

## TECHNICAL BULLETIN

## TVA2020 Temperature Compensation

**Introduction**

There have been reports of the TVA2020 drifting when monitoring temperatures are considerably different than the temperature at which the instrument was calibrated. A new menu selection in the setup menu permits you to turn on temperature compensation that may help in this condition.

**Why might this be necessary and how does this work?**

There are slight variations in electronic components from instrument to instrument and when the temperature compensation is turned on it monitors the current temperature and compares that to the temperature stored in the cal.txt file upon calibration and compensates for temperature changes. For some instruments, the compensation may help while for others (depending on the effect of temperature on the components), the impact may be minimal.

**How do I use this feature?**

Instrument firmware version 52S contains this new option and it is found under the menu tree below:

#2=Setup > #5=Other > #4=User Options > #3=More > #1=Temp Comp > #1=On and #2=Off

After installing version 52S, the temperature compensation defaults to the “ON” position. Perform a normal calibration and check performance during the day by performing a drift check. If the drift is acceptable, no actions are required. If you find that the drift is at an unacceptable level, then turning the compensation to the “OFF” position might improve performance as the impact on temperature is minimal.

Included in this new menu screen is a selection called “SN in FileName”. Activating this feature will include the instrument serial number as part of the file name when outputting logged data. Please see firmware 52S release notes “Read Me” file for more details.

Thermo Fisher Scientific  
27 Forge Parkway  
Franklin, MA 02038USA  
Telephone: (508) 520-0430  
Toll free: (866) 282-0430

[epm.techsupport@thermofisher.com](mailto:epm.techsupport@thermofisher.com)

Find out more at [thermofisher.com](http://thermofisher.com)

© 2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.