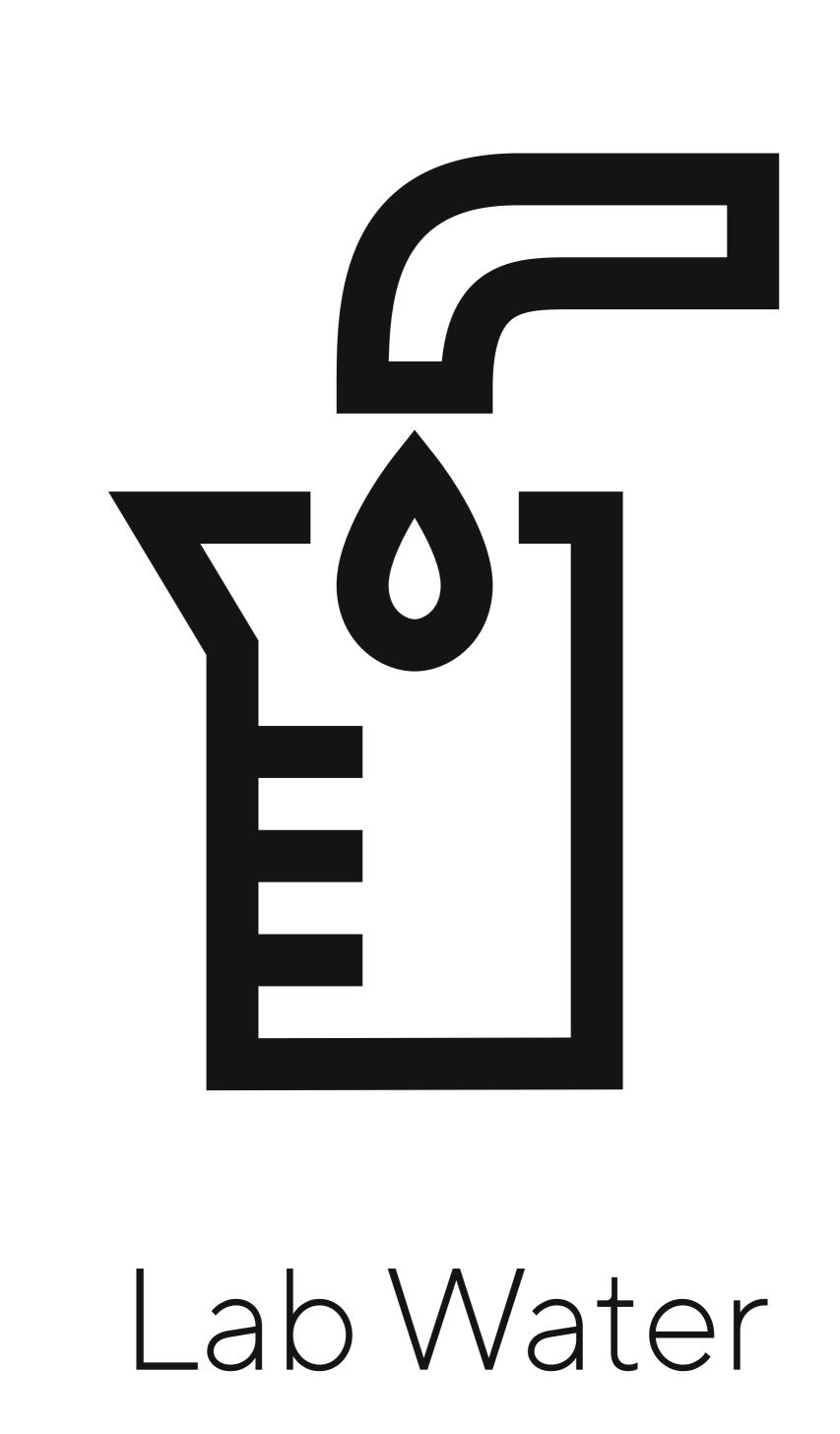
S/RICH SILS

Products and Solutions for the Laboratory

Lab Balances | Pipettes & Tips | Lab Water Systems Filtration & Microbiology Consumables | Service

Simplifying Progress







Moisture Analyzer

