

Stable  
& Fast

The Thermo Scientific Prima BT and Prima PRO mass spectrometers are the go-to-technology for online gas analysis, delivering:

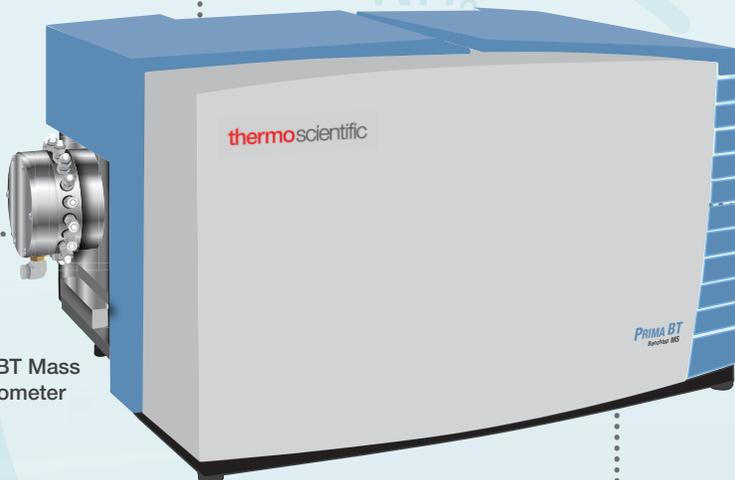
- Automatic calibration
- Low calibration frequency
- Simple-to-use software
- 99.7% up-time!

Precise

**Process optimisation**

for both research and development and manufacturing with solutions for the laboratory and production.

Prima BT Mass Spectrometer

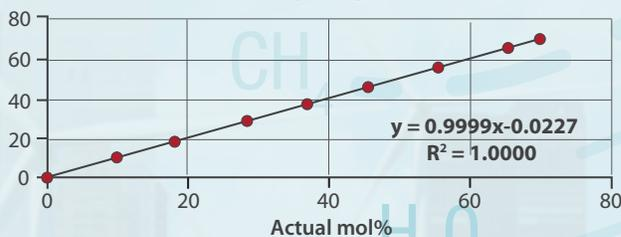


Unrivalled precision, stability and speed when measuring hydrogen when compared to other mass spectrometers and GC.

Accurate

Maximise return-on-investment by measuring up to 64 individual gas streams on one mass spectrometer.

**Hydrogen**



Dynamic measuring range from ppm to 100% of **hydrogen, oxygen, carbon dioxide** and more.

Reliable

# Real-time hydrogen measurement on the Prima Mass Spectrometer

Learn more at [thermofisher.com/PrimaBT](https://thermofisher.com/PrimaBT)

thermo scientific