

Product Name: HiLyte™ Fluor 405 acid, SE

Size: 1 mg

Catalog Number: AS-89317-1

Molecular Weight: 785.62

Solvent: DMF or DMSO

HiLyte™ Fluor 405, available in a variety of reactive forms, such as Product Description:

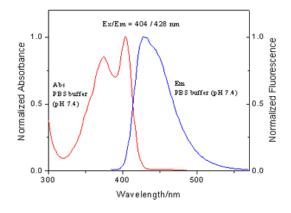
amine-reactive, thiol-reactive and amine-containing derivatives, allows for easy conjugation to biopolymers. HiLyte™ Fluor 405 has spectral characteristics similar to those of Alexa Fluor™ 405 (Invitrogen) and DyLight Fluor™ 405 (Pierce) dyes with excitation and emission wavelength at 404/428 nm in PBS buffer (pH 7.4). Conjugates of HiLyte™ Fluor 405 are well excited by the 407 nm spectral line of the Krypton laser and the 408 nm line of the violet laser diode. HiLyte™ Fluor 405 conjugates, whether conjugated to primary tag, secondary

antibodies or streptavidin, are prepared in an optimized

fluorophore/protein labeling ratio to ensure high fluorescent signal and uncompromised biological function. Fluorescence signal can be

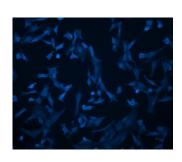
observed at Ex/Em = 404/428 nm.

Spectra: Maximum Ex/Em wavelength is 404/428 nm.



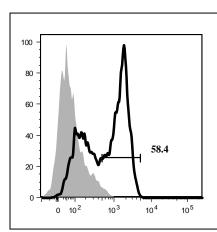
Application:

Immunofluorescence staining with HiLyte™ Fluor 405 conjugated goat Anti-mouse IgG



Alpha-tubulin in 3T3 cells probed with mouse anti -alpha tubulin antibody, visualized with HiLyte™ Fluor 405 conjugated goat anti-mouse IgG (cat# AS-28176-H405).

Flow cytometry analysis with HiLyte™ Fluor 405 labeled secondary antibody



Human PBMC cells were stained with anti-CD3 mAb (0.9 mg/ml) at 1:500 dilution. Secondary antibody goat antimouse IgG labeled with HiLyte™ Fluor 405 at 0.5 mg/ml (cat# AS-61057-H405) was used at 1:200 dilution. Filled histogram: secondary antibody staining only control.

Stock solution preparation: Make 10 mg/ml of dye stock solution (12.7 mM) in high quality anhydrous DMSO. Completely dissolve all dye content by vortexing. Dye solution should be prepared fresh immediately before an experiment. Extensive storage of the dye solution may reduce dye activity. Any solution containing the dye should be protected from light.

Shelf life and storage:

Shelf life is two years if stored at -20C, desiccated and protected from light.

Related Products:

Catalog#	Product Name
AS-89318	<u>HiLyte™ Fluor 405 amine TFA salt</u>
AS-89319	<u>HiLyte™ Fluor 405 hydrazide</u>
AS-89320	<u>HiLyte™ Fluor 405 C2 maleimide</u>
AS-72143	AnaTag™ HiLyte™ Fluor 405 Protein Labeling Kit *Ultra Convenient
AS-72142	AnaTag™ HiLyte™ Fluor 405 Microscale Protein Labeling Kit *Ultra Convenient
HiLyte™ Fluor 405 labeled Secondary Antibodies & Streptavidin	

This product is for *in vitro* research use only.