

## 651-Blue Oxazine Azide

**Catalog No:** 15417, 16417

**Format:** 1 mg, 5 x 1 mg

**Chemical Properties:** Contents: 1 mg (Catalog No. 15417) or 5 x 1 mg (Catalog No. 16417) of lyophilized 651-Blue Oxazine Azide.

Chemical name: (9-(3-azido-propoxy)-7-piperidin-1-yl-phenoxazin-3-ylidene)-dimethyl-ammonium perchlorate

Net formula:  $C_{22}H_{27}ClN_6O_6$ ; MW 506

Reagent color: dark-red

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

**Fluorescent Properties:** 651-Blue Oxazine Azide can be excited between 630 and 650 nm. Fluorescence can be detected between 660 and 680 nm.

**Molar extinction Coefficient:**  $73,000 \text{ M}^{-1}\text{cm}^{-1}$  (measured at  $A_{\text{max}}$ )

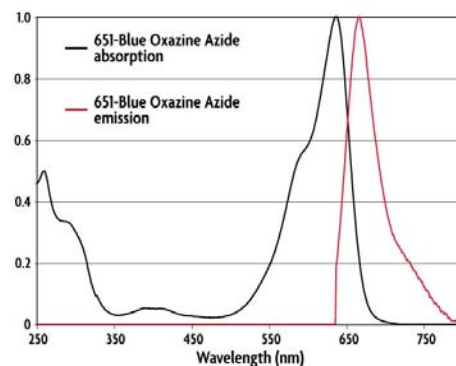
**Excitation Maximum:** 637 nm

**Emission Maximum:** 666.5 nm

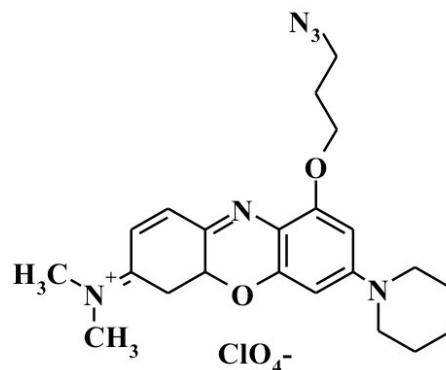
**Quality Control:** The Dye has been quality tested by TLC, spectrophotometrical evaluation, Mass Spectroscopy and analytical HPLC.

**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 651-Blue Oxazine Azide in methanol.



Structure of 651-Blue Oxazine Azide.