

# IVISense™ D-Luciferin Bioluminescent Substrate in RediJect™

## Solution

Product Number: 770504

### DESCRIPTION

IVISbrite D-Luciferin Bioluminescent Substrate in RediJect Solution is a bioluminescent *in vivo* substrate in a convenient ready-to-use, pre-formulated injectable format developed by Revvity to fit your laboratory workflow.

### CONTENTS

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

Property	Specification
Color and Form	Yellow colored solution (D-Luciferin Potassium Salt in PBS)
Concentration	30 mg/mL
Vial	Clear sterile vial
Volume per vial	850 µL of 30 mg/mL D-Luciferin
Shipping condition	Shipped on dry 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 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.

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

770504-R Rev01