



SciAps Z-9 Liquidator Specifications

- **Results in seconds**
- **No dilutions**
- **Field-portable**
- **Simple! No chemist required**
- **Simultaneously measures Li plus B, Na, Mg, K, Ca and more**

Li Brines Edition

SciAps is proud to announce the first portable LIBS powered Lithium Brines analyzer, the Z-9 Liquidator.

Combined with the Z-903 handheld LIBS, the Z-9 Liquidator provides fast and easy analysis of brines in the field or in the lab. This groundbreaking technique only needs a 1-2 mL of sample per test, with no dilution necessary. The Z-9 Liquidator samples the liquid and atomizes it into a fine mist which is then analyzed by the Z-903 LIBS. Within seconds, the concentration of Li and other key elements are displayed on nearby tablet or PC.

World's First In-Field Lithium Brines Analyzer

Fast

No more waiting days, weeks or even months for results from complicated lab techniques!

Easy

Simply feed your sample to the Z-9 Liquidator and your brines composition is reported in seconds!

Accurate

The Z-9 Liquidator is powered by the SciAps Z-903 handheld LIBS analyzer, equipped with factory built lithium brines calibration. Users can also develop a custom calibration via the Profile Builder software to further optimize performance.

Versatile

With a small footprint and light but durable construction - the SciAps Z-9 Liquidator is right at home in the lab, in the plant or even in the field! Transportation is easy for analyzing brines at the well or evaporation ponds.

LithiumBrines		
Test #370 2023-02-24 16:51:42		
	Chemistry	Uncertainty
Li		
Lithium	0.0798%	± 0.0064
Na		
Sodium	0.1533%	± 0.0069
K		
Potassium	0.0282%	± 0.0070
Ca		
Calcium	0.0307%	± 0.0027

Figure 1: Profile Builder screenshot showing lithium brines results.



LIBS
TECHNOLOGY

Have solids or rocks to analyze?

Just undock the Z-903 handheld analyzer from the Z-9 Liquidator and you're free to roam! Test outcrops, cores, pressed pellets and more!

Calibrations available for pegmatite, Nevada clays, etc.



SciAps Z-9 Liquidator Specifications

Weight	Z-9 Liquidator 15lbs + Z-903 LIBS with battery 4.93lbs
Dimensions	9.25" x 11.00" x 16.5"
Sample Volume	~ 1mL per test
Analysis Time	5s per test
Power	100-240V, 50/60Hz or optional battery
Z-903 Spectral Range	190 - 950 nm
Drift Correction	On-board automated drift correction software with factory-provided or user-provided reference materials
Regulatory	CE, RoHS, USFDA registered, Safety interlocked Class 3B laser
Calibrations	Factory built lithium brines calibration or user built calibrations via Profile Builder desktop software
Operating Software	SciAps Profile Builder for PC or Windows Tablet



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

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

SciAps