

# Correlating internalization and potency to accelerate antibody discovery and development

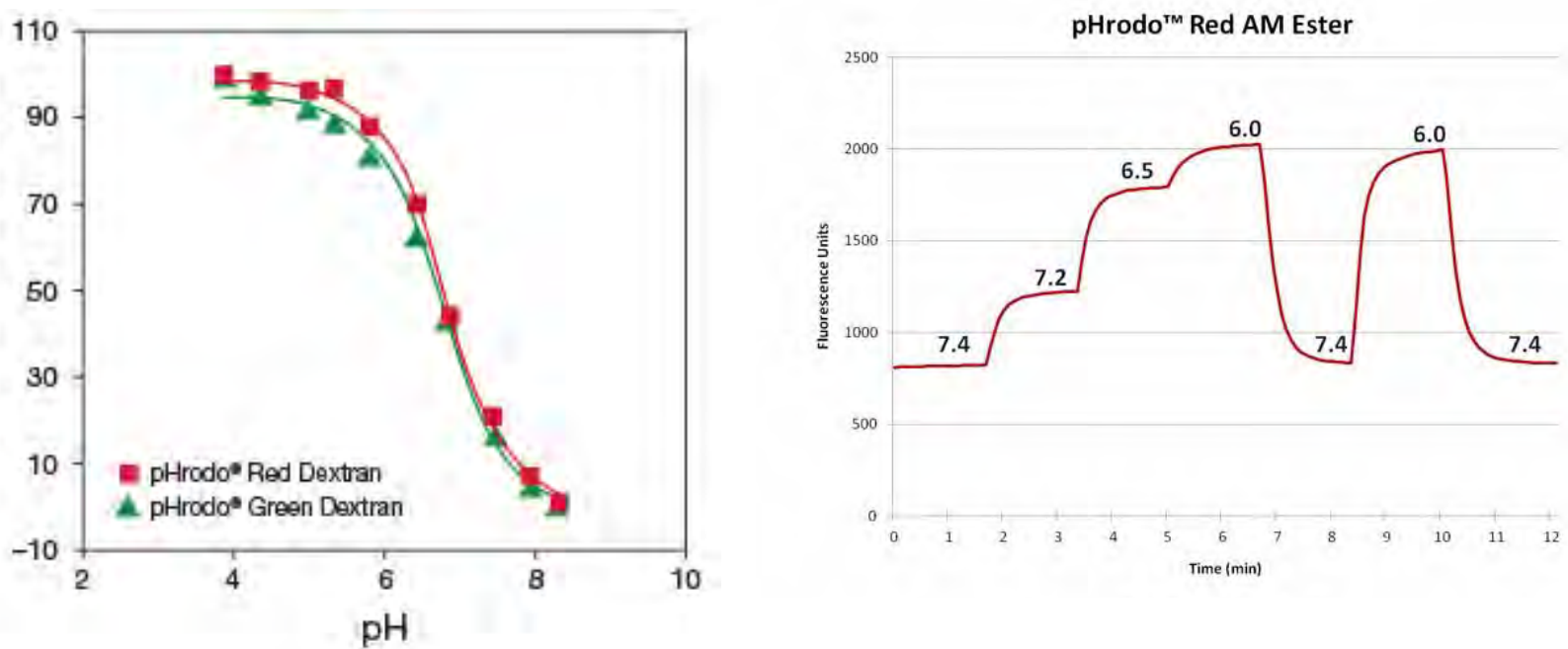
C. Langsdorf, R. Aggeler, B. Agnew, D. Beacham, J. Berlier, K. Chambers, A. Chen, A. Dix, N. Dolman, K. Gee, M. Janes, L. Smolenska, W. Zhou; Molecular Probes, part of Thermo Fisher Scientific, Eugene, OR, USA

## ABSTRACT

The basic cellular mechanisms of internalization and trafficking are important to many areas of cell biology, and especially to the proper function of therapeutic antibodies. Those antibodies intended for use as antibody drug conjugates should specifically bind to target cells and rapidly internalize into acidic compartments. Conversely, antibodies intended to kill cells via direct cell death, complement cascade, or effector cell killing should remain bound to the external surface of target cells as long as possible. However, the ability to study these internalization processes has historically been limited by the lack of tools to directly monitor the internalization and subsequent acidification of extracellular material.

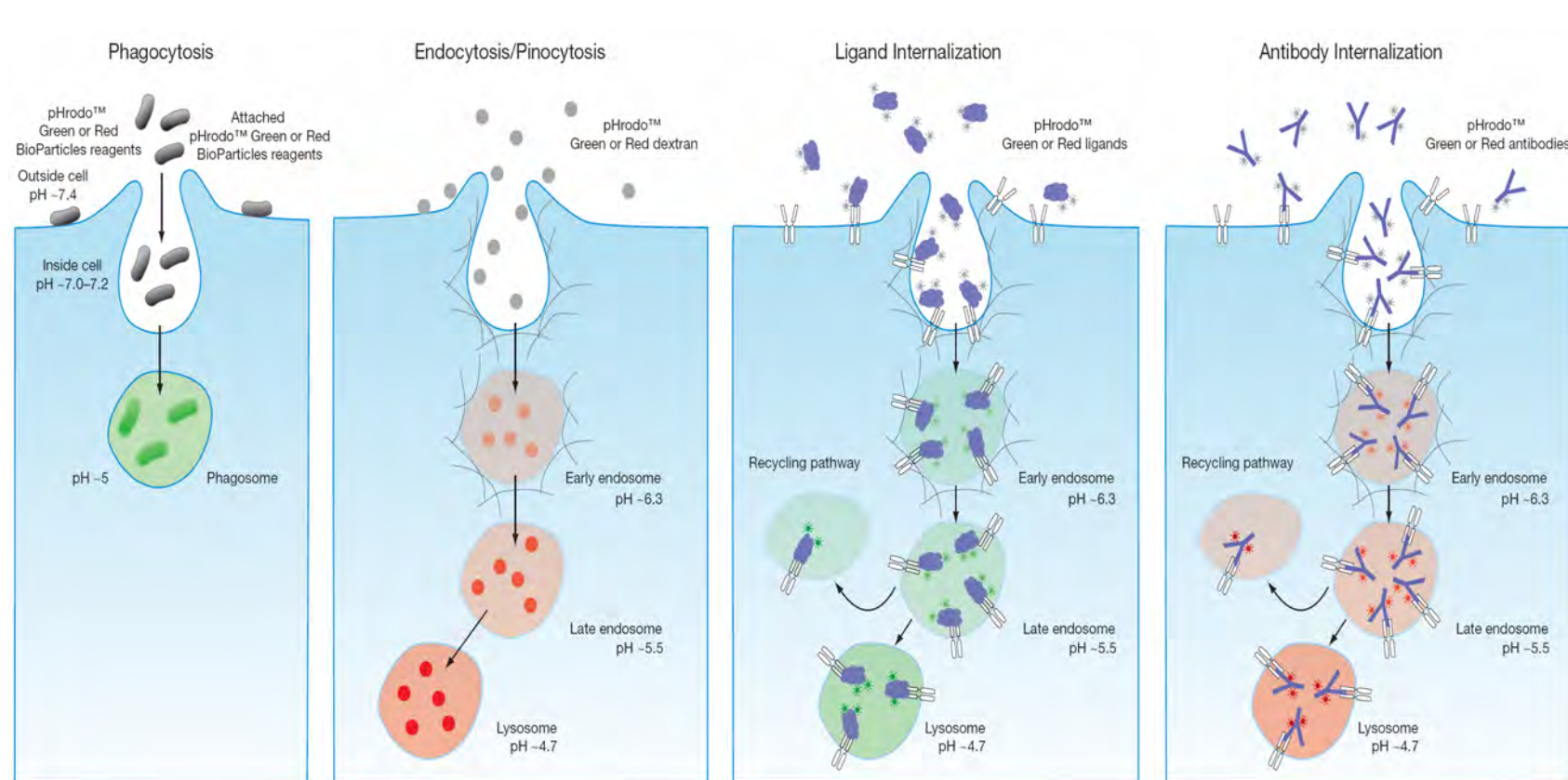
Here we present experimental data demonstrating the use of pH-sensitive fluorophores to monitor the three primary mechanisms of endocytosis in live cells. Additionally we demonstrate three approaches to directly visualize internalization of therapeutic antibodies and ADCs in live cells. Two of these techniques involve site-specific antibody labeling to screen for and further characterize internalizing clones without affecting antibody structure or function. These techniques are demonstrated using quantitative and widefield imaging to study anti-Her2 antibodies and ADCs in breast cancer models. The internalization of ADCs into Her2-positive breast cancer cells shows a strong correlation with target cell killing. We believe the insights provided by these approaches have the potential to accelerate therapeutic antibody lead generation and development.

Figure 1. pHrodo™ pH Sensor Dyes



A. pH response profile of pHrodo Red and pHrodo Green dextran conjugates. B. pHrodo fluorescence dynamically responds to pH changes. NIH-3T3 cells loaded with 5 μM pHrodo™ Red AM Intracellular pH Indicator were washed with a series of HEPES-based pH standard buffers containing 10 μM nigericin and 10 μM valinomycin to clamp intracellular pH to the indicated values.

Figure 2. Fluorogenic pH Sensing of Cellular Internalization Processes



The fluorescence intensity of pHrodo™ Red and Green dyes dramatically increases as pH drops from neutral to acidic, making them ideal tools to study endocytosis, phagocytosis, and ligand and antibody internalization.

Figure 3. Monitoring pinocytosis with fluorescently-labeled dextran

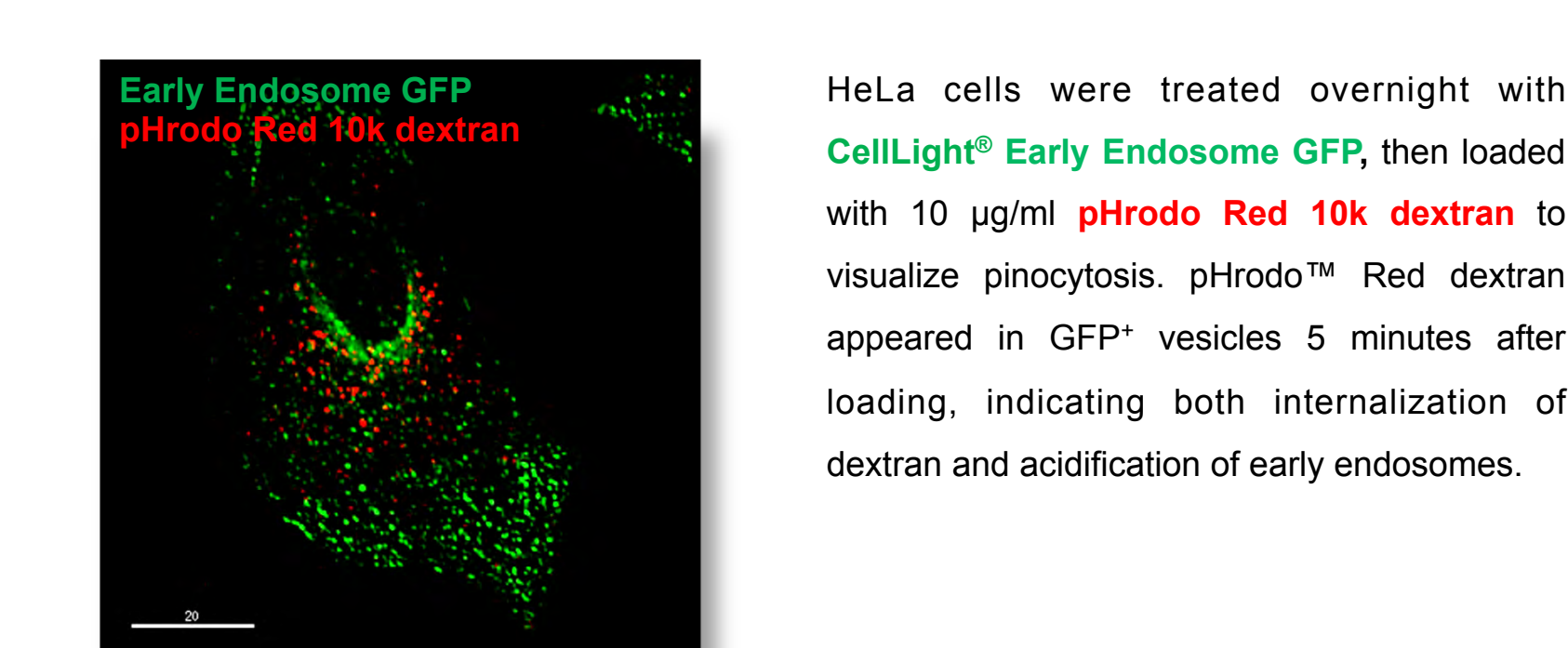


Figure 4. Fluorogenic Indication of EGF Endocytosis

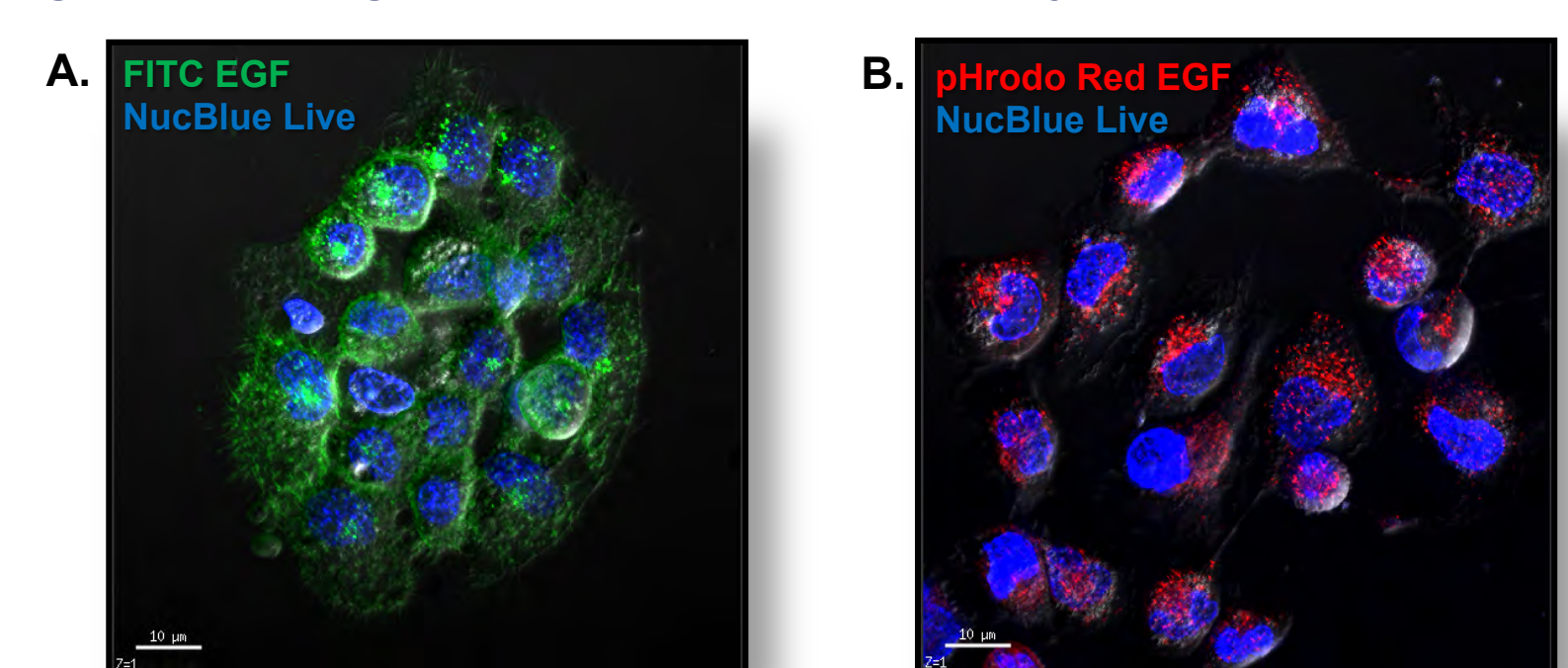
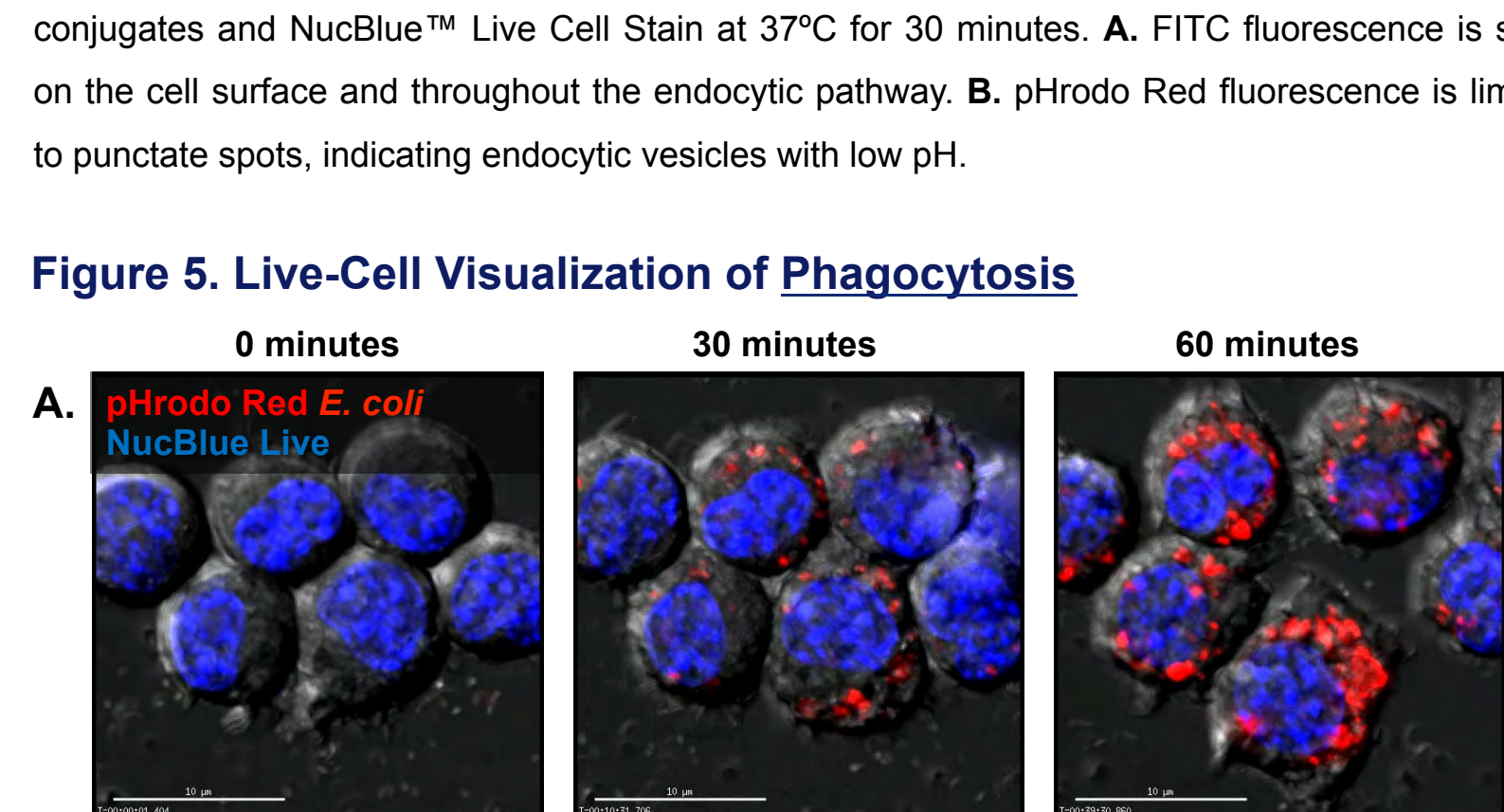
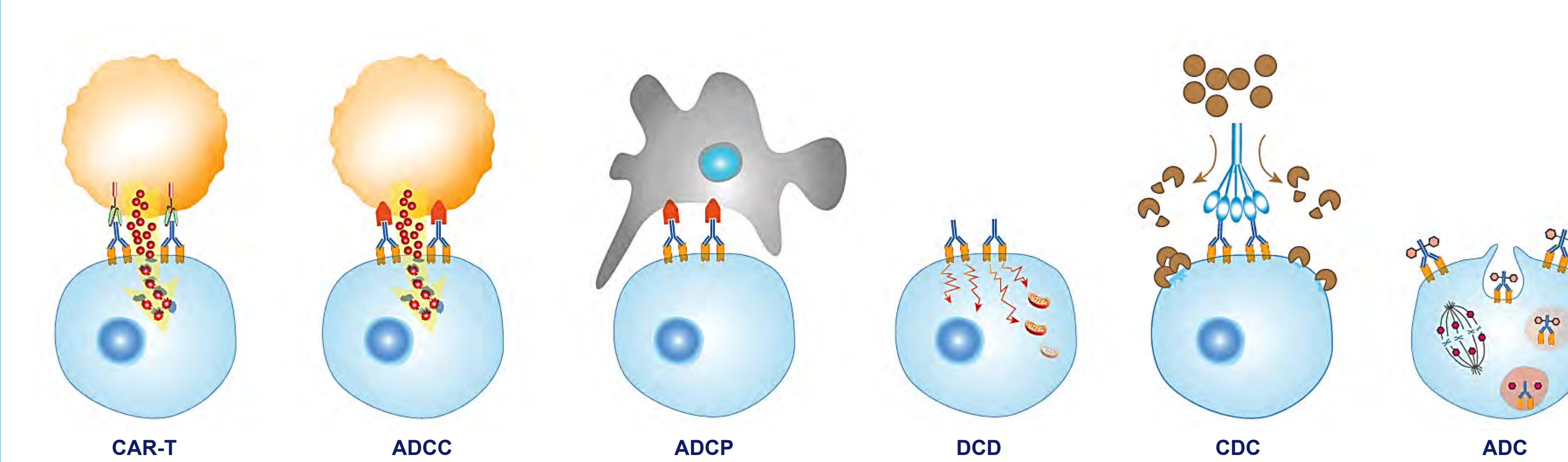


Figure 5. Live-Cell Visualization of Phagocytosis



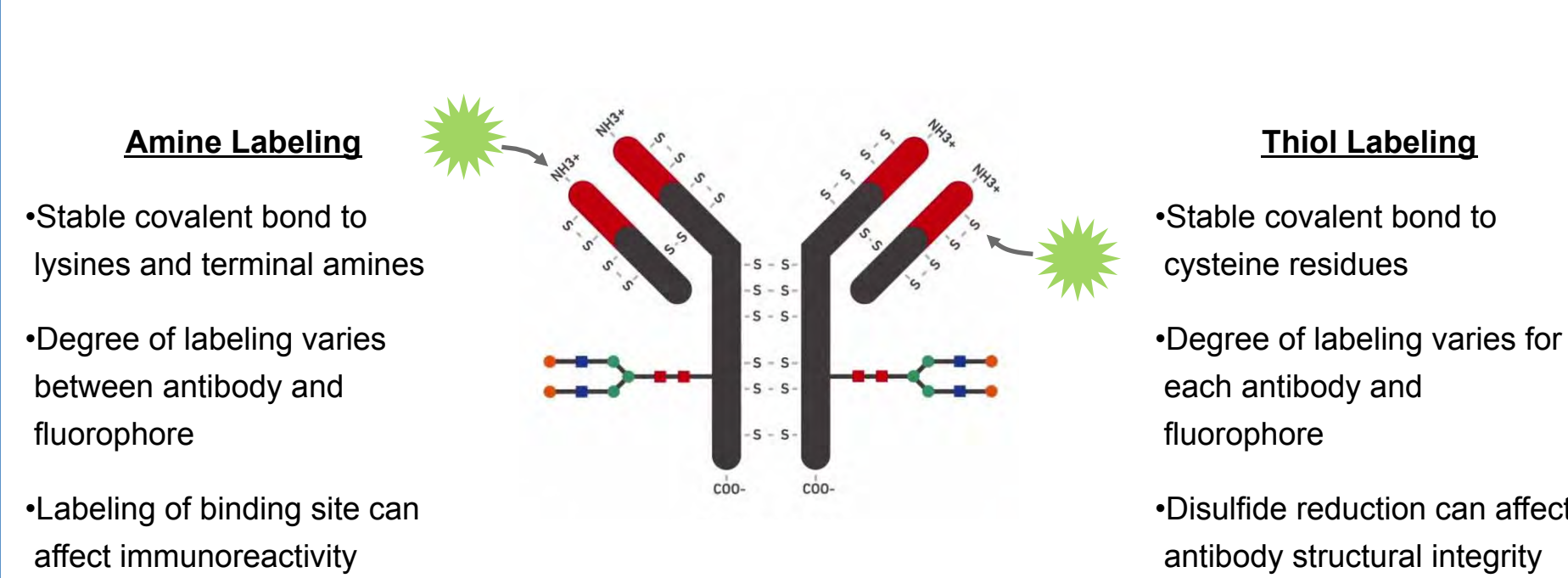
The dynamic pH-response of pHrodo dyes enable visualization of acidification processes such as the phagocytosis of bacteria by macrophages. 100 μg of pHrodo Red E. coli BioParticles® conjugates were added to MMM cells and imaged after 0, 30, and 60 minutes. The pHrodo BioParticles conjugates dramatically increase in fluorescence upon internalization and acidification.

Figure 6. Mechanisms of tumor cell killing by antibodies.



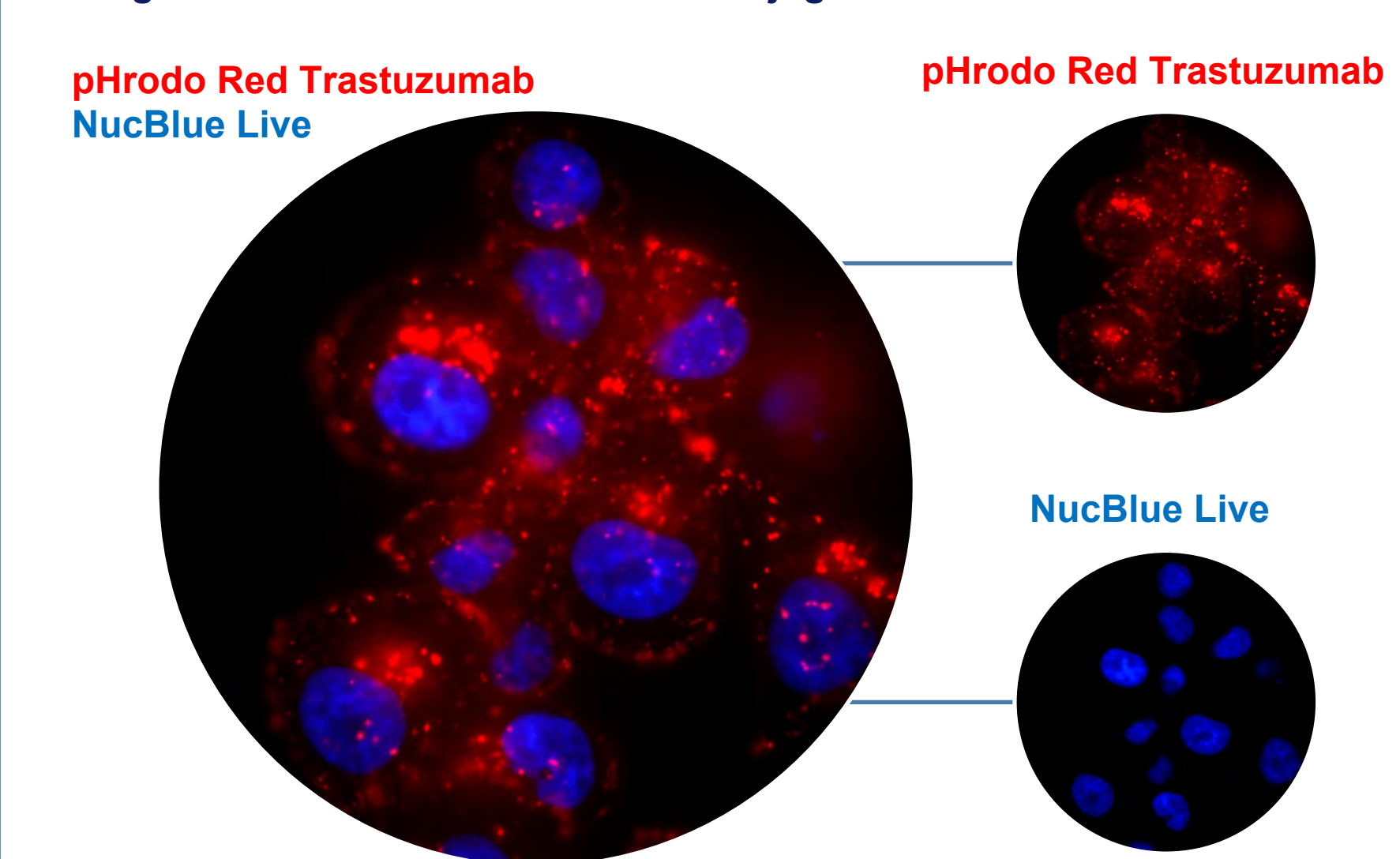
Antibodies participate in tumor cell killing by a variety of mechanisms, including: chimeric antigen receptor T cells (CAR-T), antibody-dependent cellular cytotoxicity (ADCC), antibody-dependent cellular phagocytosis (ADCP), direct cell death (DCD), complement-dependent cytotoxicity (CDC), or by the action of antibody-drug conjugates (ADC). Determination of antibody internalization kinetics is a critical step when developing therapeutic antibodies intended to participate in one of these mechanisms.

Figure 7. Conventional Antibody Labeling Strategies



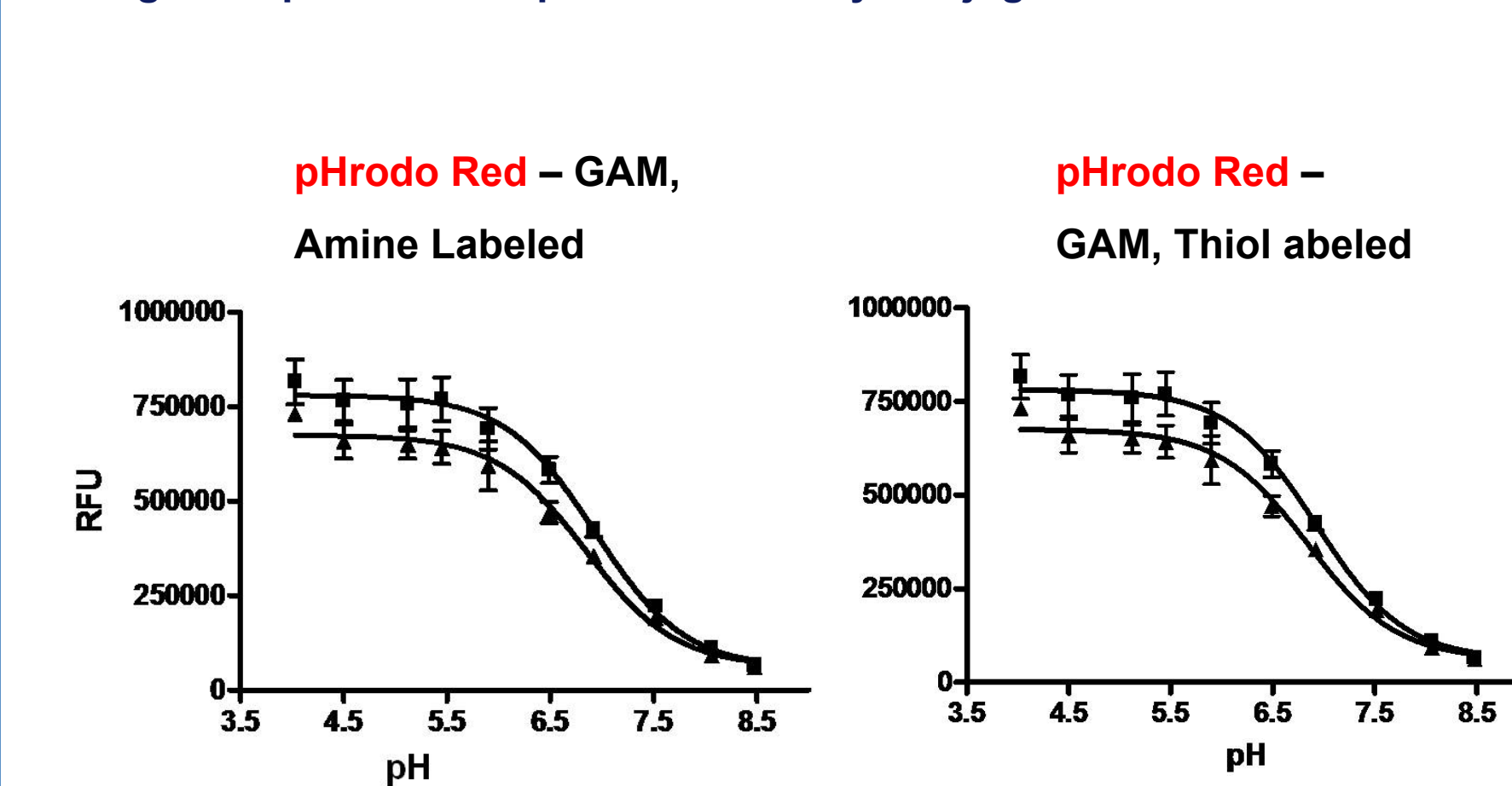
Amine- and thiol-reactive dyes are commonly used to label antibodies. However, the lack of specificity of these bioconjugation reactions can threaten immunoreactivity and lead to poorly defined constructs.

Figure 8. Internalization of amine-conjugated antibodies



Trastuzumab was labeled with amine-reactive pHrodo Red (DOL 2.8). Her2+ SK-BR-3 cells were plated in DMEM + 10% FBS in a 35 mm glass-bottom dish. Cells were treated with 10 nM pHrodo Red-labeled Trastuzumab for 16 hours at 37°C, 5% CO<sub>2</sub>. Media was removed and replaced with Live Cell Imaging Solution supplemented with 1% BSA and NucBlue Live Cell Stain for 30 minutes at 37°C. Images were captured on an EVOS FI Auto imaging system with 60x Olympus Plan Apochromat oil immersion objective and DAPI and TRITC light cubes. Red spots indicate internalization of trastuzumab into acidic vesicles.

Figure 9. pH-sensitive pHrodo-Antibody Conjugates



Antibodies were labeled with two reactive forms of pHrodo™ Red. A. First, the amine-reactive pHrodo™ Red, succinimidyl ester was used to label lysine residues on Goat anti Mouse IgG at molar ratios of 10 and 20. B. Second, the thiol-reactive pHrodo™ Red maleimide was used to specifically label cysteine residues of Goat anti Mouse IgG, with the goal of minimizing nonspecific and FC labeling. All conjugates were then tested for pH response and found to have the expected pKa of 6.8.

Figure 10. Fast, scalable screening of antibody internalization with pHrodo Fab fragments

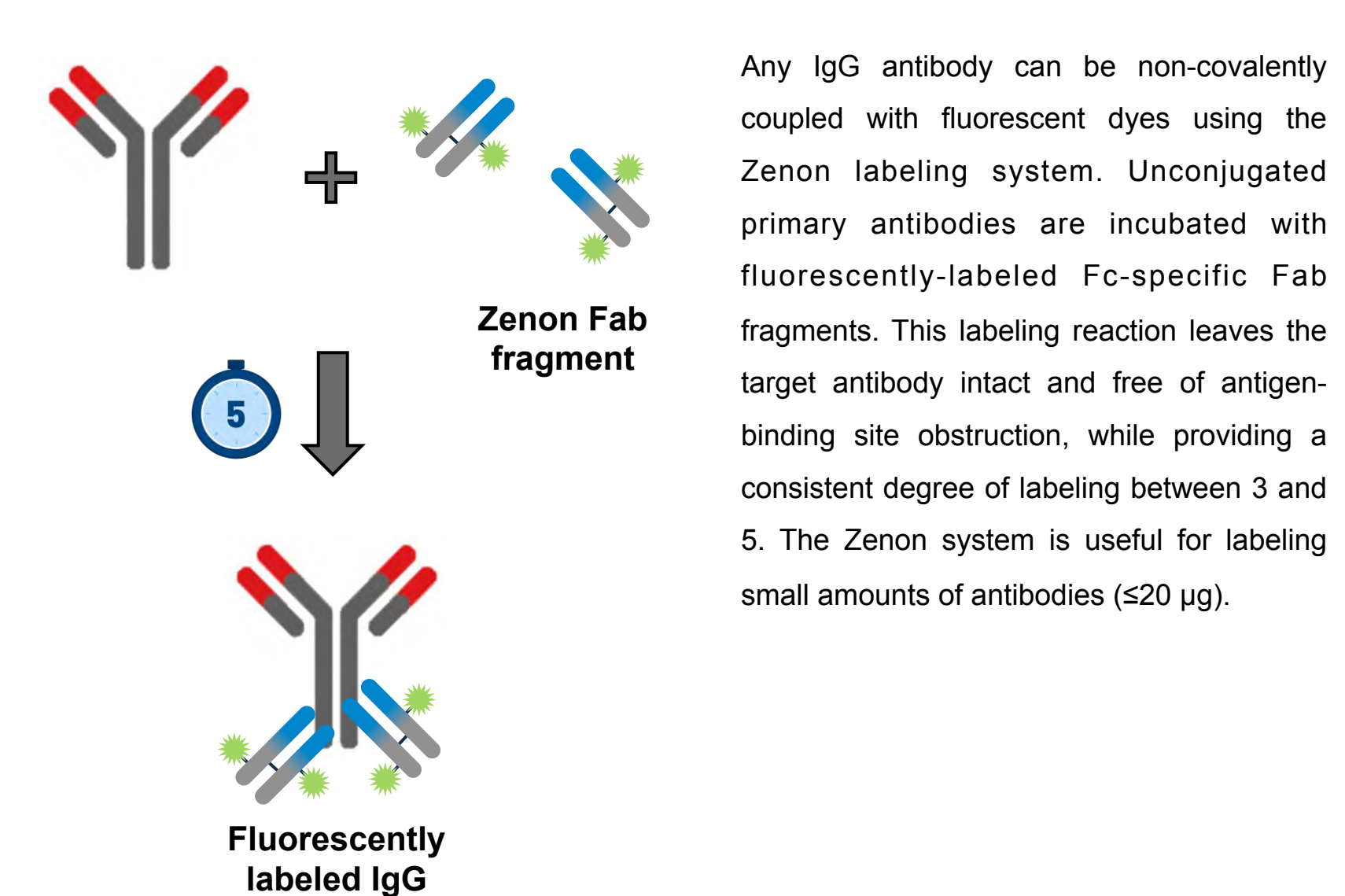


Figure 11. Trastuzumab labeled with pHrodo Red-Zenon is internalized and Trafficked to Lysosomes

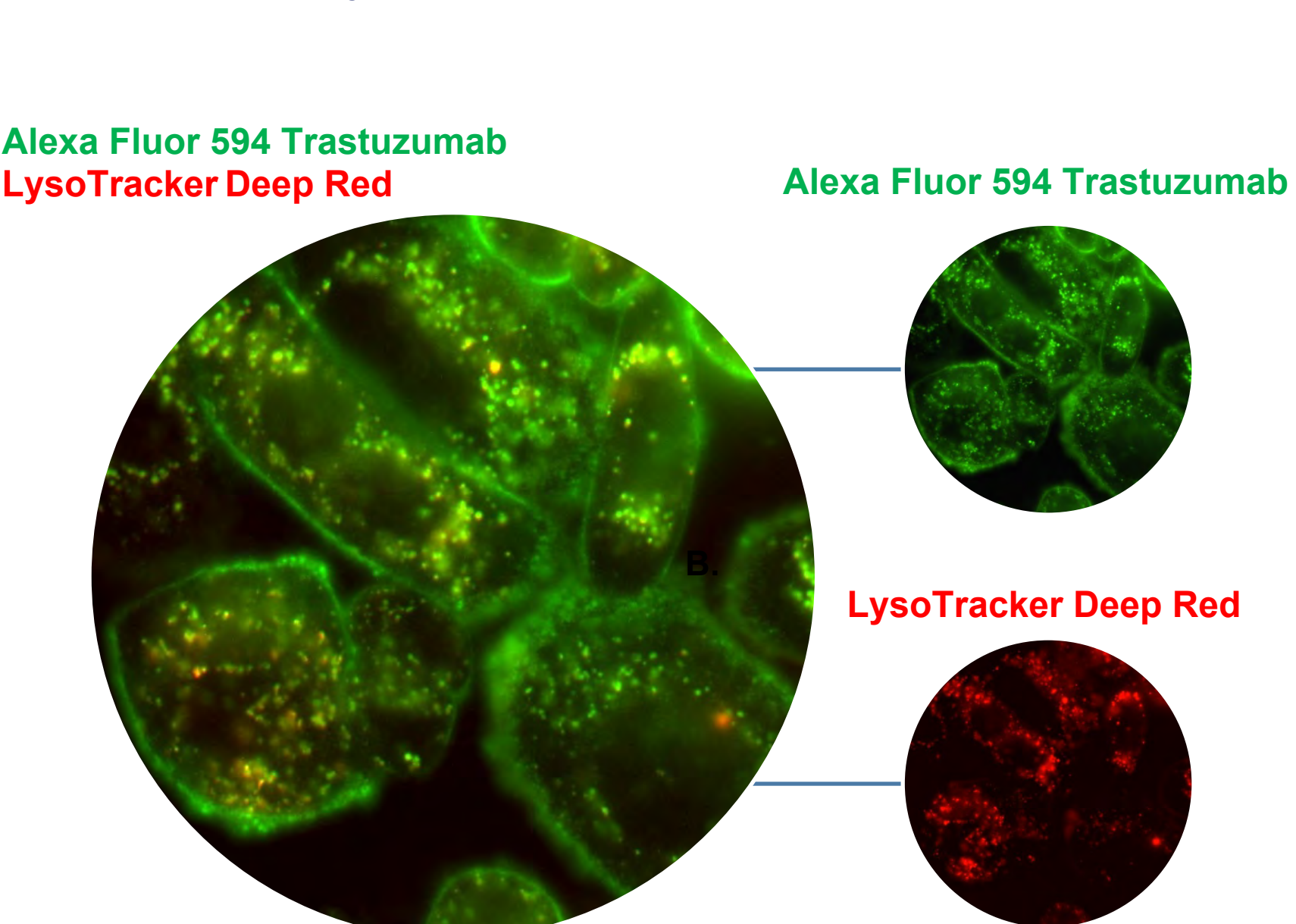
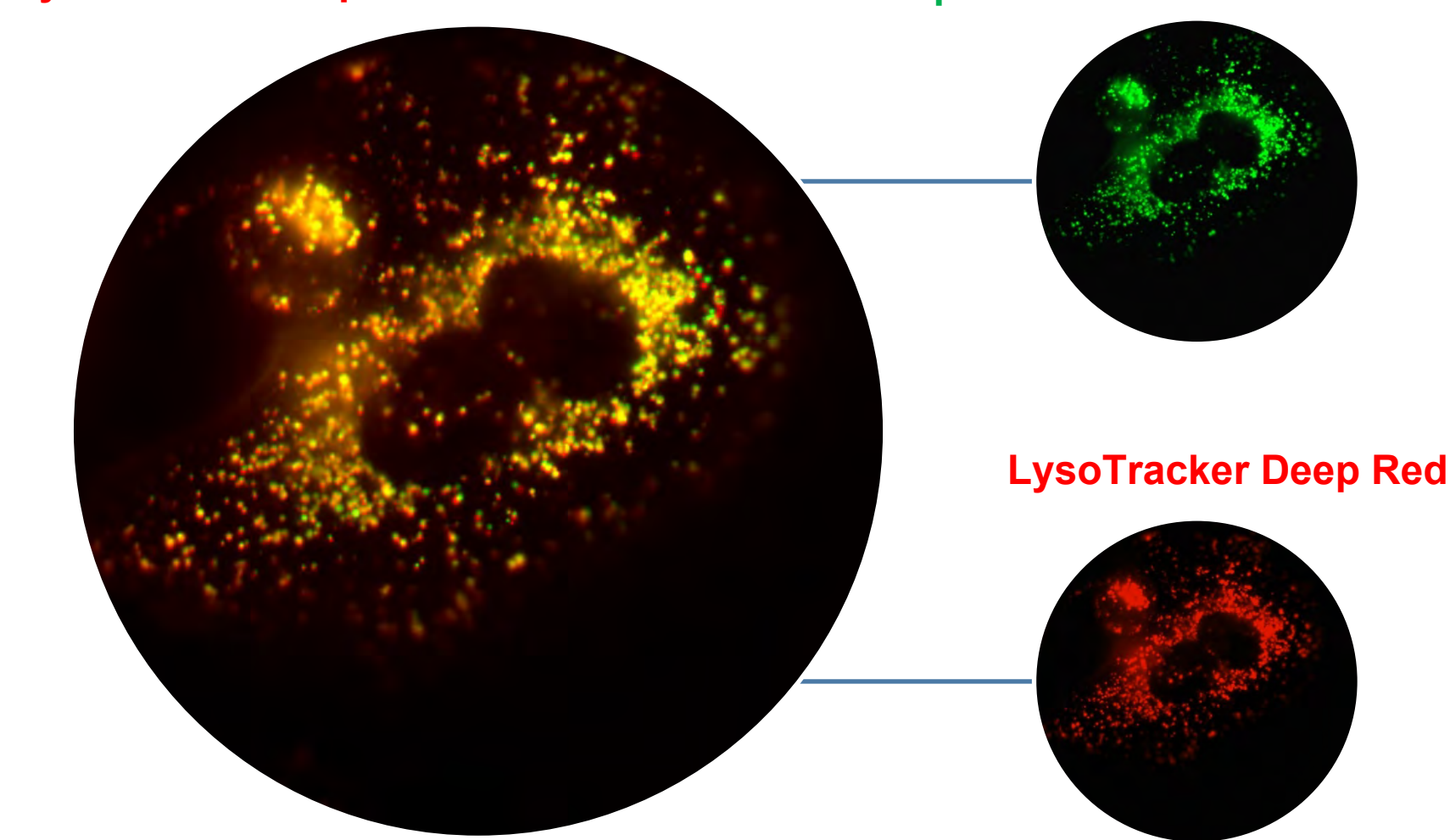


Figure 12. Site-specific labeling of heavy-chain N-linked glycans



Trastuzumab was azide-activated with the SiteClick labeling system, then click conjugated with pHrodo Red-DIBO. Cells were treated with 10 nM pHrodo Red-labeled Trastuzumab for 16 hours at 37°C, 5% CO<sub>2</sub>. Media was replaced with Live Cell Imaging Solution with 1% BSA, NucBlue Live Cell Stain, and 50 nM LysoTracker Deep Red for 30 minutes at 37°C. Red spots indicate internalization of trastuzumab into acidic intracellular vesicles that are positive for LysoTracker Deep Red.

Figure 13. SiteClick antibody labeling with pHrodo dye for site-specific labeling and antibody internalization

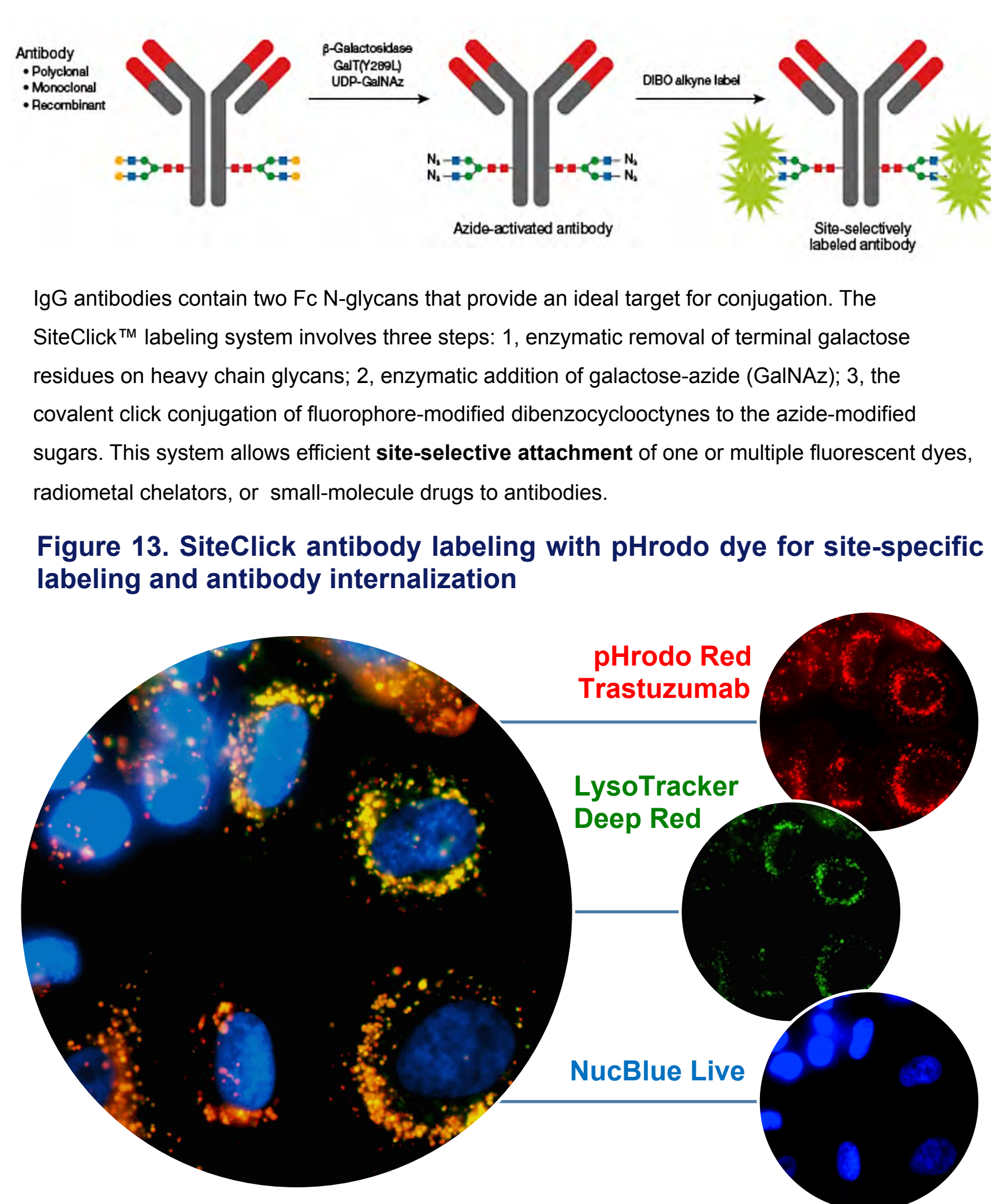


Figure 14. Direct Monitoring of ADC Internalization in Live Cells Using Dual SiteClick Alexa Fluor 647/MMAE Trastuzumab Conjugates

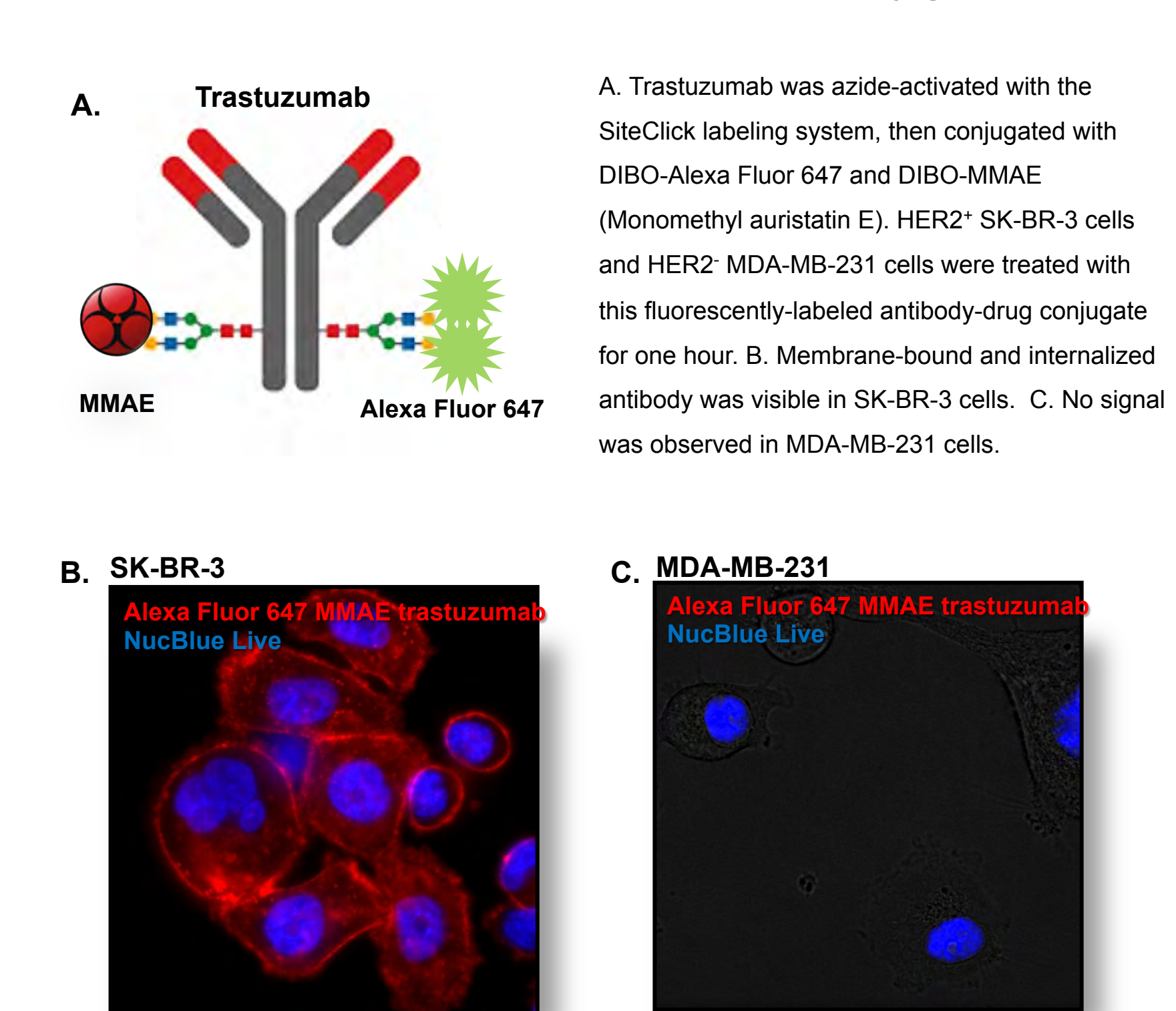
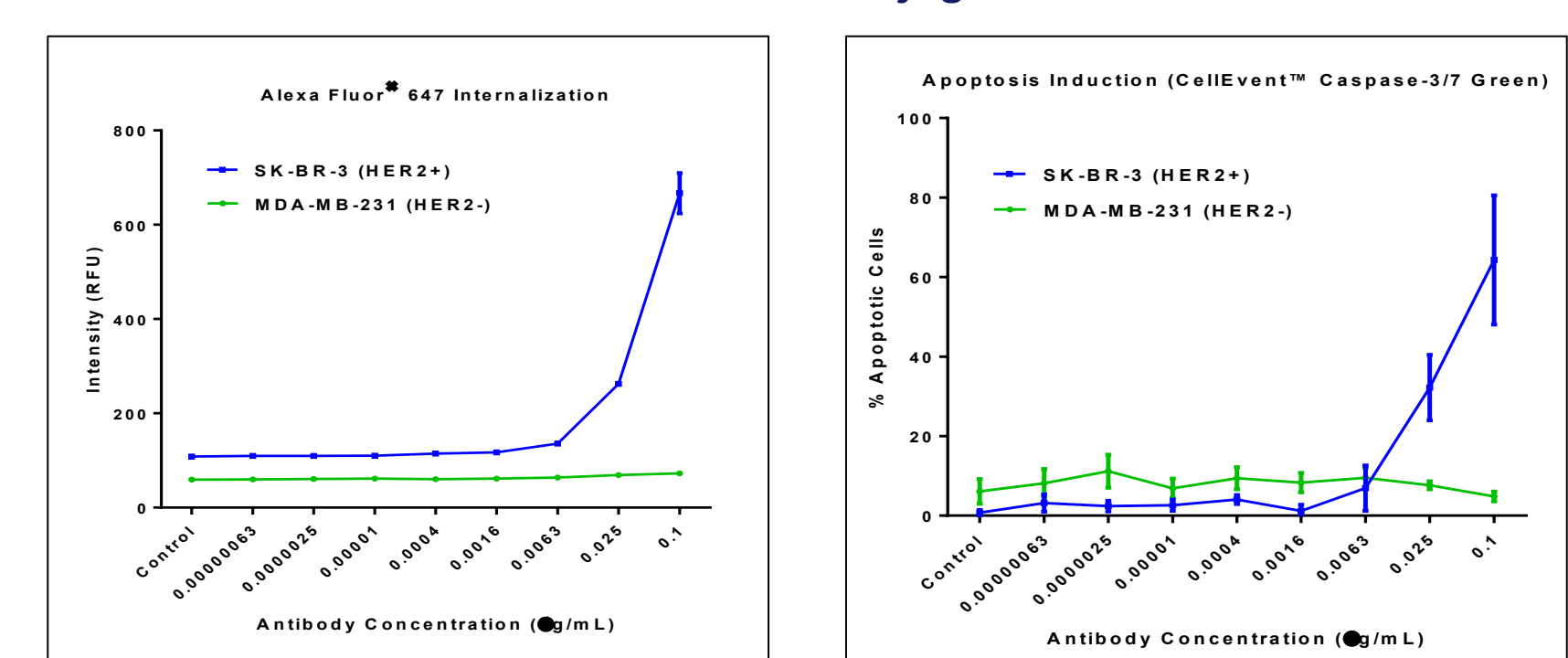


Figure 15. Specific Internalization and Cell Killing Response of Trastuzumab Alexa Fluor 647/MMAE Conjugates



## Conclusion

Studies of endocytosis, pinocytosis, and phagocytosis benefit from the ability to monitor pH changes in live cells with fluorogenic, pH-sensitive pHrodo dyes. Additionally, the internalization of antibodies can be analyzed using pHrodo dyes attached via three mechanisms.