

Flow Cytometry Compensation Beads

User Manual Version 1.2
April 30, 2024

RayBright® Flow Cytometry Universal Compensation Beads Staining Protocol

Catalog numbers: 137-00013-25 (25 tests)
137-00013-100 (100 tests)



We Provide You With Excellent
Support And Service

Tel: 770-729-2992; Fax:770-206-2393;
Web: www.raybiotech.com Email: info@raybiotech.com



RayBiotech, Inc.

RayBright® Universal Compensation Beads

TABLE OF CONTENTS

I. Introduction.....	1
II. Kit Contents.....	2
III. Assay Protocol.....	2
IV. Storage and Stability.....	3

I. INTRODUCTION

The RayBright® Universal Compensation Beads are designed to capture all isotypes of common primary antibodies to allow for easy and rapid creation of FACS compensation beads. Our beads can capture all rat, mouse, Syrian Hamster and Armenian hamster antibodies that are labeled with your fluorophores of interest. RayBright® Universal Compensation Beads are made up of two populations: negative beads and positive beads, with only the positive beads binding the fluorochrome-conjugated antibodies. It's easy and convenient to use, and a perfect tool for multi-color flow cytometry compensation.

II. KIT CONTENTS

Components	137-00013-25	137-00013-100	Part Number
	25 tests	100 tests	
RayBright® Compensation Beads (supplied in drop bottle)	1.5 mL	6 mL	Item A

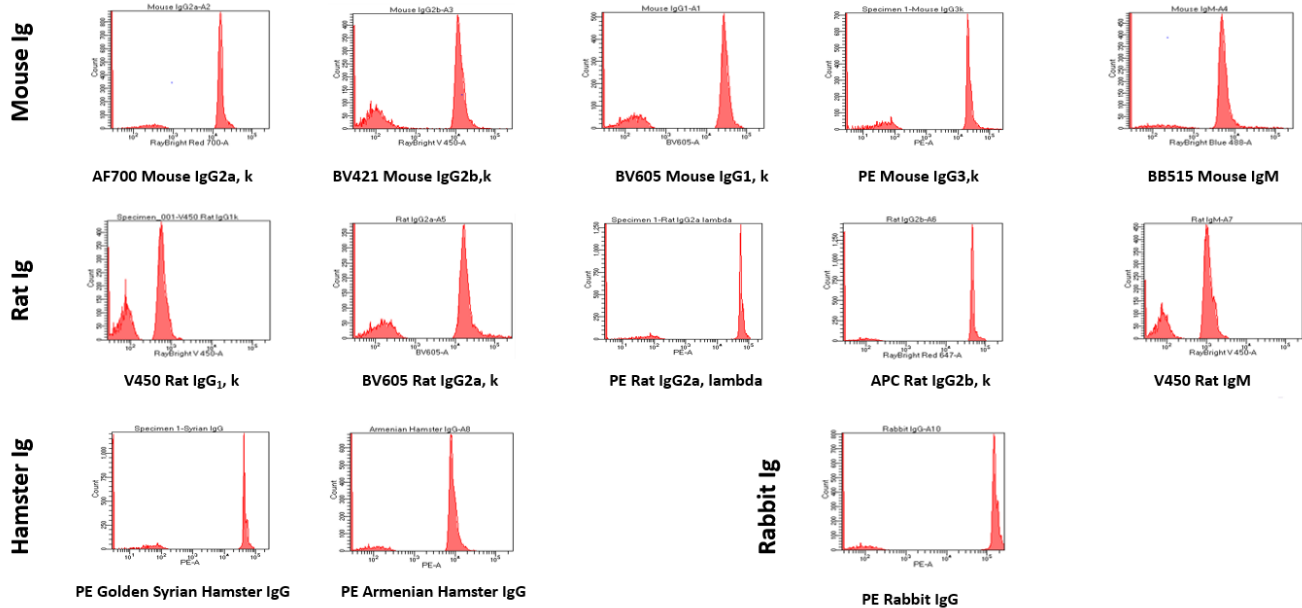
Kit can be stored at 4°C, protected from light for up to 3 months

III. Assay Protocol

1. Vortex RayBright® Compensation Beads.
2. Add 1 drop (about 50 µL) to a 5 ml FACS tube.
3. Add same amount of antibody for cell-staining and mix.
4. Incubate for 15-30 minutes on ice or 4°C.
5. Wash the beads with 1 ml of FACS buffer
6. Spin down the beads at 500g for 10 minutes at 4°C.
7. Remove the supernatant without disturbing the beads.
8. Resuspend the beads in 0.5 ml of FACS buffer. The beads are ready for compensation analysis.

Representative Data:

- 0.5 μg of RayBright[®] fluorochrome-conjugated antibodies were incubated with 50 μL of RayBright[®] Universal Compensation beads for 15 minutes at room temperature. Protocol outlined above.



IV. Storage and Stability

Store kit at 4°C, protected from light for up to three months for full functionality

This product is for research use only.