

## 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.



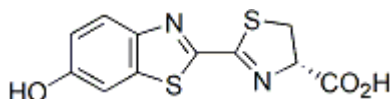
### Product attributes

CAS number	2591-17-5
------------	-----------

## 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
- $C_{11}H_8N_2O_3S_2$
- Mwt: 280
- [2591-17-5]



## References

1. BioTechniques  
7, 1116 (1989).
2. Lett. Appl. Microbiol. 1, 208 (1990).

This datasheet was generated on January 9, 2026 at 05:44:35 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/d-luciferin-free-acid/>