

IVISense™ Folate Receptor 680 Fluorescent Probe

Product Number: NEV10040

DESCRIPTION

IVISense™ Folate Receptor 680 Fluorescent Probe is a folate receptor targeted fluorescent in vivo imaging agent that can be used to image folate receptor expression in tumors.

MATERIAL

Each vial contains 24 nmol of IVISense™ Folate Receptor 680 Fluorescent Probe in dry solid form. Reconstitute IVISense™ Folate Receptor 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™ Folate Receptor 680 Fluorescent Probe per mouse.

STORAGE & HANDLING

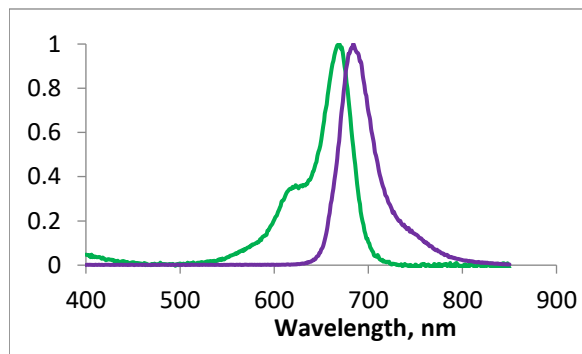
- Upon receipt, IVISense™ Folate Receptor 680 Fluorescent Probe should be IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ Folate Receptor 680 Fluorescent Probe is stable for 3 months from the date of shipment.
- Before opening the vial, check to ensure that all of the solid material is at the bottom of the vial.
- After reconstituting with PBS, gently swirl the solution to ensure that the solid is fully in solution.
- Once reconstituted with 1xPBS, the solution is stable for up to 14 days when store at 2-8 °C and protected from light.

IN VIVO IMAGING

- The recommended procedure for in vivo imaging with IVISense™ Folate Receptor 680 Fluorescent Probe is administration via intravenous injection and imaging 6-24 hours post injection.

Property	Specification
MW	1605.7 g mol ⁻¹
Fluorescence ¹	
• Excitation	670 nm
• Emission	690 nm
Absorbance ¹	670 nm
Purity ²	>90%
Appearance	Blue Solid

1. Absorbance, excitation, and fluorescence maxima in 1xPBS.



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
NEV10040-R Rev01