



**CLAISSE**  
THE FIRST AND FINEST IN FUSION®

## TAKE A WEIGHT OFF YOUR MIND!

**Have you ever thought about the amount  
of time you spend weighing flux?**

Since we have at heart your entire satisfaction and are committed to ensure efficiency in your laboratory, we now offer our same high-quality fluxes in a wide range of weights. The best fluxes on the market can now be pre-weighed to make your life easier and accelerate the sample preparation step!

**Start saving time and money today!**



# SEVEN REASONS FOR USING CLAISSÉ® PRE-WEIGHED FLUXES

## [ 1 ] SAVE TIME AND MONEY BY AVOIDING THE WEIGHING STEP

Skipping the weighing step enhances efficiency by allowing you to perform other valuable tasks in the laboratory.

## [ 2 ] HIGH PRECISION AND OUTSTANDING REPEATABILITY

Our state-of-the-art instruments automatically weigh the flux in each vial, thus avoiding variation in weight and generating accurate results.

## [ 3 ] SEALED VIALS TO AVOID CONTAMINATION

The product is therefore not altered by air, humidity or other contaminants that could be found in the laboratory.

## [ 4 ] COMPLETE TRANSFER INTO THE CRUCIBLE

Our vials are designed to prevent any waste. There is always the right amount of flux in the crucible for greater analytical reproducibility.

## [ 5 ] ANY TYPE OR QUANTITY OF CHEMICALS CAN BE PRE-WEIGHED

At Claisse, fulfilling your needs is our priority. That's why we are committed to pre-weigh any of our flux in any quantity for your convenience.

Make your life easier and try the pre-weighed fluxes offered by Claisse!

Contact your Claisse sales representative for more details.

## [ 6 ] EASY AND SAFE TO USE

The shape of our vials avoids spillage and makes them handy for everyday use.

## [ 7 ] QUICK DELIVERY

Tired of long delivery times? Our fluxes are already in stock and ready to be shipped promptly so you can maintain productivity in your laboratory.



Scan this code to obtain more information about the pre-weighed fluxes



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

