

Chromeo™ 642 Azide

Catalog No: 15288, 15289

Format: 1 mg, 5 x 1 mg

Chemical Properties: Contents: 1 mg (Catalog No. 15288) or 5 x 1 mg (Catalog No. 15289) of lyophilized Chromeo™ 642-Azide.

Net formula: C₃₈H₅₂N₆O₄P; MW 686.83

Reagent color: blue solid

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

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 M⁻¹cm⁻¹ (measured at A_{max})

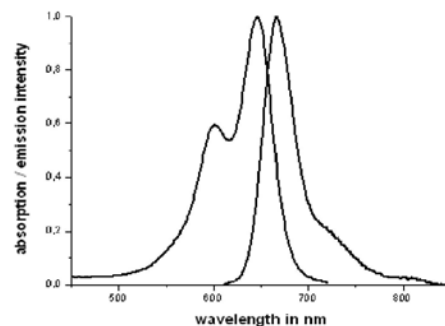
Excitation Wavelength: 646 nm

Emission Wavelength: 666.5 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™ 642 Azide in PBS.