



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP3-00499

NBP3-00499 Protocol

Flow Cytometry Protocol for Anti-Mouse Ig (H+L) Comp-Bead 3 Population (7.5 um) Kit (NBP3-00499)

Shake vigorously or vortex before use. Sonication will help to increase the number of singles. Use for routine quality control of mouse antibody conjugates used in flow cytometry. Add 1 drop from each bottle to a centrifuge tube, followed by the addition of the conjugate being evaluated. After exposure to appropriate mouse antibody conjugates, three different fluorescent peaks shall be resolved by flow cytometry.