

D-Luciferin, Sodium 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

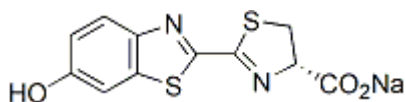
CAS number	103404-75-7
------------	-------------

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

- White solid soluble in water
- Store at -20°C and protect from light
- $\text{C}_{11}\text{H}_7\text{N}_2\text{NaO}_3\text{S}_2\cdot\text{H}_2\text{O}$
- Mwt: 302.3
- [103404-75-7]



This datasheet was generated on June 9, 2026 at 04:54:13 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/d-luciferin-sodium-salt/>