

ReadiUse™ Multiplexing Staining Buffer

Introduction

ReadiUse™ Multiplexing Staining Buffer is formulated for immunofluorescent staining applications using XV conjugates in multicolor flow cytometry experiments. The buffer is designed to reduce nonspecific staining and improve staining consistency when preparing multiplex fluorescent antibody cocktails for cell analysis. ReadiUse™ Multiplexing Staining Buffer is compatible with traditional fluorescent conjugates labeled with organic fluorochromes, Brilliant Violet dyes, and PE, APC or their tandem dye conjugates.

Specifications

Product Number: 70808 & 70809 Product Name: ReadiUse™ Multiplexing Staining Buffer

Unit Size: 100 tests & 1000 tests Volume: 25 µL/test Appearance: Aqueous buffered solution

Stabilizer: BSA and <0.1 % sodium azide

Storage Conditions: Store at 2-8°C, and protected from light.

Applications

ReadiUse™ Multiplexing Staining Buffer is used as a medium to mix fluorescent reagents with cells to decrease nonspecific staining in multicolor flow cytometry experiments. The buffer supports preparation of multiplex antibody staining cocktails for immunofluorescent cell analysis workflows.

Sample Experimental Protocol

1. Add 25 µL of ReadiUse™ Multiplexing Staining Buffer to a tube.

Note: The 25 µL ReadiUse™ Multiplexing Staining Buffer volume does not depend on the final staining volume or amount of cells used.

2. Add each fluorescent reagent at the recommended volume per test. Mix well after addition of each antibody, and then add the next one.
3. The antibody conjugate cocktail is ready to stain the aliquoted cells in tubes.

Appendix I: Product Ordering Information

Catalog Number	Product Name	Unit Size
70808	ReadiUse™ Multiplexing Staining Buffer	100 Tests
70809	ReadiUse™ Multiplexing Staining Buffer	1000 Tests