



# Dry Substance & Water Content

## Process monitoring in dairy production

### Early recognition of fluctuations

Water is a characteristic component of dairy products. It determines consistency and influences the taste as well as the quality of the final product. In most countries, the water content is subjected to severe restrictions. In Germany, for example, the water content of butter is limited to 16%.

Up to now, any deviations from the set-point could only be identified at the end of the production process. Using state-of-the-art microwave technology, it is now possible to monitor the water content/ dry substance online and during production, which provides an early recognition of fluctuations. Thus, a basis for constant high quality and cost-efficient production has been established.

## Micro-Polar LB 566

From butter to yoghurt – the measuring system Micro-Polar LB 566 provides continuous and real-time measuring data of the current water content/dry substance in a wide range of dairy products. The measurement can therefore be carried out directly on the main stream; a by-pass is not necessary.

The microwave based measurement is completely non intrusive, which means that there are no built-in components in the inner pipe. No additional time and effort is required for cleaning. The measuring performance is neither influenced by colour, viscosity, or non-homogeneity of the measured product. Due to our latest developments in the sensor technology, the Micro-Polar is a world leader for accuracy and reliability. Moreover, the instrument is extremely user-friendly arising from our customers' co-operation in the food industry.



### Online measurement for increased quality

Online-monitoring of water content or dry substance during the production process enables optimisation of the process. Deviations from the set-point can be corrected promptly, leading to a significant increase in efficiency. Furthermore, production costs can be considerably reduced. Likewise, it is ensured, that the dairy products are compliant with all statutory regulations and that the consistent high product quality meets the customer's expectations.

