

550-Red Oxazine Azide

Catalog No: 15319, 16319

Format: 1 mg, 5 x 1 mg

Chemical Properties: Contents: 1 mg (Catalog No. 15319) or 5 x 1 mg (Catalog No. 16319) of lyophilized 550-Red Oxazine Azide.

Chemical name: 1-(3-azido-propoxy)-7-dimethylamino-phenoxazine-3-one

Net formula: $C_{17}H_{17}N_5O_3$; MW 339

Reagent color: dark-blue/ black

Soluble in $CHCl_3$, MeOH and water (when pre-dissolved).

Fluorescent Properties: 550-Red Oxazine Alkyne can be excited between 520 and 580 nm. Because of the broad emission spectrum fluorescence can be detected between 610 and 650 nm.

Molar extinction Coefficient: $70,000 \text{ M}^{-1} \text{ cm}^{-1}$ (measured at A_{max})

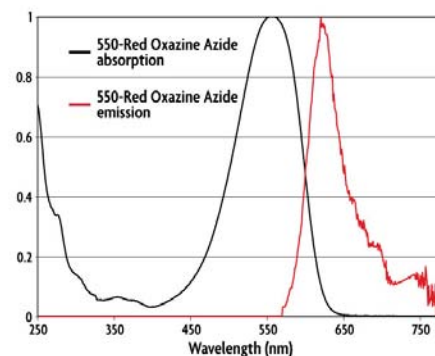
Excitation Maximum: 557 nm

Emission Maximum: 621 nm

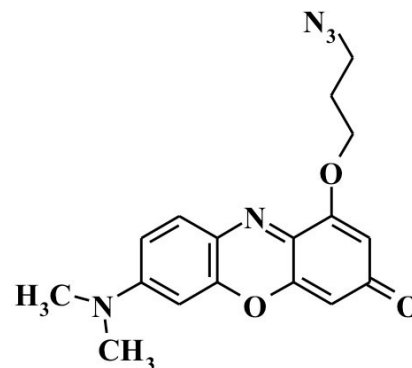
Quality Control: The Dye has been quality tested by TLC, spectrophotometrical evaluation, Mass Spectroscopy and analytical HPLC.

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 550-Red Oxazine Azide in methanol.



Structure of 550-Red Oxazine Azide.