

Chromeo™ 494 Goat anti-Rabbit IgG

Catalog No: 15042, 15062

Format: 1 mg, 100 µg

Chemical Properties:

Contents: 1 mg (Catalog No. 15042) or 0.1 mg (Catalog No. 15062) of Chromeo™ 494 conjugated Goat anti-Rabbit 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 rabbit IgG and the light chains of other rabbit immunoglobulins. No cross-reactivity with non-immunoglobulin serum proteins was observed.

Fluorescent Properties: Chromeo 494 is a unique dye with special spectral properties. It is excitable with a green laser and has a large Stokes shift of 134 nm. Chromeo 494 conjugated antibodies exhibit superior luminescent properties and stability towards photobleaching, and have absorption and emission maxima of approximately 494 and 628 nm, making them ideal partners for other 488 nm excitable dyes and conjugates.

Molar extinction Coefficient: 55,000 M⁻¹cm⁻¹ (measured at A_{max})

Quantum Yield: ~7%

Excitation Wavelength Range: 470 to 520 nm

Emission Wavelength Range: 620 to 640 nm

Applications:

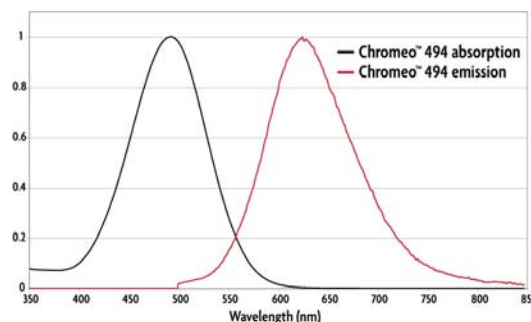
Immunofluorescence: 1:100 to 1:250 dilution

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

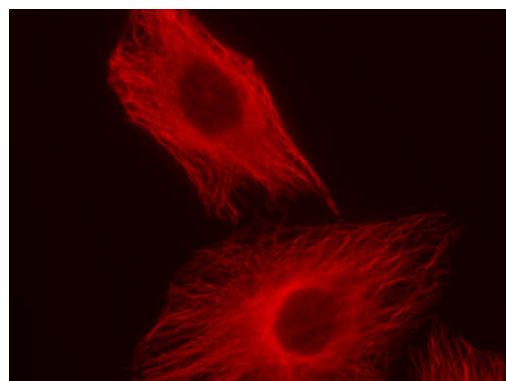
To ensure a maximum in photostability of Chromeo 494 conjugates under all experimental conditions, we recommend the use of Mowiol or TDE as mounting medium. Vectashield or MAXfluor Mounting medium might negatively influence the photostability of Chromeo 494.

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



Absorption and emission spectra of Chromeo 494 Dye.



HeLa cells stained with alpha Tubulin pAb and Chromeo 494 Goat anti-rabbit IgG.

HeLa cells were stained with alpha Tubulin rabbit pAb and Chromeo 494 Goat anti-rabbit IgG.