

ATTO 647N (STED) Goat anti-Rabbit IgG

Catalog No: 15048, 15068

Chemical Properties:

Contents: 250 μ l (Catalog No. 15048) or 35 μ l (Catalog No. 15068) of affinity purified goat anti-rabbit IgG (H+L) conjugated to ATTO 647N (STED) 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 647N (STED) belongs to a new generation of fluorescent labels for the red spectral region. Characteristic features of the label are strong absorption, excellent fluorescence quantum yield, high photo stability, excellent ozone resistance, good solubility, and very little triplet formation.

Molar extinction Coefficient: 150,000 $M^{-1}cm^{-1}$ (measured at A_{max})

Quantum Yield: 65%

Excitation Max: 644 nm

Emission Max: 669 nm

Application:

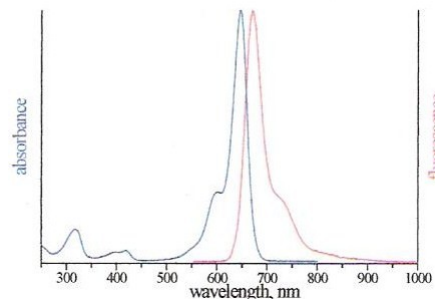
Immunofluorescence: 1:1000 to 1:2000

Quality Control:

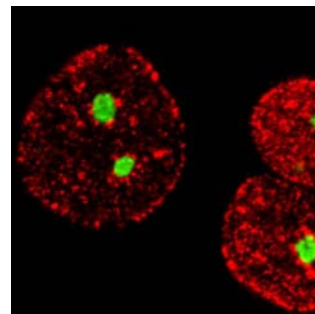
This antibody has been quality control-tested by spectra-photometrical evaluation and by immunocytochemistry.

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 647N conjugated goat anti-rabbit IgG.



Immunofluorescence: HeLa cells stained with Histone H3 trimethyl Lys27 pAb (Catalog No. 39155) at 1:500 dilution and the ATTO 647N (STED) Goat anti-rabbit secondary at 1:1,000. Green: Nucleolin antibody (Catalog No. 39541).