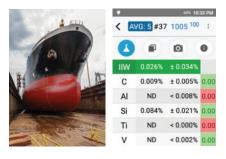


Carbon Analysis Any Place on the Planet



- Third-party inspection companies
- Pipeline inspection companies
- Chemical plants
- Refineries
- Manufacturing
- Shipbuilders
- Steel makers
- Scrap processing

SciAps Z-902 Carbon Specifications

The world's first handheld analyzer to deliver carbon content just got better. SciAps Z-902 Carbon is the key to rock-solid, in-field analysis for alloys. Confirm alloy type and grade 300 series stainless into L, H, and S series. Determine carbon equivalents for welding. Perform material verification pre- and post-fabrication.





A new standard for NDT/PMI

Identification of stainless steels and low alloy steels, including carbon analysis down to 70 ppm and instant CE (carbon equivalence) calculations using our uniquely powerful laser, onboard argon purge and high-resolution spectrometer.

- On-board argon delivers results for pennies per test.
- Powerful laser delivers 5-6 mJ/pulse on the sample, with a 50 Hz repetition rate. High energy and high repetition rate mean a clean burn.
- Laser raster can rapid-sample multiple locations in one test for fuller data.
- Analysis averaging available for multi-shot averaging and data rejection.
- Sample alignment micro-camera and LED Spotlight focus just where you want.

190 nm - 420 nm spectrometer

Includes a dedicated spectrometer for superior resolution in the 190 - 230 nm range

For more information, or to schedule a demonstration:

SciAps Inc. +1 339.927.9455



True portability

Only 4.35 lbs (1.97 kg) with battery and a narrow profile for accessing the tightest spaces; an aerospace-grade aluminum body for improved performance in high ambient temperatures; and a re-designed, Android-based user-interface all make this system the most usable platform on the market. Wifi and Bluetooth with GPS capability to print, email, and connect to virtually any information management system for efficient, real- time test data and reporting.



Weight

Security

SciAps Z-902 Carbon Specifications

4.35 lbs. (1.97 kg) with battery



SciAps Cloud Services

You use XRF for metals in alloys, because it's the best technology for most PMI. But what about carbon? Two reports? Not with SciAps Cloud. Get efficient real-time results, data merge and management, and full-featured report generation for XRF and LIBS.

10.75 x 2.875 x 8.625 inches **Dimensions** 2.7" high-brightness, color touchscreen, readable in all lighting Display conditions. Rear-facing display for easy results viewing. On-board rechargeable Li-ion battery, rechargeable inside device or with Power external charger, AC power. ARM Quad Cortex -A53 1.2 GHz Memory: 2 GB LPDDR3, 16 GB eMMC **Processing Electronics Data Storage** Results storage: 32 GB SD Built on Google's Android platform for real-time data exporting, including built-in WiFi (IEEE 802.11b/g/n), Bluetooth (BR/EDR+BLE), GPS Connectivity and USB-C to connect to virtually any information management system Integrated camera and laser target indicator for viewing sample before and during analysis for proper sample alignment. Includes second "macro Sample Viewing camera" for scanning QR or barcodes and for photo-documentation and report generation. On-board Y stage for rastering laser to discrete locations for targeted Laser Raster analysis or averaging. SciAps proprietary Opti-Purge provides an inert argon environment, improving spectral signal-to-noise ratio and improving performance in Atmosphere the UV range. 316 stainless steel standard for automated calibration and wavelength Calibration Check scale validation. On-board automated drift correction software with factory-provided or **Drift Correction** user-provided reference materials. Grade Library 500+ grades, multi-library support, libraries may be added **Grade Library** or edited. CE, RoHS, USFDA registered. Class 3b laser. Sample sensor on board, allows for operation under Class 1 conditions, subject to local LSO Regulatory approval. 190 - 420 nm **Spectral Range** Aluminum: Be, Mg, Al, Si, Ti, V, Cr, Mn, Fe, Ni, Cu, Zn, Zr, Pb, Bi, Ag, Sn Titanium Base: Al, Ti, V, Cr, Fe, Cu, Zr, Nb, Mo, Sn LAS Base: C, Al, Si, Ti, V, Cr, Mn, Fe, Cu, Ni, Nb, Mo, Pb Stainless Steels: C, Al, Si, Ti, V, Cr, Mn, Ni, Fe, Ni, Cu, Nb, Mo, W Calibrations Nickel Base: Al, Si, Ti, Cr, Mn, Fe, Co, Ni, Cu, Nb, Mo, W

Copper Base: Be, Al, Si, Cr, Mn, Fe, Ni, Cu, Zn, Ag, Sn, Pb, Bi

Specialty Bases: Mg, V, Cr, Mn, Co, Zn, Zr, Nb, Mo, Ag, Sn, Hf, Ta, W, Re,

ARM Cortex -A9 dual-core / 1.2GHz Memory: 1GB DDR2 RAM, 1GB NAND

Cobalt Base: Al, Si, Ti, Cr, Mn, Fe, Co, Ni, Cu, Nb, Mo, W

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

