

IVISense™ Hypoxia CA IX 680 Fluorescent Probe

Product Number: NEV11070

DESCRIPTION

IVISense™ Hypoxia CA IX 680 Fluorescent Probe is a Carbonic Anhydrase IX (CAIX) targeted fluorescent in vivo imaging agent that can be used to image CAIX overexpression in tumors in response to regional tumor hypoxia.

MATERIAL

Each vial contains 24 nmol of IVISense™ Hypoxia CA IX 680 Fluorescent Probe in dry solid form. Reconstitute IVISense™ Hypoxia CA IX 680 Fluorescent Probe with 1.2 mL of 1 x PBS before injecting into animals. Each vial of packaged material provides sufficient reagent for imaging approximately 10 mice (weighing ~25 grams each) when using the recommended dose of 2 nmol (100 µL) of IVISense™ Hypoxia CA IX 680 Fluorescent Probe per mouse.

STORAGE & HANDLING

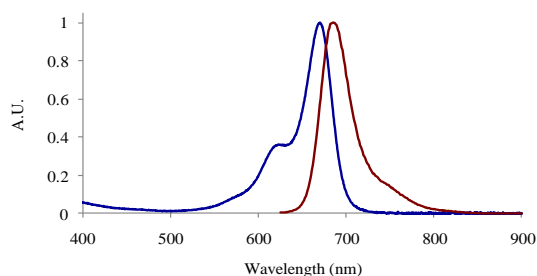
- Upon receipt, IVISense™ Hypoxia CA IX 680 Fluorescent Probe should be IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ Hypoxia CA IX 680 Fluorescent Probe is stable for six months from the date of shipment in dry solid form.
- Before opening the vial, ensure that all of the solid material is at the bottom of the vial.
- Once reconstituted, the solution is stable up to 14 days when stored at 2-8°C and protected from light.

IN VIVO IMAGING & APPLICATIONS

- The recommended procedure for in vivo imaging with IVISense™ Hypoxia CA IX 680 Fluorescent Probe is administration via intravenous injection and imaging 24-48 hours post injection. IVISense™ Hypoxia CA IX 680 Fluorescent Probe may also be injected intraperitoneally and imaged 24 hours post injection.

Property	Specification
MW	~1500 g mol ⁻¹
Fluorescence ¹	
• Excitation	670 nm
• Emission	685 nm
Absorbance ¹	670 nm
Purity ²	>90%
Appearance	Blue solid

1. Absorbance and fluorescence maxima in 1xPBS
2. As determined by RP-HPLC and measuring absorbance at 670 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
NEV11070-R Rev01