

## Rhodamine 6G (GSD) Goat anti-Mouse IgG

**Catalog No:** 15074, 15075

**Format:** 250 µl, 35 µl

### Chemical Properties:

**Contents:** 250 µl (Catalog No. 15074) or 35 µl (Catalog No. 15075) of affinity purified goat anti-mouse IgG (H+L) conjugated to Rhodamine 6G (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:** Rhodamine 6G (GSD) belongs to a class of fluorescent labels for the orange-red spectral region. Characteristic features of the label are strong absorption, excellent fluorescence quantum yield, and high photo stability.

**Molar extinction Coefficient:** 116,000 M<sup>-1</sup>cm<sup>-1</sup> (measured at A<sub>max</sub>)

**Excitation Max:** 508 nm

**Emission Max:** 558 nm

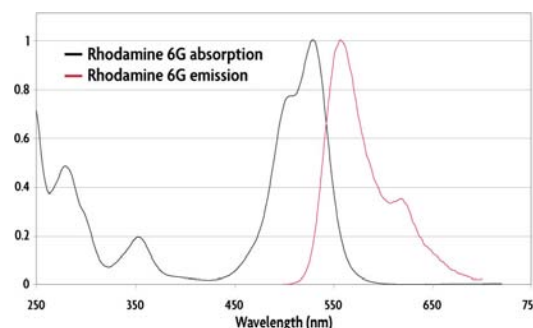
### Application:

Immunofluorescence: 1:1000 to 1:2000

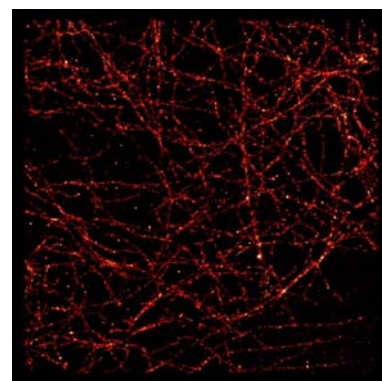
**Quality Control:** This antibody has been quality control-tested by spectrophotometrical evaluation and by immunocytochemistry.

**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 Rhodamine 6G (GSD) Dye.**



GSD fluorescence microscopy: Tubulin staining in HeLa cells using the Rhodamine 6G (GSD) Goat anti-mouse secondary. (Courtesy of Leica Microsystems, Germany.)