

---

## Streptavidin: RPE-Cy5 conjugate

(Catalog No. A16003)

---

### Description

Streptavidin: RPE-Cy5

### Purification

Streptavidin conjugated to R-Phycoerythrin - Cy5 - liquid

### Size

0.2 ml

### Formulation

Phosphate buffered saline. 0.1% Sodium Azide (NaN<sub>3</sub>) Stabilizing agent

### Specificity

Biotinylated macromolecules

### Storage

Store at 4°C. DO NOT FREEZE. This product should be stored undiluted. This product is photosensitive and should be protected from light.

### Handling Recommendations

Should this product contain a precipitate we recommend microcentrifugation before use. For maximum recovery of the products, centrifuge the vial prior to opening the cap.

### Applications & Suggested Dilutions

Flow Cytometry: Use 10ul of the suggested working dilution to label 10<sup>6</sup> cells in 100ul. Method sheets are available on request, 1/20 - 1/30; Where this product has not been tested for use in a particular technique this does not necessarily exclude its use in such procedures. Suggested working dilutions are given as a guide only. It is recommended that the user titrates the protein for use in their own system using appropriate negative/positive controls.

### Ordering Information:

| Products                        | Size   | Cat. No. |
|---------------------------------|--------|----------|
| Streptavidin: RPE-Cy5 conjugate | 0.2 ml | A16003-1 |

*This product is for research purposes only. Not intended for use in diagnostic procedures.*