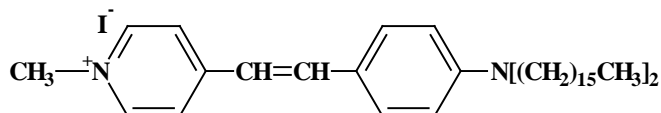


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** DiA (4-(4-dihexadecylaminostyryl)-N-methylpyridinium iodide)**CATALOG #:** 60013**MOLECULAR INFORMATION:** C<sub>46</sub>H<sub>79</sub>IN<sub>2</sub>  
Mwt: 787**PROPERTIES:**

<b>Color &amp; Form</b>	Red solid
<b>Purity</b>	≥ 95% by HPLC
<b>Solubility</b>	Soluble in DMSO and EtOH;
<b>Absorption\ Emission</b>	491 nm \ 613 nm (in MeOH)
<b>Extinction Coefficient</b>	52,000 cm <sup>-1</sup> M <sup>-1</sup> (in MeOH)

**STORAGE AND HANDLING:** Store at ≤4 °C and protect from light, especially in solution.**APPLICATION:** DiA is a green fluorescent membrane dye that diffuses much faster than DiO in cell membranes. Thus, DiA and DiI (**60010**) have been used together for two color membrane staining.Ref: 1) *Trends Neurosci.* **13**, 14(1990); 2) *J. Neurosci.* **12**, 3494(1992); 3) *J. Neurosci.* **15**, 990(1995).**TOXICITY:** Unknown

<b>FIRST AID:</b>	Potentially harmful; avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

<b>Disclaimer:</b> <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--