

Luciferase *Recombinant firefly*

Ordering Information:	Storage Conditions:
Product Number: 12500 (1mg/1mL)	Keep at -20 °C. Expiration date is one year from the date of receipt.

General Properties

CAS Number: 61970-00-1

Molecular Weight: ~60 kDa

Form: liquid (1mg/mL)

Measuring Range: ≤1 femtomole ATP sensitivity (using 0.2 µg of luciferase and suitably sensitive liquid scintillation counters or luminometers)

Biological Applications

This recombinant firefly luciferase is an enzyme expressed from a cloned gene from the North American firefly (*Photinus pyralis*) that provides the reliability and dependability needed for performing research or producing kits using bioluminescence reagents to detect ATP or luciferin substrates. This recombinant enzyme can potentially eliminate the possibility of seasonal and regional variability that may be found in luciferase purified from natural sources.

The reaction of this enzyme with luciferin, ATP, and O₂ results in the emission of light. Luciferase can be used to detect trace amounts of ATP. Firefly luciferase is also one of the most commonly utilized reporter genes for the study of gene expression. The bioluminescent reaction catalyzed by luciferase is one of the most sensitive analytical tools for measuring gene expression. Less than or equal to one femtomole of ATP can be detected by using 0.2 µg of luciferase.

References

1. McElroy, W.D. (1947) The Energy Source for Bioluminescence in an isolated System. *Proc. Natl. Acad. Sci. USA* **33**,342.
2. de Wet JR, Wood KV, Helinski DR, DeLuca M, (1985) Cloning of firefly luciferase cDNA and the expression of active luciferase in *Escherichia coli*, *Proc. Natl. Acad. Sci USA* **82**,7870-7873.
3. Khan, H.A. (2003) Bioluminometric assay of ATP in mouse brain: Determinant factors for enhanced test sensitivity, *J. Bioscience* **28**, 379-382.
4. Drew, B and C. Leeuwenburgh (2003) Method for measuring ATP production in isolated mitochondria: ATP production in brain and liver mitochondria fo Fischer-344 rats with age and caloric restriction, *Am J. Physiol. Regul. Integr. Comp. Physiol.*, **285**, R1260-R1268.
5. Hara, K. Y. and Mori, H. (2006) An efficient method for quantitative determination of cellular ATP synthetic activity, *J Biomol Screen* **11**, 310-7.
6. Sun, Y. and Chai, T. C. (2006) Augmented extracellular ATP signaling in bladder urothelial cells from patients with interstitial cystitis *Am J Physiol Cell Physiol* **290**, C27-34.
7. Stanley, P.E., A review of bioluminescent ATP techniques in rapid microbiology. *J. Biolumin. Chemilumin.* **4**, 375, (1989)
8. Rajgopal, S., and Vijayalakshmi, M.A. *Enzyme Microb. Technol.* **6**, 482-489, (1984)
9. Chappelle, E.W., et al., Determination of bacterial content in fluids., *Meth. Enzymol.* **57**, 65-72, (1978)
10. Kricka, L.J., Clinical and biochemical applications of luciferases and luciferins. *Anal. Biochem.* **175**, 14-21, (1988)

11. Liu L, Hastings JW. (2006) Two different domains of the luciferase gene in the heterotrophic dinoflagellate *Noctiluca scintillans* occur as two separate genes in photosynthetic species. *Proc Natl Acad Sci U S A*.
12. Emamzadeh AR, Hosseinkhani S, Sadeghizadeh M, Nikkhah M, Chaichi MJ, Mortazavi M. (2006) cDNA cloning, expression and homology modeling of a luciferase from the firefly *Lampyroidea maculata*. *J Biochem Mol Biol*, 39, 578.
13. Viviani VR, Ohmiya Y. (2006) Bovine serum albumin displays luciferase-like activity in presence of luciferyl adenylate: insights on the origin of protoluciferase activity and bioluminescence colours. *Luminescence*, 21, 262.
14. Schipper ML, Patel MR, Gambhir SS. (2006) Evaluation of firefly luciferase bioluminescence mediated photodynamic toxicity in cancer cells. *Mol Imaging Biol*, 8, 218.
15. Oba Y, Sato M, Inouye S. (2006) Cloning and characterization of the homologous genes of firefly luciferase in the mealworm beetle, *Tenebrio molitor*. *Insect Mol Biol*, 15, 293.
16. Palomba S, Berovic N, Palmer RE. (2006) Bioluminescence of monolayers of firefly luciferase immobilized on graphite. *Langmuir*, 22, 5451.
17. Schipper ML, Patel MR, Gambhir SS. (2006) Evaluation of Firefly Luciferase Bioluminescence Mediated Photodynamic Toxicity in Cancer Cells. *Mol Imaging Biol*.
18. Fraga H, Fernandes D, Novotny J, Fontes R, Esteves da Silva JC. (2006) Firefly luciferase produces hydrogen peroxide as a coproduct in dehydroluciferyl adenylate formation. *Chembiochem*, 7, 929.

Disclaimer: *This product is for research use only and is not intended for therapeutic or diagnostic applications. Please contact our technical service representative for more information.*