

520-Bodipy Alkyne

Catalog No: 15318, 16318

Format: 1 mg, 5 x 1 mg

Chemical Properties: Contents: 1 mg (Catalog No. 15318) or 5 x 1 mg (Catalog No. 16318) of lyophilized 520-Bodipy Alkyne.

Chemical name: 8-(4-Propynyloxyphenyl)-2,6-diethyl-1,3,5,7-tetramethyl-4,4-difluoro-4-bora-3a,4a-diaza-s-indacene

Net formula: $C_{26}H_{29}BF_2N_2O$; MW 434.2

Reagent color: red

Soluble in DMF, MeOH, DMSO and water (when pre-dissolved in DMF or DMSO).

Fluorescent Properties: 520-Bodipy Alkyne can be excited between 510 and 525 nm. Fluorescence can be detected between 530 and 550 nm.

Molar extinction Coefficient: $63,000 \text{ M}^{-1} \text{ cm}^{-1}$ (measured at A_{max})

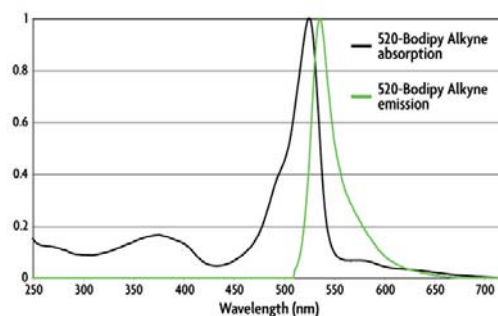
Excitation Maximum: 522 nm

Emission Maximum: 534 nm

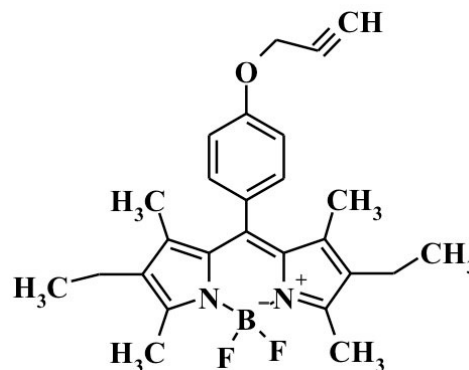
Quality Control: The Dye has been quality tested by TLC, spectrophotometrical evaluation, Mass Spectroscopy and analytical HPLC.

Storage and Guarantee: To ensure stability, the lyophilized dye should be stored at 4°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.



Absorption and Emission Spectrum of 520-Bodipy Alkyne in methanol.



Structure of 520-Bodipy Alkyne.