



NOTICE: ProZyme was purchased by Agilent in July 2018. Documents for products and product lots manufactured before August 2019 will contain references to ProZyme. For more information about these products and support, go to: www.agilent.com/en/contact-us.



PRODUCT DESCRIPTION SHEET

PRODUCT NAME: PHYCOLINK® STREPTAVIDIN-R-PHYCOERYTHRIN SAMPLER KIT

PRODUCT CODE: PJ3SX

BUFFER: Sterile-filtered solutions in 10 mM Tris-HCl, 150 mM NaCl (pH 8.2)

PRESERVATIVE: 1 µg/ml pentachlorophenol

STORAGE: Store at 2-8°C protected from light. DO NOT FREEZE.

The Sampler Kit contains six Streptavidin-R-Phycoerythrin conjugates which vary in overall size. Each conjugate was designed to address customer requirements for various applications. The applications listed are suggestions only, customers are encouraged to use the samples provided to determine the ideal conjugate for their use.

Table 1 – KIT COMPONENTS

Component Name	Quantity per Kit	Conjugate Size	Typical Application
PJRS301 Streptavidin-RPE (0.25 mg)	1 each	Large	Luminex® / Bead-Based Assays
PJRS20 Streptavidin-RPE (0.25 mg)	1 each	Large	
PJRS34 Streptavidin-RPE (0.25 mg)	1 each	Large	
PJRS25 Streptavidin-RPE (0.25 mg)	1 each	Medium	Tetramer
PJRS27 Streptavidin-RPE (0.25 mg)	1 each	Medium	
PJ39S Streptavidin-RPE (0.25 mg)	1 each	Small	Flow Cytometry

QC RELEASE

The components conform to specifications for each product – see product and lot-specific Certificates of Analysis included with the Kit for specifications and data.

APPLICATIONS

Our range of Streptavidin-R-Phycoerythrin (SA-RPE) conjugates differ in overall size. We recommend that you test the conjugates to find out which is the best for your system.

PJRS301 is our newest conjugate, developed for long-term signal consistency in Luminex assays.

We do not have specific dilutions protocols for our SA-RPE conjugates, as this may depend on your system & application. Below are some suggestions based on application:

Luminex

Dilution of SA-RPE to 2-20 µg/ml for addition to wells is a good starting point. For our SA-RPE Luminex QC assay we dilute to 10 µg/ml and add 25 µl (0.25 µg) per well.

Flow cytometry

The final conjugate concentration needed in your system should be determined by titration to optimize signal & background. However, dilutions covering the following concentration range might be a decent starting point:

0.01 µg/ml, 0.03 µg/ml, 0.1 µg/ml, 0.3 µg/ml

MHC tetramers

We suggest using the NIH Emory Tetramer Core Facility protocol as a starting point, which uses PJRS25 (SA-RPE). PJRS27 may also be appropriate. PJ27S (SA-APC) is also used in the Emory protocol.

<http://tetramer.yerkes.emory.edu/support/protocols>