



SciAps Z-901 CSi Specifications

A dedicated handheld LIBS for carbon and silicon measurements

- The perfect complement to handheld XRF
- Measures C + Si in steels, L-grade & H-grade stainless
- Ultra compact at 3.87 lbs., handles like XRF

X-ray analysis is great on transition and heavy metals, making it the preferred tool for most PMI operators. But additional data is required for users of carbon steels and L- and H-grade stainless. Z-901 CSi provides dedicated information on carbon and silicon, making it fast and economical to meet critical performance standards with a high- performance handheld LIBS analyzer.



Now widely accepted for carbon testing:

SciAps breakthrough LIBS analyzer is included in API 578 2nd Edition, accepted at every major refinery and in use worldwide by fabricators, pipeline integrity management programs, power plants and other users of carbon steels and L- and H-grade stainless.

Add carbon to complement XRF

An alternative to spark OES, handheld LIBS analyzers use a miniature laser to measure what X-ray can't: Li, Be, B, C, F, Na and more. For industries that focus on carbon content in steels and stainless, the CSi offers a dedicated solution. Features include advanced spectrometer design for high resolution and wide range, 190-260 nm, and microanalysis capability with 100 um laser spot size. Meets multi-test averaging protocols.

Fully featured analyzer

CSi has an internal camera for precise targeting of analysis locations, especially welds; a macro camera for photo-documentation of samples, reading bar codes and QR codes; a patented "sample sensor" that allows Class 1 operation; intuitive Android operating system; high-resolution rear-facing display; rugged metal body; narrow snout for welds or difficult-to-access test locations; and user-replaceable argon that provides hundreds of tests at pennies per test.

Easy reports

XRF users who add CSi to their PMI toolkit can efficiently merge tests using SciAps cloud data management and report generation, with results displayed for specific parameters like carbon equivalents and residuals. Using another company's XRF? No problem. Merge our LIBS data with any XRF to get the job done.



SciAps X-550 and Z-901 CSi are the best combination for widest variety of PMI and highest throughput. Get them both in OneBox.

For more information, or to schedule a demonstration:

SciAps Inc.
+1 339.927.9455

SciAps



**A dedicated
handheld LIBS for
carbon and silicon
measurements**

SciAps Z-901 CSI Specifications

Weight	3.87 lbs. (1.76 kg) with battery
Dimensions	10.75" x 2.38" x 8.63" (273 x 60 x 219 mm)
Display	2.7" high brightness, color touchscreen, readable in all lighting conditions. Rear facing display for easy results viewing.
Power	On-board rechargeable Li-ion battery, rechargeable inside device or with external charger, AC power.
Processing Electronics	1.2GHz ARM Cortex-A53 quad-core. Memory: 2GB LPDDR3, 16GB eMMC
Data Storage	Results Storage: 32 GB SD card
Connectivity	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 and USB-C to connect to virtually any information management system.
Sample Viewing	Integrated camera and laser target indicator for viewing sample before and during analysis for proper sample alignment. Includes second "macro camera" for scanning QR or barcodes and for photo-documentation and report generation.
Laser Raster	On-board Y stage for rastering laser to discrete locations for targeted analysis or averaging. Typically, raster in 3 locations.
Atmosphere	SciAps proprietary Opti-Purge provides an inert argon environment improving spectral signal-to-noise ratio and improving performance in the UV range.
Calibration Check	316 stainless steel standard for automated calibration and wavelength scale validation.
Drift Correction	On-board automated drift correction software with factory-provided or user-provided reference materials.
Regulatory	CE, RoHS, USFDA registered. Class 3b laser. Sample sensor on board, allows for operation under Class 1 conditions, subject to local LSO approval.
Spectral Range	188 - 260 nm
Calibrations	Low Alloy Steel: C (0.008 to 1.25%), Si Stainless Steel: C (0.0075 to 0.25%)
Security	Password protected; Multi-user support with configurable access settings