



SciAps X-200 Specifications

Totally re-engineered

- Weighs only 3.1 lbs. with battery
- Redesigned heat sinking can operate continuously up to 130 °F (54 °C) ambient temperature
- New internal circuit boards
- New housing and upgraded metal components
- Up-to-date software, user interface, and calibration

SciAps X-200 XRF analyzer is the perfect blend of high performance and attractive price. It's a top-performing SDD with highly optimized X-ray tube and detector geometry, plus improved heat dissipation, reduced weight, and improved durability. Operates on Android OS using Wi-Fi and software with the ease of a smart phone for revolutionary speed, connectivity, and test data management.



High performance and value

The X-200 Series are the lowest priced XRF guns on the market, yet still offer best-in-class analytical and speed performance, with all the accuracy and throughput of SciAps flagship X-550 and X-505.

X-ray tube

40 kV Rh anode for alloys
50 kV Au anode for
geochem, soil, RoHS



Alloys

Fast on every alloy family, SciAps X-200 is a great choice for scrap processing, non-destructive testing and manufacturing quality control.



Geochemistry/Soils

Industry-comparable precision, LODs and elemental range for environmental, pathfinder, exploration and mining. Analyzers maybe factory calibrated with fundamental parameters, Compton Normalization (EPA Method 6200), or user-defined derivative calibrations.



RoHS

Fast, precise analysis for Pb, Hg, Cd, Br and Cr, plus other restricted materials and Halogen Free requirements. Operates at three beam conditions for optimized LODs across the element suite. Automated sample type recognition (polymer, alloy, or mixed type).

For more information , or to
schedule a demonstration:

SciAps Inc.
+1 339.927.9455

SciAps

All internal electronics now use less power and operate
at higher temperatures - no delays due to drift or shutdown.



SciAps X-200

Specifications

Weight	3.1 lbs. with battery
Dimensions	9.38in (238mm) x 11.15in (283mm) x 3.34in (84mm)
Power	On-board rechargeable Li-ion battery, rechargeable inside device or with external charger, and AC power.
Display	3.5" color touch screen, smartphone-type display – PowerVR SGX540 3D graphic
Sample Viewing	Internal camera for viewing sample before and during analysis for proper sample alignment. Second macro-camera for scanning QR or barcodes and for photo-documentation and report generation.
Comms/Data Transfer	Wi-Fi, Bluetooth, USB. Connectivity to most devices, including SciAps Profile Builder PC software.
Excitation Source	5 W X-ray tube. 50 kV max, 200 uA max, Rh or Au anode depending on application.
Detector	20mm ² silicon drift detector (active area), <140eV resolution FWHM at 5.95Mn Kalpha line.
X-Ray Filtering	6-position filter wheel for beam optimization
Processor	ARM Cortex -A9 dual-core / 1.2GHz Memory: 1GB DDR2 RAM, 1GB NAND
Pulse Processor	14-bit ADC with digitization rate of 80 MSPS 8K channel MCA USB 2.0 for high speed data transfer to host processor; digital filtering implemented in FPGA for high throughput pulse processing 50nS – 24uS peaking time
Calibration	Fundamental Parameters, Compton Normalization, and/or Compton Normalization for Alloy. Geochem, Environmental and Soil calibrations.
Calibration Check	External 316 stainless check standard for calibration verification and energy scale validation
Environmental Temp. Range	10°F to 130°F at 25% duty cycle.
Security	Password protected usage (user level) and internal settings (admin)
Regulatory	CE, RoHS, USFDA registered, Canada RED Act

SciAps Inc.
sales@sciaps.com
SciAps.com
+1 339.927.9455

 [YouTube.com/SciAps](https://www.youtube.com/SciAps)

SciAps