

FRACTURING SYSTEM

Nuclear density meter for fracturing and cementing applications

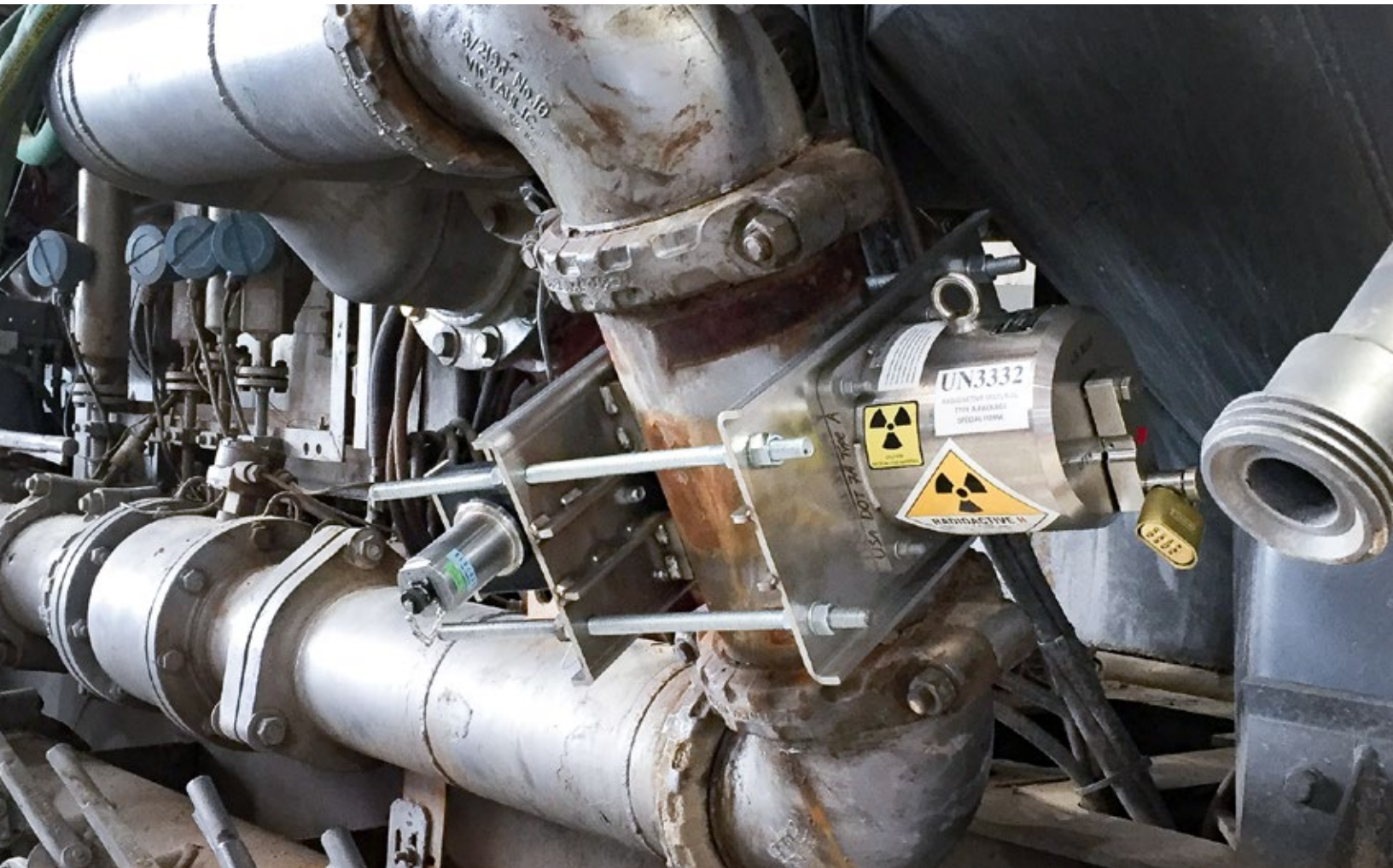


 **BERTHOLD**

DENSITY METER FOR FRACTURING APPLICATIONS

Berthold offers radiometric density systems that are specifically designed for use in hydraulic fracturing and cementing applications. The systems are installed on blender trucks, high pressure lines, and cementing lines; providing an online measurement of PPA, PPG, density, specific gravity, or percent solids. In addition to their high repeatability, our fracturing systems stand out with their ruggedness in the face of mechanical shock and vibration, and for the low source activities required.

Berthold has supplied nuclear density meters to the fracturing industry for more than 10 years. During that time, we have learned that the fracturing environment is tougher than any other industry – we have learned how to address these special needs. Therefore, we offer a complete line of specialized products that exactly meet our customers’ demands in the fracturing industry – no matter if a new system or a retrofitting solution is required. With more than 1,000 density systems installed on fracturing applications, we have proven our capability to provide the products, support, and service that this industry deserves.



The easy-to-use measurement system

Berthold’s fracturing system stands out with its excellent performance including high accuracy and repeatability. The transmitter DuoXpert LB 475 easily connects to the FracSENS LB 6770 detector and links directly to the customer control system. No calculations must be done in the control system and the transmitter converts the measured count rate into the output values specified by the customer for low pressure or high-pressure applications.

DuoXpert LB 475 Transmitter

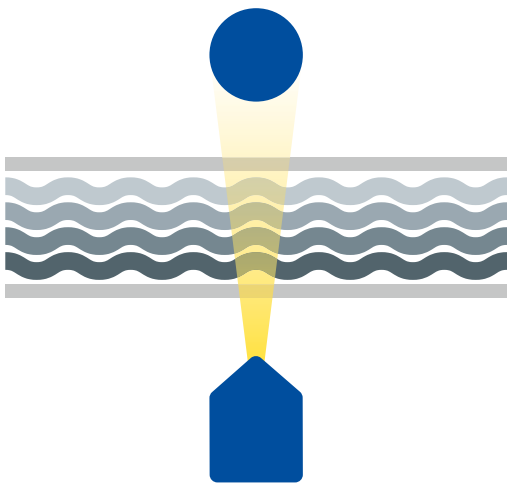
Operators benefit from the new simplified user-friendly interface where calibration is as simple as possible. Only two-points – a low and a high density, e.g. air and water – have to be acquired by the customer. Advanced self-diagnostics and monitoring features ensure safe function of the system. Operators also benefit from the new service functionalities like data logging, which can be accessed either remotely through ethernet connection or through the USB port at the front panel of the system.

Technology

- Backlit 3.5” color touchscreen
- Dual Core DSP/ARM controller
- Ethernet connection for remote access
- USB port for USB storage media or mouse
- Optional in-field user-interface for FracSENS LB 6770

Features

- Easy to use with advanced touch HMI interface
- Specially designed for hydraulic fracturing and cementing applications
- Measures density, outputting PPA, PPG, percent solids, specific gravity, or density
- Simplified calibration procedure
- Advanced self-diagnostics and monitoring features



Clamp-on density measurement on blender trucks, high pressure lines, and cementing lines



THE EXPERTS IN MEASUREMENT TECHNOLOGY

Berthold Technologies stands for excellent know-how, high quality and superior reliability. The customer is always the focus of our solution. We know our business!

Using our varied product portfolio, our enormous specialized knowledge and extensive experience, we develop suitable solutions together with our customers for new, individual measurement tasks in a wide variety of industries and applications. Berthold Technologies is specialized in radiometric process measurements for 70 years. This is our core competence with state-of-the-art and innovative products and solutions covering a vast range of industries and applications.

We are here for you – worldwide!

The engineers and service technicians from Berthold Technologies are wherever you need them. Our global network assures you fast and above all competent and skilled assistance in case of need. No matter where you are, our highly qualified experts are willing and ready to assist you in a timely manner with the best possible solution for you, even in the most challenging applications..

Berthold Technologies U.S.A., LLC

99 Midway Lane · Oak Ridge · TN 37830
+1 (865) 483 1488 · berthold-us@berthold.com · www.berthold.com/us

