

## D-Luciferin, Potassium Salt

D-Luciferin is a substrate for the enzyme firefly luciferase. Chemiluminescence at 560 nm is generated when luciferin is oxidized by an ATP-dependent process, catalyzed by luciferase.



### Product attributes

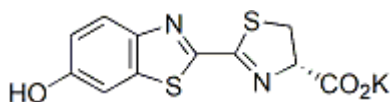
CAS number	115144-35-9
------------	-------------

## Product Description

D-Luciferin is a substrate for the enzyme firefly luciferase. Chemiluminescence at 560 nm is generated when luciferin is oxidized by an ATP-dependent process, catalyzed by luciferase. Luciferase is encoded by the *luc* gene, which is widely used as a reporter gene in a variety of cells. Because of the intrinsic low background of chemiluminescence technique, detection of the *luc* gene expression can be made at very low level. In addition, luciferin/luciferase has been used to measure  $10^{-15}$  molar quantity of ATP.

D-luciferin, potassium salt is a water-soluble form of D-luciferin.

- Light yellow solid soluble in water
- Store at  $-20^{\circ}\text{C}$  and protect from light
- $\text{C}_{11}\text{H}_7\text{KN}_2\text{O}_3\text{S}_2$
- Mwt: 318.4
- [115144-35-9]



## References

1. Appl Opt. 50(10), 1389 (2011).
2. PLoS ONE 5(10), e13545 (2010).
3. Antioxidants and Redox Signaling (2010) doi: 10.1089/ars.2010.3203
4. Biomaterials 31, 1429 (2010).
5. Methods in Molecular Biology (2010) DOI 10.1007/978-1-60761-901-7\_11
6. Plant Physiology 139, 1557 (2005).
7. Current Biology 15, 1352 (2005).
8. Cancer Research 64, 6684 (2004).
9. The Plant Cell 16, 769 (2004).
10. PNAS (2003) doi:10.1073/pnas.2435285100

This datasheet was generated on June 5, 2026 at 10:49:40 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/d-luciferin-potassium-salt/>