



# UNLOCK

THROUGHPUT. COMPLIANCE.  
DATA INTEGRITY.

## TECHNICAL NOTE

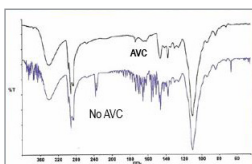
All at one go with **PerkinElmer Spectrum Two™ FT-IR**



**Nitrogen purging and desiccant change for maintenance can be time consuming and cost intensive**

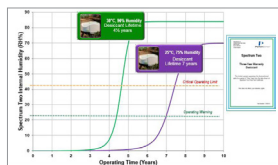
### OpticsGuard™ Technology- a differentiator in itself!

- Unique solution to FT-IR's nemesis – humidity
- Innovative design to protect the optics providing a barrier to the external environment
- No need of Nitrogen purging/ regular desiccant replacement
- Typical >5 years desiccant lifetime (90% RH)



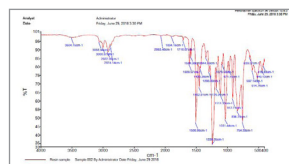
**Sample spectra with interference of atmospheric H<sub>2</sub>O and CO<sub>2</sub> peaks is a big concern**

Atmospheric Vapor Compensation™ (AVC) - No purge and no reference curves required to address interference of H<sub>2</sub>O and CO<sub>2</sub>. Highest accuracy of results



**High power consumption is too much to handle**

Spectrum Two™ unique power management system and **OpticsGuard™** Technology can be switched in and out of standby without risk to sensitive optical components, significantly reducing power consumption and long term cost of ownership.



**Noisy baseline poses a challenge always**

### PerkinElmer Spectrum Two™ with UATR sample spectra is free of any noise in baseline

PerkinElmer Spectrum Two™ Spectral data is free of baseline noise in both Transmittance and reflectance mode due to high signal to noise ratio achieved over the complete spectral region

**More no. of scans in FT-IR sample analysis is often a challenging prospect**

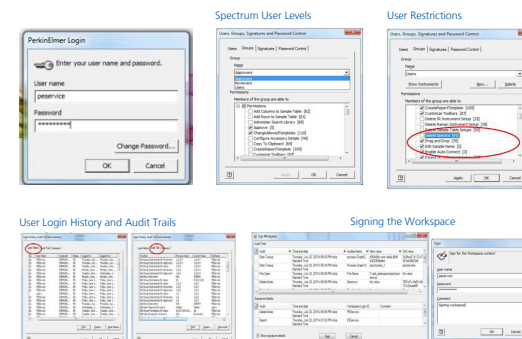
Provided the molecule is pure, good spectral quality can be achieved with lesser number of scans which ultimately reduces sample analysis time and increases the sample throughput



**FT-IR 21CFR part 11 compliance software should meet current FDA Regulatory Demands**

PerkinElmer's solution assures validation of all computerized system with incorporation of:

- User access controls (password protected User account)
- User rights
- Audit trails
- Electronic data security & Electronic signature



**PerkinElmer (India) Pvt. Ltd.**

G-Corp Tech Park, 8<sup>th</sup> Floor, Near Hypercity Mall, Ghodbunder Road, Kasarvadavali, Thane (West) – 400 615, Maharashtra, India.

Toll free: 1800 221 660 | E-mail: marketing.india@perkinelmer.com

For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright ©2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

**PerkinElmer**  
For the Better