

# AnaTag™ Protein LABELING KITS

Performance + Convenience

Bovine pulmonary artery endothelial cells probed with biotin-conjugated phalloidin and mouse anti-oxphos complex V; visualized with HiLyte Fluor™ 555 labeled Streptavidin (Cat# 60666), and HiLyte Fluor™ 488 labeled goat anti-mouse (Cat# 61057-H555). Nucleus stained with Hoechst 33342 (Cat# 83218).

Incorporating some of the industry's brightest and most photostable dyes in an ultra-fast, ultra-convenient kit, AnaTag™ Protein Labeling Kits are the perfect marriage of performance and convenience.

- Highest performance
- Easy to use
- Cost effective

## Superior Performance & Convenience

*Fast labeling - as short as 30 minutes*

*Available in two convenient reaction sizes:*

- Microscale kit: up to 200 ug protein x 3 labelings
- Larger scale kit: up to 5 mg protein x 3 labelings

*High Performance HiLytePlus™ & HiLyte Fluor™ Dyes*

- Protein conjugates brighter than Cy3 & Cy5 protein conjugates

*Classic Dyes*

- AMCA, FAM, FITC, ROX, TAMRA



ISO 9001:2000 Company

AnaTag™ Labeling Kits			
Number of Reactions			
Fluorophore	Abs / Em (nm / nm)	3 x 5 mg	3 x 200 ug
AMCA-X	353 / 442	72055	72056
5-FITC	494 / 519	72059	72060
5-FAM	495 / 520	72053	72054
HiLyte Fluor™ 488	499 / 523	72047	72048
5 TAMRA	547 / 574	72063	72064
HiLytePlus™ 555	552 / 567	72091	72092
HiLyte Fluor™ 555	553 / 568	72045	72046
5-ROX	573 / 602	72061	72062
HiLyte Fluor™ TR	591 / 622	72051 (3x1mg)	72052
HiLytePlus™ 647	649 / 669	72093	72094
HiLyte Fluor™ 647	652 / 669	72049	72050
HiLytePlus™ 750	751 / 779	72089	72090
HiLyte Fluor™ 750	754 / 778	72043	72044

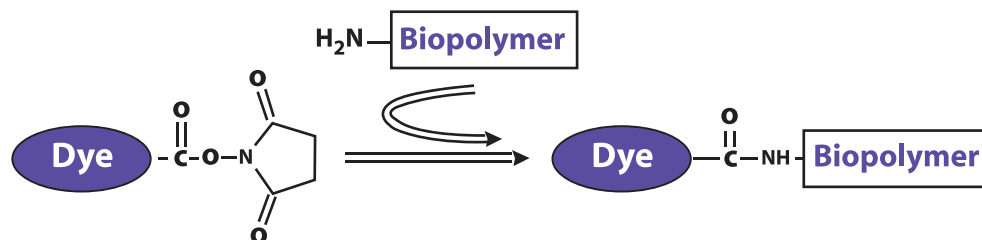


Figure 1. Labeling of an amino group (for instance, a lysine) on a biopolymer with a succinimidyl ester of a dye.

