



Optimized D-Luciferin

IVISbrite™ D-Luciferin Potassium Salt Bioluminescent Substrate for *in vivo* imaging

Catalog number: 122799

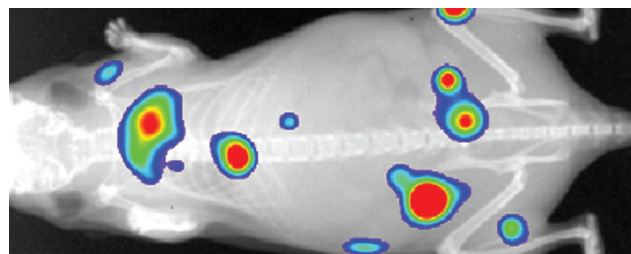
Molecular information: C₁₁H₇KN₂O₃S₂ (MW: 318.4)

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 Revvity, a 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* imaging due to properties of its emission spectra.

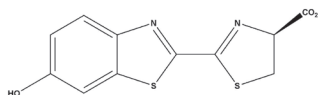
IVISbrite D-Luciferin can be used in a number of ways. It can be used in a variety of *in vitro* assays, where the production of light can be monitored with either a luminometer or a scintillation counter. It can also be used to monitor light production *in vivo*, and can be monitored with Revvity's 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 Revvity's IVISbrite™ luciferase expressing tumor cell lines and our IVISbrite Red F-luc lentiviral particles.

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



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!



Learn more at www.revvity.com