

Claisse<sup>®</sup> is a Malvern Panalytical brand

# EAGON 2<sup>®</sup> FUSION INSTRUMENT

Keeping ahead through expertise in sample preparation by fusion



# SAFE, SIMPLE, HIGH PERFORMANCE

The Eagon 2 is a fully automatic fusion instrument that prepares glass disks for XRF analysis. Its innovative patented design ensures performance, operator safety and ease of use. The concept of the Eagon 2 instrument makes fusion and the consequent benefits of accurate XRF analysis easily reachable.



#### The Eagon 2's value to sample preparation by fusion

- Ultimate safety
- Optional exhaust adapter for minimum infrastructure requirements
- Optimized method development



The Eagon 2 is a great quality control tool leading to very high analytical performance and allowing

The versatility of the Eagon 2 is convenient when it comes to analyzing cosmetics samples.

With this simple and low maintenance instrument, you can quickly switch from producing glass beads for XRF to producing solutions for ICP analysis. It then facilitates your experiments.

The safety door of the instrument prevents spills and allows a clean and safe preparation of food

The Eagon 2 is a great quality control tool leading to very high analytical performance and allowing

The Eagon 2 is a versatile fusion instrument that is easy to use and that requires low maintenance.

In addition to providing excellent reproducibility, the Eagon 2 instrument is very helpful in analyzing

The Eagon 2 instrument is useful for the preparation of soils and sediments since it leads to high-

### WHAT IS FUSION?

Fusion is a sample preparation method developed in the mid 50s. It consists of dissolving at high temperatures a fully oxidized sample in a suitable solvent (a flux) in a platinum, zirconium or graphite crucible. The melted mixture is agitated and poured into a mold to create a glass disk for XRF analysis. It can be poured into a beaker to create a solution for AA or ICP analysis.



\*\*\*The Eagon 2 instrument can't be used to make solutions for AA or ICP analysis.

#### Why should I use fusion in my laboratory?

This universal technique has numerous benefits when you compare it with other sample preparation methods such as pressed pellets or acid digestion.

	Fusion	Pressed pellets
Affected by mineralogy	No	Yes
Affected by particle size	No	Yes
Desirable size of powder (microns)	50-100 (easy)	5-30 (difficult)
Accuracy	≤1%	≤10%
Easy calibration with synthetic standards	Yes	No
Application of matrix correction	Yes	No

# WHY INVEST IN THE EAGON 2 **FUSION INSTRUMENT**?

#### High analytical performance

- Reproducible
- Fully automatic operation that guarantees a
- perfect repetition of fusion cycles
- Superior quality ceramic holders for crucibles and
  - molds to ensure the lowest contamination
  - · Optimized fusion conditions for all materials.

#### Programmable fusion parameters

- Temperature setting
- Duration
- Oxidation steps
- · Non-wetting agent injection
- Agitation · Cooling.

### QUICK RETURN ON INVESTMENT (ROI)

#### Low cost of ownership

- Casting dish sensors
- Ensure safe and reliable operation no possible damage of the instrument related to pouring without the platinum mold
- · Crucible and mold holders can easily be dismounted for cleaning
- Platinumware friendly, no damaging temperature gradient.



### **Optimized method development**

 Pause and inspection function to visualize the fusion process during the fusion cycle.

#### Ultimate safety

- Cold-to-cold operation
- The operator is protected against hot material and surfaces by interlocked doors during the entire fusion cycle.

#### Easy to use

- One-touch operation
- Library of predefined fusion methods
- · Definition of customizable fusion method.

#### Minimal infrastructure required

- Simple electrical connection
- No O<sub>2</sub>, compressed air or water cooling system needed
- The optional exhaust adapter limits the infrastructural requirements.



#### Tailored solutions through expertise

Malvern Panalytical experts are the reference for all steps of the fusion process or development of new applications. Their personalised solutions allow customers to properly manage risks, reach a quick ROI, and benefit from a large database of preparation methods for a broad range of samples.

# TECHNICAL SPECIFICATIONS

#### Productivity

Fully automatic fusion instrument that prepares glass disks for XRF ar
 Two fusion positions with independent synchronous or asynchronous

#### Heating

- Temperature operation range: 300 1200°C
- Temperature monitored by a type R thermocouple inside the heating Resistance-based heating system

#### Electrical

- Voltage: 220-230 V
  Current: 32 A
  Frequency: 50-60 Hz Electrical

#### Dimensions

XINV

Alex.

Dimensions • Width: 64.5 cm (25.4 in.) • Depth: 68.5 cm (27.0 in.) • Height: 56.5 cm (22.2 in.) • Weight: 78 kg (172 lb.)

#### Programmable Fusion Parameters

- Ramp-dwell steps (temperature /time)
- Full temperature range solid phase oxidation
- Crucible angle and agitation speed
- Non-wetting agent injection
  Crucible pouring angle
  Passive or forced cooling (2 speeds)

#### **Control and Operation**

- One-touch operation
- Precise temperature control (± 5°C)
- Full microprocessor control
  Pause and inspection function to visualize the fusion process during Alarm when the cycle is completed

#### Software and Communication

- Front panel operation
- Optional alpha /numeric programming via a PC for additional conver Up to 32 user-defined methods

#### Safety

- Certified CE and machinery directive 98/37/EC compliant
- Outer doors are interlocked during fusion cycle to enhance safety
- Maximum external temperature of 70°C
   Casting dish sensors ensure safe and reliable operation
   User operation levels are protected by password

nalysis <b>us</b> operation
chamber
g the fusion cycle
ience



# WHY CHOOSE MALVERN PANALYTICAL?

We are global leaders in materials characterization, creating superior, customerfocused solutions and services which supply tangible economic impact through chemical, physical and structural analysis.

Our aim is to help you develop better quality products and get them to market faster. Our solutions support excellence in research, and help maximize productivity and process efficiency.

Malvern Panalytical is part of Spectris, the productivity-enhancing instrumentation and controls company.

www.spectris.com

### SERVICE & SUPPORT

Malvern Panalytical provides the global training, service and support you need to continuously drive your analytical processes at the highest level. We help you increase the return on your investment with us, and ensure that as your laboratory and analytical needs grow, we are there to support you.

Our worldwide team of specialists adds value to your business processes by ensuring applications expertise, rapid response and maximum instrument uptime.

- · Local and remote support
- · Full and flexible range of support agreements
- · Compliance and validation support
- Onsite or classroom-based training courses
- e-Learning training courses and web seminars
- Sample and application consultancy



### MALVERN PANALYTICAL

Grovewood Road, Malvern, Worcestershire, WR14 1XZ, United Kingdom

Tel. +44 1684 892456 Fax. +44 1684 892789 Lelyweg 1, 7602 EA Almelo, The Netherlands

Tel. +31 546 534 444 Fax. +31 546 534 598

info@malvernpanalytical.com www.malvernpanalytical.com

www.malvernpanalytical.com/claisse