



SciAps reveNIR Specifications

Why choose the SciAps reveNIR?

- Rapid decision-making with real-time data
- Lightweight and compact for extreme field conditions
- Detects minerals critical for ore vectoring and exploration
- A reliable, indispensable tool for modern geotechnical teams

The SciAps reveNIR (Near-infrared spectroscopy) handheld analyzer is a cutting-edge tool designed for mineral exploration and reconnaissance. By combining advanced vis-NIR technology with portability and efficiency, the reveNIR delivers real-time mineralogical analysis in remote and challenging environments. Whether you're analyzing geological formations or identifying pathfinder minerals, this device provides unparalleled performance and reliability.

Filling the Gap in the Market

The discontinuation of the ASD Halo left a significant gap in the market—one that existing solutions couldn't fill. SciAps responded by developing the reveNIR, a compact, powerful handheld analyzer designed to meet this critical need. Built with advanced technology, the reveNIR offers precision and portability for accurate field analysis, redefining standards for NIR spectrometry in mineral exploration.

Power Packed Details

Instrument Type:

Handheld vis-NIR Spectrometer

Primary Application:

Mineral identification and reconnaissance

Target Users:

Geologists, exploration teams, and mining professionals

Portability:

Lightweight, compact design suitable for remote locations

The reveNIR's lightweight construction and robust design ensure optimal functionality in rugged terrains. At under 4.5 lbs (with battery), it's 2 lbs lighter and 50% smaller than the ASD Halo. Its ergonomic design makes it ideal for use in remote locations, whether accessed by helicopter or small plane.

Reflectance Spectra Analysis

The reveNIR records reflectance spectra to predict the mineral composition of samples, making it a critical tool for understanding geological formations and identifying pathfinder minerals.

- Spectral range - 350 to 2510 nm.
- A basic mineral spectral library of 50 important exploration minerals is included to get started immediately.
- Capability for user spectral additions and subtractions and converting publicly available USGS SpecLib 7a downloads from the internet.
- Real time Spectral Scalars calculations of Al-OH, ISM (Illite Spectral Maturity), Mg-OH, and Fe3t which indicate crystallinity, composition and/or geothermal conditions that serve as additional vectors to potential mineralization.

Transform Exploration with the SciAps reveNIR

The SciAps reveNIR is redefining material identification for mineral exploration professionals. With its advanced capabilities, customizable features, and field-ready design, it's the ideal tool for modern geological analysis. Contact us today to learn how the reveNIR can enhance your exploration efforts.



NEW!
reveNIR

For more information, or to
schedule a demonstration:

SciAps Inc.

+1 339.927.9455


SciAps



SciAps reveNIR Specifications

**The Essential Tool for
Mineral Exploration**

Weight	4.5lbs with battery
Dimensions	10.74in x 10.64in x 3.15in
Excitation Source	Halogen Bulb
Detector	VIS1: CCD array (>2000 pixels) NIR1: InGaAs array (256 pixels) NIR2: Extended InGaAs (256 pixels)
Available Apps	Mining
Environmental Temperature Range	5°C to 40°C
Analytical Range	350nm to 2500nm
Processing Electronics and Host Processing	Powerful quad Arm® Cortex®-A53 processor with a Neural Processing Unit (NPU) operating at up to 2.3 TOPS
Power	On-board rechargeable Li-ion battery, rechargeable inside device or with external charger, AC power, hot-swap capability (60 s max swap time)
Display	3.5" color touchscreen Smartphone type display - PowerVR SGX540 3D graphic
Comms/Data Transfer	USB-C, Bluetooth, Wi-Fi, GPS
Calibration	Built-in mineral library
Calibration Check	Unit is wavelength calibrated. During regular operation, a white external reference calibration is needed every 30-60 minutes when the unit is warmed up. The unit's calibration is automatically checked against its internal references to correct fast drift
Security	Password protected usage (user level) and internal settings (admin)
Dual Cameras	Two 5-megapixel cameras
Regulatory	CE, RoHS

 [YouTube.com/SciAps](https://www.youtube.com/SciAps)
sales@sciaps.com
SciAps.com
 +1 339.927.9455



**THE LIGHTEST & FASTEST
HANDHELD VIS-NIR
ON THE MARKET**