

OPTIMIZED D-LUCIFERIN

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

Catalog Number: 122799

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.

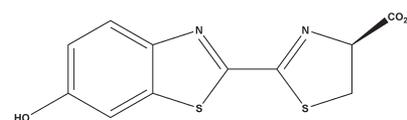
Luciferin is a chemical substance found in the cells of various bioluminescent organisms. When Luciferin is oxidized under the catalytic effects of luciferase and ATP, light is produced. Because the reaction is dependent on ATP, it allows researchers to determine the presence of energy or life. Firefly 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 PerkinElmer's IVIS® imaging platform. Because luciferin can penetrate cell membranes, it allows transformed cells to be monitored for luciferase activity. Luciferin can also be used with PerkinElmer's Bioware® Brite luciferase expressing oncology cell lines and our Red-FLuc lentiviral particles.

There are many considerations when choosing a luciferin substrate and it is important to know that you are using the highest quality Luciferin for your experiments.

Use the luciferin proven in countless peer reviewed publications!



For research use only. Not for use in diagnostic procedures.

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 ©2012-2016, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.