



FCS Express Software for ZE5 Cell Analyzer Users

To provide a more complete flow cytometry workflow, Bio-Rad is partnering with De Novo Software to make FCS Express available for ZE5 Cell Analyzer users. FCS Express Software is a fully integrated analysis, statistics, graphing, and reporting tool that utilizes integrated spreadsheets so your graphs and statistics update as your gates change. It is optimized to support and display ZE5 Cell Analyzer data, and the analysis can be exported to PDF, PowerPoint, and six image formats at publication-quality resolution.

- All new ZE5 Cell Analyzer purchases will ship with a free 1-year FCS Express Software Internet license for noncommercial usage
- Existing ZE5 customers have the option to purchase a stand-alone 1-year license from Bio-Rad

Visit bio-rad.com/ZE5 for more information.

Ordering Information

Catalog #	Description
12013092	FCS Express Software , 1-year Internet license for academic customers
12013095	FCS Express Software , 1-year Internet license for nonacademic customers

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

