

Epsilon 1 Wells for faster well interventions



Contact us for more
information

Production chemistry data within minutes

On-asset solution for Production Chemistry and Flow Assurance challenges.

The Epsilon 1 wells is an on-asset monitoring solution that gives you critical production chemistry and flow assurance data within minutes. It maximizes uptime and production by giving you the ability to make data driven decisions about production parameters for scale-squeezes and other interruptions in real time.



Fast on-site
measurements



Easy to
operate and
maintain



Real-time
data



No sample
alterations

With the Epsilon 1 Wells you can predict with confidence when a scale squeeze needs to take place, saving millions of dollars. Shell has been successfully using this technology for several years.

www.malvernpanalytical.com/epsilon1



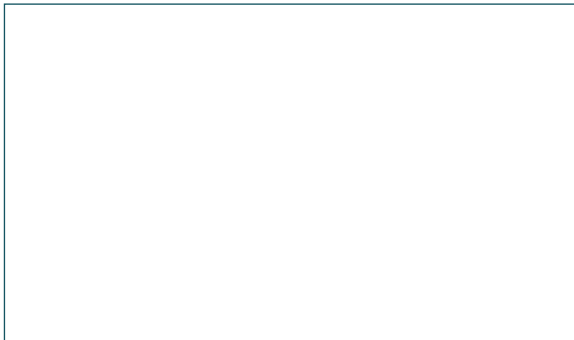
**Malvern
Panalytical**
a spectris company

Applications (all sample types except gas)

- Water composition: produced water, injection water, flowback water, aqueous treatment chemicals, scale formation, proppants, corrosion inhibitors.
- Scale composition: barium, calcium, strontium, sulfur, phosphorus.
- Oil composition: sulfur, nickel, vanadium and other contaminants.
- Sand/solid/metal composition, emulsion/sludge estimation.
- Production contaminants monitoring (e.g. mercury and others), condensate impurities.

Solution capabilities

- Fast on-site measurements (lab quality data in 20 mins).
- Better insight by minimal sample alterations due to long transportation times.
- Offshore optimized Benchtop XRF analyzer – DNV-GL certification pending.
- Easy to operate and maintain – minimal consumables (10mL sampling cup).



Malvern Panalytical

Groveswood 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