

Chromeo™ 505 Carboxylic Acid

Catalog No: 15610, 16610

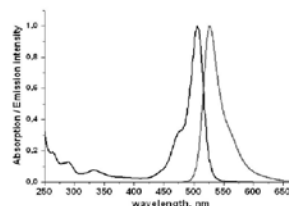
Format: 1 mg, 5 x 1 mg

Chemical Properties: Contents: 1 mg (Catalog No. 15610) or 5 x 1 mg (Catalog No. 16610) of lyophilized Chromeo™ 505 Carboxylic Acid

Net formula: $C_{21}H_{12}Cl_2O_5$; MW 415.23

Reagent color: yellow-orange

Soluble in DMF, ethanol and methanol.



Spectrum of Chromeo™ 505 in PBS.

Fluorescent Properties: Chromeo 505 can be excited in a range from 480 -505 nm which includes excitation by a green laser. Even if the emission spectrum is shifted towards longer wavelength by some nm, Filters designed for fluorescein can be used with Chromeo 505 and its conjugates. In contrast to fluorescein, Chromeo 505 exhibits a large tolerance to a broad pH range and high photostability.

Molar extinction Coefficient: $70,000 \text{ M}^{-1}\text{cm}^{-1}$ (measured at A_{max})

Excitation Wavelength: 505 nm

Emission Wavelength: 526 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.