

IVISense™ 645 MAL Fluorescent Dye (5mg)

Product Number: NEV11274

DESCRIPTION

IVISense™ 645 MAL Fluorescent Dye is a thiol (-SH, sulfhydryl) reactive maleimide-containing red fluorochrome for molecular labeling applications. IVISense™ 645 MAL Fluorescent Dye enables optimized and integrated imaging applications in both in vivo and in vitro applications. IVISense™ 645 MAL Fluorescent Dye's superior brightness makes it particularly well suited for the labeling of thiol and cysteine containing molecules for a wide range of imaging applications. The 645 nm fluorescence wavelength is ideal for biological imaging applications on in vitro and in vivo fluorescence systems, including advanced microscopes, high content imaging and flow cytometry systems.

MATERIAL

Available as a solid in 1 mg or 5 mg vials.

STORAGE & HANDLING

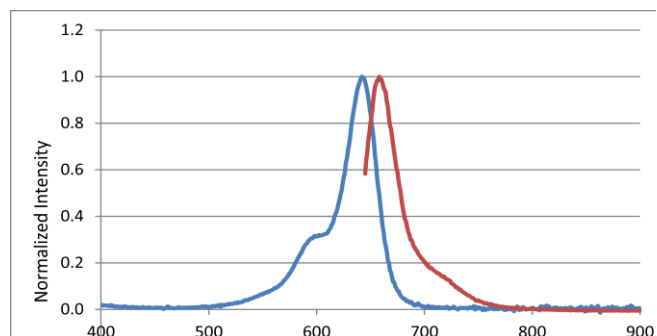
- Upon receipt, IVISense™ 645 MAL Fluorescent Dye should be IMMEDIATELY STORED AT -20 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ 645 MAL Fluorescent Dye is stable for up to 6 months from the date of shipment.
- Allow to equilibrate to room temperature before opening.

IN VIVO IMAGING

- Red fluorophore optimized for labeling target molecules including peptides, small molecules, nanoparticles, and macromolecules such as proteins and antibodies
- Highly hydrophilic and suitable for single or multiple dye loading.

Property	Specification
MW	1418 g mol ⁻¹
Fluorescence ¹ • Emission	660 ± 3nm
Absorbance ¹	643 ± 3 nm
Extinction	~210,000 M ⁻¹ cm ⁻¹
Purity ²	>90%
Appearance	Royal blue Solid

1. Absorbance and fluorescence maxima in 1xPBS.
2. As determined by RP-HPLC, measuring absorbance at 645nm



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS, OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF
NEV11274-R Rev01