



CLAISSE
THE FIRST AND FINEST IN FUSION®

CERTIFIED REFERENCE MATERIALS READY FOR CALIBRATION

Claisse has a new turnkey solution for your XRF applications. Our fused borate disks are made with **Certified Reference Materials (CRMs)** or **customized mixtures of synthetic oxides**. They are **custom-made** to fulfill your personal analytical needs.

Fused borate disks:

- Help you build the best XRF calibration curves for your application
- Offer a great alternative to a XRF drift monitor when a more specific element needs to be monitored
- Can replace silica based drift monitors when unavailable



FIVE REASONS FOR USING CLAISSÉ FUSED BORATE DISKS

1. HIGH-QUALITY DISK PREPARATION

Claisse produces borate disks that **exceed the industry's quality standards**. The **high accuracy** and **precision** of our fused borate disks can help you build your XRF application calibration curves and obtain accurate results for your samples.

2. TIME SAVING

Claisse will save you a significant amount of time by **finding the appropriate CRMs** for your calibration needs and **preparing fused disks for you**.

3. QUICK APPLICATION START-UP

With your fused disks already **prepared for the calibration curve**, you can integrate the new application rapidly in your day-to-day routine and start analyzing your samples right away.

4. LOW COST

The cost of a fused disk is **less than 50% of the cost of CRMs** when bought as a powder.

5. DIVERSITY

Claisse can produce **almost any mix of oxides**, thus extending the calibration possibilities. For **synthetic standards**, we can perform the tedious task of **preparing perfectly homogeneous mixes to cover the calibration range needed** for the elements.

Did you know that Claisse produces many fused disks that are used as drift monitor every year?

HERE ARE A FEW EXAMPLES OF OUR AVAILABLE PRODUCTS

- PbZn calibration (25 disks) with synthetic oxide
- NIST srm 1880 to 1889
- JCA cement standard kit (JCA 601a XRF1 to XRF15)
- Cement and raw material calibration curves (31 disks)
- Iron ore exploration (28 disks)

Scan this code to obtain more information about our fused borate disks made with CRMs



CLAISSÉ
THE FIRST AND FINEST IN FUSION®

claisse.com

Claisse Headquarters
350 Franquet St., suite 45
Québec City, Quebec
G1P 4P3 CANADA

Phone: +1 418. 656. 6453
Fax: +1 418. 656. 1169

