

LB 350

ONLINE MOISTURE ANALYSER

Radiometric moisture analyser
for representative and robust
measurements



 **BERTHOLD**

ONLINE MOISTURE MEASUREMENT ON BULK MATERIAL

The LB 350 is a moisture measurement system for real-time determination of the moisture content of a wide variety of different bulk materials such as:

- Coke
- Sinter mixture
- Ore
- Sand
- Building material

and is used on many different types of vessels.

Neutron moisture measurement is very robust and a proven technology with a long track record of success for moisture measurement on bulk materials; it functions where most other moisture measurement technologies fail for various reasons. The large measuring volume of up to one meter in diameter ensures a representative measured value. Once the system has been installed and calibrated, it can be left alone for several years providing reliable and maintenance-free moisture monitoring of bulk products.

The LB 350 can additionally be equipped with bulk density compensation. Bulk density compensation increases the accuracy of the moisture measurement if the bulk density of the product fluctuates.



Moisture measurement on coke
with surface probe

LB 350

Online moisture analyser

Possible applications

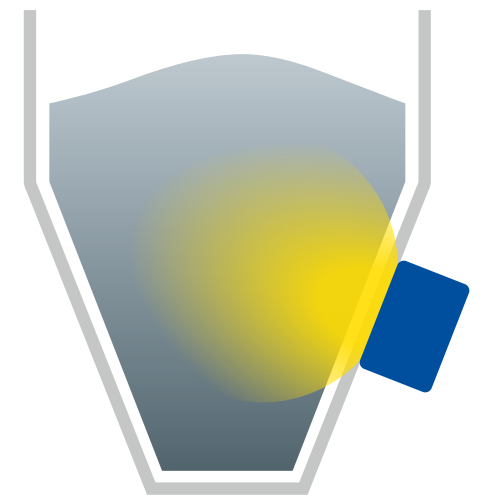
- Bunkers
- Silos
- Weighing tanks
- Feeding tanks

Customer benefits

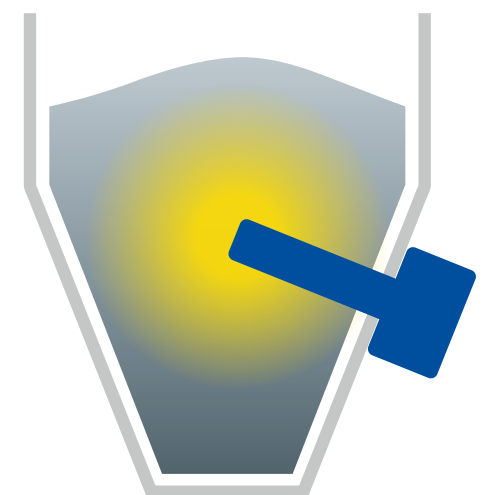
- Representative measured values due to large measuring volume
- Accurate moisture measurement
- Not effected by e.g. temperature, conductivity, pressure, pH-value etc.
- Different sensor probes for measuring in or through the vessel.
- Can be supplied with density compensation

Technical Properties

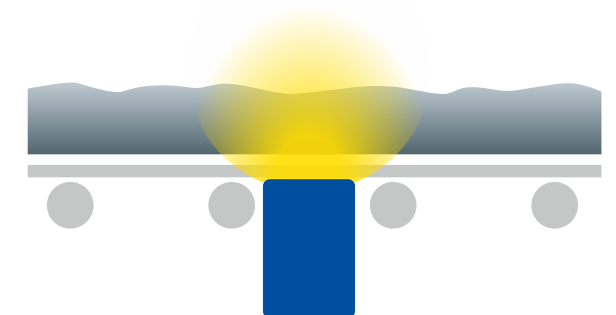
- Proven neutron moisture measurement technology
- Attainable accuracy ± 0.1 % depending on application
- Analog output from sensor (0/4...20 mA)
- Digital input for "hold signal"
- Digital outputs for system status and min and max values
- Sensor operating temperature: -20...+50 °C
- Density compensation enabled



Moisture measurement with surface probe



Moisture measurement with bunker probe



Moisture measurement with surface probe on conveyor



THE EXPERTS IN MEASUREMENT TECHNOLOGY

Berthold Technologies stands for excellent know-how, high quality and 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 specialised in radiometric process measurements for 70 years. This is our core competence with state-of-the-art and cutting edge 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 and specialists are ready and waiting and will be with you in no time at all with the ideal solution for even the most difficult measurement task.

Berthold Technologies GmbH & Co. KG

Calmbacher Straße 22 · 75323 Bad Wildbad · Germany
+49 7081 1770 · industry@berthold.com · www.berthold.com

