (BD™ OBS) is one whose fluorescence is quenched by molecular oxygen in a predictable fashion. For many applications, knowing the relative concentration of oxygen is sufficient. For certain applications, it may be advantageous to be able to calculate actual oxygen concentrations from the fluorescence. Understanding the theory which relates oxygen concentration to signal will allow the user to make such calculations.

## Stern-Volmer Theory

For the fluor used in the BD OBS, and others like it, fluorescence is inversely proportional to the concentration of oxygen in the vicinity of the fluor. The behavior of fluorophores quenched this way follows a mathematical model called the Stern-Volmer equation. At its simplest, the Stern-Volmer equation takes the following form:<sup>1</sup>

$$[1] \quad I_0 / I = 1 + K_{SV} * [O_2]$$

where  $[O_2]$  is the oxygen concentration,  $K_{SV}$  is the Stern-Volmer constant,  $I$  is the fluorescence intensity at  $[O_2]$  and  $I_0$  is the fluorescence intensity at the reference condition (zero oxygen).

## Definitions

### Normalized Relative Fluorescence (NRF)

The ratio of fluorescence at the oxygen concentration of interest to that at the ambient reference condition.  $NRF = I / I_A$ , where  $I$  indicates fluorescence intensity, and the  $A$  subscript represents "ambient".

### Dynamic Range (DR)

The maximum possible signal occurs at zero oxygen, so the DR is the ratio of fluorescence at zero oxygen to that at ambient oxygen, or  $DR = I_0 / I_A$ . For a given set of experimental conditions (temperature, wavelengths, etc.), DR will be a constant. To convert fluorescence to actual oxygen concentration, it is necessary to know DR.

## Use of Ambient as the Reference Condition

Zero oxygen has traditionally been chosen as the reference condition because it yields the simplest form of *Equation 1*, although the choice is entirely arbitrary. Because it is impractical to collect fluorescence intensities in a zero-oxygen environment, "ambient" oxygen is used as the reference condition. Thus, we normalize fluorescence intensities to those at ambient oxygen concentration instead of zero oxygen (see definition of "NRF" above).

The use of an alternate reference condition changes the form of *Equation 1* slightly, but not its fundamental essence, as will be shown in the following discussion. At ambient  $[O_2]$ , the Stern-Volmer equation takes the form:

$$[2] \quad I_0 / I_A = 1 + K_{SV} * [O_2]_A$$

where the subscript "A" designates ambient.

$I_0 / I_A = DR$ , so *Equation 2* can be rearranged to yield:

$$[3] \quad K_{SV} = (DR - 1) / [O_2]_A$$

Thus, by measuring DR, it is possible to calculate  $K_{SV}$ . It can be seen that DR and  $K_{SV}$  are related constants.

Furthermore, it can be shown that the ratio of DR to NRF is equivalent to  $I_0 / I$ . Substituting this into *Equation 1* yields:

$$[4] \quad DR / NRF = 1 + K_{SV} * [O_2]$$

*Equation 4* is equivalent to *Equation 1*. Solving for NRF yields:

$$[5] \quad NRF = DR / (1 + K_{SV} * [O_2])$$

*Equation 5* is the expression of the Stern-Volmer law pertinent to typical applications of the BD OBS. It relates the NRF signal to oxygen concentration. It can be seen from *Equation 5* that fluorescence (NRF) varies inversely with  $[O_2]$ , and that  $[O_2]$  may be calculated from NRF if  $K_{SV}$  and DR are known.

*continued*

## BD Biosciences

Clontech  
Discovery Labware  
Immunocytometry Systems  
Pharminggen



## How to Determine DR and $K_{SV}$

DR may be determined by a simple experiment, or from the sodium sulfite positive control well (100 mM sodium sulfite in PBS buffer) we recommend for each plate. Dynamic range will be the ratio between the signal once it has reached its maximum (typically 30-60 minutes) and that same well's fluorescence at ambient conditions. Stated another way, DR is the NRF of the sulfite well, once the sulfite has removed all the oxygen from the well. DR may then be used to calculate  $K_{SV}$ , per *Equation 3*.

DR is a characteristic property of the sensor and how it is formulated, and should thus be constant for all wells. Hence, it is possible to calculate DR for a plate from a single well. DR does vary slightly with temperature and from reader to reader, however, so it should be calculated on the same reader, by the same method and under the same experimental conditions as the experimental data it will be used to analyze.

A typical value of DR is 6, which yields a  $K_{SV}$  of 23.9 atm<sup>-1</sup> (i.e., for  $[O_2]_A = 0.209$  atm) or 0.031 mmHg<sup>-1</sup> (for  $[O_2]_A = 159$  mmHg).

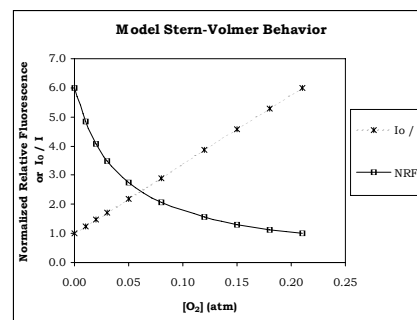
## Calculating Oxygen Concentrations

Using *Equation 5*, it is possible to convert fluorescence values into actual oxygen concentrations as shown in *Figure 1*. To solve for  $[O_2]$ , simply rearrange *Equation 5* to yield:

$$[6] \quad [O_2] = (DR / NRF - 1) / K_{SV}$$

Thus, it is possible to calculate  $[O_2]$  from NRF, provided one knows the values of DR and  $K_{SV}$ . These two constants (defined above) may be determined experimentally from a single experiment.

Because BD™ OBS is in equilibrium with the atmosphere, it should be noted that the concentration of oxygen in the well depends on the rate at which it is being consumed by the content of the well (e.g., the cells), as well as the rate at which oxygen diffuses in from the atmosphere. We are still developing a simple method that will allow one to factor out the diffusion term and thereby determine the rate of oxygen consumption in the well.



**Figure 1:** Theoretical response of a Stern-Volmer fluorophore (such as that utilized in the BD Oxygen Biosensor System) to varying concentrations of oxygen. It can be seen that signal varies inversely with oxygen concentration. This example was calculated using the equations outlined in the text, assuming DR = 6.

## Reference

1. Stern, O. and Volmer, M., Phys. Z., **20:183** (1919).

For additional information on the BD Oxygen Biosensor System, visit our website at:

[www.bdbiosciences.com/discovery\\_labware/Products/drug\\_discovery/oxygen\\_biosensor\\_system](http://www.bdbiosciences.com/discovery_labware/Products/drug_discovery/oxygen_biosensor_system)

**BD Biosciences**  
Two Oak Park  
Bedford, MA 01730 USA  
tel: 800.343.2035  
fax: 800.743.6200

**BD**  
2771 Bristol Circle  
Oakville, Ontario  
Canada L6H 6R5  
tel: 905.855.5550  
fax: 905.829.5405

**Nippon BD**  
Akasaka DS Bldg.  
5-26 Akasaka 8-chome  
Minato-ku, Tokyo 107 Japan  
tel: (81) 24 593 5405  
fax: (81) 24 593 5761

**BD Biosciences**  
Singapore Branch  
30 Tuas Avenue 2  
Singapore 639461  
tel: (65) 6861 0633  
fax: (65) 6860 1590

**BD Biosciences**  
Discovery Labware Europe  
11 rue Aristide Bergès BP 4  
38800 Le Pont De Claix, France  
tel: (33) 04 76 68 32 98  
fax: (33) 04 76 68 55 10

**BD Biosciences**  
4 Research Park Drive  
Macquarie University Research Park  
North Ryde NSW 2113 Australia  
tel: (612) 8875 5239  
fax: (612) 8875 7200

