

Protonex™ Green 500-Latex Bead Conjugate

Catalog number: 21223

Unit size: 1 mL

Component	Storage	Amount (Cat No. 21223)
Protonex™ Green 500-Latex Bead Conjugate	Refrigerated (2-8 °C), Minimize light exposure	1 mL

OVERVIEW

The Protonex™ Green 500-Latex Bead Conjugate is a ready-to-use reagent designed to study phagocytosis and phagosome acidification in live cells. This conjugate combines synthetic latex beads with Protonex™ Green 500, a novel pH-sensitive fluorophore that remains non-fluorescent at neutral pH and becomes highly fluorescent upon entering acidic environments such as maturing phagosomes and phagolysosomes.

As a standalone reagent, it enables users to integrate phagocytic detection into their own custom assays. The FITC-like excitation/emission properties of Protonex™ Green 500 make it compatible with a wide range of fluorescence imaging and detection systems. These conjugates can be used in combination with red fluorescent dyes like RFP, Calbryte™ 630 calcium dye, calcein red, or Cy5-labeled antibodies for multiplexed cell functional analysis. It is ideal for immunological research, drug discovery, and mechanistic studies of innate immune function, autophagy, or particle uptake.

AT A GLANCE

Chemical and Physical Properties

Solvent:	Water
Solids Content:	1% in PBS
Number of Microspheres per mL:	~4e+10
Ex/Em:	445/503 nm
Mean Diameter:	0.72 µm

KEY PARAMETERS

Fluorescence microscope

Emission	FITC filter set
Excitation	FITC 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 10^7 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™ Green 500-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 FITC filter set.

Note: For fluorescence microplate readers, monitor the fluorescence intensity at Ex/Em = 450/510 nm (Cutoff = 490 nm).

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 if you have any questions.