



CLAISSSE
THE FIRST AND FINEST IN FUSION®

KEEPING AHEAD THROUGH CLAISSSE EXPERTISE IN SAMPLE PREPARATION BY FUSION

Scale up your analytical precision



LE DOSER™

HOW TO REACH EFFICIENCY WITH CLAISSE EXPERTISE?



LE **DOSER**

Claisse offers a global solution in sample preparation by fusion to improve efficiency in the laboratory. Our knowledge and experience combined with the reputation of PANalytical allow us to constantly innovate to fulfill our customers' needs as well as to help them obtain accurate and precise analytical results.



Consumables



Services



Expertise





CLAISSE

THE FIRST AND FINEST IN FUSION®

LeDoser™ instrument is more than a weighing device: it is an essential instrument to achieve constant, on target results and efficient laboratory management. It is specifically designed to weigh and dispense borate flux with high precision for the preparation of glass disks for XRF and solutions for ICP. This instrument minimizes repetitive strain injuries (RSI) arising from weighing related tasks, hence improving the operator's safety. Three different weighing modes and smart flux flow software control makes it fully adaptable to your needs, without compromise. LeDoser's industrial grade balance incorporates the most cutting-edge technology weighing system with unrivalled robustness. LeDoser instrument is synonymous with peace of mind, quick return on investment and outstanding analytical performance.





Scale up your analytical precision



WHY INVEST IN LeDOSER AUTOMATIC WEIGHER

EFFICIENT LABORATORY MANAGEMENT

- Automated and reproducible weighing
- Leads to a more efficient use of the operator's time during the weighing process
- Eliminates calculation errors and fatigue for increased analytical accuracy and reproducibility
- Promotes a safer work environment, decreasing the risk of operator injuries caused by repetitive tasks

PEACE OF MIND

- 3 different weighing modes: Sample-to-flux ratio – Catch weight – Absolute weight
- Seamless integration in laboratory Standard Operation Procedure (SOP)
- Tank stirrer option to ensure constant performance for high and low throughput
- LIMS ready – simple and straightforward data transfer

HIGH PERFORMANCE

- Cutting-edge technology weighing system for high productivity
- High precision (up to 0.0001 g)
- Smart flux flow software control for constant yield and maximized weighing speed
- Runs autonomously for up to 250 weighings*

* According to the requested quantity of flux per fusion

EASY TO USE

- Intuitive touchscreen interface
- One-touch operation
- No training required for operation
- No calculations required to determine flux weight in relation to sample quantity

QUICK RETURN ON INVESTMENT (ROI)

Low cost of ownership

- Save up to 80% weighing time
- Simple and efficient mechanism made of durable parts
- Robust industrial grade balance for unrivalled robustness
- Protective equipment prevents infiltration of flux in the balance mechanism
- Low maintenance

Minimal infrastructure required

- Small and compact (benchtop instrument)
- Easy installation



TECHNICAL SPECIFICATIONS



PRODUCTIVITY

- Weighs and dispenses flux for XRF, AA and ICP analysis
- Runs autonomously for up to 250 weighings*
- Maximized weighing speed depends on the requested tolerance, flux quantity and flux type**

* According to the requested quantity of flux per fusion

** Weighing speed will also be influenced by the environmental condition (installation)

ELECTRICAL

- Voltage: 100-240 VAC
- Current: 1 A
- Frequency: 50-60 Hz
- Power: 100 VA
- Phase: 1 phase

DIMENSIONS

- Height: 50.5 cm (20 in.)
- Depth: 50.5 cm (20 in.)
- Width: 30.5 (12 in.)

WEIGHT (WITH BALANCE)

- 19 kg (42 lb.)

CONTROL AND OPERATION

- One-touch operation
- Touchscreen user interface
- Detailed weighing parameters and results display
- Automated sliding doors
- Can be used as a balance
- Bar code reader can be plugged directly in the instrument
- Crucible detector
- Low volume detector located in the flux tank

PROGRAMMABLE WEIGHING PARAMETERS

- Sample-to-flux ratio
- Number of dosing steps
- Target weight and tolerance for each dosing step
- Calculation method of the tolerance for each dosing step

SOFTWARE AND COMMUNICATION

- 2 USB ports
- Limitless weighing program storage (4G Compact Flash)
- Smart flux flow software control
- Dosing results transferable through Ethernet connection
- Management tool for different flux composition
- Balance calibration function available from the interface

SAFETY

- User operation levels are protected by password
- Certified CE and CSA
- Conformal coated PCB for high corrosion resistance
- Meets UL 94 flammability standard
- Meets RoHS requirements



Scan this QR code to obtain more information about LeDoser instrument.



CLAISSSE
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

