

Chromeo™ 488 Azide

Catalog No: 15280, 15281

Format: 1 mg, 5 x 1 mg

Chemical Properties: Contents: 1 mg (Catalog No. 15280) or 5 x 1 mg (Catalog No. 15281) of lyophilized Chromeo™ 488 Azide.

Net formula: $C_{33}H_{37}N_5O_6$; MW 599.69

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})

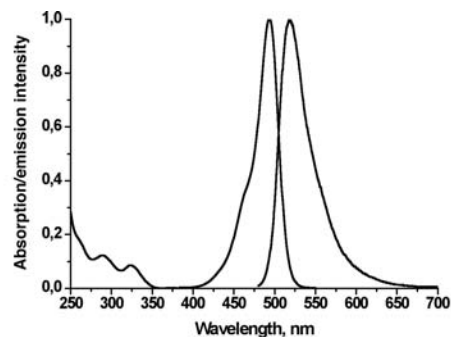
Excitation Wavelength: 488 nm

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. In the absence of an express written agreement to the contrary, all products are sold by Active Motif for research purposes only and the exclusive use of the original purchaser, and are not to be sold.



Spectrum of Chromeo™ 488 in PBS.