

IVISense™ Osteo 680 Fluorescent Probe

Product Number: NEV10020EX

DESCRIPTION

IVISense™ Osteo 680 Fluorescent Probe is a fluorescent in vivo bisphosphonate imaging agent. IVISense™ Osteo 680 Fluorescent Probe images areas of microcalcifications and bone remodeling and enables imaging of bone growth and resorption.

MATERIAL

Each vial contains 24 nmol of IVISense™ Osteo 680 Fluorescent Probe as a lyophilized solid. The solution has been filtered (0.2µ) prior to lyophilization. Upon dilution with 1.2 mL of 1 x PBS, this material provides sufficient reagent for imaging approximately 10 mice (weighing ~25 grams each) when using the recommended dose of 2 nmol/100 µL 1xPBS of IVISense™ Osteo 680 Fluorescent Probe per mouse.

STORAGE & HANDLING

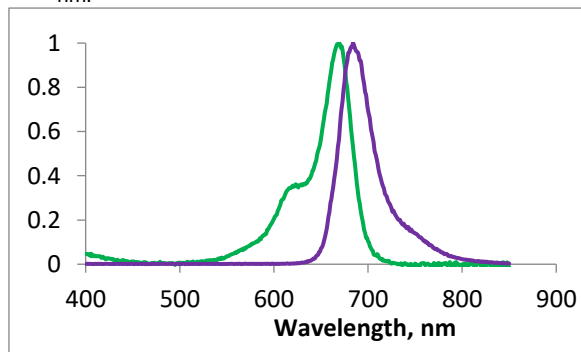
- Upon receipt, IVISense™ Osteo 680 Fluorescent Probe should be IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ Osteo 680 Fluorescent Probe is stable for up to six months from the date of shipment.
- Once reconstituted, the PBS solution is stable up to 14 days when stored at 2-8 °C.

IN VIVO IMAGING & APPLICATIONS

- The recommended procedure for in vivo imaging with IVISense™ Osteo 680 Fluorescent Probe is administration via tail vein injection and imaging 24 hours post injection.
- Imaging Bone Growth: IVISense™ Osteo 680 Fluorescent Probe can be used to measure the effects of therapeutic stimulation of bone growth.
- Imaging Bone Remodeling: IVISense™ Osteo 680 Fluorescent Probe can be used to characterize bone remodeling associated with animal models of arthritis.

Property	Specification
MW	1470.5 g mol ⁻¹
Fluorescence ¹	
• Excitation	668 ± 10 nm
• Emission	687 ± 10 nm
Absorbance ¹	668 ± 5 nm
Purity ²	>95%
Appearance	Blue solid

1. Absorbance, excitation, and fluorescence maxima of IVISense™ Osteo 680 Fluorescent Probe in 1xPBS.
2. Based on concentration resulting in absorbance of 0.3 to 0.5 AU.
3. As determined by RP-HPLC and measuring absorbance at 680 nm.



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
NEV10020EX-R Rev01