

Chromeo™ 494 Azide

Catalog No: 15282, 15283

Format: 1 mg, 5 x 1 mg

Chemical Properties: Contents: 1 mg (Catalog No. 15282) or 5 x 1 mg (Catalog No. 15283) of lyophilized Chromeo™ 494 Azide.

Net formula: $C_{29}H_{37}BrN_6O_3$ MW 597.56

Reagent color: light-red

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

Fluorescent Properties: Chromeo 494 can be excited with a green laser. As the dye shows a large Stokes shift of 134 nm, the emission should be measured at a wavelength of 630 nm. Its fluorescent properties make Chromeo 494 an ideal partner for multiplexing with other 488-excitable dyes and conjugates.

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

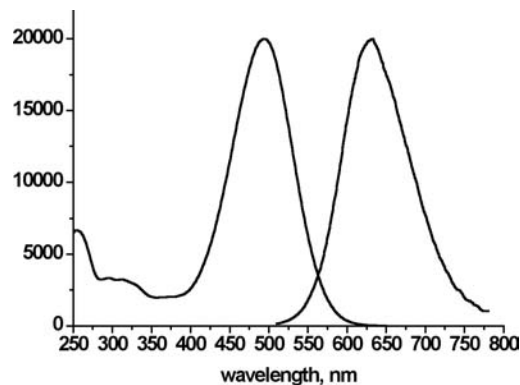
Excitation Wavelength: 494 nm

Emission Wavelength: 628 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™ 494 in PBS.