

revvity

IVISense™ fluorescent imaging panels

Fluorescent panel user guide

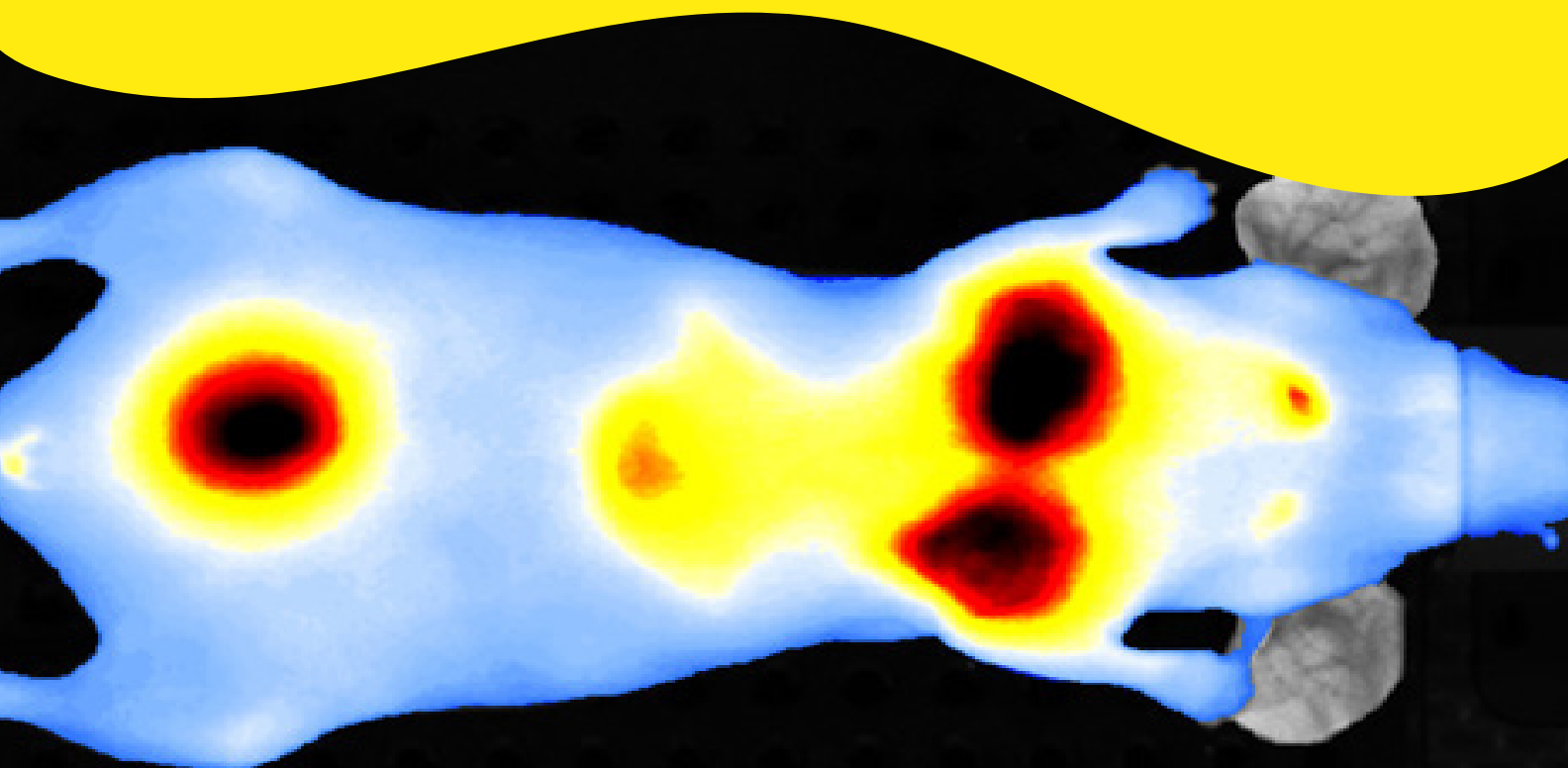


Table of contents

Protocol For Monoplex, Multiplex, and Longitudinal Imaging with <i>In Vivo</i> Probe Panels	3
Study Preparation	3
Single Probe Imaging	3
Two Probe Imaging	4
Longitudinal Imaging	4
Toxicology Probe Cocktail	4
Probe Dose and Clearance Table	5

Fluorescence molecular imaging

Revvity's imaging probes are developed through an extensive R&D process and designed to incorporate drug-like biodistribution properties for optimal target delivery and performance. The table summarizes proper dosages, imaging time points, routes of metabolism, and probe clearance kinetics. Probes can be used either singly or in pairs by combining appropriate pairs of 680 nm and 750 nm probes. These probes can also be used for longitudinal studies by working within the parameters of the tissue pharmacokinetics and only reinjecting upon complete clearance.

Protocol for monoplex, multiplex, and longitudinal imaging with *in vivo* probe panels

Study preparation

1. Two weeks before the imaging study, switch mice to low fluorescence chow. Regular mouse chow contains chlorophyll that auto fluoresces around 700 nm which can interfere with imaging.
2. On the study day, it is essential to prepare ahead of time for optimal results. Group and number the mice to be injected and imaged. Appropriate study design should include both positive control and negative control (i.e. un-diseased) mice injected with probe(s).
3. Mouse hair removal is essential for sensitive, high-quality fluorescence imaging. Either genetically hairless mice (SKH-1E) or normal, haired mice (BALB/c, C57BL/6, etc.) with depilation, must be used for optimal fluorescence tomographic imaging. This can be performed under injectable or inhaled anesthesia.
 - To minimize light scattering and absorption in fluorescence imaging, hair is removed from the appropriate body region of all mice; depilatory cream (Nair lotion, Church and Dwight Co., Inc., Princeton, NJ) is applied thickly on hair over the imaging region of each mouse, rinsed off thoroughly with warm water, and reapplied until all hair has been removed. [Rinsing must be done carefully and thoroughly to minimize any introduction of skin lesions that can cause imaging artifacts.]
 - Care should be taken to remove hair from an area larger than just the region of focus to assure that you can capture target and surrounding background fluorescence. For tomographic imaging, hair must be removed from front, sides, and back for the region of focus.
4. Establish the readiness of the imaging system by checking the anesthesia chamber and connections to the system. Activate the anesthesia, setting evaporator to the appropriate settings for your particular set-up.
5. Make sure you know ahead of time the proper positioning of the mouse you will use to facilitate acquisition of the best quality data for your particular animal model.

Single probe imaging

1. Prepare the imaging probe according to included instructions. All probes must be injected systemically either through the retro-orbital plexus or the tail vein. For most of the probes intraperitoneal (IP) or subcutaneous (SC) injection will either not work at all or will be extremely variable and with high injection site signal.
2. Place a heating pad beneath the anesthesia induction chamber to keep the body temperature of the mice constant. Be careful not to overheat. Anesthetize the first mouse by placing it in a gas anesthesia induction chamber.
3. Remove the mouse from the induction chamber when it appears completely anesthetized, and confirm the depth of anesthesia through unresponsiveness to toe pinch.
4. Inject the appropriate volume of probe (100 - 150 μ L, as per each probe's specific instructions) via the retro-orbital plexus (or tail vein) of the anesthetized mouse. Record the injection time.
5. Return the mouse to the cage for recovery and go to the next mouse for injection.

- Repeat steps 1 - 6 until all mice are injected.
- Imaging is performed at the suggested time(s) using a single excitation/emission filter pair optimal for the wavelength of the probe to be imaged (see table). Anesthetize mice using inhaled anesthesia and place them carefully in the appropriate orientation in the imaging system. Multiple images can be acquired with little or no concern for photobleaching of the probes.

Two probe imaging

- Prepare the first imaging probe as described in that probe's instructions. Either use the Probe 1 solution to solubilize Probe 2 in order to minimize injection volume or make each probe at half-volumes for mixing. [Note: Bear in mind that three of the probes (IVISense™ Annexin-V 750, IVISense MMP 680, and IVISense Pan Cathepsin 680) come in 10X liquid form]. Multiple specific strategies for preparation are possible, but it is ideal to keep mouse injection volumes under 250 μ L.
- Prepare and inject mice as described above.
- Imaging should be performed as described above, but for both 680 nm and 750 nm using the appropriate excitation and emission filter pairs. Correct times for acquisition should be noted; although most of the probes are optimal for 24 h imaging, some can be imaged earlier, and some should be imaged earlier (see table). For example, IVISense Annexin-V 750 imaging is optimal for most applications at 2 h, whereas IVISense Pan Cathepsin 680 is optimal at 24 h. You can either image both wavelengths at both 2 and 24 h, or you can image 750 nm at 2 h and 680 nm at 24 h.
- Multiple repeat acquisitions can be performed with little or no concern for photobleaching of the probes.

Longitudinal imaging

- Revvity probes are well characterized with respect to tissue clearance kinetics, providing guidance for longitudinal imaging strategies. Depending on the probe, reinjection generally can be performed three to seven days following the first image acquisition.

- IVISense Osteo 680 bone turnover imaging, however, requires a different strategy due to the very long tissue clearance kinetics. Secondary imaging time points must be performed using a pre-imaging strategy; briefly, mice should be imaged immediately prior to each additional probe injection to allow subtraction correction of additional imaging datasets.

Toxicology probe cocktail

- Revvity recently published research showing the utility of an IVISense Annexin-V/MMP/Transferrin Receptor cocktail of 750 nm probes (AMT-750) in the imaging of drug-induced liver injury (Vasquez & Peterson, J Pharmacol Exp Ther 2017; 361:87-98). The Toxicology Panel of probes provides an extra vial of IVISense MMP 750 FAST to allow the preparation of AMT-750.
- Prepare the AMT-750 as below:

IVISense probe	Prepared stock concentration	Volume
Annexin-V 750	As supplied in liquid form	0.5 mL
MMP 750 FAST	Prepare as 80 μ M stock in PBS	0.5 mL
Transferrin Receptor 750	Prepare as 10 μ M stock in PBS	0.5 mL
		Total 1.5 mL Mouse Dose 150 μL

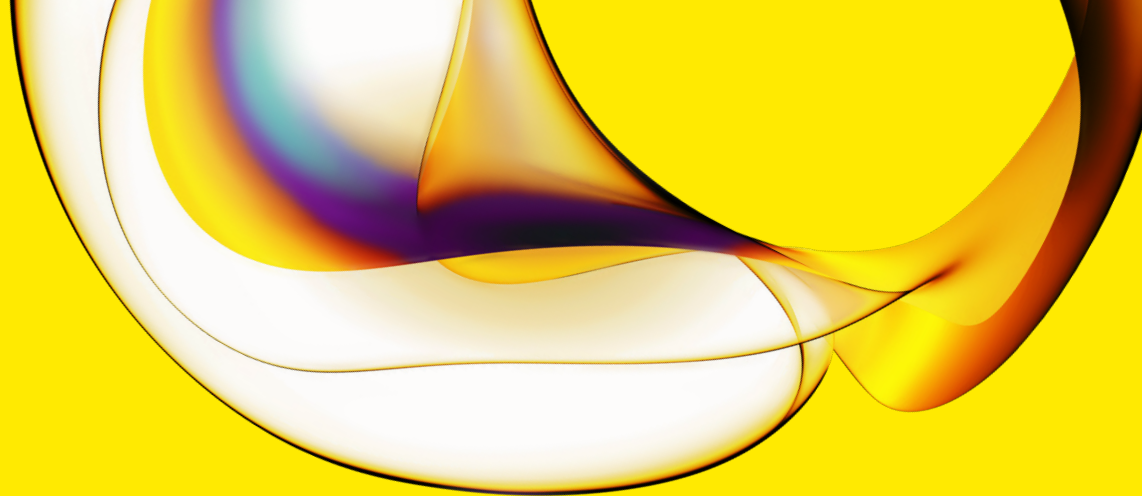
The ratio of the cocktail is specifically designed to optimize liver injury signal while minimizing normal background signal in liver and kidneys.

- AMT-750 should be injected retro-orbitally or via tail vein at 2 h or 24 h post-drug treatment. Some drugs (given as a single IP bolus of 100 - 300 mg/kg) will induce early biological changes predictive of tissue injury, whereas others require 24 h to manifest tissue biological changes. Imaging time is optimized for 24h post-AMT-750 injection (Vasquez & Peterson, 2017).

Probe dose and clearance table

Part number	IVISense probe	Packaged amount	Mouse dose (25 g)	Optimal imaging time	Probe clearance	Route of metabolism/background tissue(s)	ex/em wavelengths (IVIS® Spectrum)
NEV10054EX	Vascular 680	24 nmol	2 nmol	24 h	6-7 d	Low Liver, Lung	675/720
NEV10723EX	Vascular 680 sample size	8 nmol	2 nmol	24 h	6-7 d	Low Liver, Lung	675/720
NEV10011EX	Vascular 750	24 nmol	2 nmol	24 h	6-7 d	Low Liver, Lung	745/800
NEV11053	Annexin-V 750	1 mL	100 µL	2 h	3 d	Kidneys (High), Liver	745/800
NEV10090	Bombesin Receptor 680	24 nmol	2 nmol	24 h	6-7 d	Pancreas, Kidney	675/720
NEV11000	CAT K 680 FAST	24 nmol	2 nmol	6-24 h	3 d	Kidney > Liver	675/720
NEV11112	CAT B 680 FAST	24 nmol	2 nmol	6-24 h	3 d	Salivary Glands > Liver, Kidneys	675/720
NEV11098	CAT B 750 FAST	48 nmol	4 nmol	6-24 h	3 d	Salivary Glands > Liver, Kidneys	745/800
NEV10040	Folate Receptor 680	24 nmol	2 nmol	6 h (6-24)	6-7 d	Kidneys	675/720
NEV10645	Integrin Receptor 680	24 nmol	2 nmol	24 h	14 d	Kidneys	675/720
NEV10873	Integrin Receptor 750	24 nmol	2 nmol	24 h	4-6 d	Kidneys	745/800
NEV10878	Integrin Receptor 750 Sample Size	7 nmol	2 nmol	24 h	4-6 d	Kidneys	745/800
NEV11169	Neutrophil Elastase 680 FAST	48 nmol	4 nmol	3-6 h	2 d	Bladder > Liver, Intestines	675/720
NEV10126	MMP 680	20 nmol	2 nmol	24 h (24-36)	6-7 d	Liver	675/720
NEV10168	MMP 750 FAST	24 nmol	2 nmol	24 h (12-24)	6-7 d	Liver > Kidneys	745/800
NEV10932	MMP 750 FAST sample size	7 nmol	2 nmol	24 h (12-24)	6-7 d	Liver > Kidneys	745/800
NEV10020EX	Osteo 680	24 nmol	2 nmol	3-24 h	4 weeks	Bladder	675/720
NEV10003	Pan Cathepsin 680	20 nmol	2 nmol	24 h (24-48)	6-7 d	Liver	675/720
NEV10001EX	Pan Cathepsin 750	24 nmol	2 nmol	24 h	6-7 d	Low Liver, Intestine	745/800
NEV10972EX	Pan Cathepsin 750 Sample size	8 nmol	2 nmol	24 h	6-7 d	Low Liver, Intestine	745/800
NEV11171	Pan Cathepsin 750 FAST	48 nmol	4 nmol	6-24 h	3 d	Low Liver, Bladder	745/800
NEV11079	Renin 680 FAST	24 nmol	2 nmol	24 (12-24) h	4 d	Kidney, lung, liver	675/720
NEV10091	Transferrin Receptor 750	24 nmol	2 nmol	24 (6-24) h	4 d	Liver, kidney	745/800
NEV11118	680 NHS Fluorescent Labeling Kit	Kit: 2 x 0.25 mg	Labeling reaction	Dependent on the protein that is conjugated	Dependent on the protein that is conjugated	Dependent on the protein that is conjugated	675/720
770504	IVISbrite™ D-Luciferin Substrate in RediJect™ Solution	10 x 850 µL of 30 mg/mL	125 µL per 25 g mouse (or 150 µL per 30 g mouse)	10-20 min or as determined by kinetic curve*	NA	NA	NA

* See "Determining the Luciferin Kinetic Curve for Your Model" available on our website or by contacting Global Technical Support at global.techsupport@revvity.com



www.revivity.com

revvity

Revvity, Inc.
940 Winter Street, Waltham, MA 02451 USA
(800) 762-4000 | www.revivity.com

For a complete listing of our global offices, visit www.revivity.com
Copyright ©2023, Revvity, Inc. All rights reserved.

471627