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

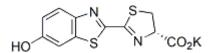
115144-35-9

Call us: 800-304-5357 Email: btinfo@biotium.com

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°C and protect from light
- C₁₁H₇KN₂O₃S₂
- 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 November 4, 2025 at 12:15:38 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/d-luciferin-potassium-salt/