

Analysis of Glass by X-Ray Fluorescence

ARL OPTIM'X WD-XRF Spectrometer

Key Words

- ARL OPTIM'X
- Glass
- WDXRF
- X-Ray Fluorescence

Introduction

The simplest form of glass is the single component fused silica (SiO_2). However it is both difficult to process and expensive. To reduce these difficulties, some other oxides are added imparting specific properties to the resultant glass.

Most of glasses are composed of about 70 % silica, which is a glass former, soda as a flux in the form of carbonate and sulfate (about 14 %), and lime as a stabilizer in the form of limestone (about 10 %). Other types of oxides like alumina or magnesia improve the physical characteristics of glass, particularly the resistance to atmospheric conditions.

In-depth coloring is obtained by incorporation of various metallic oxides: oxides of chromium, iron, manganese or copper.

Instrumentation



An ARL OPTIM'X XRF spectrometer from Thermo Electron Corporation has been used to derive limits of detection and precision for the analysis of glasses. The ARL OPTIM'X is a wavelength dispersive system which provides superior resolution and light elements capability. It is fitted with an Air-cooled Rh End-Window Tube with thin Be window (0.075 mm) and has a

maximum power of 50 Watts. Thanks to close coupling between the X-ray tube anode and the sample the performance of the ARL OPTIM'X is equivalent to a 200W conventional WD-XRF instrument. The instrument can be equipped with the unique SmartGonio™, a series of Multichromators™ or both.

Table 1 shows limits of detection for various elements in soda-lime glasses prepared as pressed powders.



Calibration and limits of detection

A series of pressed glass samples have been measured on an ARL OPTIM'X. Calibration curves have been derived by relating intensities for each oxide (or element) to concentrations in the standard samples. X-ray fluorescence measures elements, but the results can be related directly to the oxide forms of these elements when only one single form is present in the sample. Using the calibration curves, limits of detection have been derived using the SmartGonio for the most common oxides/elements found in soda-lime glasses (Table 1).

OXIDE/ ELEMENT	LINE	CRYSTAL	DETECTOR	LOD [PPM]
Na_2O	$\text{K}\alpha_{1,2}$	AX-06	FPC	100
MgO	$\text{K}\alpha_{1,2}$	AX-06	FPC	60
Al_2O_3	$\text{K}\alpha_{1,2}$	PET	FPC	47
SiO_2	$\text{K}\alpha_{1,2}$	PET	FPC	N.R.
P_2O_5	$\text{K}\alpha_{1,2}$	PET	FPC	48
SO_3	$\text{K}\alpha_{1,2}$	PET	FPC	23
Cl	$\text{K}\alpha_{1,2}$	PET	FPC	24
K_2O	$\text{K}\alpha_{1,2}$	LiF 200	FPC	14
CaO	$\text{K}\alpha_{1,2}$	LiF 200	FPC	12
TiO_2	$\text{K}\alpha_{1,2}$	LiF 200	FPC	12
Cr_2O_3	$\text{K}\alpha_{1,2}$	LiF 200	FPC	9
MnO	$\text{K}\alpha_{1,2}$	LiF 200	FPC	9
Fe_2O_3	$\text{K}\alpha_{1,2}$	LiF 200	FPC	9
ZnO	$\text{K}\alpha_{1,2}$	LiF 200	SC	3.6
SrO	$\text{K}\alpha_{1,2}$	LiF 200	SC	2.4
ZrO_2	$\text{K}\alpha_{1,2}$	LiF 200	SC	1.8
BaO	$\text{L}\beta_1$	LiF 200	FPC	51
PbO	$\text{L}\beta_1$	LiF 200	SC	9

Table 1: Analytical parameters and limits of detection for various oxides/element in soda-lime glass (100 sec. counting time)

N.R. = LoD is not relevant for major elements

FPC = flow proportional counter

SC = scintillation counter

Excitation conditions: 40 kV / 1.25 mA

Collimator: 0.29°

Precision tests

Precision tests have been carried out by analyzing repeatedly the same pressed pellet sample for eleven consecutive analyses. Eighteen elements/oxides are determined using a counting time of 36 seconds per analytical line. The results are summarized below for two different glass samples (Tables 2 and 3). In the case when precision should be improved for some elements this counting time could be increased. Doubling the counting time would improve the precision by a factor of about 1.4 (square root of 2).

Sample A

RUN	Na ₂ O %	MgO %	Al ₂ O ₃ %	SiO ₂ %	K ₂ O %	CaO %	Fe ₂ O ₃ %	SO ₃ PPM	TiO ₂ PPM
Time (s)	36	36	36	36	36	36	36	36	36
Run 1	13.98	0.185	1.79	72.59	0.588	10.85	0.330	582	579
Run 2	13.93	0.193	1.81	72.60	0.582	10.82	0.333	640	563
Run 3	13.97	0.177	1.80	72.64	0.588	10.82	0.330	608	563
Run 4	14.01	0.178	1.80	72.64	0.582	10.87	0.330	645	581
Run 5	13.95	0.182	1.80	72.60	0.588	10.83	0.329	576	564
Run 6	13.94	0.177	1.81	72.61	0.589	10.82	0.329	573	569
Run 7	13.86	0.185	1.80	72.64	0.588	10.83	0.330	658	569
Run 8	13.92	0.186	1.81	72.59	0.585	10.84	0.331	652	566
Run 9	13.94	0.184	1.81	72.63	0.591	10.82	0.334	651	579
Run 10	13.98	0.183	1.80	72.63	0.586	10.87	0.332	617	526
Run 11	13.95	0.188	1.78	72.62	0.588	10.83	0.330	619	561
Avg.	13.95	0.183	1.80	72.62	0.587	10.84	0.331	620	565
Std.Dev.	0.04	0.005	0.01	0.02	0.003	0.02	0.0015	32	15

Table 2: Repeatability for the analysis of the major and minor oxides in Sample A

Sample B

RUN	Na ₂ O %	MgO %	Al ₂ O ₃ %	SiO ₂ %	K ₂ O %	CaO %	Fe ₂ O ₃ PPM	SO ₃ %	TiO ₂ PPM
Time (s)	36	36	36	36	36	36	36	36	36
Run 1	13.35	0.180	1.67	73.07	0.556	10.77	773.3	0.177	556.0
Run 2	13.33	0.180	1.68	73.08	0.564	10.76	757.9	0.181	568.0
Run 3	13.28	0.186	1.67	73.08	0.554	10.81	789.6	0.180	555.1
Run 4	13.28	0.185	1.66	73.11	0.559	10.83	768.2	0.186	587.2
Run 5	13.35	0.181	1.67	73.05	0.554	10.79	763.9	0.181	594.7
Run 6	13.32	0.172	1.67	73.11	0.566	10.80	767.3	0.186	541.4
Run 7	13.33	0.185	1.67	73.06	0.554	10.79	758.9	0.180	570.3
Run 8	13.26	0.185	1.69	73.04	0.555	10.78	771.7	0.185	565.2
Run 9	13.33	0.180	1.64	73.11	0.561	10.82	775.7	0.183	553.7
Run 10	13.30	0.193	1.68	73.08	0.556	10.80	764.1	0.188	572.6
Run 11	13.31	0.184	1.66	73.06	0.561	10.78	785.3	0.186	566.0
Avg.	13.31	0.183	1.67	73.08	0.558	10.80	770.5	0.183	566.4
Std.Dev.	0.03	0.01	0.01	0.03	0.004	0.02	10	0.003	15

Table 3: Repeatability for the analysis of the major and minor oxides in sample B

Conclusion

All limits of detection obtained show that the ARL OPTIM'X can deliver adequate analysis results, notably for bottle glass application. Repeatability of analysis is excellent for major and minor elements even for Na₂O and MgO. Longer counting time may be used in case elements present below 100 ppm need to be controlled precisely. These results show that the ARL OPTIM'X spectrometer is well suited to produce precision results for the determination of the main oxides and the coloring agents in glasses.

In addition to these offices, Thermo Fisher Scientific maintains a network of representative organizations throughout the world.

RUN	P ₂ O ₅ PPM	Cl PPM	Cr ₂ O ₃ PPM	MnO PPM	As ₂ O ₃ PPM	SrO PPM	ZrO ₂ PPM	BaO PPM	PbO PPM
Time (s)	36	36	36	36	36	36	36	36	36
Run 1	166.9	113.4	93.6	48.4	101.6	127.4	209.1	454.3	228.2
Run 2	146.3	129.5	91.4	44.7	101.5	124.8	204.6	392.6	218.9
Run 3	193.3	111.2	91.1	43.8	95.9	127.1	207.0	361.8	197.8
Run 4	199.2	104.6	96.2	29.9	103.8	127.0	205.4	375.7	234.6
Run 5	158.1	111.8	94.6	41.8	103.7	122.7	204.7	385.2	228.5
Run 6	171.3	107.9	85.2	49.5	95.0	126.3	203.8	355.4	194.8
Run 7	203.6	113.4	88.9	40.3	96.0	125.0	205.4	434.1	234.4
Run 8	190.4	135.6	94.5	44.7	96.8	125.9	203.4	315.1	207.3
Run 9	150.7	110.1	88.6	43.1	114.9	127.1	206.6	401.2	220.8
Run 10	255.0	104.0	83.6	41.9	99.2	125.8	206.2	402.2	214.6
Run 11	218.3	97.9	80.6	38.9	97.7	126.8	203.1	429.8	197.8
Avg.	186.6	112.7	89.9	42.4	100.5	126.0	205.4	391.6	216.1
Std.Dev.	32.5	11	5	5.2	5.7	1.4	1.8	39.7	15

RUN	P ₂ O ₅ PPM	CL PPM	CR ₂ O ₃ PPM	MNO PPM	AS ₂ O ₃ PPM	SRO PPM	ZRO ₂ PPM	BAO PPM	PBO PPM
Time (s)	36	36	36	36	36	36	36	36	36
Run 1	200.6	100.7	63.4	9.3	118.2	122.4	227.2	883.9	895.6
Run 2	159.5	111.2	65.9	3.9	112.0	122.2	226.1	960.6	914.5
Run 3	193.3	115.7	64.7	18.2	115.6	119.3	225.3	925.4	911.5
Run 4	156.6	103.5	74.7	8.7	105.6	126.5	225.1	891.3	900.0
Run 5	187.4	97.9	63.6	12.3	106.8	126.3	226.2	948.8	904.8
Run 6	183.0	114.6	59.5	14.7	115.2	125.5	226.3	960.5	904.6
Run 7	193.3	101.8	67.9	6.1	113.5	124.9	226.9	960.6	910.1
Run 8	191.8	95.2	66.9	9.3	101.6	124.7	227.4	918.0	919.7
Run 9	219.7	113.4	61.5	9.3	109.6	124.6	222.9	916.9	913.4
Run 10	243.2	101.3	64.4	1.0	103.4	123.0	226.0	980.8	912.7
Run 11	191.8	109.0	70.8	13.8	106.9	122.1	227.7	950.9	875.6
Avg.	192.8	105.8	65.8	9.7	109.9	123.8	226.1	936.2	905.7
Std.Dev.	24.2	7.2	4.3	4.9	5.4	2.2	1.4	31.2	12

Australia
+61 2 8844 9500

Austria
+43 1 333 50340

Belgium
+32 2 482 30 30

Canada
+1 800 532 4752

China
+86 10 5850 3588

Denmark
+45 70 23 62 60

France
+33 1 60 92 48 00

Germany
+49 6103 408 1014

India
+91 22 6742 9434

Italy
+39 02 950 591

Japan
+81 45 453 9100

Latin America
+1 608 276 5659

Netherlands
+31 76 587 98 88

South Africa
+27 11 570 1840

Spain
+34 91 657 4930

Sweden/Norway/Finland
+46 8 556 468 00

Switzerland
+41 21 694 71 11

UK
+44 1442 233555

USA
+1 800 532 4752

www.thermo.com



Thermo Electron SA, Ecublens, Switzerland is ISO certified.

©2007 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

AN41708_E 06/07C