

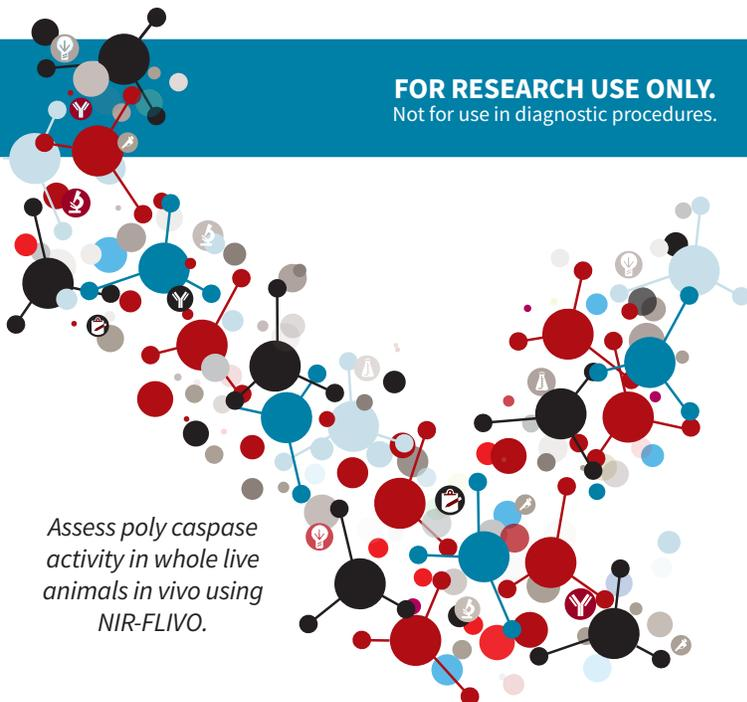
## 1. INTRODUCTION

FLIVO® (FLuorescence in vivo) is a powerful method for assessing caspase activity *in vivo*. Like our FLICA® probes<sup>1,2</sup>, but optimized for whole live animal imaging, FLIVO Tracers are non-cytotoxic fluorescent inhibitors of caspases. ICT's near-infrared (NIR)-FLIVO poly caspase tracers contains the preferred binding sequence for most caspases, Val-Ala-Asp (VAD). This preferred poly caspase tripeptide binding sequence is labeled at the amino terminus end with \*Dylight® 690 or 747 NIR label and linked at the carboxyl end to a fluoromethyl ketone (FMK) reactive entity.

Apoptosis is an evolutionarily conserved process of programmed cell suicide. It is centered on a cascade of proteolytic enzymes called caspases that are triggered in response to pro-apoptotic signals. Like most other proteases, caspases are synthesized as pro-form precursors that undergo proteolytic maturation, either autocatalytically or in a cascade by enzymes with similar specificity<sup>3</sup>. Active caspase enzymes consist of two large (~20 kD) and two small (~10 kD) subunits that non-covalently associate to form a two heterodimer, tetrameric active caspase<sup>4-6</sup>. Once activated, caspases cleave protein substrates leading to the eventual disassembly of the cell. Caspases have been identified in organisms ranging from *C. elegans* to humans. Mammalian caspases play distinct roles in both apoptosis and inflammation.

FLIVO kits provide a simple yet accurate method to detect caspase activity *in vivo*. To label cells containing elevated levels of active caspases, inject the FLIVO Tracer intravenously and let it circulate. Because the FLIVO Tracer is cell-permeant, it readily diffuses in and out of all cells it encounters as it circulates throughout the body. If there are active caspase enzymes inside a cell, FLIVO will form an irreversible covalent bond with a reactive cysteine on the large subunit of the caspase heterodimer, thereby inhibiting further enzymatic activity. The bound FLIVO Tracer will remain inside the cell if the cell membrane is intact. Any unbound FLIVO is removed from the circulation of the animal in about an hour. The remaining NIR fluorescent signal in the tissue is a direct measure of caspase activity that occurred at the time the reagent was injected. Apoptotic cells will retain a higher concentration of FLIVO Tracer and fluoresce brighter than non-apoptotic cells. There is no interference from pro-caspases or inactive forms of the enzyme. If the treatment is causing cell death via apoptosis, apoptotic cells will have an elevated level of caspase activity relative to non-apoptotic or negative control cells and fluoresce near-infrared with FLIVO.

NIR-FLIVO Tracers are used in conjunction with NIR-FLIVO Free Dye Control Assays (sold separately). The NIR-FLIVO Free Dye Control Assay uses the NIR fluorescent dye molecule. When injected, both



*Assess poly caspase activity in whole live animals in vivo using NIR-FLIVO.*

the Free Dye Control and the Tracer reagents will generate a fluorescent signal, but only the Tracer should bind to active caspases and remain inside an apoptotic cell. The fluorescent signal generated from the Free Dye Control reveals where the fluorescent reagent may have become trapped inside the cell, tissues, or body of the animal without specifically binding with an active caspase. In the context of the overall experiment, this base level of fluorescence is considered background noise compared with the signal generated in animals that were injected with the NIR-FLIVO Tracer.

An initial experiment may be necessary to determine when and how much NIR-FLIVO to inject based on the size of the animal, tissue type, experimental conditions, rate of apoptosis, and method of analysis. Generally, the longer FLIVO circulates, the lower the non-specific background signal; however, some apoptotic cells may be lost over time. After 60 minutes, most of the unbound NIR-FLIVO Tracer or Free Dye Control will have cleared the bloodstream. The bound NIR-FLIVO Tracer will remain inside an apoptotic cell and generate a positive signal if the cell membrane is still intact.

Once the animals have been injected with FLIVO, they are ready for analysis and no further staining is necessary. Because FLIVO is a direct stain, it eliminates any false positives that may arise from manipulation of the tissue. This gives a true representation of the induction of apoptosis *in vivo* as a result of the experimental condition. Live animals may be analyzed in a whole animal imager, and /or tissues may be prepared and further analyzed by histological methods. Tissues labeled with FLIVO can be counter-stained with other reagents such as DAPI, and fixed or frozen for future analysis. The fluorescence intensity can be quantified by excising the tissue and analyzing cells with a flow cytometer. NIR-FLIVO 690 Tracer optimally excites at 690 nm and has a peak emission at 709 nm. NIR-FLIVO 747 Tracer optimally excites at 747 nm and has a peak emission at 776 nm.

## 2. KIT CONTENTS

### Catalog #9112 contains:

- 2 vials of NIR-FLIVO 690 Tracer (\*Dylight® 690-VAD-FMK), 10 Tests per vial, 71.4 µg per vial, #6306
- 1 bottle of 10X Injection Buffer, 5 mL, #6220

### Catalog #9114 contains:

- 2 vials of NIR-FLIVO 747 Tracer (\*Dylight® 747-VAD-FMK), 10 Tests per vial, 77.3 µg per vial, #6308
- 1 bottle of 10X Injection Buffer, 5 mL, #6220

## 3. STORAGE

Store the unopened kit and each unopened component at -20°C until the expiration date. Once reconstituted with DMSO, use NIR-FLIVO Tracer immediately, or store at ≤-20°C for 6 months protected from light and thawed no more than twice during that time.

## 4. SAFETY DATA SHEETS (SDS)

SDS are available at online at [www.immunochemistry.com](http://www.immunochemistry.com) or by calling 1-800-829-3194 or 952-888-8788.

## 5. RECOMMENDED MATERIALS

- DMSO, 50 µL per vial to reconstitute NIR-FLIVO Tracer
- DiH<sub>2</sub>O, 45 mL to dilute 10X Injection Buffer

- 0.2 µm syringe filter to sterilize Injection Buffer
- Injection materials such as a syringe and needle
- Experimental and control animals ready to be assessed
- Optional: tools to dissect, extract, and examine labeled tissues
- Optional: NIR-FLIVO® 690 / 747 Free Dye Control Assays (kits #9113 and #9115, respectively)

## 6. DETECTION EQUIPMENT

Detection equipment such as instrumentation designed for non-invasive live whole animal imaging, alternatively, excised tissues can be analyzed by fluorescence microscope or flow cytometer.

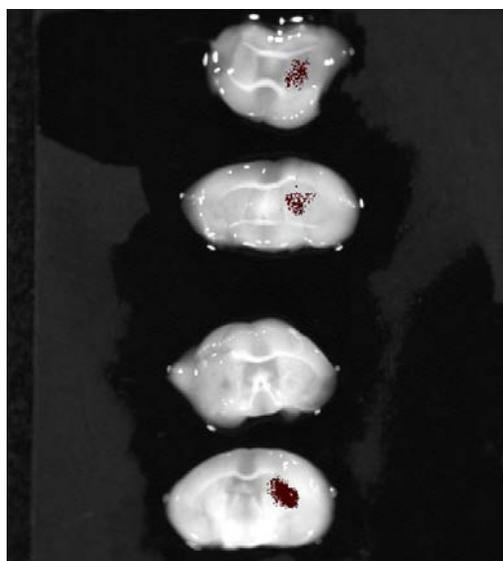
- NIR-FLIVO 690 Tracer optimally excites at 690 nm and has a peak emission at 709 nm.
- NIR-FLIVO 747 Tracer optimally excites at 747 nm and has a peak emission at 776 nm.

## 7. EXPERIMENTAL PREPARATION

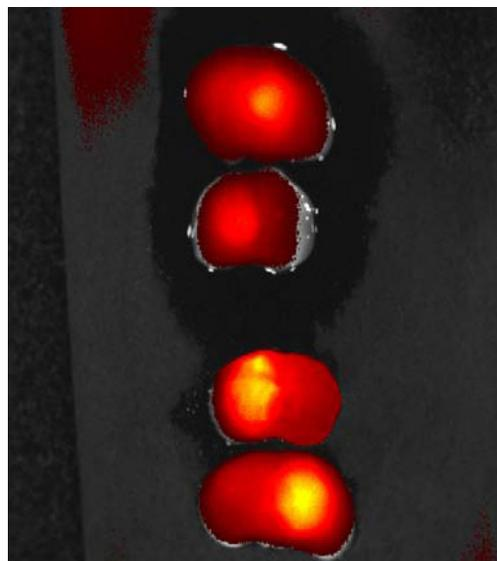
Plan your experiment so that NIR-FLIVO Tracer can be reconstituted, diluted and then injected at the time when caspase activity is expected to be occurring in the animal. It may be necessary to set up an initial experiment to determine when and how much NIR-FLIVO Tracer to inject as the resulting positive fluorescent signal is a direct measure of caspase activity that occurred at the time of injection. The amount of NIR-FLIVO Tracer may need to be adjusted to

### FIGURE 1. DETECT APOPTOSIS IN MURINE BRAIN ABSCESSSES USING NIR-FLIVO 690 TRACER

Brain abscesses were induced in mice following the intracerebral inoculation of live *Staphylococcus aureus*. Animals received intravenous injections of NIR-FLIVO 690 (\*DyLight® 690-VAD-FMK, kit #9112) or NIR-FLIVO 690 Free Dye Control (kit #9113) at 17 hours post-infection, whereupon signals were acquired 1 hour later from brain tissues *ex vivo* using an IVIS® Spectrum™ (Caliper Life Sciences). Strong caspase activity was associated with brain abscesses following administration of the NIR-FLIVO 690 tracer (right image), whereas minimal signal was detected in animals injected with the NIR-FLIVO 690 Free Dye Control (left image).



FREE DYE CONTROL



NIR-FLIVO 690 TRACER

accommodate the experimental model and research conditions being investigated.

1. Expose the test animals to your experimental conditions to assess caspase activity in the target tissue(s).
2. Prepare the animals for intravenous NIR-FLIVO injection.

## 8. PREPARATION OF 1X INJECTION BUFFER

ICT's Injection Buffer is an isotonic solution used for diluting and injecting NIR-FLIVO Tracers.

1. 10X Injection Buffer may form precipitates during cold storage. If this happens, gently warm it until all crystals have dissolved. Do not boil.
2. Dilute 10X Injection Buffer 1:10 in diH<sub>2</sub>O. For example, add 1 mL 10X Injection Buffer to 9 mL diH<sub>2</sub>O for a total of 10 mL.
3. Sterilize the 1X Injection Buffer by filtering through a 0.2 μm syringe filter or equivalent.
  - 1X Injection Buffer may be stored at 2-8°C and used within 1 week or frozen and used within 6 months.

## 9. PREPARATION OF NIR-FLIVO TRACER

NIR-FLIVO Tracers are supplied as a lyophilized powder that may be slightly visible as an iridescent sheen inside the vial. Protect from light and use gloves and eye protection when handling. Each vial

**FOR RESEARCH USE ONLY.**  
Not for use in diagnostic procedures.

of NIR-FLIVO 690 Tracer contains 71.4 μg. Each vial of NIR-FLIVO 747 Tracer contains 77.3 μg. This amount is typically enough to inject 10 mice or 2 young rats at approximately 300 nanomoles/Kg.

1. Reconstitute each vial of NIR-FLIVO Tracer with 50 μL DMSO to form the 10X stock solution. The stock solution may have a blue-green color. Once reconstituted, it may be stored at ≤-20°C for 6 months protected from light and thawed no more than twice during that time. Wear gloves and eye protection.
2. Immediately prior to injection into the animal, further dilute the 10X stock NIR-FLIVO Tracer 1:10 by adding 450 μL sterile 1X Injection Buffer to each vial to form 500 μL of the 1X NIR-FLIVO Tracer solution. Inject 1X NIR-FLIVO Tracer within 1 hour of dilution into aqueous buffer; protect from light during handling.

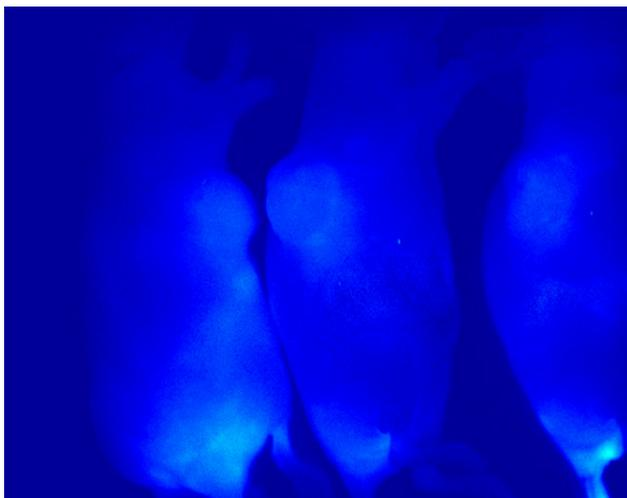
## 10. INTRAVENOUS INJECTION

The recommended dose of NIR-FLIVO Tracer is 300 nanomoles/Kg per injection, but the necessary amount will depend on the extent of apoptosis in the animal. Each vial of reagent will test 10 tumor-bearing mice or 2 tumor-bearing young rats at approximately 300 nanomoles/Kg. Animal models that do not involve tumors may require less reagent. An initial experiment may be necessary to determine how much of the NIR-FLIVO Tracer to inject.

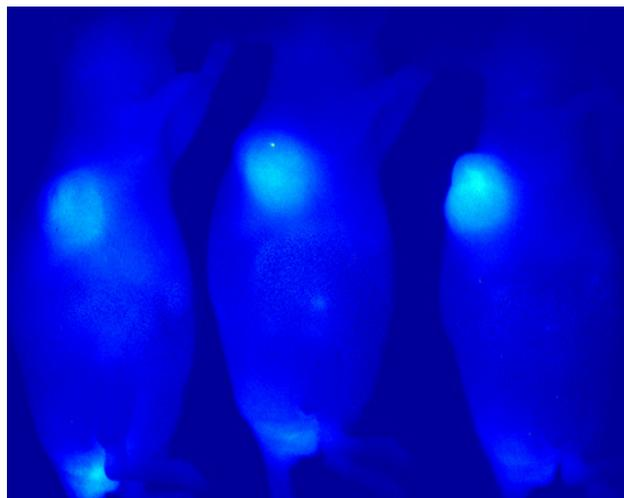
1. Inject 50 μL 1X NIR-FLIVO Tracer into the tail vein or other large vein. Intraperitoneal or intravitreal injection is not advised. The exact IV location and amount of NIR-FLIVO Tracer injected may vary depending on the size of the animal, the target tissue, and the experimental conditions.

### FIGURE 2. *IN VIVO* MONITORING OF APOPTOSIS DURING EXPERIMENTAL ANTI-CANCER TREATMENT

Apoptotic tissues in murine tumor models were imaged *in vivo* with NIR-FLIVO™ 747 (\*DyLight® 747-VAD-FMK, kit #9114) using the CRi Maestro™ imaging system. NIR-FLIVO 747 was used to non-invasively monitor the efficacy and time kinetics of an experimental anti-cancer treatment. After receiving a placebo or an experimental treatment, negative control mice (left image) and experimentally treated mice (right image) were injected intravenously with NIR-FLIVO 747 Tracer and imaged non-invasively at various time points with the CRi Maestro™ imaging system. Apoptotic tumor tissues fluoresce bright blue in the experimental group subjects, indicating that the experimental treatment had an apoptotic effect.



CONTROL GROUP



EXPERIMENTAL GROUP



- Optional: If a larger injection volume is preferred, dilute NIR-FLIVO Tracer with a greater amount of 1X Injection Buffer and inject 1/10th of the final volume into the animal.
- Allow NIR-FLIVO Tracer to circulate within the animal for at least 30 minutes. After 60 minutes, most of the unbound FLIVO will have cleared from the bloodstream. Generally, the longer the reagent circulates, the lower the background signal; however, some positive cells (apoptotic or pyroptotic) may be lost over time. NIR-FLIVO Tracer will remain inside a cell containing active caspases as long as the cell membrane is intact.

## 11. CIRCULATION

When determining the circulation time, it is important to consider both the reagent clearance rate as well as the rate of cellular apoptosis in the model system. NIR-FLIVO Tracers will clear from the circulating bloodstream within an hour and clear the liver in 3-4 hours. In general, non-specific background signal will decrease with increased circulation time, but caspase-positive cells may be lost over time via natural degradation. The NIR-FLIVO Tracer will be retained inside an apoptotic cell as long as the cell membrane is intact; it will generate the specific signal. The NIR-FLIVO Free Dye Control should not remain in the body; any remaining signal is background fluorescence. The Free Dye Control should circulate for the same length of time as the Tracer. An initial experiment may be necessary to determine the optimal circulation time, which will vary by the method of analysis and the tissue of interest:

- Analysis by *in vivo* imaging: Allow the NIR-FLIVO Tracer to circulate at least 4 hours before imaging. Depending on the rate of cellular apoptosis in the animals receiving the NIR-FLIVO Tracer, the animals should be able to be reinjected with another dose of the NIR-FLIVO Tracer and reanalyzed within 1-4 days if desired.
- Analysis of non-hepatic histological samples *ex vivo*: Allow NIR-FLIVO Tracer to circulate *in vivo* at least 30-60 minutes before preparing samples. Protect samples from light.
- Analysis of hepatic histological samples *ex vivo*: Allow NIR-FLIVO Tracer to circulate *in vivo* at least 4 hours before preparing samples. Protect samples from light.

## 12. PREPARATION OF ANIMALS FOR LIVE IMAGING

Anesthetize animals according to the experimental protocol and place in optical imaging machine. Consult instrument manufacturer for assistance with equipment set up and data collection.

## 13. PREPARATION OF HISTOLOGY SAMPLES

Prepare the animal tissues according to your desired protocol. We have listed several methods here:

- Excise the tissue, freeze, and make thin tissue sections for histology.
- Perfuse with formalin or a non-methanol, non-ethanol fixative or embedding agent.
- Extract the tissue and prepare cells for flow cytometry.

## 14. REFERENCES

- Amstad, P. A. et al. Detection of caspase activation in situ by fluorochrome-labeled caspase inhibitors. *Biotechniques* 31, 608-610, 612, 614, passim (2001).
- Bedner, E., Smolewski, P., Amstad, P. & Darzynkiewicz, Z. Activation of caspases measured in situ by binding of fluorochrome-labeled inhibitors of caspases (FLICA): correlation with DNA fragmentation. *Exp Cell Res* 259, 308-313, doi:10.1006/excr.2000.4955 (2000).
- Kumar, S. Mechanisms mediating caspase activation in cell death. *Cell Death Differ* 6, 1060-1066, doi:10.1038/sj.cdd.4400600 (1999).
- Wilson, K. P. et al. Structure and mechanism of interleukin-1 beta converting enzyme. *Nature* 370, 270-275, doi:10.1038/370270a0 (1994).
- Walker, N. P. et al. Crystal structure of the cysteine protease interleukin-1 beta-converting enzyme: a (p20/p10)<sub>2</sub> homodimer. *Cell* 78, 343-352 (1994).
- Rotonda, J. et al. The three-dimensional structure of apopain/ CPP32, a key mediator of apoptosis. *Nat Struct Biol* 3, 619-625 (1996).

**FOR RESEARCH USE ONLY.**

Not for use in diagnostic procedures.



**BRIGHT MINDS, BRIGHT SOLUTIONS.**

ImmunoChemistry Technologies, LLC gratefully acknowledges the significant contributions made by one of its founders, Brian W. Lee, Ph.D in the development of this product, including the creation and illustration of its strategy and protocol.

**ImmunoChemistry Technologies, LLC**

9401 James Ave. S. #155, Bloomington, MN 55431 USA 800-829-3194

Copyright ©2019 ImmunoChemistry Technologies, LLC. FLICA® and FLIVO® are registered trademarks of ImmunoChemistry Technologies, LLC. \*Dylight® is a registered trademark of Thermo Fisher Scientific.

immunochemistry.com