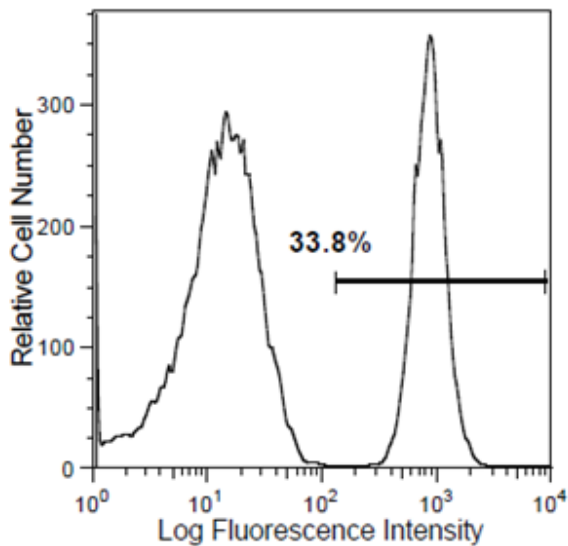
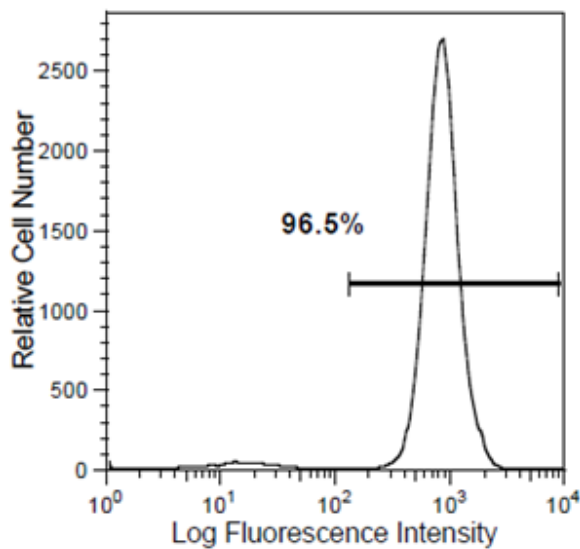


MojoSort™ Mouse Pan B Cell Isolation Kit

Before Isolation



After MojoSort™ Isolation



A single cell suspension from C57BL/6 mouse lymph nodes cells was prepared to isolate B cells using the MojoSort™ Mouse Pan B cell Isolation Kit. Cells were stained with CD19 (clone 6D5) APC. Dead cells were excluded by 7-AAD.