D-Luciferin, Free Acid

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

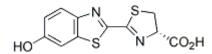
2591-17-5

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.

- White solid soluble in DMSO and slightly soluble in water
- Store at -20°C and protect from light
- C₁₁H₈N₂O₃S₂
- Mwt: 280
- [2591-17-5]



References

1. BioTechniques

7, 1116 (1989).

2. Lett. Appl. Microbiol. 1, 208 (1990).

This datasheet was generated on November 8, 2025 at 12:15:01 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/d-luciferin-free-acid/