

625-Nile Blue Azide

Catalog No: 15415, 16415

Format: 1 mg, 5 x 1 mg

Chemical Properties: Contents: 1 mg (Catalog No. 15415) or 5 x 1 mg (Catalog No. 16415) of lyophilized 625-Nile Blue Azide.

Chemical name: N-(5-(3-azidopropylamino)-9H-benzo(a)-phenoxazin-9-ylidene)-N-methyl-methanaminium chloride

Net formula: $C_{21}H_{22}ClN_6O$; MW 409.6

Reagent color: dark-blue

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

Fluorescent Properties: 625-Nile Blue Azide can be excited between 600 and 635 nm. Fluorescence can be detected between 650 and 680 nm.

Molar extinction Coefficient: $18,000 \text{ M}^{-1} \text{ cm}^{-1}$ (measured at A_{max})

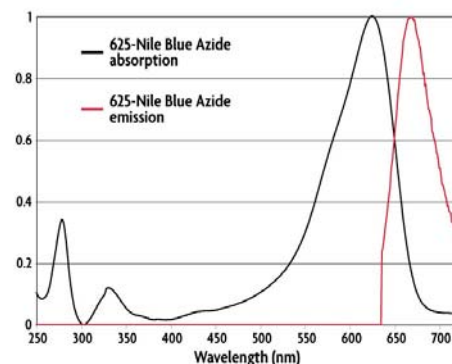
Excitation Maximum: 625 nm

Emission Maximum: 669.5 nm

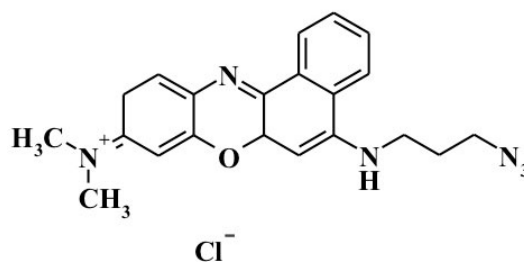
Quality Control: The Dye has been quality tested by TLC, spectrophotometrical evaluation, Mass Spectroscopy and analytical HPLC.

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 625-Nile Blue Azide in methanol.



Structure of 625-Nile Blue Azide.