



## QualitySpec® TREK

### Making field spectroscopy easy

Finally! A truly portable full-range near-infrared (NIR) contact spectrometer that is fast, accurate, and easy to use. From the makers of the FieldSpec®, LabSpec®, and TerraSpec® product lines comes QualitySpec® Trek, redefining true instrument field-portability while maintaining the highest standards of spectral data quality you expect from ASD Inc., a PANalytical company.

QualitySpec Trek delivers precision full-range spectral measurements through a hand-held system designed around a radically streamlined cable-free workflow. With one pull of the trigger this revolutionary hand-held full-range spectrometer delivers accurate spectral results in seconds. By optimizing spectral collection workflow, portability, and ease of use, QualitySpec Trek saves time in the field and maximizes your research budget investment. There simply is no easier way to collect accurate full-range spectra.

QualitySpec Trek is a general purpose, all-in-one, full-range portable spectrometer measuring the visible and shortwave infrared regions (350-2500 nm).



#### Contact NIR spectroscopy

Spectral researchers and scientists, regardless of application, often rely on non-destructive contact measurement techniques to acquire spectral signatures. In most cases these measurements are obtained with an instrument configuration involving an optical accessory with an internal light source that is tethered to a portable spectrometer through a fiber optic cable and power cable. In turn, the spectrometer is typically controlled through a portable laptop or PDA.

By eliminating the need for cables and controller accessories, QualitySpec Trek offers a turnkey solution that simplifies your spectral collection work flow and is well-suited for a variety of applications including:

- Spectral profiling of soil horizons
- Spectral characteristics of man-made materials
- Plant physiology studies
- Spectral characterization of snow/ice profiles
- Characterization of industrial raw and finished materials

## QualitySpec Trek features and benefits

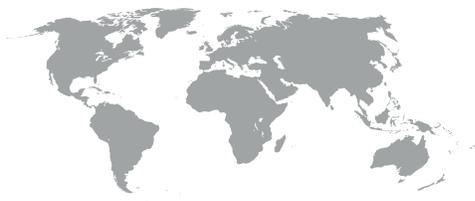
- **Portability:** Lightweight, rugged design allows unprecedented portability in a full-range highly accurate NIR spectrometer.
- **Fast and accurate:** Enhanced spectrometer configuration builds on the trusted legacy of ASD for unparalleled accuracy and performance.
- **Cable free:** Remove cable management from your data collection workflow with a truly cable-free integrated system.
- **Integrated computer and LCD:** On board computing and LCD display eliminates the need for third-party laptops or hand held devices. Simplify your workflow with this all-in-one instrument.
- **Audio voice recorder:** Automate your field notes with the convenience of voice memos. QualitySpec Trek's audio voice recorder feature can record useful descriptive information for each data point measured.
- **Internal wavelength and white reference:** On-board internal referencing features ensure instrument performance, high data quality, and liberates the operator from the necessity of frequent white reference measurements.
- **On-board GPS:** Automatically populate accurate coordinate and elevation data for every measurement.
- **Sample clip:** Each QualitySpec Trek is shipped with a magnetic sample clip for easy measurements of thin samples like parchment, foliage, or fabric.
- **Data and project management:** Review spectra on the instrument or sync to your laptop for advanced project management functionality with the Trek PC Manager. Laptop software facilitates communication with the device and allows for easy spectrum, file and project management with user friendly file naming, spectrum review and tagging, and easy import/export capability.
- **Multiple languages:** QualitySpec Trek is ready to address global markets with fully integrated English, Mandarin Chinese, and Spanish language options.

## QualitySpec Trek specifications

Spectral range	350 - 2500 nm
Spectral Resolution	3 nm @ 700 nm 9.8 nm @ 1400 nm 8.1 nm @ 2100 nm
Height x width x depth	cm: 31 x 10 x 30 in: 12.3 x 4.0 x 11.7
Weight with battery	2.5 kg 5.5 lbs
Weight without battery	2.0 kg 4.3 lbs
Languages	English, Spanish, Mandarin Chinese



### Global and near



ASD Inc.,  
a PANalytical company  
2555 55th Street, Suite 100  
Boulder, CO 80301  
United States of America  
T +1 (303) 444-6522  
F +1 (303) 444-6825  
NIR.info@panalytical.com  
www.asdi.com

PANalytical B.V.  
Lelyweg 1, 7602 EA Almelo  
P.O. Box 13, 7600 AA Almelo  
The Netherlands  
T +31 (0) 546 534 444  
F +31 (0) 546 534 598  
info@panalytical.com  
www.panalytical.com