

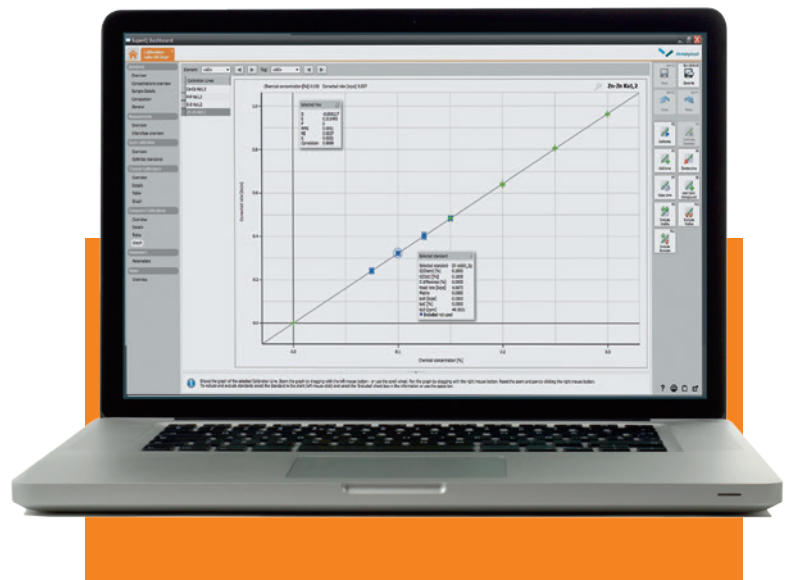


PANalytical
get insight



OIL-TRACE

Accurate elemental analysis for
fuels, oils and wear metals





Accurate results, easy operation

Savings on setup time, standards and maintenance

One calibration for many different oils, fuels and petrochemicals

| Your challenges | Our solution |
|---|---|
| Accurate analyses using multiple calibrations for a wide range of petrochemical products | Single calibration enabled by: - advanced FP-Compton matrix corrections - multi-scatter corrections |
| Maintenance of multiple calibrations | Number of calibrations dramatically reduced |
| Reproducible sample preparation due to infinite thickness problems when dealing with X-rays | Largely independent of sample mass/volume. Advanced Fluorescent Volume Geometry modeling for infinite thickness corrections |
| Getting best accuracy, regardless of varying oxygenates and hydrocarbon composition | Automatic oxygen and hydrocarbon composition corrections |
| International norm compliance | Application templates to simplify setup and analysis according to international norms and test methods |
| Sourcing standards of varying composition | One generic set of oil standards |



Superior accuracy

Single calibration - many applications

Traditionally, separate calibrations must be made for quantifying important elements in different types of fuels, oils and other petrochemicals. A single Oil-Trace calibration removes the time-consuming maintenance of many calibrations.

ASTM

D2622, D6334, D4927, D6443, D6376, D5059, D4294, D6481

ISO

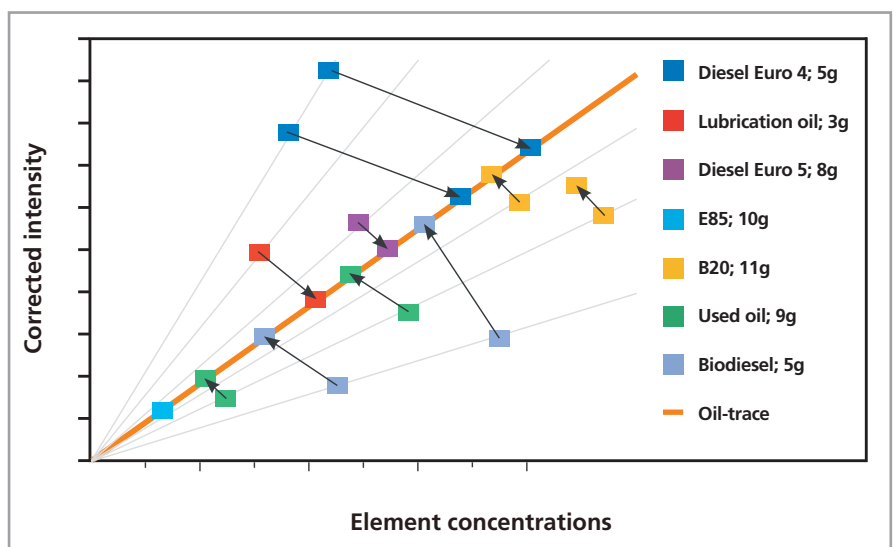
14596, 14597, 20884, 20847, 8754, 13032

IP

336, 496

JIS K2541-4

Differences in oxygen content and sample weights/volumes are compensated by Oil-Trace



Calibrate with a single fuel, oil or petrochemical standards set of your choice and Oil-Trace allows for accurate analysis of any fuel, oil or other petrochemical.

Completing the solution

Integrated in SuperQ and Epsilon 3 software

Oil-Trace is closely integrated in the Zetium - Petro edition SuperQ and Epsilon 3 software environments giving access to the many useful features present within these mature packages:

- Routine analysis and sample handling
- Calibration update and instrument monitoring
- Statistical process control (SPC)
- Data storage and reporting
- Data transmission

Oil-Trace comprises everything required for successful oil analysis:

- Oil-Trace software
- Liquid standards for C, H and O characterization
- Compact weighing balance
- Volumetric flask
- Comprehensive user setup documentation (CD-ROM)



Oil-Trace:
accurate elemental
analysis using fewer
calibrations

Tailored system solutions

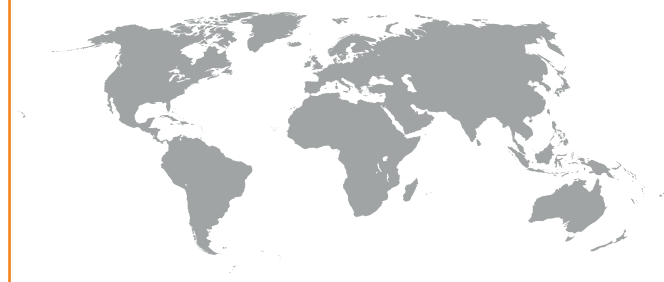
Epsilon 3^X and Epsilon 3^{XLE} spectrometers –
Compliant and compact



Petro edition of Zetium spectrometer -
Compliant and complete



Global and near



PANalytical B.V.
Lelyweg 1, 7602 EA Almelo
P.O. Box 13, 7600 AA Almelo
The Netherlands

T +31 546 534 444
F +31 546 534 598

info@panalytical.com
www.panalytical.com

Regional sales offices

Americas
T +1 508 647 1100
F +1 508 647 1115

Europe, Middle East, Africa
T +31 546 834 444
F +31 546 834 969

Asia Pacific
T +65 6741 2868
F +65 6741 2166