

Examining the Kinetics of Neutrophil Phagocytosis by Flow Cytometry

Author

Agilent Technologies, Inc.

Abstract

Studying neutrophil phagocytosis is essential to understanding the initial host response to pathogens. This application note demonstrates the analysis of neutrophil phagocytosis under various conditions using the Agilent NovoCyte flow cytometer, and quantifies the effect of a potent actin-remodeling inhibitor, cytochalasin D. The data generated show that measuring phagocytosis with the NovoCyte flow cytometer enables the rapid analysis of millions of cells, making results easier to achieve.

Examining the kinetics of neutrophil phagocytosis by flow cytometry

The immune system has evolved specialized cells that engulf and kill invading pathogens in a process known as phagocytosis. These cells, known as phagocytes, are essential for elimination of dangerous micro-organisms, as well as apoptotic bodies, from the host. Phagocytosis can be divided into two key stages: the uptake of particles by ingestion, followed by clearance of the particle by degradation.

Initiation of phagocytosis begins with the interaction of receptors on the phagocyte with ligands on the target particle. These ligands can be components of the target particle, such as lipopolysaccharide (LPS), or host proteins that coat a target particle and are recognized by phagocytes, also known as opsonization. Proteins used for opsonization include immunoglobulins, components of the complement cascade, and other serum proteins.

There are several types of phagocytes including neutrophils, macrophages, dendritic cells, and others. Neutrophils account for 50 to 70% of circulating leukocytes and are the first line of defense against many pathogens. During an inflammatory response, neutrophils are rapidly recruited to the site of infection and fight pathogens by both phagocytosis and release of antimicrobial molecules. Neutrophils are extremely efficient phagocytes that can internalize IgG-opsonized targets in less than 20 seconds. Therefore, studying neutrophil phagocytosis is essential to understanding the initial host response to pathogens.

Two methods for measuring phagocytosis are routinely used: microscopy and flow cytometry. Both techniques measure the capability of phagocytic cells to ingest a fluorescent particle (labeled micro-organisms or beads). There are several advantages to using flow cytometry to measure phagocytosis. Microscopy requires time-consuming manual counting of ingested particles, which limits the total number of cells analyzed. Conversely, flow cytometry can analyze thousands of cells per second, facilitating statistical analysis while offering multiparametric measurements for simultaneous evaluation of neutrophil features.

Finally, neutrophils are easy to prime and activate by manipulation in the lab, which is a necessary process for microscopic analysis of phagocytosis. Flow cytometry allows the measurement of neutrophil phagocytosis in white blood cells (WBCs) without creating artifacts, and may be more representative of phagocytosis *in vivo*. This application note measured neutrophil phagocytosis on the NovoCyte flow cytometer under various conditions, and quantified the effect of a potent actin-remodeling inhibitor, cytochalasin D.

Neutrophil phagocytosis increases with incubation time and particle concentration

Cell phagocytosis can be measured using flow cytometry by quantifying the engulfment of fluorescent particles. These studies often use cultured bacteria, yeast polysaccharides, or latex/polystyrene beads as the target particle for phagocytosis. Particles

opsonized by immunoglobulins are recognized by Fcγ receptors that specifically interact with the Fc portion of IgG, resulting in receptor activation and the initiation of phagocytosis.

The kinetics of neutrophil phagocytosis were measured by adding IgG-coated, FITC-labeled latex beads to leukocytes, and quantifying the percentage of cells that had engulfed the beads (Figure 1). Under these conditions, phagocytosis of opsonized beads steadily increases over time, reaching ~70% at 160 minutes. Increased phagocytosis also results in an increase in cell side scatter (SSC) of neutrophils because the cells have ingested solid beads (Figure 1A). The phagocytosis of control beads did not change over time, indicating that these results are due to specific IgG-mediated phagocytosis.

Next, we examined how an increased ratio of target beads to neutrophils affects the rate of phagocytosis (Figure 2). Phagocytosis directly correlates with the increased ratio of opsonized beads to neutrophils. Almost 90% of cells had undergone phagocytosis at a ratio of 50:1, where an apparent increase in SSC is observed (Figure 2A). However, when the beads-to-neutrophils ratio is greater than 10:1, a noticeable increase in phagocytosis occurs in the control beads, reaching 20% phagocytic cells at 50:1. These data indicate that the ideal ratio of beads to neutrophils that should be used for these assay conditions is 10:1, to avoid an increase in nonspecific phagocytosis of target beads.

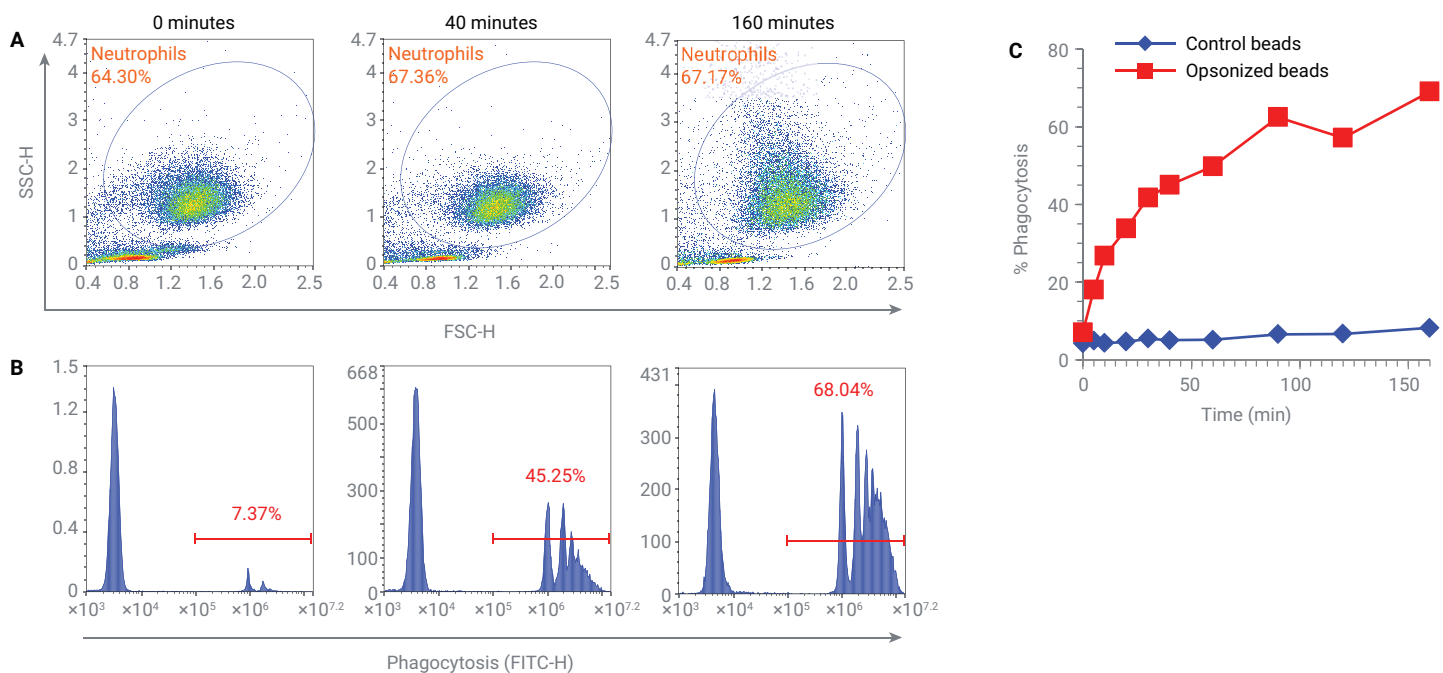


Figure 1. Neutrophil phagocytosis increased with incubation time. FITC latex beads (2 μ m) coated with IgG were mixed with isolated WBCs at a ratio of 10:1. Cells were incubated for the indicated times (0 to 160 minutes) then washed with cold PBS and 0.02% EDTA, followed by the addition of Trypan blue to quench the extracellular fluorescent signal. Monocytes adhered to the plate, leaving only granulocytes and lymphocytes. Lymphocytes are eliminated by FSC (A). Representative plots of phagocytosis of opsonized beads at 0, 40, and 160 minutes are shown for FSC/SSC (A) and FITC (B). Kinetic data for frequency of phagocytic cells are shown for both opsonized (red) and control FITC beads (blue) (C).

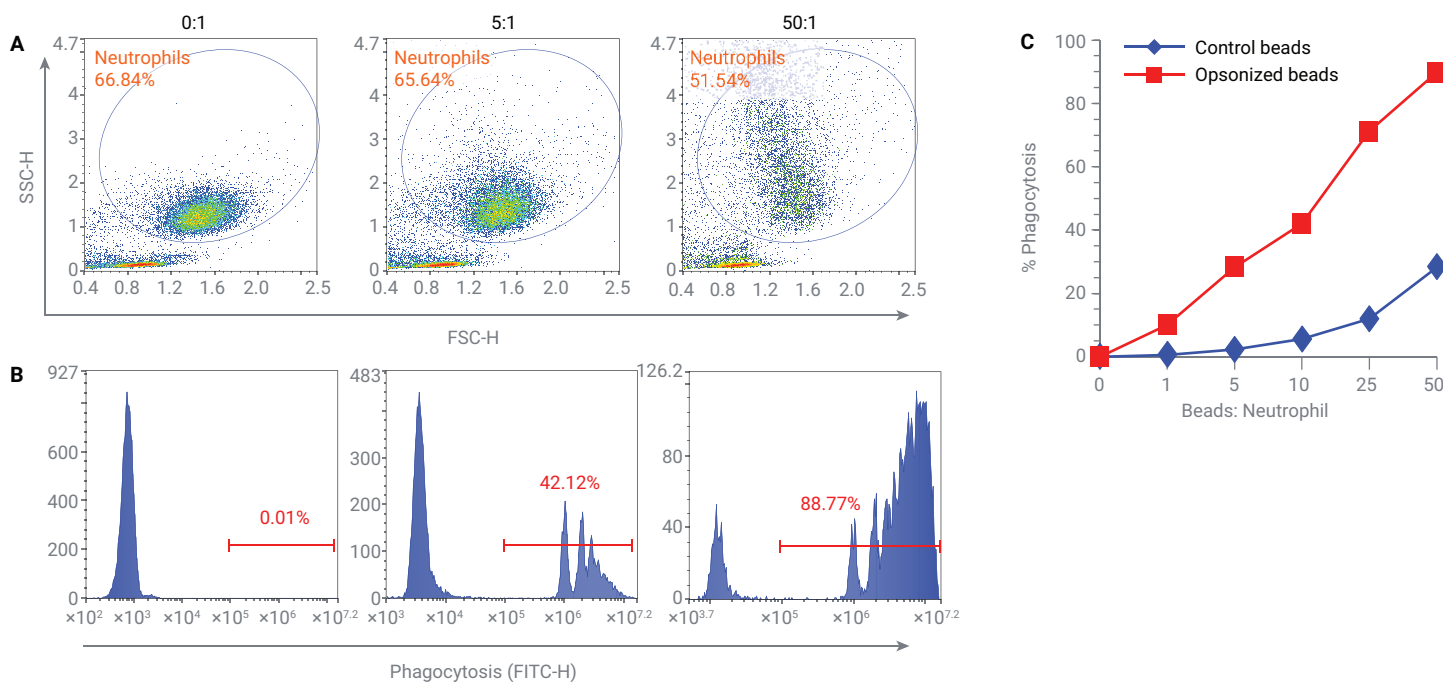


Figure 2. Neutrophil phagocytosis increases with increased bead to cell ratio. FITC latex beads (2 μ m) coated with IgG were mixed with isolated WBCs at the indicated ratio of beads to neutrophils for 40 minutes, then washed with cold PBS and 0.02% EDTA, followed by addition of Trypan blue to quench the extracellular fluorescent signal. Representative plots for ratios of 0:1 (no beads), 5:1, and 50:1 are shown for FSC/SSC (A) and FITC (B). Frequency of phagocytic cells are quantified for both opsonized (red) and control FITC beads (blue) (C).

Inhibition of phagocytosis with cytochalasin D

Many conditions and agents have been reported to impair phagocytosis, including inhibition of actin rearrangement, which is necessary during the uptake of large particles. Cytochalasins are antagonists of actin

remodeling, and commonly used to inhibit phagocytosis. Therefore, the effect of cytochalasin D on neutrophil phagocytosis was examined. Increasing concentrations of cytochalasin D dramatically inhibited neutrophil phagocytosis, which was nearly halved at 370 nM as the IC₅₀ was calculated to be 0.439 μ M (Figure 3).

Complete ablation of phagocytosis was achieved with 10 μ M cytochalasin D, and a dose-dependent effect was observed. These data demonstrate that cytochalasin D has a strong concentration-dependent inhibitory effect on phagocytosis, and measurements by flow cytometry are specific and sensitive enough for these types of assays.

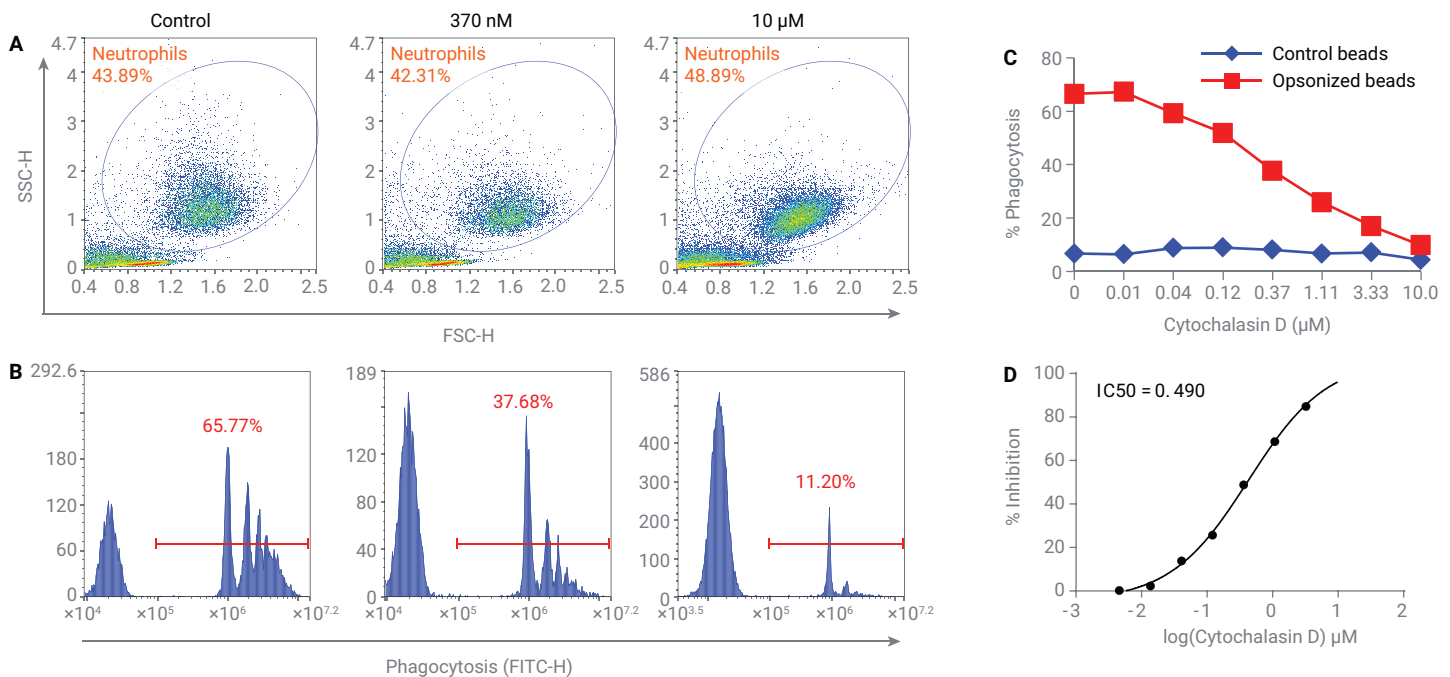


Figure 3. Neutrophil phagocytosis decreases with the addition of the actin-remodeling inhibitor, cytochalasin D. FITC latex beads (2 μ m) coated with IgG were mixed with isolated WBCs at a ratio of 20:1 with indicated concentrations of cytochalasin D. Cells were incubated for 30 minutes, then washed with cold PBS and 0.02% EDTA, followed by addition of Trypan blue to quench the extracellular fluorescent signal. Representative plots of control (no inhibitor), 370 nM, and 10 μ M cytochalasin D are shown for FSC/SSC (A) and FITC (B). The frequency of phagocytic cells is shown for both opsonized (red) and control FITC beads (blue) (C). The cytochalasin D concentration-dependent inhibitory effect on phagocytosis IC₅₀ is 0.439 μ M (D).

Conclusion

This application note investigated the effect of time and bead-to-cell ratio on neutrophil phagocytosis. Increasing both the time and bead concentrations enhanced IgG-induced neutrophil phagocytosis. It was also seen that cytochalasin D is a potent inhibitor, capable of complete abrogation of neutrophil phagocytosis. Measuring phagocytosis with the NovoCyte flow cytometer allows rapid analysis of millions of cells, making results easier to achieve.

References

1. Bassøe, C. F. Assessment of Phagocyte Functions by Flow Cytometry. *Curr. Protoc. Cytom.* **2002** Aug, Chapter 9, Unit 9.19.
2. Greenberg, S.; Grinstein, S. Phagocytosis and Innate Immunity. *Curr. Opin. Immunol.* **2002**, 14, 136–45.
3. Kolaczowska, E.; Kubes, P. Neutrophil Recruitment and Function in Health and Inflammation. *Nature Reviews Immunology* **2013**, 13, 159–175.
4. Chow, C. W.; Downey, G. P.; Grinstein, S. Measurements of phagocytosis and phagosomal maturation. *Curr. Protoc. Cell Biol.* **2004** May, Chapter 15, Unit 15.7
5. Heyman, B. Regulation of Antibody Responses Via Antibodies, Complement, and Fc Receptors. *Annu. Rev. Immunol.* **2000**, 18, 709–737.

www.agilent.com/chem

For Research Use Only. Not for use in diagnostic procedures.

This information is subject to change without notice.