

## Protonex™ Red 600-Latex Bead Conjugate

Catalog number: 21209

Unit size: 1 mL

Component	Storage	Amount (Cat No. 21209)
Protonex™ Red 600-Latex Bead Conjugate	Refrigerated (2-8 °C), Minimize light exposure	1 vial (1 mL)

### OVERVIEW

Protonex™ Red-latex bead conjugate demonstrated pH-dependent fluorescence. Unlike most of the existing fluorescent dyes that are more fluorescent at higher pH, acidic conditions enhance the fluorescence of Protonex™ Red-latex bead conjugate. The fluorescence of Protonex™ Red-latex bead conjugate dramatically increases as pH decreases from neutral to the acidic, making it a robust tool to study phagocytosis and its regulation by drugs and/or environmental factors. The lack of fluorescence outside the cell eliminates the wash steps. Protonex™ Red-latex bead conjugate provides a powerful tool to study phagocytosis. Protonex™ Red-latex bead conjugate is low fluorescent outside the cells, but fluoresce brightly red in acidic compartments (such as phagosomes, lysosomes and endosomes). This Protonex™ Red-latex bead conjugate can be also used for multiplexing cell functional analysis with green dyes such as GFP, Fluo-8, calcein, or FITC-labeled antibodies. Protonex™ Red has the spectral properties similar to those of Texas Red, making the common filter set of Texas Red readily available to the assays of Protonex™ Red.

### AT A GLANCE

#### Chemical and Physical Properties

<b>Solvent:</b>	Water
<b>Solids Content:</b>	1% in PBS
<b>Number of Microspheres per mL:</b>	~4e+10
<b>Ex/Em:</b>	575/597 nm
<b>Mean Diameter:</b>	0.72 µm

### KEY PARAMETERS

#### Fluorescence microscope

Emission	Texas Red filter set
Excitation	Texas Red filter set
Recommended plate	Black wall/clear bottom

### SAMPLE EXPERIMENTAL PROTOCOL

#### Important

The following is a recommended protocol for granulocytes. This protocol only provides a guideline and should be modified according to your specific experimental conditions.

#### Protocol

1. Prepare cells as desired. For example, prepare the granulocytes at 10<sup>7</sup> cells/mL with Hanks and 20 mM Hepes buffer (HHBS), and

add 100 µL to a polypropylene tube.

**Note:** Each cell line should be evaluated on an individual basis to determine the optimal cell density.

2. Add 1-10 µL of the Protonex™ Red 600-Latex Bead Conjugate to the tube and incubate with gentle shaking for 30 minutes at 37 °C.

**Note:** Each cell line should be evaluated on an individual basis to determine the optimal incubation time.

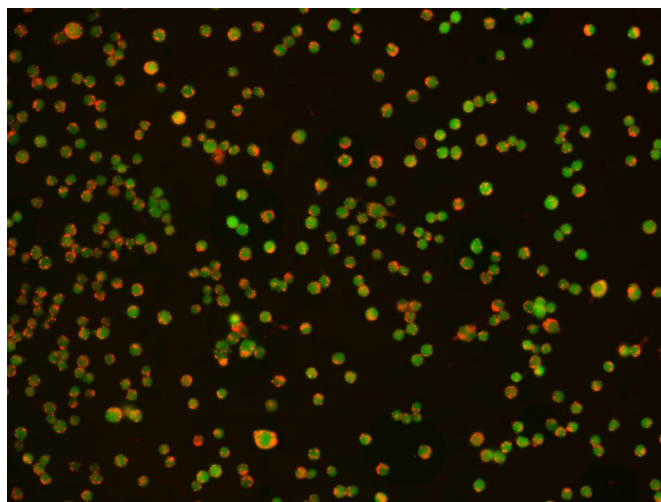
3. Prepare an identical sample that is incubated at 4 °C and label it as a control.
4. At the end of the 30-minute incubation, stop phagocytosis by adding 2mL of ice-cold HHBS and mix well.

5. Wash the cells 2 times with cold HBSS.

6. Resuspend the cells in 500 µL of cold HBSS, keep the samples at 4 °C, and analyze immediately using a fluorescence microscope equipped with a Texas Red® filter set.

**Note:** For fluorescence microplate readers, monitor the fluorescence intensity at Ex/Em = 570/600 nm (Cutoff = 585 nm).

### EXAMPLE DATA ANALYSIS AND FIGURES



**Figure 1.** Phagocytosis was examined in RAW 264.7 cells by Protonex™ Red 600-Latex Bead Conjugate (Cat # 21209). The cells were incubated with Protonex™ 600 Latex Beads in a growth medium for 4 hours. CytoTrace™ Green CMFDA (Cat # 22017) was used to track live cells. The image (20X) was taken using Keyence

Fluorescence Microscopy.

**DISCLAIMER**

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email [info@aatbio.com](mailto:info@aatbio.com) if you have any questions.