

Chromeo™ 642 Carboxylic Acid

Catalog No: 15310, 16310

Format: 1 mg, 5 x 1 mg

Chemical Properties: Contents: 1 mg (Catalog No. 15310) or 5 x 1 mg (Catalog No. 16310) of lyophilized Chromeo™ 642 carboxylic acid

Net formula: $C_{35}H_{45}N_2O_5P$; MW 604,73; m.p. >300 °C;

Reagent color: blue solid

Soluble in DMF, ethanol and methanol.

Fluorescent Properties: Chromeo 642 shows similar fluorescent properties as the red-fluorescent Cy5 dye. Filters designed for Cy5 can be used with Chromeo 642 and its conjugates. Chromeo 642 exhibits a large tolerance to pH and unmatched photostability, which provides more time for image capture.

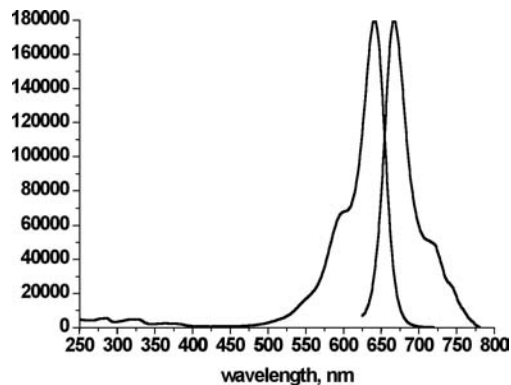
Molar extinction Coefficient: $180,000 \text{ M}^{-1}\text{cm}^{-1}$ (measured at A_{max})

Excitation Wavelength: 642 nm

Emission Wavelength: 660 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™ 642 in PBS.