



**Malvern
Panalytical**



**SAVE ADDITIONAL
€265.000/YEAR
FOR REFINING CATALYST
PROTECTION***

www.panalytical.com/Xflow

EPSILON XFLOW

Save additional €265.000/year by using Epsilon Xflow, a low-maintenance and highly reproducible sulfur, iron, nickel and vanadium cat-killer analyzer compared to any other analyzer.

- Extend catalyst life time by simultaneous on-line analysis of S, Fe, Ni and V: **savings € 100.000/year**
- Reduce conflicts between lab and on-line results (ASTM D2622, ISO 13032) saving precious time for product release, reducing temporarily higher give-away and costly re-analysis: **savings €100.000/year.**
- Quickly switch between process streams: **savings € 50.000/year**
- Reduce instrument- and calibration maintenance: **savings €15.000/year and 5 days/year**

www.panalytical.com/Xflow

MORE INFO:

EMEA

juergen.wess@panalytical.com

APAC

xue.shilei@panalytical.com

AMEC

daniel.raymond@panalytical.com

**Monetary savings are based on 400.000 bbl/day refining capacity and typical operating costs. No warranties or guarantees are given.*



EASY ON-LINE PRODUCTION CONTROL OF LIQUID PROCESS STREAMS



Customized integration of the Epsilon Xflow into your production process guarantees results right from the start. Malvern Panalytical's specialists offer a complete solution for integration, including required sample conditioning systems, probes, fast loop pumps and more.

The Epsilon Xflow on-line liquid analyzer, is specifically designed for continuous elemental composition analysis of any liquid. The on-line solution provides real-time elemental insight in your processes, thereby leading to more efficient plant operations and reduced costs.

- Multi-element analyzer with low detection limits
- Monitor and control your processes via Plant Control System
- Chemical resistance for a wide range of liquids
- Designed according to ATEX requirements
- Low maintenance
- Excellent long-term repeatability

WHY CHOOSE MALVERN PANALYTICAL?

Malvern Panalytical is a global leader in materials characterization, creating expert solutions for the challenges associated with maximizing productivity, developing better quality products and getting them to market faster. We provide superior, customer-focused solutions and services which deliver tangible economic impact through chemical, physical and structural analysis of materials.

www.malvernpanalytical.com

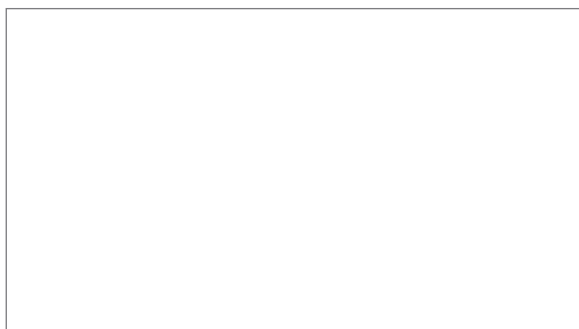
Malvern Panalytical is part of Spectris, the productivity-enhancing instruments and controls company.
www.spectris.com

SUPPORT AND SERVICE

Malvern Panalytical helps you increase your laboratory's productivity and the return on your investment with our global training, service and support network.

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
- Validation support
- Onsite or classroom-based training courses
- e-Learning courses and web seminars
- Sample and application consultancy



MALVERN PANALYTICAL

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

Tel. +44 1684 892456
Fax. +44 1684 892789

Lelyweg 1, 7602 EA Almelo,
Netherlands
P.O.Box 13, 7600 AA Almelo,
Netherlands

Tel. +31 546 534 444
Fax. +31 54 534 598

info@panalytical.com
www.malvernpanalytical.com

Disclaimer: Although diligent care has been used to ensure that the information in this material is accurate, nothing herein can be construed to imply any representation or warranty as to the accuracy, correctness or completeness of this information and we shall not be liable for errors contained herein or for damages in connection with the use of this material. Malvern Panalytical reserves the right to change the content in this material at any time without notice. Copyright © Malvern Panalytical 2017. This publication or any portion thereof may not be copied or transmitted without our express written permission. 9498-707-65411-PN11368