



## PODS+ Alarms/Notifications explanations

### Instrument Settings (2) Menu

- **Auto off timeout:** The period of time in minutes that elapse before the instrument automatically shuts down completely.

### Notifications Setup Menu

- **Calibration Due Warning:** Warns the operator that the calibration is due during power up of the instrument.
- **Calibration Due Alert:** Alerts the operator that the calibration is due at the beginning of every sample.
- **Coincidence Alarm:** This alarm is activated when the internally set coincidence value has been exceeded. The sample will continue to run normally until completion.
- **Particle Difference limit:** This is a percentage value that is used if the Particle Difference Alarm is enabled.
- **Particle Difference Alarm:** This alarm is activated when the Particle Difference (percentage) limit has been exceeded between the minimum and maximum values on a given channel(s) between runs of an individual sample.

**Example:** Particle Difference limit is 70, the Sample recipe calls for 3 runs. For run 1 the 4 $\mu$ m channel reports 8000 counts, the 2<sup>nd</sup> run shows 4000 counts, and the 3<sup>rd</sup> run reports 1200 counts. The count percentage difference between 8000 and 4000 is only a 50% difference so this would not exceed the Particle Difference limit (70). However, 8000 counts (run 1) compared to 1200 counts (run 3) is a difference of 85% which would exceed the Particle Difference limit of 70 and this will cause an alarm to be generated. The purpose for this feature is to detect a dramatic count change between multiple runs of the same sample that could indicate heavy particle settling quickly or some other phenomenon.

### I/O Setup Menu (Located under the Notifications Setup Menu)

**It is important to note that these alarms when enabled ONLY affect control of the I/O port. I.e. if any one of these alarms are triggered AND have been enabled within the I/O Setup menu then the I/O port would be electronically disabled immediately.**

- System Alarm
- Count Alarm
- Count Difference Alarm
- Coincidence Alarm