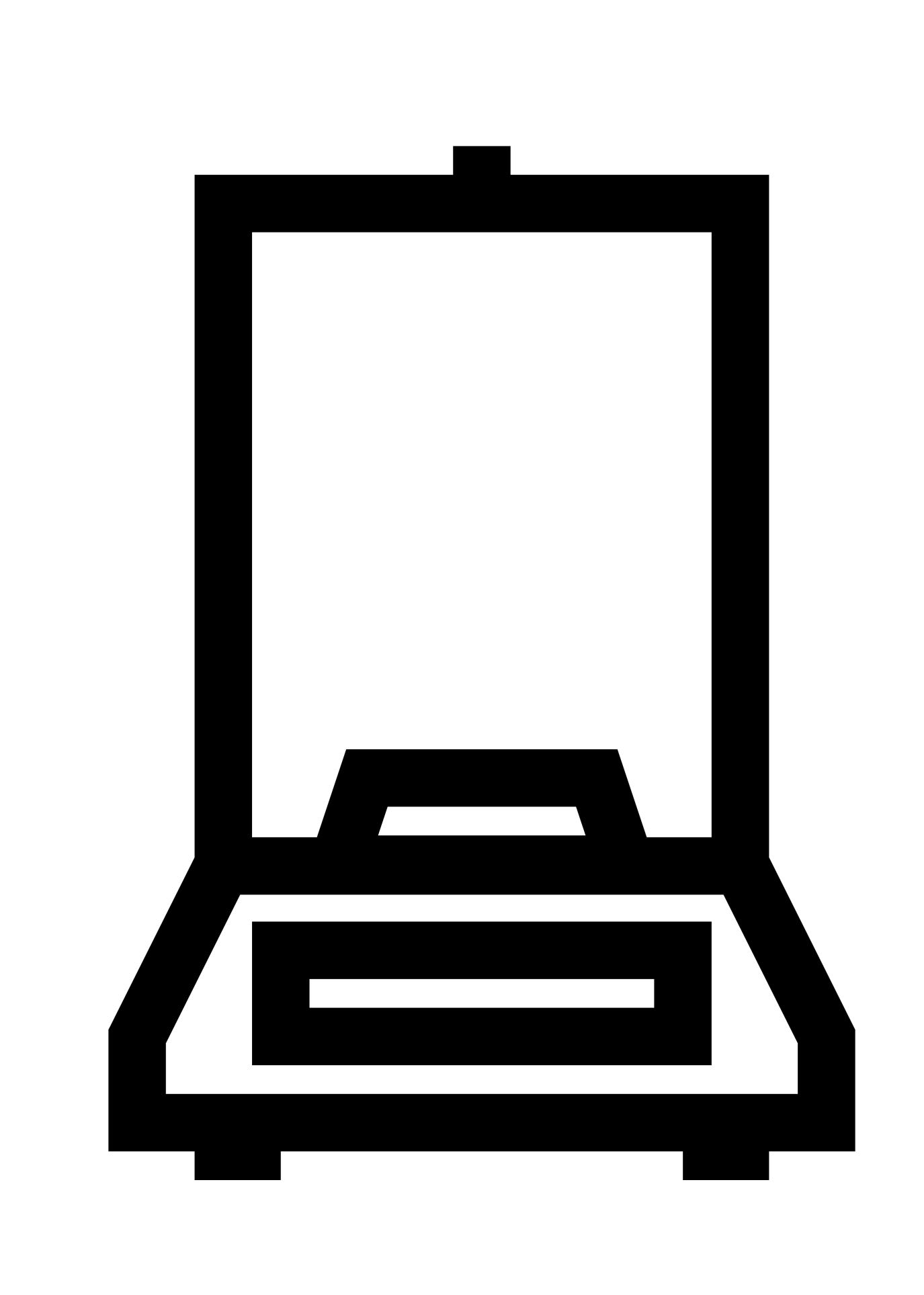
SARTORIUS

Simplifying Progress

Products and Solutions for the Laboratory

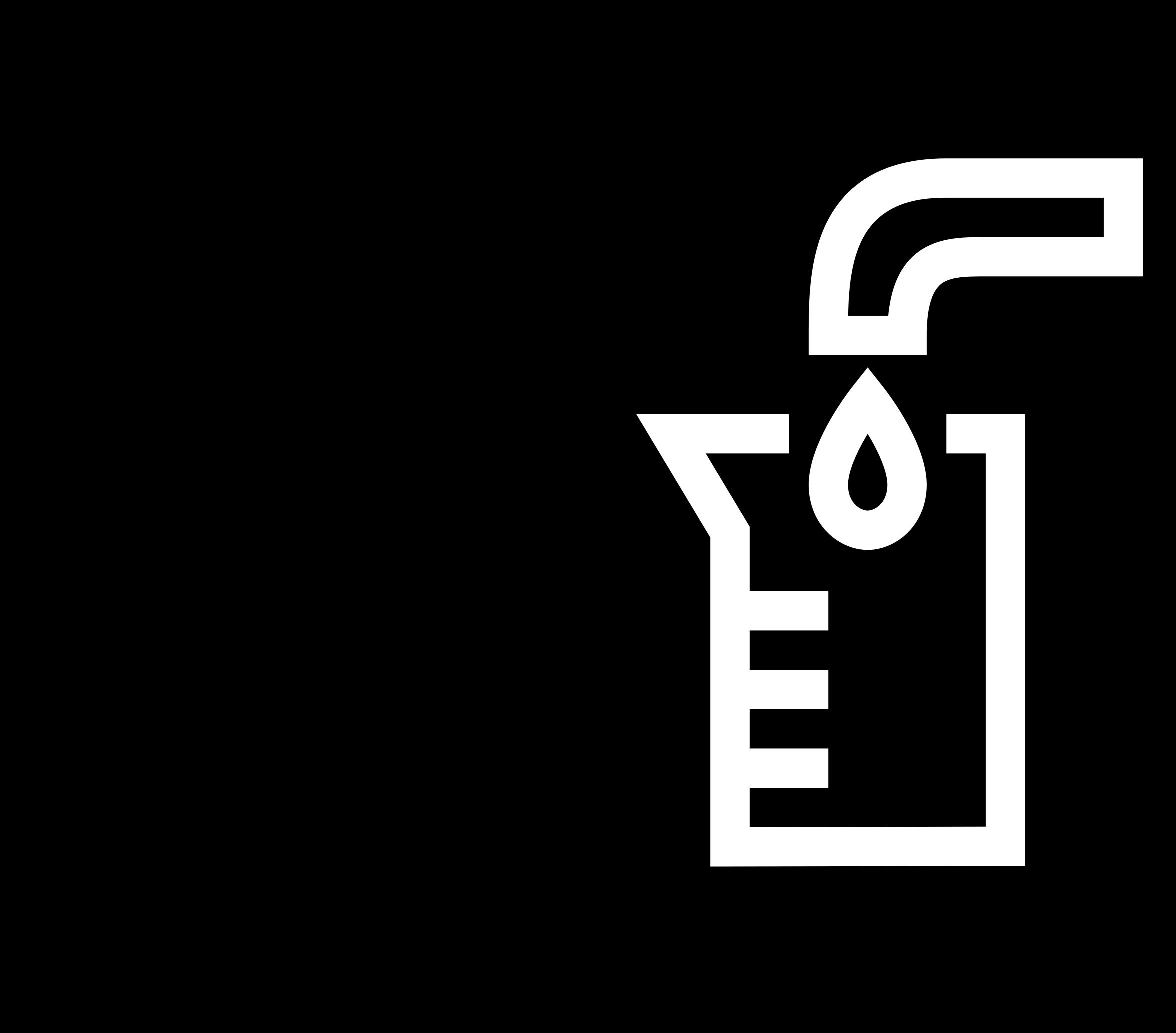
Lab Balances | Pipettes & Tips | Lab Water Systems Filtration & Microbiology Consumables | Service



Lab Weighing



Liquid Handling



Lab Water

