

# In Vivo imaging Protocol with IVISbrite D-Luciferin RediJect Solution.

## IVISbrite™ D-Luciferin Bioluminescent Substrate - RediJect™ Solution

<b>Part number</b>	770504
<b>Properties</b>	Yellow colored solution (D-Luciferin Potassium salt in PBS)
<b>Concentration</b>	10 Sterile Vials each containing 850 µL of 30 mg/ml D-Luciferin
<b>Storage and handling</b>	Store below -70 °C. Repeated freeze Thaw is not recommended.

- Just before your experiment, remove a vial from the kit and place it in a 37 °C water bath for five minutes. Vortex the tube for one minute and it is ready to use.
- For *in vivo* imaging studies, we recommend injection of IVISbrite™ D-Luciferin - RediJect™ Solution at 150 mg/kg (150 µL/mouse injection\*) using a 25 gauge needle, usually with 1 cc syringe. Injections can be performed intraperitoneally, subcutaneously or intravenously.
- A Luciferin kinetic curve should be performed for each new animal model to determine signal plateau duration. **Please see our 'Determining the Luciferin kinetic curve for your model' instruction sheet available for download on our website.**
- Once plateau is determined, allow D-Luciferin to distribute in animals under conditions consistent with those the animals were under during kinetic curve generation, i.e. under anesthesia and warmed to 37 °C.
- Place fully anesthetized animals in the IVIS imaging system and perform bioluminescence imaging.

\*Calculations based on a 30 g mouse

