



CLAISSSE
THE FIRST AND FINEST IN FUSION®

PLATINUM AND LABWARE

Platinum that meets the highest
quality standards



HOW TO REACH EFFICIENCY WITH CLAISSE EXPERTISE?



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.



Instruments



Services



Expertise





CLAISSE

THE FIRST AND FINEST IN FUSION®

Claisse is the one-stop shop for any special needs in platinum or labware for sample preparation by fusion!

Claisse offers high quality molds and crucibles made of 95% platinum and 5% gold. Zirconium stabilized molds are also available for special needs in sample preparation by fusion. All of our molds and crucibles are sold in a wide range of sizes and thicknesses to help you obtain the best analytical results. The heavy molds are highly recommended since they are well known for their durability and because they can be polished several times.

We provide our customers with other precious metal accessories used in laboratories such as dishes, electrodes, tongs and tweezers.

Our experts know that peroxide fusion requires labware made of different materials, that's why nickel or zirconium crucibles are available at Claisse.

**Platinum that meets the highest
quality standards**



WHY INVEST IN OUR PLATINUMWARE?

- High quality
- Available from stock for immediate delivery
- Wide choice of platinumware for sample preparation by fusion
- The 95% platinum and 5% gold accessories prevent sticking
- Can be shipped everywhere in the world
- Precious metal scrap policy

POLISHING YOUR PLATINUMWARE

It is essential to keep the surface of your platinum molds and crucibles flat and smooth to maintain a high quality of analysis. We offer a specialized



polishing service to help you quickly recover the mirror aspect of your platinum accessories. We sell a polishing kit, mold reshapers and a training program for customers who would like to polish their labware on their own.

PRECIOUS METAL SCRAP POLICY

Did you know that Claisse can buy your precious metal accessories at a competitive market rate? A credit is given as a partial payment on any of our products or on new platinumware.



CATALOG

MOLDS

95% Pt – 5% Au

M03030-B-0A	30 mm Claisse mold, 0.8 mm thick, 21 g
M03130-B-0A	30 mm Claisse heavy mold, 1 mm thick, 27 g
M03032-B-0A	32 mm Claisse mold, 0.8 mm thick, 23 g
M03132-B-0A	32 mm Claisse heavy mold, 1.2 mm thick, 36 g * Available from stock
M03001-0A	32 mm Claisse shallow mold, 1 mm thick, 31 g
M03009-B-0A	32 mm Claisse Eagon 2 mold, 0.8 mm thick, 40 g
M03010-B-0A	32 mm Claisse heavy Eagon 2 mold, 1.2 mm thick, 60 g
M03035-B-0A	35 mm Claisse mold, 0.8 mm thick, 26 g
M03135-B-0A	35 mm Claisse heavy mold, 1.2 mm thick, 40 g * Available from stock
M03003-0A	35 mm Claisse shallow mold, 1 mm thick, 36 g
M03038-B-0A	38 mm Claisse mold, 0.8 mm thick, 34 g
M03040-B-0A	40 mm Claisse mold, 0.8 mm thick, 35 g
M03140-B-0A	40 mm Claisse heavy mold, 1,1 mm thick, 48 g * Available from stock
M03002-0A	40 mm Claisse shallow mold, 1 mm thick, 44 g
M03011-B-0A	40 mm Claisse Eagon 2 mold, 1 mm thick, 58 g
M03012-B-0A	40 mm Claisse heavy Eagon 2 mold, 1.2 mm thick, 69 g

95% Pt – 5% Au - DPHzr (Zr Grain Stabilized)

M03018-B-0A	32 mm Claisse Pt-DPH Eagon 2 mold, 0.8 mm thick, 40 g
M03019-B-0A	32 mm Claisse heavy Pt-DPH Eagon 2 mold, 1.2 mm thick, 60 g
M03020-B-0A	40 mm Claisse Pt-DPH Eagon 2 mold, 1 mm thick, 58 g
M03021-B-0A	40 mm Claisse heavy Pt-DPH Eagon 2 mold, 1.2 mm thick, 69 g

All weights are nominal.

Claisse offers a mold polishing service to ensure a better durability of your platinumware and to maintain the high quality of your analysis.



Scan this QR code to obtain more information about the platinum and labware offered by Claisse.

CRUCIBLES AND LIDS

95% Pt – 5% Au

CR3133-B-0A	Claisse Optimix™ crucible (with three humps), 30 g * Available from stock
CR3325-B-0A	Claisse Bis!™ crucible, 26 g
CR3330-B-0A	Claisse heavy Bis! crucible, 30 g * Available from stock
CR3120-B-0A	Claisse flat bottom crucible, 20 mL, 24 g
CR3125-B-0A	Claisse flat bottom crucible, 25 mL, 26 g
CR3130-B-0A	Claisse flat bottom crucible, 30 g * Available from stock
CR3010-B-0A	Claisse Eagon 2 flat bottom crucible, 40 g
CR3011-B-0A	Claisse Eagon 2 crucible with dimple, 40 g

95% Pt – 5% Au - DPHzr (Zr Grain Stabilized)

CR3008-B-0A	Claisse Pt-DPH Eagon 2 flat bottom crucible, 40 g
CR3009-B-0A	Claisse Pt-DPH Eagon 2 crucible with dimple, 40 g

100% Zr

CR1900-0B	Zirconium crucible, 25 mL
AC3101-0A	Zirconium lid for crucible Cr1900
P-0430-02	Zirconium crucible, 50 mL
P-0431-02	Zirconium lid for crucible P-0431-02

100% Ni

CR1902-0A	Nickel crucible, 25 mL
AC3102-0A	Nickel lid for crucible Cr1902
P-0450-02	Nickel crucible, 50 mL
P-0451-02	Nickel lid for crucible P-0450-02

All weights are nominal.

ACCESSORIES

PRK-RS010	Bis!™ crucible reshaper
PRK-RS012	Flat bottom crucible reshaper
PRK-RS002	30 mm mold reshaper
PRK-RS004	32 mm mold reshaper
PRK-RS005	35 mm mold reshaper
PRK-RS007	38 mm mold reshaper
PRK-RS008	40 mm mold reshaper
PRK-PK001	Claisse mold polishing kit (120 V)
PRK-PK002	Claisse mold polishing kit (240 V)
P-0600-20	Polishing recharge 1
P-0600-30	Polishing recharge 2
PRK-VM001	Claisse VortexMixer™ (110 V)
PRK-VM002	Claisse VortexMixer (220 V)



CLAISSE
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

