

## Chromeo<sup>™</sup> 488 Carboxylic Acid

## **Catalog No: 15510, 16510 Format:** 1 mg, 5 x 1 mg

## **Chemical Properties:**

Contents: 1 mg (Catalog No. 15510) or 5 x 1 mg (Catalog No. 16510) of lyophilized Chromeo<sup>™</sup> 488 Carboxylic Acid and 1 ml MAXfluor<sup>™</sup> Mounting Medium.

## Chromeo™ 488 Carboxylic Acid

Net formula:  $C_{30}H_{31}NO_7$ ; MW 517.58 Reagent color: yellow-orange Soluble in DMF, ethanol and methanol.

**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. Chromeo 488 exhibits a large tolerance to pH and high photostability, which provides more time for image capture.

Molar extinction Coefficient: 73,000 M<sup>-1</sup>cm<sup>-1</sup> (measured at A<sub>max</sub>)

Excitation Wavelength: 488 nm

Emission Wavelength: 520 nm

To ensure a maximum in photostability of Chromeo 488 conjugates under all experimental conditions, we recommend MAXfluor Mounting medium in fluorescence microscopy experiments. MAXfluor Mounting Medium is a non-hardening, glycerol-based, aqueous mounting medium which provides optimal fluorescent stability and inhibits photobleaching. The effectiveness of MAXfluor Mounting Medium may depend on which fluorescent dye it is used with. MAXfluor Mounting Medium is designed to be dispersed over an entire coverslip. Recommended usage is 15-20 µl per 22 mm square coverslip. The coverslips may be sealed with nail polish or other sealants for long-term storage.

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

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



Spectrum of Chromeo<sup>™</sup> 488 in PBS.