

Cut analysis time in half

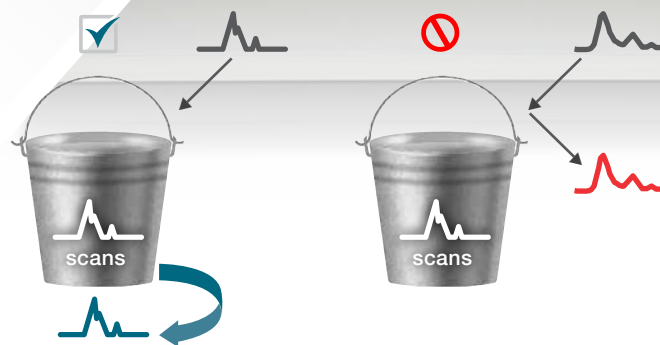
Rely on Smart Background

We are in the third decade of the 21st Century and you're doing important work. So why use an old instrument that has less computing power than your cell phone? It's time to upgrade and leave the complicated analysis to us – so you can focus on what's important, your results. The Thermo Scientific™ Nicolet™ Summit™ FTIR Spectrometers are simple, yet advanced instruments. The Smart Background feature is just one of the many intelligent ways we save you time in the lab.

What would it be like if you could literally cut your analysis time in half? Smart Background is a proprietary technology that completely removes the background collection step from your workflow. We automatically make sure you have the most recent, accurate background from your instrument, even when you aren't using it.

When you turn on Smart Background, the instrument will first collect a fresh set of backgrounds to establish a "reference bucket" of spectra. Once completed, the Lightbar (see guide to the right) will pulse blue to indicate that the instrument is constantly taking new background scans. Each new background scan is instantly compared to the reference bucket. If the new scan is similar to the reference bucket, then the oldest scan is removed from the bucket and the new one is added in. This rolling process ensures the background is up to date and accurate. If the new background scan is not similar to the reference bucket scans (i.e., you bump the instrument or place a sample on the ATR crystal), then it is automatically rejected. See the figure to the right for an illustration of the process.

The caliber of your work should match the caliber of your instruments. Make the upgrade with a Nicolet Summit FTIR Spectrometer. With features like Smart Background, you will be more efficient than ever before. Available models include the Summit LITE, Summit X, Summit PRO, and Summit OA.



Smart Background Lightbar Status Guide

-  A blue light moves back and forth as the initial reference scans are taken.
-  The Lightbar pulses blue to signify that new background scans are being taken.
-  Two solid segments in the center of the Lightbar indicate that spectra are being rejected.

Learn more at thermofisher.com/summit