

Beetle Luciferin, Potassium Salt (D-luciferin)

来自北美萤火虫 (*Photinus pyralis*) 和其它甲虫的萤光素酶基因, 通常作为报告基因用于研究真核细胞瞬时转染分析中的转录调控, 以及作为标记物研究真核细胞稳定转染。甲虫萤光素 (也称为 D- 萤光素) 作为单钾盐来合成, 是众多甲虫萤光素酶报告基因系统中的底物。D- 萤光素为那些选择自行配制检测试剂进行体外或体内萤光素酶活性检测的研究者而准备。

应用方向	产品引文
神经炎症 免疫调节	Jana Van Broeckhoven, <i>et al.</i> (2022) Macrophage-based delivery of interleukin-13 improves functional and histopathological outcomes following spinal cord injury. <i>J Neuroinflammation</i> . 19(1):102. PMID: 35488301
全细胞生物传感器	Maria Maddalena Calabretta, <i>et al.</i> (2022) A Luciferase Mutant with Improved Brightness and Stability for Whole-Cell Bioluminescent Biosensors and In Vitro Biosensing. <i>Biosensors (Basel)</i> . 12(9):742. PMID: 36140127
体外翻译和抑制	Dov Borovsky, <i>et al.</i> (2022) The Ribosome Is the Ultimate Receptor for Trypsin Modulating Oostatic Factor (TMOF). <i>Biomolecules</i> . 12(4):577. PMID: 35454167
非病毒性基因治疗	Haifei Gao, <i>et al.</i> (2021) CEBA: A new heterobifunctional reagent for plasmid DNA functionalization by click chemistry. <i>Int J Pharm</i> . 601:120566. PMID: 33812974
信号通路	P Thamarasseri Dhanagovind, <i>et al.</i> (2021) IL-6 Signaling Protects Zebrafish Larvae during Staphylococcus epidermidis Infection in a Bath Immersion Model. <i>J Immunol</i> . 207(8): 2129-2142. PMID: 34544800
植物育种	Mengjia Zhuang, <i>et al.</i> (2021) The wheat SHORT ROOT LENGTH 1 gene TaSRL1 controls root length in an auxin-dependent pathway. <i>J Exp Bot</i> . 72(20):6977-6989. PMID: 34328188
小鼠肾上腺诱导物 Dkk3/REIC	Hirofumi Fujita, <i>et al.</i> (2020) Dkk3/REIC, an N-glycosylated Protein, Is a Physiological Endoplasmic Reticulum Stress Inducer in the Mouse Adrenal Gland. <i>Acta Med Okayama</i> . 74(3):199-208. PMID: 32577017
嗜盐微生物基因调控	Chris R. Davis, <i>et al.</i> (2020) A bioluminescent reporter for the halophilic archaeon Haloferax volcanii. <i>Extremophiles</i> . 24(5): 773–785. PMID: 32749548
MAPK 信号通路	Charles S Goldsmith, <i>et al.</i> (2019) Inhibition of p38 MAPK activity leads to cell type-specific effects on the molecular circadian clock and time-dependent reduction of glioma cell invasiveness. <i>BMC Cancer</i> . 19(1). PMID: 30674294

Beetle Luciferin 使用引文

应用方向	产品引文
酿酒酵母代谢节律	Archana Krishnamoorthy, <i>et al.</i> (2015) Dual-Color Monitoring Overcomes the Limitations of Single Bioluminescent Reporters in Fast-Growing Microbes and Reveals Phase-Dependent Protein Productivity during the Metabolic Rhythms of <i>Saccharomyces cerevisiae</i>. <i>Appl Environ Microbiol.</i> 81(18): 6484–6495. PMID: 26162874
生物节律	J. Brian Robertson, <i>et al.</i> (2008) Real-time luminescence monitoring of cell-cycle and respiratory oscillations in yeast. <i>Proc Natl Acad Sci U S A.</i> 105(46): 17988–17993. PMID: 19004762
光动力疗法 (PDT) 与癌症治疗	Theodossiou, T., <i>et al.</i> (2003) Firefly luciferin-activated Rose Bengal: In vitro photodynamic therapy by intracellular chemiluminescence in transgenic NIH 3T3 cells. <i>Cancer Res.</i> 63, 1818-1821. PMID: 12702568
植物生长调节剂的调控机制研究	Bouquin, T., <i>et al.</i> (2001) Control of specific gene expression by gibberellin and brassinosteroid. <i>Plant Physiol.</i> 127, 450-458. PMID: 11598220
癌基因 Myc 与视网膜母细胞瘤	Santoni-Rugiu, E., <i>et al.</i> (2000) Involvement of Myc activity in a G(1)/S-promoting mechanism parallel to the pRb/E2F pathway. <i>Mol. Cell. Biol.</i> 20(10), 3497-3509. PMID: 10779339

如需了解更多萤光素酶底物，请点击下方链接：

wechat.promega.com.cn

产品	规格	目录号
Beetle Luciferin, Potassium Salt	5mg	E1601
Beetle Luciferin, Potassium Salt	50mg	E1602
Beetle Luciferin, Potassium Salt	250mg	E1603
Beetle Luciferin, Potassium Salt	1g	E1605



欢迎关注 Promega 生命科学