



# CELL ANALYSIS

## ZE5 Cell Analyzer Data Scaling Settings in Third-Party Analysis Software

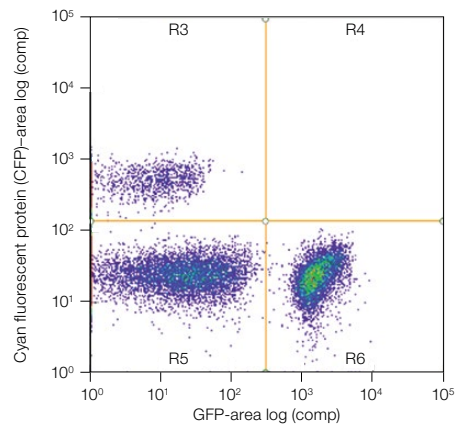
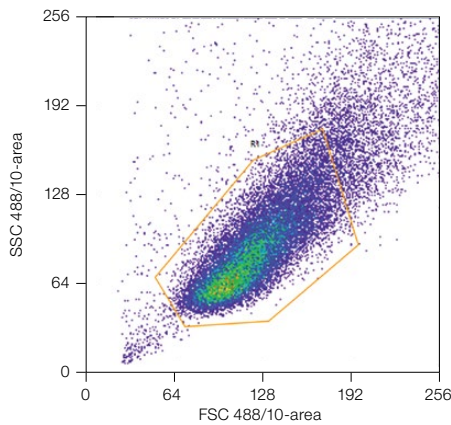
### Quick Guide

## FCS Express Software

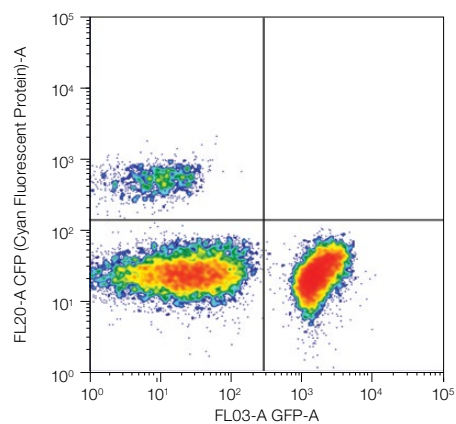
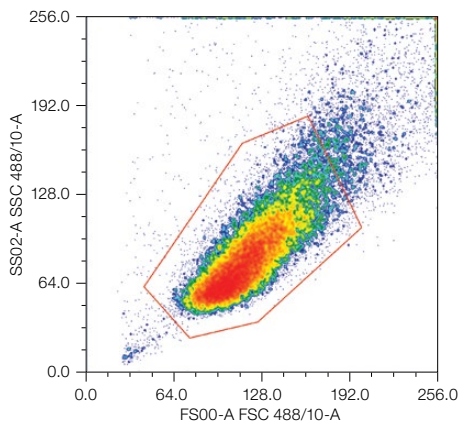
Flow cytometry data files generated on the ZE5 Cell Analyzer are exported with either “ZE5” (Everest Software v2.0 and later) or “YETI” (Everest Software v1.4 and earlier) as the keyword. The latest versions of FCS Express Software (v7.04 and greater) are able to identify both “ZE5” and “YETI” as keywords and apply ZE5 Cell Analyzer-specific settings automatically after data loading. As a result, ZE5 Cell Analyzer data display exactly the same as shown in Everest Software, and there is no need to manually adjust settings when using FCS Express Software.

Examples of data plots displayed identically in Everest and FCS Express Software:

### Everest Software




### FCS Express Software

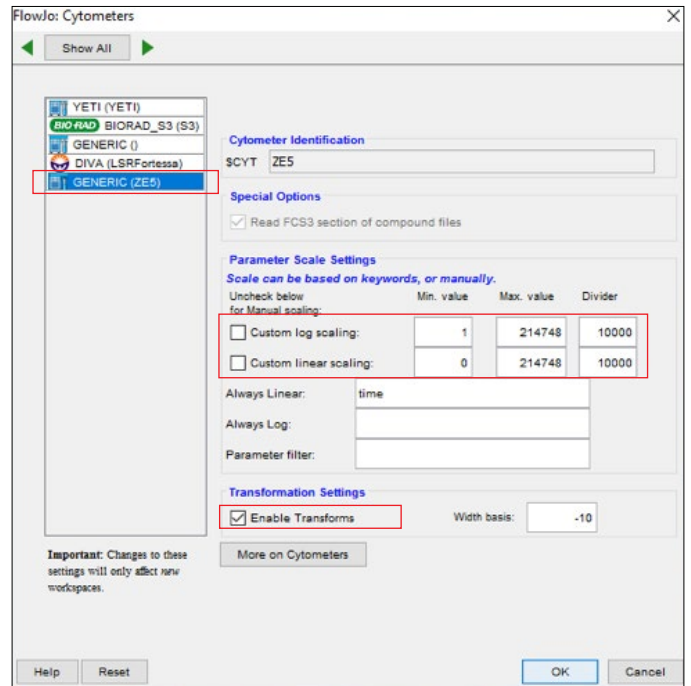


# FlowJo Software

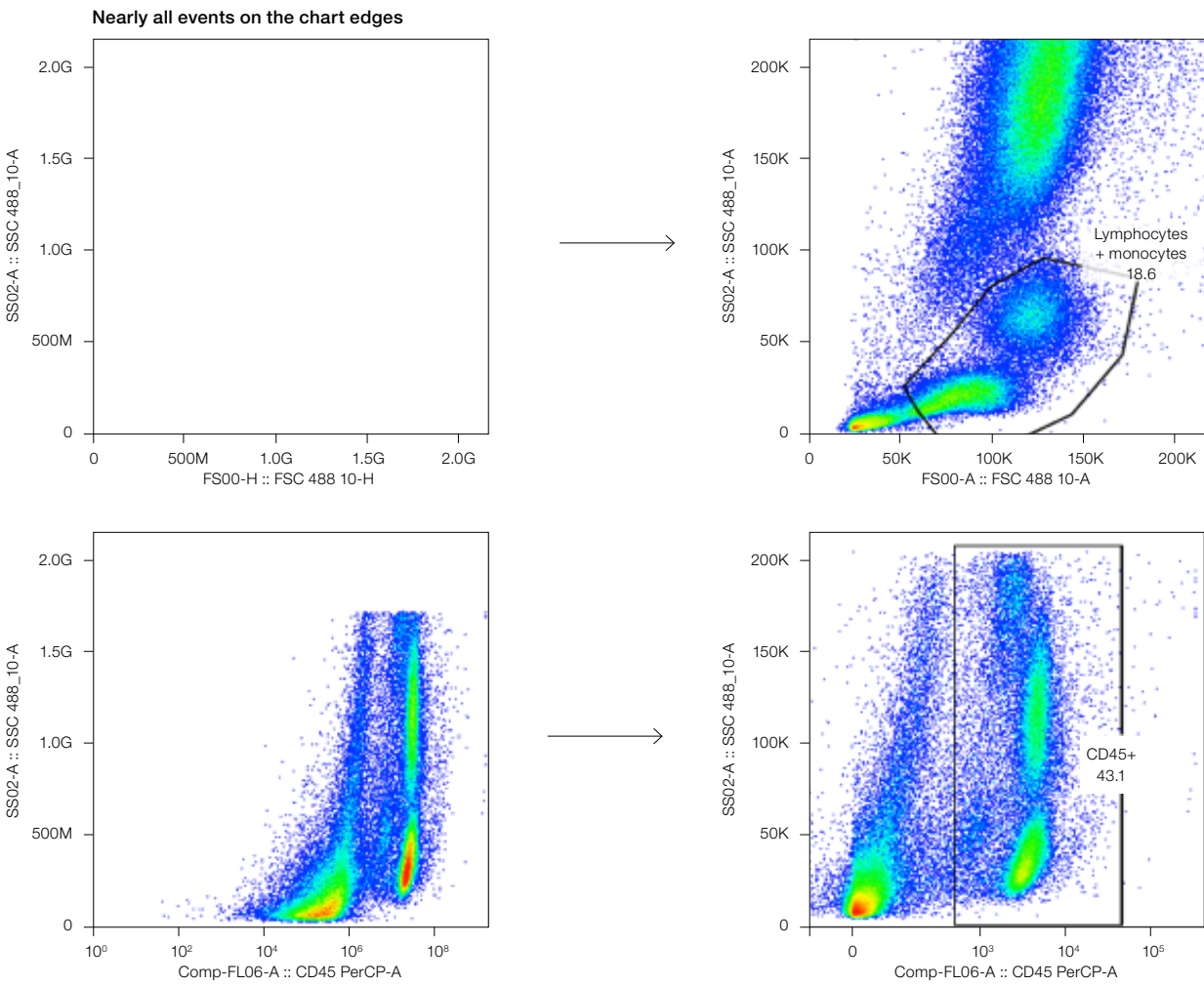
In certain versions of FlowJo Software (particularly v10.5 or earlier), it has been observed that neither “ZE5” nor “YET1” are included in the keyword library. As a result, users need to manually adjust cytometer-specific settings to properly display ZE5 data.

## Instructions

1. Click the  icon in the toolbar in FlowJo Software to access Preferences.
2. Select **Cytometers**.
3. Click the **GENERIC (ZE5)** option in the cytometer list.
4. Uncheck the **Custom log scaling** and **Custom linear scaling** boxes.
5. Set the fields in **Custom log scaling** as follows:
  - **Min. value:** 1
  - **Max. value:** 214748
  - **Divider:** 10000
6. Set the fields in **Custom linear scaling** as follows:
  - **Min. value:** 0
  - **Max. value:** 214748
  - **Divider:** 10000
7. Check the **Enable Transforms** box.
8. Click **OK** and close and relaunch FlowJo Software to apply the modified preferences.



Examples of FlowJo Software data plots before and after the manual scaling adjustment:



Visit [bio-rad.com/ZE5](https://www.bio-rad.com/ZE5) for more information.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. in certain jurisdictions.  
All trademarks used herein are the property of their respective owner.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Website** [bio-rad.com](http://bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399  
**Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23  
**France** 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300  
**Israel** 0 3 9636050 **Italy** 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Luxembourg** 00 800 00 24 67 23  
**Mexico** 52 555 488 7670 **The Netherlands** 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23  
**Portugal** 00 800 00 24 67 23 **Russian Federation** 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23  
**Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23 **Switzerland** 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311  
**United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

