

Raman process analysis

## Five things you didn't know about the Ramina Process Analyzer and the MarqMetrix BallProbe

### 1. If you can touch a sample, you can take a measurement

The Thermo Scientific™ Ramina™ Process Analyzer is designed to natively use a variety of MarqMetrix® probes. TouchRaman® technology allows you to simply touch the end of your BallProbe® to the sample and take a Raman measurement with your Ramina system. The BallProbe's short focal length can accurately analyze liquids, solids, slurries, powders, and heterogeneous mixtures, while eliminating user to user variation.



Bioreactor BallProbe with TouchRaman immersion technology

### 2. BallProbe is available in multiple shapes and sizes

The BallProbe comes in various sizes for different applications. Whether you need a probe that can withstand harsh environments, moderate chemical exposure, or demanding bioprocessing settings, we offer a range of BallProbes that will work with your application.



Process BallProbe

### 3. BallProbe is built to last

Ultra-resilient, BallProbe is designed to withstand temperatures from -20c to 150c and pressures up to 6,000 psi. No matter what you want to measure, BallProbe can take the heat and pressure.

### 4. BallProbe is compatible with multiple Raman systems

While BallProbe is designed for use with the Ramina Process Analyzer, it is also compatible with a number of other Raman systems from other vendors. All Ramina probes are incredibly easy to swap, giving our customers additional flexibility. BallProbe self-aligns, eliminating annoying downtime associated with probe swaps with other vendors.

### 5. You can customize your BallProbe

We know you have unique applications that require powerful solutions. In partnership with MarqMetrix, BallProbes can be manufactured to custom specs and from the most suitable materials for your application upon request.



FlowCell with BallProbe technology

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



**thermo** scientific