

# IVISense™ D-Luciferin Ultra Bioluminescent Substrate in RediJect™ Solution

Product Number: 770505

## DESCRIPTION

IVISbrite D-Luciferin Ultra Bioluminescent Substrate in RediJect Solution is preformulated with a rapidly clearing fluorescent dye to validate the integrity of your substrate injection. With a quick fluorescent image on your IVIS system, you know immediately if an injection failed and you can normalize your bioluminescence results.

## CONTENTS

- 10 vials – IVISbrite D-Luciferin Ultra Bioluminescent Substrate in RediJect Solution format (1 vial images 5 mice)

Property	Specification
Color and Form	Green colored solution (D-Luciferin Potassium Salt in PBS with fluorescent tracer)
Concentration	30 mg/mL
Vial	Amber sterile vial
Volume per vial	850 µL of 30 mg/mL D-Luciferin
Shipping condition	Shipped on blue ice

## STORAGE & HANDLING

- Upon receipt, IVISbrite D-Luciferin should be IMMEDIATELY STORED AT  $\leq -20^{\circ}\text{C}$  and protected from air and light.
- When stored and handled properly, contents are stable until expiration date printed on vial label.
- Just before use, thaw required number of vials in a  $37^{\circ}\text{C}$  water bath and vortex.
- IVISbrite D-Luciferin Ultra supports intravenous, subcutaneous, or intraperitoneal injection.
- Repeated freeze thaw is not recommended.

## PROTOCOL

For *in vivo* imaging studies, we recommend intraperitoneal injection of 150 mg/kg (150 µL/mouse\*). Load a 1 mL syringe directly from the vial and inject using a 25 gauge needle. \*Calculations based on a 30g mouse.

- For fluorescence imaging, select exposure time between 1 and 5 seconds and image using 745 nm excitation and 800 nm emission filter set (ICG filter set for IVIS instruments). Select a route of injection (ROI) on the scruff of the neck on the dorsal side to evaluate the accuracy of your intraperitoneal injection. For ventral imaging, ROI should be drawn around the thoracic region.

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

770505-R Rev01