

ATTO 532 (GSD) Goat anti-Mouse IgG

Catalog No: 15070, 15071

Format: 250 µl, 35 µl

Chemical Properties:

Contents: 250 µl (Catalog No. 15070) or 35 µl (Catalog No. 15071) of affinity purified goat anti-mouse IgG (H+L) conjugated to ATTO 532 (GSD) 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 mouse IgG and shows minimum cross-reactivity with bovine, chicken, goat, guinea pig, horse, human, rabbit, rat and sheep serum proteins. No cross-reactivity was observed with non-immunoglobulin serum proteins.

Fluorescent Properties: ATTO 532 (GSD) belongs to a class of fluorescent labels for the orange-red spectral region, like Rhodamine 6G. The dye is very photostable, exhibits a strong absorption and an excellent fluorescence quantum yield. ATTO 532 (GSD) is very soluble in water, allowing conjugations with high dye-to-protein ratios.

Molar extinction Coefficient: 115,000 M⁻¹cm⁻¹ (measured at A_{max})

Excitation Max: 534 nm

Emission Max: 560 nm

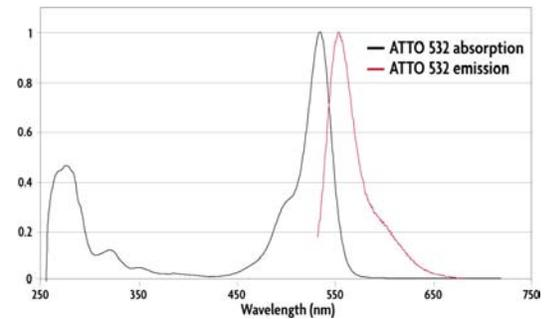
Application:

Immunofluorescence: 1:1000 to 1:2000

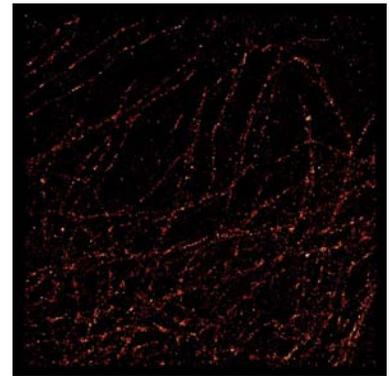
Quality Control: This antibody has been quality control-tested by spectrophotometrical evaluation and by immunocytochemistry.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Upon receipt, unconjugated antibodies may be stored at -20°C for up to 2 years. Fluorophore- & enzyme-conjugated antibodies should be stored at 4°C. Fluorophore-conjugated antibodies should be protected from light. Keep reagents on ice when not in storage; to avoid repeated freeze/thaw cycles, we recommend aliquoting items that will be stored frozen into single-use fractions prior to freezing. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



Absorption and emission spectra of ATTO 532 (GSD) Dye.



GSD fluorescence microscopy: Tubulin staining in HeLa cells using the ATTO 532 (GSD) Goat anti-mouse secondary. (Courtesy of Leica Microsystems, Germany.)