



# Fast and accurate automated method for free sulfite analysis in wine

## Authors

Eric Naigeon,<sup>1</sup> Marco Rastetter,<sup>2</sup>  
Mari Klemm,<sup>3</sup> and Annu  
Suoniemi-Kähärä<sup>3</sup>

<sup>1</sup>Thermo Fisher Scientific,  
Paris, France

<sup>2</sup>Thermo Fisher Scientific,  
Dreieich, Germany

<sup>3</sup>Thermo Fisher Scientific,  
Vantaa, Finland

## Keywords

Discrete Analyzer, Photometric  
Analyzer, Gallery Analyzer, Arena

## Goal

To develop a method for free sulfite  
analysis in wine using automated  
discrete photometry

## Introduction

Sulfur dioxide (SO<sub>2</sub>) is added to control the process of wine making. It serves many useful functions, for example, it acts as an enzyme inhibitor in musts preventing juice from browning. As a microbiological control agent, SO<sub>2</sub> is added to the winemaking process to prevent oxidation in the finished product.

Sulfur dioxide can be found in wine in its free forms, SO<sub>2</sub> (gas) and bisulfate ion (HSO<sub>3</sub><sup>-</sup>), or bound to compounds that incorporate a carbonyl group, such as acetaldehyde. Free forms of SO<sub>2</sub> are pH and temperature dependent and because of the acidic nature of wines, SO<sub>2</sub> is usually present and measured as bisulfate ion (HSO<sub>3</sub><sup>-</sup>). Results are reported as SO<sub>2</sub>.

The presence of total SO<sub>2</sub>, both free and bound, is regulated and, as a result, a warning statement is required on wine labels because sulfite is considered an allergen. The European Union established a maximum permitted level of total SO<sub>2</sub> in wine varying from 150 to 500 mg/L which is dependent upon the sugar level of the product.

In the USA, the maximum level of total SO<sub>2</sub> permitted is 350 mg/L.<sup>1</sup> The measurement of both total and free SO<sub>2</sub> can be automated using Thermo Scientific™ system reagents and Thermo Scientific™ Gallery™ discrete analyzers<sup>2</sup>.

In this study, an automated method to measure free SO<sub>2</sub> in wine samples is presented. The method is based on the reaction between sulfur dioxide, *p*-rosaniline hydrochloride, and formaldehyde. This method is designed to use optimal reagent concentrations and volumes to provide accurate results. The concentration of free SO<sub>2</sub> in the sample is calculated automatically from a calibration curve. This method enables a laboratory to fully automate SO<sub>2</sub> determinations and replace traditional time consuming reference and distillation methods.

As a result of using bar-coded system reagents, this new automated method to measure free SO<sub>2</sub> is very quick and easy to use. Analysis of 60 samples takes 35 min with only an additional 10 min for daily calibration and analyzer start-up. From the same samples, various sugars, acids, color, and total SO<sub>2</sub> can be run automatically.<sup>3</sup> Compared to Flow Injection Analysis (FIA), the photometric method requires only small volumes of reagents, thus becoming the more economically and environmentally friendly choice.<sup>4</sup>

## Experimental

### Materials and methods

#### Instruments

A Gallery discrete analyzer was used for this study.

Reference values were obtained from a commercially available FIA analyzer and the reference method was based on *p*-rosaniline hydrochloride and formaldehyde reactions.

#### Method principle to analyze free SO<sub>2</sub>

The method is based on the reaction between SO<sub>2</sub>, *p*-rosaniline hydrochloride, and formaldehyde and is performed at 37 °C, using a 575 nm filter and a 700/750 nm filter for side wavelength. The purpose of measuring a side wavelength is to remove the effect of bubbles which may appear in cuvettes. The side wavelength is determined from the spectrum area where no reaction occurs. See Figure 1 for the reaction details.

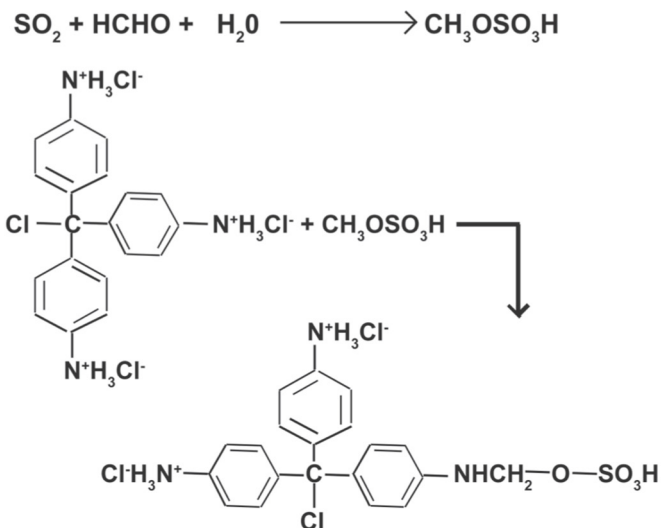


Figure 1. Method principle to analyze free SO<sub>2</sub>

#### Application to wine samples

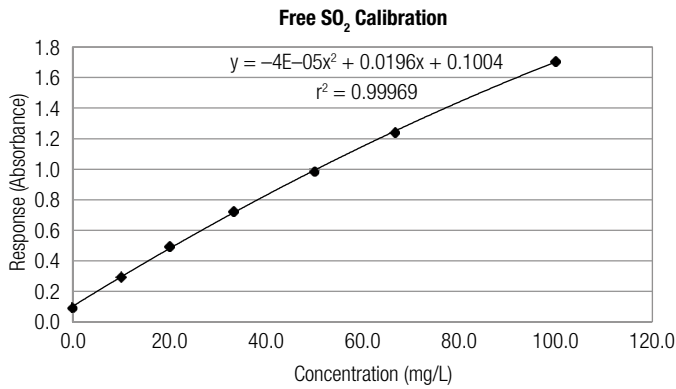
The automated application to measure free SO<sub>2</sub> on Gallery discrete analyzers consists of two reagents, an end-point measurement with a sample blank, and a second order calibration curve used for result calculation. The measuring range of the method is from 2 to 100 mg/L and can be extended by automated dilution of the sample.

First, 100 µL of SO<sub>2</sub> Free R1 reagent and 10 µL of sample with 10 µL of water are incubated for 30 seconds and the sample color is blanked. Then, 100 µL of R2 reagent is added. After 75 seconds of incubation the reaction is measured at 575 nm with a side wavelength measurement at 700/750 nm.

#### Reagents and calibrator

Thermo Scientific system reagents for Gallery discrete analyzers were used. The SO<sub>2</sub> Free kit is available using the P/N 984634.

Calibration was performed with a self-made standard prepared from sodium metabisulfite (Na<sub>2</sub>S<sub>2</sub>O<sub>5</sub> MW = 190.11 g/mol). The standard concentration was 200 mg/L and calibration points were automatically diluted by the analyzer (Figure 2).



**Figure 2. Method calibration for free SO<sub>2</sub> using an automated dilution of the calibration stock solution**

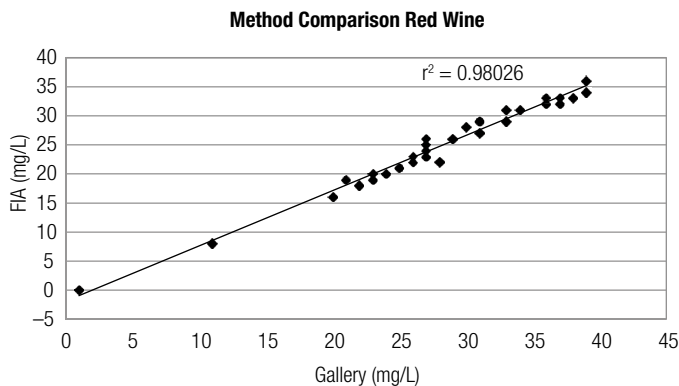
### Samples

Fifty-seven (57) wine samples were analyzed. Sample types included both red (N = 35) and white wines (N = 22). Reference values were determined using the *p*-rosaniline method on a FIA analyzer. No sample pretreatment was done prior to the analysis.

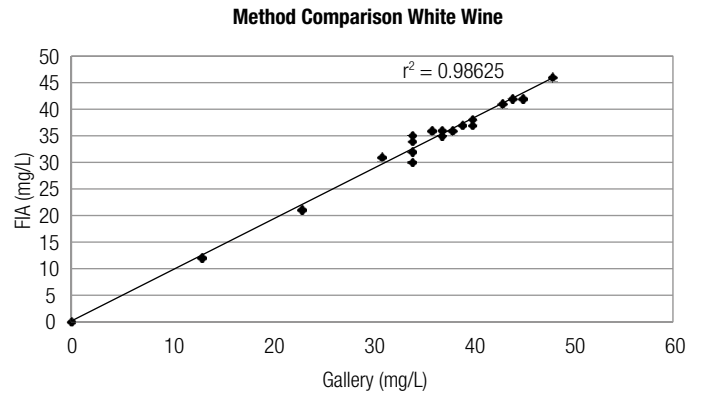
### Results and discussion

The results of free SO<sub>2</sub> in 35 red wines and 22 white wines using the Gallery discrete analyzer and FIA are compared in Table 1 (page 4). The concentration of free SO<sub>2</sub> in samples varied from 0 to 46 mg/L with an average of 28 mg/L. The average bias was +3 mg/L when calculated using red wine samples and +1 mg/L when calculated using white wine samples.

The method correlation study is shown graphically in Figures 3 and 4. Concentrations of red wine samples varied from 20 to 40 mg/L and concentrations of white wine samples from 35 to 50 mg/L.



**Figure 3. Red wine sample comparison between Gallery and FIA methods**



**Figure 4. White wine sample comparison between Gallery and FIA methods**

Very good correlation between the Gallery discrete analyzer and the FIA analyzer was shown by both sample types.

The coefficient of determination was slightly better for white wines than for red wines, respectively  $r^2 = 0.9863$  and  $0.9803$ . When the data from all samples were combined into a single plot, the correlation coefficient ( $r^2$ ) was  $0.9755$ .

In addition, precision for one of the red wine samples (N = 40) was calculated (data not shown) using the Gallery method and was 1.8% within run, and 2.1% between runs, which correlates with the allowed EBC/ASBC standard ranges.

### Conclusions

The automated photometric Thermo Scientific SO<sub>2</sub> Free method correlates well with the FIA method with a coefficient of determination ( $r^2$ ) of  $0.9755$ . The concentration of samples varied from 0 to 46 mg/L with an average of 28 mg/L. Samples (N = 57) including different types of wines were tested. Precision for the red wine samples (N = 40) was 1.8% within run, and 2.1% between runs.

This new automated SO<sub>2</sub> Free method is very quick and easy to use. Analysis of 60 samples takes only 35 min and allows simultaneous analysis of various sugars and acids, color, and total SO<sub>2</sub>. Compared to the FIA method, the photometric method requires only small volumes of reagents, thus being the more economically and environmentally friendly choice.

**Table 1. Bias of the two Free SO<sub>2</sub> methods**

Red Wine Sample	Gallery (mg/L)	FIA (mg/L)	Bias (mg/L)	White Wine Sample	Gallery (mg/L)	FIA (mg/L)	Bias (mg/L)
Red Wine 1	38	33	5	White Wine 1	40	37	3
Red Wine 2	1	0	1	White Wine 2	43	41	2
Red Wine 3	11	8	3	White Wine 3	39	37	2
Red Wine 4	20	16	4	White Wine 4	34	34	0
Red Wine 5	22	18	4	White Wine 5	37	35	2
Red Wine 6	26	22	4	White Wine 6	34	30	4
Red Wine 7	23	20	3	White Wine 7	38	36	2
Red Wine 8	23	19	4	White Wine 8	34	32	2
Red Wine 9	24	20	4	White Wine 9	40	38	2
Red Wine 10	31	29	2	White Wine 10	37	35	2
Red Wine 11	29	26	3	White Wine 11	36	36	0
Red Wine 12	27	26	1	White Wine 12	34	35	-1
Red Wine 13	37	32	5	White Wine 13	34	35	-1
Red Wine 14	39	36	3	White Wine 14	13	12	1
Red Wine 15	39	34	5	White Wine 15	23	21	2
Red Wine 16	36	33	3	White Wine 16	0	0	0
Red Wine 17	34	31	3	White Wine 17	31	31	0
Red Wine 18	36	33	3	White Wine 18	48	46	2
Red Wine 19	33	29	4	White Wine 19	37	36	1
Red Wine 20	31	29	2	White Wine 20	39	37	2
Red Wine 21	25	21	4	White Wine 21	45	42	3
Red Wine 22	21	19	2	White Wine 22	44	42	2
Red Wine 23	21	19	2				
Red Wine 24	36	32	4				
Red Wine 25	28	22	6				
Red Wine 26	30	28	2				
Red Wine 27	26	23	3				
Red Wine 28	27	23	4				
Red Wine 29	27	24	3				
Red Wine 30	27	25	2				
Red Wine 31	33	31	2				
Red Wine 32	36	33	3				
Red Wine 33	37	33	4				
Red Wine 34	31	27	4				
Red Wine 35	25	21	4				

## References

1. Jackowetz, N; Li, E; Mira de Orduña, R. *Sulphur Dioxide Content of Wines: the Role of Winemaking and Carbonyl Compounds*. Research Focus, Cornell Viticulture and Enology Program, 2011–3.
2. Automated Discrete Photometry. [Online] <http://www.thermoscientific.com/en/products/environmental-industrial-testing-kits.html> (accessed November 3, 2014).
3. Cudiamat, G. Continuous Flow Analysis & Discrete Analyzers. LC/GC [Online] 2010, <http://www.chromatographyonline.com/lcgc/Research/Continuous-Flow-Analysis-amp-Discrete-Analyzers/ArticleStandard/Article/detail/678395> (accessed Dec 10, 2014).
4. Hicks, M.; Schenken, J.; Steinrauf, M. *Laboratory Instrumentation*, 3rd Edition; J. B. Lippincott & Co.: Pennsylvania, 1987; pp 284–292.

Find out more at [thermofisher.com/discreteanalysis](http://thermofisher.com/discreteanalysis)

**ThermoFisher**  
SCIENTIFIC