

X-ray fluorescence

Analysis of glass by X-ray fluorescence with the ARL OPTIM'X WD-XRF Spectrometer

Keywords

ARL OPTIM'X, glass, WDXRF, X-ray fluorescence



ARL OPTIM'X XRF spectrometer

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°

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

Introduction

The simplest form of glass is the single component fused silica (SiO₂). 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 Fisher Scientific is used to derive limits of detection and precision for the analysis of glasses. The ARL OPTIM'X is a wavelength dispersive system that 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 200 W conventional WD-XRF instrument. The instrument can be equipped with the unique SmartGonio™, a series of monochromators or both.

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

| Oxide element | Line | Crystal | Detector | LoD (ppm) |
|--------------------------------|-------|---------|----------|-----------|
| Na ₂ O | Ka1,2 | AX06 | FPC | 100 |
| MgO | Ka1,2 | AX06 | FPC | 60 |
| Al ₂ O ₃ | Ka1,2 | PET | FPC | 47 |
| SiO ₂ | Ka1,2 | PET | FPC | N.R. |
| P ₂ O ₅ | Ka1,2 | PET | FPC | 48 |
| SO ₃ | Ka1,2 | PET | FPC | 23 |
| Cl | Ka1,2 | PET | FPC | 24 |
| K ₂ O | Ka1,2 | LiF200 | FPC | 14 |
| CaO | Ka1,2 | LiF220 | FPC | 12 |
| TiO ₂ | Ka1,2 | LiF200 | FPC | 12 |
| Cr ₂ O ₃ | Ka1,2 | LiF200 | FPC | 9 |
| MnO | Ka1,2 | LiF200 | FPC | 9 |
| Fe ₂ O ₃ | Ka1,2 | LiF200 | FPC | 9 |
| ZnO | Ka1,2 | LiF200 | FPC | 3.6 |
| SrO | Ka1,2 | LiF200 | FPC | 2.4 |
| ZrO ₂ | Ka1,2 | LiF220 | FPC | 1.8 |
| BaO | Lβ1 | LiF200 | FPC | 51 |
| PbO | Lβ1 | LiF220 | FPC | 9 |

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 are determined using the SmartGonio™ for the most common oxides/elements found in soda-lime glasses (Table 1). The recommended crystals for glass application are AX06, PET and LiF200.

Precision tests

Precision tests have been carried out by analyzing repeatedly the same pressed pellet sample for eleven consecutive analyses. Eighteen oxide/elements 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).

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.

Sample A

| RUN | Na ₂ O % | MgO % | Al ₂ O ₃ % | SO ₂ % | K ₂ O % | CaO % | Fe ₂ O ₃ % | So ₂ ppm | TiO ₂ ppm | P ₂ O ₅ % | Cl % | Cr ₂ O ₃ % | MnO % | As ₂ O ₃ % | SrO % | ZrO ₂ % | BaO ppm | PbO ppm |
|----------|---------------------|-------|----------------------------------|-------------------|--------------------|-------|----------------------------------|---------------------|----------------------|---------------------------------|-------|----------------------------------|-------|----------------------------------|-------|--------------------|---------|---------|
| Run 1 | 13.98 | 0.185 | 1.79 | 72.59 | 0.588 | 10.85 | 0.330 | 582 | 579 | 166.9 | 113.4 | 93.6 | 48.4 | 101.6 | 127.4 | 209.1 | 454.3 | 228.2 |
| Run 2 | 13.93 | 0.193 | 1.81 | 72.60 | 0.582 | 10.82 | 0.333 | 64 | 563 | 146.3 | 129.5 | 91.4 | 44.7 | 101.5 | 124.8 | 204.6 | 392.6 | 218.9 |
| Run 3 | 13.97 | 0.177 | 1.80 | 72.64 | 0.588 | 10.82 | 0.330 | 608 | 563 | 193.3 | 111.2 | 91.1 | 43.8 | 95.9 | 127.1 | 207.0 | 361.8 | 197.8 |
| Run 4 | 14.01 | 0.178 | 1.80 | 72.64 | 0.582 | 10.87 | 0.330 | 645 | 581 | 199.2 | 104.6 | 96.2 | 29.9 | 103.8 | 127.0 | 205.4 | 375.7 | 234.6 |
| Run 5 | 13.95 | 0.182 | 1.80 | 72.60 | 0.588 | 10.83 | 0.329 | 576 | 564 | 158.1 | 111.8 | 94.6 | 41.8 | 103.7 | 122.7 | 204.7 | 385.2 | 228.5 |
| Run 6 | 13.94 | 0.177 | 1.81 | 72.61 | 0.589 | 10.82 | 0.329 | 573 | 569 | 171.3 | 107.9 | 85.2 | 49.5 | 95.0 | 126.3 | 203.8 | 355.4 | 194.8 |
| Run 7 | 13.86 | 0.185 | 1.80 | 72.64 | 0.588 | 10.83 | 0.330 | 658 | 569 | 203.6 | 113.4 | 88.9 | 40.3 | 96.0 | 125.0 | 205.4 | 434.1 | 234.4 |
| Run 8 | 13.92 | 0.186 | 1.81 | 72.59 | 0.585 | 10.84 | 0.331 | 652 | 566 | 190.4 | 135.6 | 94.5 | 44.7 | 96.8 | 125.9 | 203.4 | 315.1 | 207.3 |
| Run 9 | 13.94 | 0.184 | 1.81 | 72.63 | 0.591 | 10.82 | 0.334 | 651 | 579 | 150.7 | 110.1 | 88.6 | 43.1 | 114.9 | 127.1 | 206.6 | 401.2 | 220.8 |
| Run 10 | 13.98 | 0.183 | 1.80 | 72.63 | 0.586 | 10.87 | 0.332 | 617 | 526 | 255.0 | 104.0 | 83.6 | 41.9 | 99.2 | 125.8 | 206.2 | 402.2 | 214.6 |
| Run 11 | 13.95 | 0.188 | 1.78 | 72.62 | 0.588 | 10.83 | 0.330 | 619 | 561 | 218.3 | 97.9 | 80.6 | 38.9 | 97.7 | 126.8 | 203.1 | 429.8 | 197.86 |
| Avg. | 13.95 | 0.183 | 1.80 | 72.62 | 0.587 | 10.84 | 0.331 | 620 | 565 | 186.6 | 112.7 | 89.9 | 42.4 | 100.5 | 126.0 | 205.4 | 391.6 | 216.1 |
| Std.Dev. | 0.04 | 0.005 | 0.01 | 0.02 | 0.003 | 0.02 | 0.0015 | 32 | 15 | 32.5 | 11 | 5 | 5.2 | 5.7 | 1.4 | 1.8 | 39.7 | 15 |

Table 2. Repeatability for the analysis of the major and minor oxides in Sample A at 36 seconds per oxide/element.

Sample B

| RUN | Na ₂ O % | MgO % | Al ₂ O ₃ % | SiO ₂ % | K ₂ O % | CaO % | Fe ₂ O ₃ % | SO ₂ ppm | TiO ₂ ppm | P ₂ O ₅ % | Cl % | Cr ₂ O ₃ % | MnO % | As ₂ O ₃ % | SrO % | ZrO ₂ % | BaO ppm | PbO ppm |
|----------|---------------------|-------|----------------------------------|--------------------|--------------------|-------|----------------------------------|---------------------|----------------------|---------------------------------|-------|----------------------------------|-------|----------------------------------|-------|--------------------|---------|---------|
| Run 1 | 13.35 | 0.180 | 1.67 | 73.07 | 0.556 | 10.77 | 773.3 | 0.177 | 556.0 | 200.6 | 100.7 | 63.4 | 9.3 | 118.2 | 122.4 | 227.2 | 883.9 | 895.6 |
| Run 2 | 13.33 | 0.180 | 1.68 | 73.08 | 0.564 | 10.76 | 757.9 | 0.181 | 568.0 | 159.5 | 111.2 | 65.9 | 3.9 | 112.0 | 122.2 | 226.1 | 960.6 | 914.5 |
| Run 3 | 13.28 | 0.186 | 1.67 | 73.08 | 0.554 | 10.81 | 789.6 | 0.180 | 555.1 | 193.3 | 115.7 | 64.7 | 18.2 | 115.6 | 119.3 | 225.3 | 925.4 | 911.5 |
| Run 4 | 13.28 | 0.185 | 1.66 | 73.11 | 0.559 | 10.83 | 768.2 | 0.186 | 587.2 | 156.6 | 103.5 | 74.7 | 8.7 | 105.6 | 126.5 | 225.1 | 891.3 | 900.0 |
| Run 5 | 13.35 | 0.181 | 1.67 | 73.05 | 0.554 | 10.79 | 763.9 | 0.181 | 594.7 | 187.4 | 97.9 | 63.6 | 12.3 | 106.8 | 126.3 | 226.2 | 948.8 | 904.8 |
| Run 6 | 13.32 | 0.172 | 1.67 | 73.11 | 0.566 | 10.80 | 767.3 | 0.186 | 541.4 | 183.0 | 114.6 | 59.5 | 14.7 | 115.2 | 125.5 | 226.3 | 960.5 | 904.6 |
| Run 7 | 13.33 | 0.185 | 1.67 | 73.06 | 0.554 | 10.79 | 758.9 | 0.180 | 570.3 | 193.3 | 101.8 | 67.9 | 6.1 | 113.5 | 124.9 | 226.9 | 960.6 | 910.1 |
| Run 8 | 13.26 | 0.185 | 1.69 | 73.04 | 0.555 | 10.78 | 771.7 | 0.185 | 565.2 | 191.8 | 95.2 | 66.9 | 9.3 | 101.6 | 124.7 | 227.4 | 918.0 | 919.7 |
| Run 9 | 13.33 | 0.180 | 1.64 | 73.11 | 0.561 | 10.82 | 775.7 | 0.183 | 553.7 | 219.7 | 113.4 | 61.5 | 9.3 | 109.6 | 124.6 | 222.9 | 916.9 | 913.4 |
| Run 10 | 13.30 | 0.193 | 1.68 | 73.08 | 0.556 | 10.80 | 764.1 | 0.188 | 572.6 | 243.2 | 101.3 | 64.4 | 1.0 | 103.4 | 123.0 | 226.0 | 980.8 | 912.7 |
| Run 11 | 13.31 | 0.184 | 1.66 | 73.06 | 0.561 | 10.78 | 785.3 | 0.186 | 566.0 | 191.8 | 109.0 | 70.8 | 13.8 | 106.9 | 122.1 | 227.7 | 950.9 | 875.6 |
| Avg. | 13.31 | 0.183 | 1.67 | 73.08 | 0.558 | 10.80 | 770.5 | 0.183 | 566.4 | 192.8 | 105.8 | 65.8 | 9.7 | 109.9 | 123.8 | 226.1 | 936.2 | 905.7 |
| Std.Dev. | 0.03 | 0.01 | 0.01 | 0.03 | 0.004 | 0.02 | 10 | 0.003 | 15 | 24.2 | 7.2 | 4.3 | 4.9 | 5.4 | 2.2 | 1.4 | 31.2 | 12 |

Table 3. Repeatability for the analysis of the major and minor oxides in Sample B at 36 seconds per oxide/element.