

Chromeo™ 488 Alkyne

Catalog No: 15290, 15291

Format: 1 mg, 5 x 1 mg

Chemical Properties: Contents: 1 mg (Catalog No. 15290) or 5 x 1 mg (Catalog No. 15291) of lyophilized Chromeo™ 488 Alkyne.

Net formula: $C_{33}H_{34}N_2O_6$; MW 554.64

Reagent color: yellow-orange

Soluble in DMF, MeOH, chloroform, DMSO and H₂O (when pre-dissolved in a small volume of DMF or DMSO).

Fluorescent Properties: Chromeo 488 can be excited with a green laser and shows similar fluorescent properties as fluorescein or other 488-excitable dyes. Filters designed for fluorescein can be used with Chromeo 488 and its conjugates.

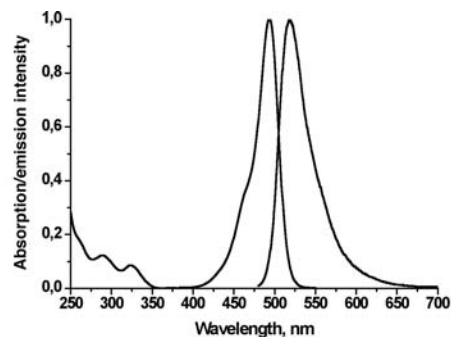
Molar extinction Coefficient: 73,000 M⁻¹cm⁻¹ (measured at A_{max})

Excitation Wavelength: 488 nm

Emission Wavelength: 517 nm

Quality Control: The Dye has been quality tested by TL, spectrophotometrical evaluation, Mass Spectroscopy and analytical HPLC.

Storage and Guarantee: To ensure stability, the lyophilized dye should be stored at -20°C in the dark. As the dye is moisture sensitive, it should be stored in the original foil pouch with desiccant. This product is guaranteed for 6 months from the date of arrival.



Spectrum of Chromeo™ 488 in PBS.