



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

# Luciferase (Firefly, Recombinant)

(Cat. # 786-1308, 786-1309)



think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)

## INTRODUCTION

Firefly luciferase is a unique photoprotein used in bioluminescence for high sensitive detection and quantification of ATP and as a reporter for studying gene function and regulation.

Luciferase (Firefly, Recombinant) is recombinant 62 kDa protein expressed in *E. coli*.

## ITEM(S) SUPPLIED

Cat. #	Description	Size
786-1308	Luciferase (Firefly, Recombinant)	100 µg
786-1309	Luciferase (Firefly, Recombinant)	0.5 mg

## STORAGE CONDITIONS

Luciferase (Firefly, Recombinant) is shipped with dry ice. Store at -20°C in small one time use aliquots. When stored properly as advised the enzyme is stable for 1 year. For long term storage store aliquots at -70°C

## SPECIFICATIONS

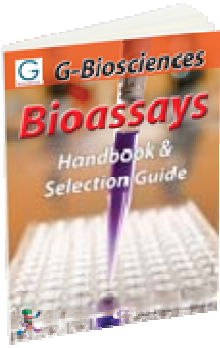
- **Single band purity**, >99 % by SDS-PAGE
- **High specific activity**:  $\geq 1 \times 10^{10}$  units/mg.
- **Molecular weight**: 62kDa
- **Protein amount**: determined based on absorbance at 280 nm and molecular extinction co-efficient of Luciferase
- **Source**: *E. coli*.

## PREPARATION BEFORE USE

Luciferase (Firefly, Recombinant) is a temperature sensitive enzyme. Thaw the enzyme on ice and keep it on ice when using it immediately. Avoid repeated thawing. For making enzyme dilutions it is recommended to add 1 % carrier protein such as BSA to the dilution buffer. Do not vortex enzyme any time. Mixing should be done gently using a pipette.

## RELATED PRODUCTS

Download our Bioassay Handbook.



<http://info2.gbiosciences.com/complete-bioassay-handbook>

For other related products, visit our website at [www.GBiosciences.com](http://www.GBiosciences.com) or contact us.



[www.GBiosciences.com](http://www.GBiosciences.com)