

OPTIMIZED D-LUCIFERIN

D-Luciferin K⁺ Salt Optimized for *In Vivo* Imaging

Catalog Number: 122796

Molecular Information: C₁₁H₇KN₂O₃S₂ (MW: 318.4)

Luciferin is essential in performing your bioluminescent assay – and the quality of your research will depend on the quality of your luciferin. That's why PerkinElmer, the world leader in the area of *in vivo* preclinical imaging, offers high quality Luciferin at an affordable price.

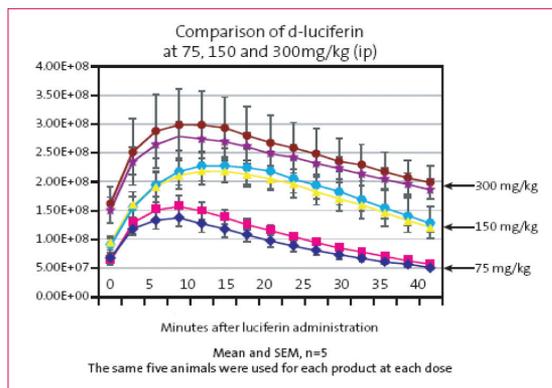
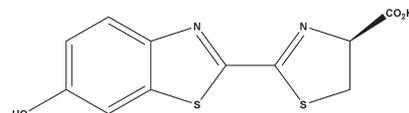
Luciferin is a chemical substance found in the cells of various bioluminescent organisms. When Luciferin is oxidized under the catalytic effects of luciferase and ATP, a bluish-green light is produced. Because the reaction is dependent on ATP, it allows researchers to determine the presence of energy or life. Firefly luciferin is a particularly good reporter for *in vitro* biophotonic imaging due to properties of its emission spectra.

Luciferin can be used in a number of ways. It can be used in a variety of *in vitro* assays, where the production of light can be monitored with either a luminometer or a scintillation counter.

It can also be used to monitor light production *in vivo*, and can be monitored with a PerkinElmer IVIS® Imaging System. Because luciferin can penetrate cell membranes, it allows transformed cells to be monitored for luciferase activity.

There are many considerations when choosing a luciferin substrate such as dosing and toxicity. It is important to know that you are using the highest quality Luciferin for your experiments. You might want to ask:

- Has your Luciferin been validated by PerkinElmer scientists?
- Is your Luciferin used exclusively by PerkinElmer physicists to calibrate IVIS Imaging Systems?
- Has your Luciferin been used in countless publications?



PerkinElmer *in vivo* imaging reagents are intended for animal research and not for use in humans.

Learn more at www.perkinelmer.com/invivoreagents

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2012, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.