

Chromeo™ 494 Carboxylic Acid

Catalog No: 15110, 16110

Format: 1 mg, 5 x 1 mg

Chemical Properties: Contents: 1 mg (Catalog No. 15110) or 5 x 1 mg (Catalog No. 16110) of lyophilized Chromeo™ 494 carboxylic acid

Net formula: $C_{26}H_{32}BrN_2O_4$; MW 516.45

Reagent color: light red.

Soluble in DMF, ethanol and methanol.

Fluorescent Properties: Chromeo 494 can be excited with a green laser. As the dye exhibits 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.

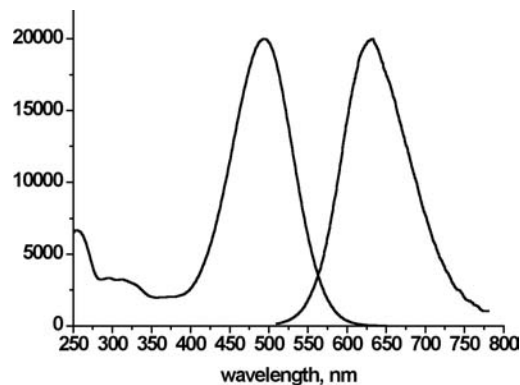
Molar extinction Coefficient: $55,000 \text{ M}^{-1} \text{ cm}^{-1}$ (measured at A_{max})

Excitation Wavelength: 494 nm

Emission Wavelength: 628 nm

Quality Control: The Dye has been quality tested by TLC and spectrophotometrical evaluation.

Storage and Guarantee: To ensure stability, the lyophilized dye should be stored at 4°C in the dark. This product is guaranteed for 6 months from the date of arrival.



Spectrum of Chromeo™ 494 in PBS.