

Import Dialog Change (version 1.0.2)

Version 1.0.2 adds a checkbox to the Import Dialog that enables applying automatically computed correction factors to the imported records. This is meant for the following scenario: You have an above-canopy data set, and a below canopy data set. The below canopy data set contains a few above-canopy readings obtained by placing the below sensor next to the above sensor and logging (on the below unit) one or two A readings. These are logged at the beginning (or end, or both) of the below canopy data file(s).

Thus, the below canopy file data section might look like this:

```
:
### Contributing Sensors
Sensor W1      PCH-3004      4072    1477    1142    1110    961
### Observations
A      64      20100413 14:18:50    W1      141.70 229.80 247.80 125.50 29.55
B      65      20100413 14:21:52    W1      3.825  6.566  2.599  3.519  0.709
B      66      20100413 14:22:10    W1      13.52  3.187  4.179  2.189  0.739
B      67      20100413 14:22:37    W1      2.614  2.141  5.129  5.542  0.621
:
B      123     20100413 14:40:24    W1      140.50 181.60 135.40 88.48  58.30
A      124     20100413 14:40:29    W1      142.50 182.70 136.60 88.56  59.24
```

Meanwhile, the above unit was logging at regular intervals, but was never matched to the below unit.

```
:
A      341     14:17:02      1.689 11.32 7.852 4.581 2.590
A      342     14:17:17      1.598 10.57 8.211 4.585 2.607
A      343     14:17:32      1.519 9.771 8.438 4.642 2.613
A      344     14:17:46      1.489 9.337 8.482 4.762 2.615
A      345     14:18:02      1.495 9.019 8.355 4.911 2.635
A      346     14:18:17      1.067 0.709 2.919 0.101 0.065
A      347     14:18:32      61.99 50.01 49.23 20.84 4.550
A      348     14:18:48      25.40 36.08 42.53 23.68 7.634
A      349     14:19:02      25.30 34.80 42.79 23.49 7.693
:
```

If the above unit were an LAI-2000 (as was the case here), and the below an LAI-2200, then they really couldn't be easily matched anyway. This is not a problem when the below canopy file has some A readings that correspond to some readings in the above file.

To import A records for each of the B records in the below file, and “recalibrate” them to match the A readings that are already there, set the Import Dialog as follows (note the **Adjust...** checkbox on the third line): The ‘below’ file is the selected file, and the ‘above’ file (1) is checked.

Import Records into Day05.txt

For every record of type **B** found in **selected** files in Day05.txt

Insert a(n) **A** record from the files below, using **closest in time**

☒ Adjust imported values to match any existing A records

Source	File	Date	From	To	#A	#B	#?	#1	#2
▼ Day05.txt									
☐	T3	20100413 14:11:50	12:55:12	14:11:09	4	59	0	0	0
☒	1	20100413 12:50:14	12:52:02	14:45:46	456	0	0	0	0
☐	T4	20100413 14:17:29	12:55:18	14:16:59	4	59	0	0	0
☐	T6	20100413 14:49:41	14:15:10	14:47:10	3	58	0	0	0

☒ Show only source files with the same date as target files

Cancel **OK**

After the import, the “Contributing Sensors” section will show the sensor information for the imported records (if it was an LAI-2000, no serial number is shown, as is the case below). The calibration values in this case are the multipliers the program calculated to adjust the sensor’s readings.

```

:
### Contributing Sensors
Sensor W1      PCH-3004      4072   1477   1142   1110   961
Sensor W2                5.644  6.122  5.570  4.925  6.456  <- The multipliers for adjusting each ring
### Observations
A      64      20100413 14:18:50   W1      141.70 229.80 247.80 125.50 29.55
A      360     20100413 14:21:46   W2      251.46 258.16 190.11 102.79 53.99  <- Added and adjusted
B      65      20100413 14:21:52   W1       3.825 6.566 2.599 3.519 0.709
A      362     20100413 14:22:17   W2      234.25 260.74 194.73 102.35 53.31  <- Added and adjusted
B      66      20100413 14:22:10   W1       13.52 3.187 4.179 2.189 0.739
A      363     20100413 14:22:32   W2      237.75 259.51 194.73 102.74 52.72  <- Added and adjusted
B      67      20100413 14:22:37   W1       2.614 2.141 5.129 5.542 0.621
:
B      122     20100413 14:40:15   W1      113.70 169.60 110.30 79.36 49.49
A      434     20100413 14:40:17   W2      159.85 188.07 153.24 94.13 39.63  <- Added and adjusted
B      123     20100413 14:40:24   W1      140.50 181.60 135.40 88.48 58.30
A      124     20100413 14:40:29   W1      142.50 182.70 136.60 88.56 59.24

```

In practice, multiple A readings at the start and end of the below file are a good idea, especially when using the smaller view caps, where slight variations in direction or level can make a significant difference in the reading.

