

EN

SO₂ Free

REF 984634

● 3 x 20 ml Reagent 1
● 3 x 20 ml Reagent 2

INTENDED USE

Reagent for photometric determination of free Sulfur Dioxide in homogenous liquid samples using automated Thermo Scientific Arena or Gallery analyzer.

METHOD

The method is based on the reaction between sulfur dioxide, pararosaniline and formaldehyde. Method is performed at 37 °C, using 575 nm filter and for side wavelength 700 or 750 nm filter.

REAGENT INFORMATION

Ready-to-use reagent		Barcode id
Reagent 1	3 x 20 ml	A40
Reagent 2	3 x 20 ml	A41

Note: Labels of reagent vials have two barcodes. For Arena analyzers, turn the short barcode to the barcode reader. For Gallery analyzers, turn the long barcode to the barcode reader.

Concentrations

Pararosaniline ≤ 0.1 %
Formaldehyde ≤ 0.2 %

Precautions

SO₂ Free R2 is hazardous. See separate sheet inside the kit for Hazardous- and Precautions-phrases: H314, P280, P303 + P361 + P353, P305 + P351 + P338. EUH208: Contains formaldehyde. May produce an allergic reaction.

Exercise the normal precautions required for handling all laboratory reagents.

The product has to be disposed of as laboratory chemical in accordance with local regulations.

Preparation

The reagents are ready-to-use.

Note: Check that there are no bubbles on the surface of the reagent when you insert vials into the analyzer.

Storage and Stability

Reagents in unopened vials are stable at 2...8 °C until the expiry date printed on the label. Do not freeze the reagents.

Refer to the Application Notes of your analyzer for the on board stability of reagents.

SAMPLES**Sample Type**

Food and other sample material e.g wine.

Sample concentration and Arena/Gallery application

All method related details are in the separate application note.

Sample preparation

If the sample has substances interfering the measurement, please handle it according to the following suitable preparation procedure:

Note: Polyvinylpyrrolidone (PVP) cannot be used with this reagent because of the inhibition of color forming.

- Use clear, colorless and practically neutral liquid samples directly.
- Filter or centrifuge turbid solutions.

TEST PROCEDURE

See the separate Arena or Gallery System Application note for an automated procedure. Due to the differences in sample matrixes, all performance should be evaluated by the user.

Materials required but not provided

Distilled water (aseptic and free of heavy metals) and general laboratory equipment.

Calibration

A fresh solution is used for the calibration. Weigh precisely 0.0303 g of sodium metabisulfite (Na₂S₂O₅ e.g. S9000/ Sigma-Aldrich, MW = 190.11 g/mol, purity ≥ 98 %) into a 250 ml volumetric flask and fill up to

mark with distilled water. The solution has a sulfur dioxide concentration of 80 mg/l. The standard must be used fresh.

Quality Control

Use quality control samples at least once a day and after each calibration and every time a new bottle of reagent is used. It is recommended to use two level of controls. The control intervals and limits must be adapted to the individual laboratory requirements. The results of the quality control sample(s) should fall within the limits preset by the laboratory.

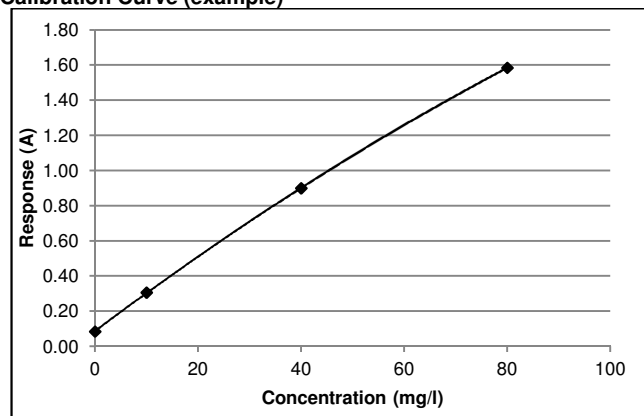
CALCULATION OF RESULTS

The results are calculated automatically by the analyzer using a calibration curve.

Conversion factors:

$$\text{mmol/l} \times 64.054 = \text{mg/l}$$

$$\text{mg/l} \times 0.0156 = \text{mmol/l}$$

Calibration Curve (example)

Calibrator	Response (A)	Conc. (mg/l)
Water	0.084	0
SO ₂ F 10	0.305	10
SO ₂ F 40	0.899	40
SO ₂ F 80	1.584	80

Note that the calibration curve is lot dependent. This calibration curve is performed by Gallery analyzer.

LIMITATIONS OF THE PROCEDURE**Interference**

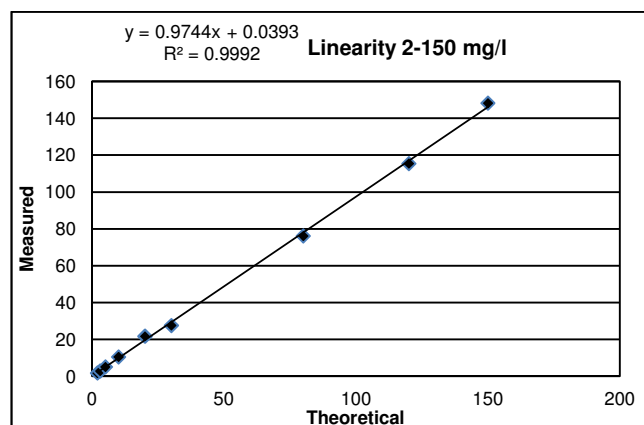
The determination is specific for free sulfur dioxide. No interference was observed.

MEASURING RANGE

The test has been developed to determine free sulfur dioxide concentrations within a measuring range from 2 to 150 mg/l.

PERFORMANCE CHARACTERISTICS

The results obtained in individual laboratories may differ from the performance data given. Linearity testing has been performed with water based solutions. Different matrixes may change the linearity limits of the test.



Determination limit (=Test limit low)

The determination limit is the lowest concentration that can be measured quantitatively. The determination limit for this method is 2 mg/l.

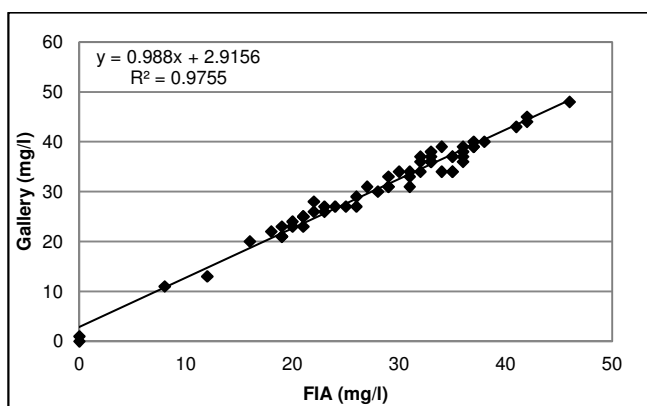
Precision**Gallery analyzer**

	White wine		Rosé wine		Red wine	
	N	39	N	40	N	40
	Mean	31	Mean	5	Mean	24
	SD	SD	SD	CV %	SD	CV %
Within run	0.480	1.5 %	0.168	3.4 %	0.442	1.8 %
Between run	0.752	2.4 %	0.156	3.2 %	0.502	2.1 %
Total	0.892	2.9 %	0.230	4.6 %	0.669	2.8 %

A precision study was performed using the Gallery analyzer, with the number of measurements being n = 40.

Method Comparison

Graph below shows the correlation of wine samples measured with Gallery system SO2 Free method and similar type of FIA pararosaniline method.

**OTHER REMARKS**

Note that part of the application performance has been verified with pure chemicals dissolved in deionized water. The results obtained in individual laboratories may differ from the given performance data due to e.g. sample matrix, concentrations or analysis environment. Each laboratory is responsible to verify the method to prove the analysis performance.

WASTE MANAGEMENT

Please refer to local legal requirements. It is recommended to empty the analyzer cuvette waste bin and waste water daily. Emptying should be done immediately after the analysis when using hazardous reagents/solutions.

Note: If using reagents/solutions that react with each other, cuvette waste bin and waste water should be emptied and washed between use of these reagents.

ADDITIONAL MATERIAL

Certificate of analysis and SDS are available at www.e-labeling.eu/TSF

Applications for Gallery and Arena automated analyzers are available upon request from the local sales representative. Information in the Application note can change without prior notice.

MANUFACTURER

Thermo Fisher Scientific Oy
Ratastie 2, P.O. Box 100, FI-01621 Vantaa, Finland
Tel. +358 10 329200

CONTACT INFORMATION

www.thermoscientific.com
e-mail: system.support.fi@thermofisher.com

Date of revision

2015-05-22

Changes from previous version

Precautions updated.

General updates.