



Flow Cytometry Experiment Planner

Use our helpful, customizable Experiment Planner as an aid in the design of your flow cytometry experiments, taking you step by step through all the aspects you need to consider for successful results. Simply fill in the information about your experiment setup in the spaces provided to help you keep track of the details and allow smoother execution of your flow cytometry analysis.

Experiment name: _____

Date: _____

Basic Details

Hypothesis: _____

What will be measured by flow cytometry? _____

Sample type: _____

Target species: _____

Number of samples to be tested: _____

Total number of cells or volume of blood required: _____

Staining format (for example, FACS tubes or 96-well plate): _____

Controls (fill in all that apply)

Unstained control: _____

Positive control: _____

Negative control: _____

Secondary only control: _____

Isotype control: _____

Compensation controls on cells: _____

Compensation controls on beads: _____

Fluorescence minus one: _____

Flow cytometer information

Flow cytometer: _____

Lasers: _____

Band pass/long pass filters: _____

Speed of acquisition: _____

Number of cells to acquire: _____

See appendix for

- Plate layout
- Expanded antibody selection table

Visit [bio-rad-antibodies.com/flow-cytometry](https://www.bio-rad-antibodies.com/flow-cytometry) for further advice on flow cytometry.

Appendix
96-Well Plate Layout



