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

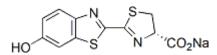
Call us: 800-304-5357 Email: btinfo@biotium.com

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
- C₁₁H₇N₂NaO₃S₂.H₂O
- Mwt: 302.3
- [103404-75-7]



This datasheet was generated on November 3, 2025 at 09:32:39 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/d-luciferin-sodium-salt/