

## IVISbrite™ D-Luciferin Potassium Salt Bioluminescent Substrate

**Product Number: 122799**

**Lot Number: K9973PE**

**Date of Manufacture: 22<sup>nd</sup> November 2023**

**Storage temperature: -18°C**

**Expiry date: 21<sup>st</sup> November 2028**

**Approval date: 18<sup>th</sup> January 2024**

**Vacuum dried, packed under argon in amber glass vial with crimp cap**

### DESCRIPTION

D-Luciferin is a chemical substance found in the cells of organisms that generate bioluminescence. When luciferin is oxidized under the catalytic effects of firefly luciferase and ATP, light is produced. Luciferin is able to penetrate cell membranes and can be used to monitor activity of cells of interest in vivo that have been transformed to express luciferase. Because the reaction with luciferase is ATP dependent, cellular viability can also be determined. Use IVISbrite™ D-Luciferin Potassium Salt Bioluminescent Substrate with confidence in your in vivo and in vitro studies. \*

Property	Specification
Color and Form	Nearly white powder
Molecular Formula	C <sub>11</sub> H <sub>7</sub> KN <sub>2</sub> O <sub>3</sub> S <sub>2</sub>
Molecular Weight	318.41
CAS#	115144-35-9
Solubility	up to 40 mg/ml in H <sub>2</sub> O and Phosphate Buffered Saline
Chemical Purity	>99% by HPLC
Biochemical Performance	>98.5 % +/- 1.2 with ATP/ Luciferase
Optical Purity	>98% by HPLC
Vial	Amber vial filled under argon gas
Volume per vial	1 g D-Luciferin
Shipping condition	Shipped on blue ice

## STORAGE & HANDLING

- Upon receipt, IVISbrite™ D-Luciferin should be stored at -18°C and protected from air and light.
- When stored and handled properly, an unopened vial of lyophilized D-Luciferin is stable for up to five (5) years.
- Luciferin is a light sensitive reagent and should be kept out of direct light as much as possible. We recommend that the luciferin be protected from light (e.g., covered with a light blocking material such as tin foil) during use, from the stock solution preparation until completion of the procedure.
- Repeated freeze thaw is not recommended.

\* For imaging protocols, D-Luciferin solution preparation, and solution storage, please see Preparation of IVISbrite™ D-Luciferin for In Vitro and In Vivo Bioluminescent Assays available on the Revvity website.

**This document has been generated electronically and does not require a signature**

**Limited Warranty:** This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. 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.

The Revvity logo, consisting of the word "revvity" in a lowercase, sans-serif font.

Revvity  
940 Winter Street  
Waltham, MA 02451 USA

(800) 762-4000  
www.revvity.com

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.