

## Chromeo™ 505 Goat anti-Mouse IgG

**Catalog No:** 15030, 15050

**Format:** 1 mg, 100 µg

### Chemical Properties:

**Contents:** 1 mg (Catalog No. 15030) or 0.1 mg (Catalog No. 15050) of Chromeo™ 505 conjugated Goat anti-Mouse IgG (H+L). The antibody concentration is 2 mg/ml in 0.01 M potassium phosphate, 0.15 M sodium chloride pH 7.4, containing 2 mM sodium azide.

**Specificity:** This antibody was purified by immunoaffinity chromatography. It reacts with whole molecular mouse IgG and the light chains of other mouse immunoglobulins. No cross-reactivity with non-immunoglobulin serum proteins was observed.

**Fluorescent Properties:** Chromeo 505 can be excited in a range from 480-515 nm, which enables excitation by a green laser. Even if the emission spectrum is shifted towards longer wavelength by some nm, filters designed for fluorescein can be used with Chromeo 505 conjugated antibodies. In contrast to fluorescein, Chromeo 505 exhibits a large tolerance to a broad pH range and high photostability. The Chromeo 505 conjugated antibodies have absorption and emission maxima of approximately 514 and 530 nm, which are compatible with common excitation sources and filter sets.

**Molar extinction Coefficient:** 70,000 M<sup>-1</sup>cm<sup>-1</sup> (measured at A<sub>max</sub>)

**Quantum Yield:** ~30%

**Excitation Wavelength Range:** 485 to 515 nm

**Emission Wavelength Range:** 525 to 540 nm

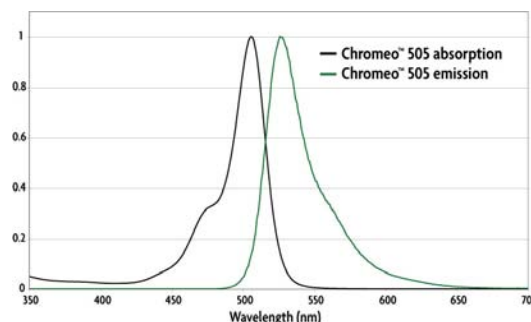
### Applications:

Immunofluorescence: 1:1000 to 1:2000 dilution

Plate-based assays: 1:800 to 1:1000 dilution

**Quality Control:** This antibody has been quality control-tested by spectrophotometrical evaluation, by immunohistochemistry (IHC) and by plate-based assays.

**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



**Absorption and emission spectra of Chromeo 505 Dye.**



**Chromeo 505 staining in HeLa cells.**

HeLa cells were stained with alpha Tubulin mouse mAb (Clone 5-B-1-2; Cat. No. 39527) and Chromeo 505 Goat anti-mouse IgG.