

ATTO 594 Goat anti-Rabbit IgG

Catalog No: 15047, 15067

Chemical Properties:

Contents: 250 μ l (Catalog No. 15047) or 35 μ l (Catalog No. 15067) of affinity purified goat anti-rabbit IgG (H+L) conjugated to ATTO 594 in PBS, pH 7.4, 0.01% sodium azide.

Specificity: This antibody was purified by immunoaffinity chromatography and cross-adsorbed. It reacts with whole molecule rabbit IgG and shows minimum cross-reactivity with bovine, chicken, goat, guinea pig, horse, human, mouse, rat and sheep serum proteins. No cross-reactivity was observed with non-immunoglobulin serum proteins.

Fluorescent Properties:

ATTO 594 is a new fluorescent label belonging to the class of Rhodamine dyes. Characteristic features of the label are strong absorption, high fluorescence quantum yield, high thermal and photo stability, excellent water solubility, and very little triplet formation.

Molar extinction Coefficient: 120,000 $M^{-1}cm^{-1}$ (measured at A_{max})

Quantum Yield: 85%

Excitation Max: 601 nm

Emission Max: 627 nm

Application:

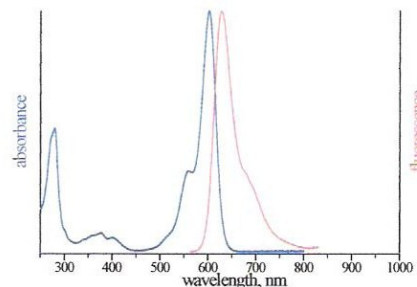
Immunofluorescence: 1:500 to 1:2000

Quality Control:

This antibody has been quality control-tested by spectra-photometrical evaluation and by immunocytochemistry.

Storage:

For short-term storage, the conjugated antibody should be stored at 4°C protected from light. This product is guaranteed for 6 months from the date of arrival.



Spectrum of ATTO 594 conjugated goat anti-rabbit IgG.



Immunofluorescence: HeLa cells stained with Histone H3 trimethyl Lys27 pAb (Catalog No. 39155) at 1:500 and the ATTO 594 Goat anti-rabbit secondary at 1:1,000. DAPI staining shown in blue.