

# Fluorescent Imaging Agent

Caution: For Laboratory Use. A product for research purposes only.

## HypoxiSense 680

Product Number: NEV11070

**DESCRIPTION:** *HypoxiSense 680* 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:** (Needs to be reconstituted)

**CONTENTS:** Each vial contains 24 nmol of *HypoxiSense 680* in dry solid form. Reconstitute *HypoxiSense 680* 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 *HypoxiSense 680* per mouse.

**PROPERTIES:** The physical characteristics of *HypoxiSense 680* can be found in **Table 1** and **Figure 1**.

### STORAGE & HANDLING:

- Upon receipt, *HypoxiSense 680* should be **IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT**.
- When stored and handled properly, *HypoxiSense 680* is stable for up to 6 months in dry solid form 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 1 x PBS, the solution is stable up to 10 days when stored at 2-8 °C and protected from light.**

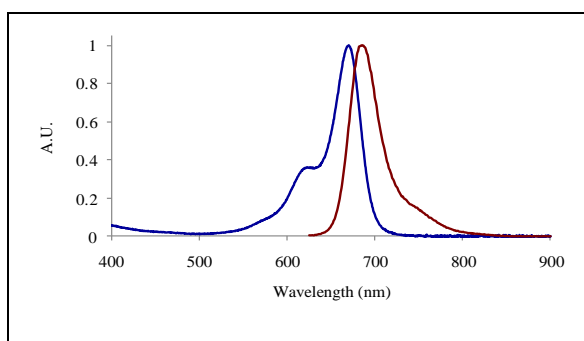
### IN VIVO IMAGING AND APPLICATIONS:

- The generally recommended procedure for *in vivo* imaging with *HypoxiSense 680* is administration via intravenous injection and imaging **24-48 hours post injection**. *HypoxiSense 680* may also be injected intraperitoneally and imaged 24 hours post injection.

**Table 1.** *HypoxiSense 680* Characteristics

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

1. Absorbance and fluorescence maxima of *HypoxiSense 680* in 1x PBS.  
2. As determined by RP-HPLC and measuring absorbance at 670 nm.



**Figure 1.** Normalized absorbance (blue) and fluorescence emission (red) spectra of *HypoxiSense 680* in 1x PBS.

## NOTES:

- *HypoxiSense 680* is intended for research purposes only and is not for human use. It must be used by or directly under the supervision of a technically qualified individual experienced in handling potentially hazardous materials. Please read the Material Safety Data Sheet (MSDS) provided for this product.
- Several of *PerkinElmer's* products and product applications are covered by U.S. and foreign patents and patents pending. Our products are not available for resale or other commercial uses without a specific agreement with *PerkinElmer*.

PerkinElmer, Inc.  
549 Albany Street  
Boston, MA 02118 USA  
P: (800) 762-4000 or (+1) 203-925-4602  
[www.perkinelmer.com/nenradiochemicals](http://www.perkinelmer.com/nenradiochemicals)

**For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)**

Copyright ©2010, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc.  
All other trademarks are the property of their respective owners.