

Chromeo™ 546 Alkyne

Catalog No: 15294, 15295

Format: 1 mg, 5 x 1 mg

Chemical Properties: Contents: 1 mg (Catalog No. 15294) or 5 x 1 mg (Catalog No. 15295) of lyophilized Chromeo™ 546 Alkyne.

Net formula: $C_{38}H_{51}N_3O_4P$ MW 644.81

Reagent color: purple solid.

Soluble in DMF, MeOH, acetonitrile, ethanol and H_2O (when pre-dissolved in a small volume of DMF or acetonitrile).

Fluorescent Properties: Chromeo 546 shows similar fluorescent properties as the yellow-orange Cy3 dye. Filters designed for Cy3 can be used with Chromeo 546 and its conjugates. Chromeo 546 exhibits a large tolerance to pH and unmatched photostability, which provides more time for image capture.

Molar extinction Coefficient: $96,800 M^{-1} cm^{-1}$ (measured at A_{max})

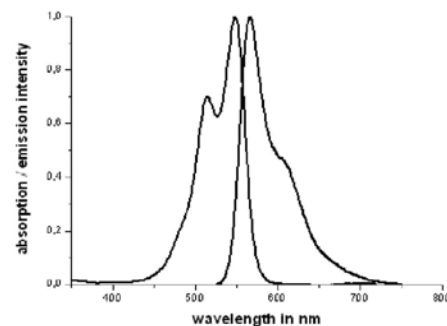
Excitation Wavelength: 550 nm

Emission Wavelength: 567 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 $-20^{\circ}C$ 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™ 546 Azide in PBS.