

# STOP

Before using this product, please read the Limited Use License statement below:

## **Important Limited Use License information for pSELECT-zeo-Lucia**

The purchase of the pSELECT-zeo-Lucia vector conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes.

The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes.

Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic, or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research.

If the purchaser is unwilling to accept the limitations of this limited use statement, InvivoGen is willing to accept return of the product with a full refund. The product must be returned in resaleable condition. For information on purchasing a license to this product for purposes other than research, contact 10515 Vista Sorrento Parkway San Diego, CA 92121 USA. Tel: 858-457-5873 Fax: 858-457-5843.

---

### TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)



# pSELECT-zeo-Lucia

A secreted luciferase reporter gene system selectable with Zeocin®

Catalog code: psetz-lucia  
<https://www.invivogen.com/lucia-gene>

For research use only

Version 23A19-MM

## PRODUCT INFORMATION

### Contents:

- 20 µg of pSELECT-zeo-Lucia provided as lyophilized DNA
- 1 ml of Zeocin® (100 mg/ml)
- 1 tube of QUANTI-Luc™ 4 Reagent, a Lucia luciferase detection reagent (sufficient to prepare 25 ml)

### Storage and Stability:

- Product is shipped at room temperature.
- Lyophilized DNA should be stored at -20°C.
- Resuspended DNA should be stored at -20°C and is stable for up to 1 year.
- Store Zeocin® at 4°C or at -20°C. The expiry date is specified on the product label.
- Store QUANTI-Luc™ 4 Reagent at -20°C. After preparation, the working solution is stable for 48 hours at 4°C and for 1 month at -20°C. Prepare aliquots to avoid repeated freeze-thaw cycles. Note: This product is photosensitive and should be protected from light.

### Quality Control:

- Plasmid construct has been confirmed by restriction analysis and full-length ORF sequencing.
- Plasmid DNA was purified by ion exchange chromatography.

## GENERAL PRODUCT USE

InvivoGen provides the Lucia luciferase reporter gene in the pSELECT-zeo plasmid. The plasmid pSELECT-zeo-Lucia can be used *in vivo* and *in vitro* to transfect mammalian cells stably or transiently. Lucia luciferase gene expression is driven by the EF-1α/HTLV composite promoter that combines the elongation factor 1 alpha core promoter and the 5' untranslated region of the Human T-cell Leukemia Virus. The pSELECT-zeo-Lucia plasmid contains the Zeocin® resistance marker for selection in both mammalian cells and bacteria.

Lucia luciferase is a completely novel and optimized luciferase with strong bioluminescent activity. It is expressed by a synthetic gene designed on natural secreted luciferase genes from marine copepods. Lucia luciferase is a secreted coelenterazine-utilizing luciferase that generates 1000-fold higher bioluminescent signal compared to the commonly used Firefly and *Renilla* luciferases. Lucia luciferase is designed for high and prolonged expression in mammalian cells.

Lucia luciferase activity can be determined using QUANTI-Luc™ 4 Lucia/Gaussia, a Lucia and Gaussia luciferase detection reagent.

## PLASMID FEATURES

### First expression cassette

- hEF1-HTLV prom is a composite promoter comprising the Elongation Factor-1α (EF-1α) core promoter<sup>1</sup> and the R segment and part of the U5 sequence (R-U5') of the Human T-Cell Leukemia Virus (HTLV) Type 1 Long Terminal Repeat<sup>2</sup>. The EF-1α promoter exhibits a strong activity and yields long lasting expression of a transgene *in vivo*. The R-U5' has been coupled to the EF-1α core promoter to enhance stability of RNA.

- Lucia luciferase is a synthetic CpG-free gene that codes for a secreted coelenterazine-utilizing luciferase. ORF size (from the ATG to the stop codon): 634 bp.

## PLASMID FEATURES

- **SV40 pAn:** The Simian Virus 40 late polyadenylation signal enables efficient cleavage and polyadenylation reactions resulting in high levels of steady-state mRNA<sup>3</sup>.

- **pMB1 Ori** is a minimal *E. coli* origin of replication with the same activity as the longer Ori.

### Second expression cassette

- **CMV enh/prom:** The human cytomegalovirus immediate-early gene 1 promoter/enhancer was originally isolated from the Towne strain and was found to be stronger than any other viral promoters

- **EM7** is a bacterial promoter that enables the constitutive expression of the antibiotic resistance gene in *E. coli*.

- **Sh ble:** Resistance to Zeocin® is conferred by the *Sh ble* gene from *Streptoalloteichus hindustanus*. The *Sh ble* gene is driven by the CMV promoter/enhancer in tandem with the bacterial EM7 promoter. Therefore, Zeocin® can be used to select stable mammalian cells transfectants and *E. coli* transformants.

- **βGlo pan:** The human beta-globin 3'UTR and polyadenylation sequence allows efficient arrest of the transgene transcription<sup>4</sup>.

1. Kim DW. *et al.*, 1990. Use of the human elongation factor 1α promoter as a versatile and efficient expression system. *Gene* 91(2):217-23. 2. Takebe Y. *et al.*, 1988. SR alphapromoter: an efficient and versatile mammalian cDNA expression system composed of the simian virus 40 early promoter and the R-U5 segment of human T-cell leukemia virus type 1 long terminal repeat. *Mol Cell Biol.* 8(1):466-72. 3. Carswell S. & Alwine JC., 1989. Efficiency of utilization of the simian virus 40 late polyadenylation site: effects of upstream sequences. *Mol Cell Biol.* 9(10):4248-58. 4. Yu J. & Russell JE., 2001. Structural and functional analysis of an mRNP complex that mediates the high stability of human β-globin mRNA. *Mol Cell Biol.* 21(17):5879-88.

## METHODS

### Plasmid resuspension

Quickly spin the tube containing the lyophilized plasmid to pellet the DNA. To obtain a plasmid solution at 1 µg/µl, resuspend the DNA in 20 µl of sterile H<sub>2</sub>O. Store resuspended plasmid at -20°C.

### Plasmid amplification and cloning

Plasmid amplification and cloning can be performed in commonly used laboratory *E. coli* strains, such as DH5α.

### Zeocin® usage

This antibiotic can be used for *E. coli* at 25-50 µg/ml in liquid or solid media and at 50-200 µg/ml to select Zeocin®-resistant mammalian cells.

### QUANTI-Luc™ 4 Reagent

1. Dilute the total volume of the 20X tube (1.25 ml) of QUANTI-Luc™ 4 Reagent into 23.75 ml sterile water to obtain 25 ml of working solution.
2. Vortex very briefly (a few seconds).
3. Use the working solution immediately or store until required for use.

## RELATED PRODUCTS

Product	Description	Cat. Code
QUANTI-Luc™ 4 Lucia/Gaussia Zeocin®	Luminescence detection kit Selection antibiotic	rep-qlc4lg1 ant-zn-1

## TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)



# QUANTI-Luc™ 4 Reagent

A coelenterazine-based luminescence assay reagent

<https://www.invivogen.com/ quanti-luc>

For research use only

Version 23A16-MM

## PRODUCT INFORMATION

### Contents

- 1 tube of QUANTI-Luc™ 4 Reagent (20X)

One tube of QUANTI-Luc™ 4 Reagent is sufficient for 5 x 96-well plates (25 ml standard Flash/end-point detection).

**Note:** This sample cannot be sold separately from the QUANTI-Luc™ 4 Lucia/Gaussia kit.

QUANTI-Luc™ 4 Lucia/Gaussia comprises two liquid components:

- QUANTI-Luc™ 4 Reagent 20X (coelenterazine substrate)
- QUANTI-Luc™ 4 Stabilizer 25X (optimized Glow assay reagent)

Find more information at <https://www.invivogen.com/ quanti-luc>.

### Storage and Stability

- Store QUANTI-Luc™ 4 Reagent at -20°C for up to 12 months.
- After preparation, the working solution is stable for 48 hours at 4°C and 1 month at -20°C. Prepare aliquots to avoid repeated freeze-thaw cycles.

**Note:** This product is photosensitive and should be protected from light.

### Quality Control

Each lot is thoroughly tested to ensure the absence of lot-to-lot variation.

- Physicochemical characterization (pH, appearance).
- Functional assays using recombinant Lucia protein or reporter cells.

## DESCRIPTION

QUANTI-Luc™ 4 Reagent is a component of the QUANTI-Luc™ 4 Lucia/Gaussia kit. It contains the coelenterazine substrate for the detection of secreted Lucia or Gaussia activity in live-cell supernatants, and of intracellular Renilla after cell lysis. The light signal produced correlates to the amount of luciferase protein expressed. It is quantified using a luminometer and expressed as relative light units (RLUs).

## METHODS

### Preparation of QUANTI-Luc™ 4 Reagent working solution (1X):

1. Dilute the total volume of the 20X tube (1.25 ml) of Reagent into 23.75 ml of sterile water to obtain 25 ml of working solution.
2. Vortex **very briefly** (a few seconds).
3. Use the working solution immediately or store until required for use. QUANTI-Luc™ 4 Reagent working solution can be stored for 48 hours at 4°C or 1 month at -20°C.

### Flash detection of luciferase activity from cell culture medium:

To obtain **end-point readings** using a luminometer **with an injector**.

1. Set the luminometer with the following parameters: 50 µl of injection, end-point measurement with a 4 second start time and 0.1 second reading time.
2. Pipet 20 µl of sample per well into a 96-well white (opaque) or black plate, or a luminometer tube.
3. Prime the injector with QUANTI-Luc™ 4 Reagent 1X and proceed **immediately** with the measurement.

To obtain **end-point readings** using a luminometer **without injectors**.

1. Set the luminometer with a 0.1 second reading time.
2. Pipet 20 µl of sample per well into a 96-well white (opaque) or black plate, or a luminometer tube.
3. Add 50 µl of QUANTI-Luc™ 4 Reagent 1X to each well or tube.
4. Gently tap the plate several times to mix (do **not** vortex).
5. Proceed **immediately** with the measurement.

## RELATED PRODUCTS

Product	Cat. Code
QUANTI-Luc™ 4 Lucia/Gaussia Kit	
500 tests	rep-qlc4lg1
2 x 500 tests	rep-qlc4lg2
5 x 500 tests	rep-qlc4lg5

### TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

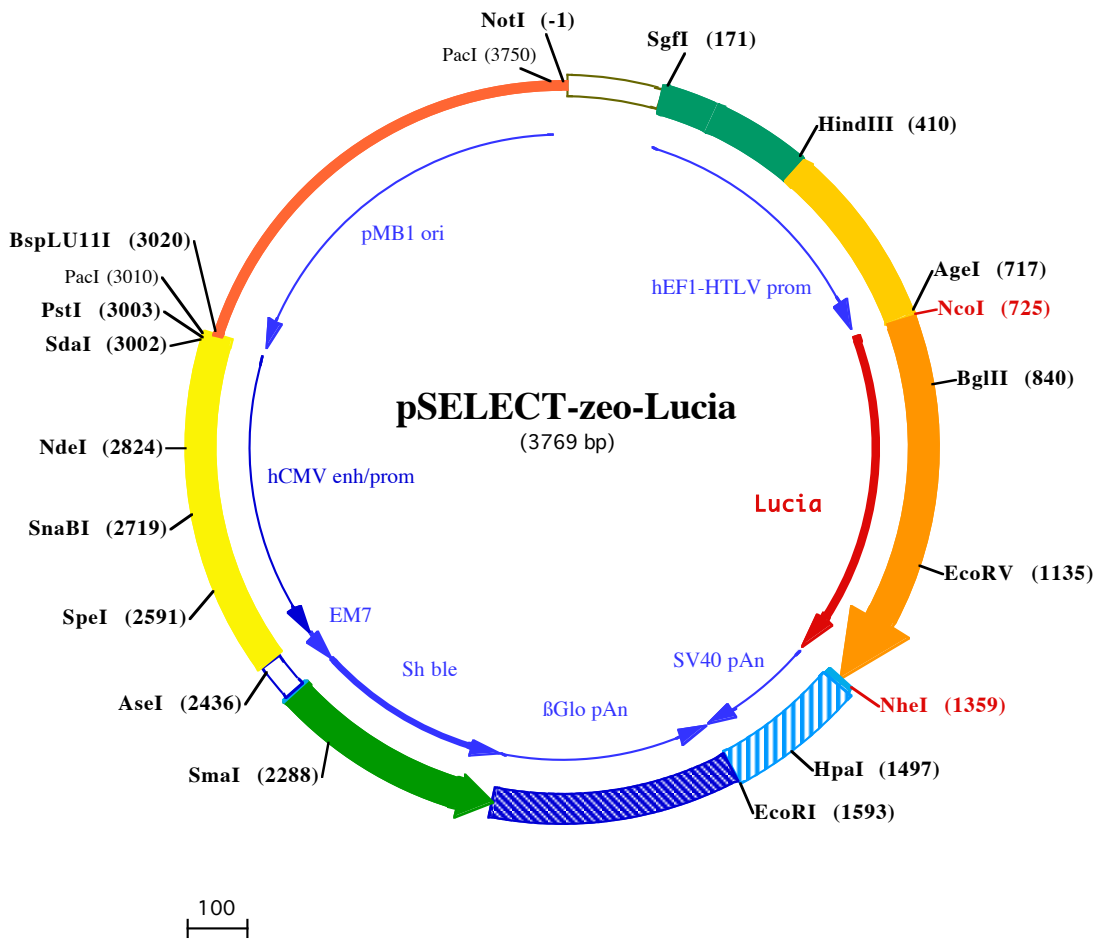
InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)





**NotI (-1)**  
1 CGCGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACA  
**SgfI (171)**  
101 AAACAACTAGCAAATAGGCTGTCCCGAGTGAAGTGCAGGTGCCAGAACATTTCTCTATCGAAGGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGCA  
201 GAGCGCACATCGCCACAGTCCCGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGTGCCTAGAGAAGTGGCGGGGTAACCTGGGAAAGTGATG  
301 TCGTGTACTGGCTCCGCTTTTTCCCGAGGTGGGGGAGAACCCTATATAAGTCAGTAGTCGCGCGTGAACGTTCTTTTTCCCAACGGGTTTCCGCCAC  
**HindIII (410)**  
401 AACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCCTTACCGCGCCGCCCTACCTGAGGCGCCATCCACGCCGTTGAGTGCCTTCTGCCGCT  
501 CCCGCTGTGGTGCCTCCTGAATGCGTCCCGCTAGTAAGTTAAAGCTCAGTGCAGACCGGCTTTGTCCGGCGCTCCCTTGAGCCTACCTA  
601 GACTCAGCGGCTCCACGCTTTCCTGACCTGCTTGTCTCAACTCTACGCTTTGTTCGTTTTCTGTTCTGCGCGTTACAGATCCAAGCTGTGACC  
**AgeI (717)**  
701 GCGCCTACCTGAGATCACCGGTACCATGGAATCAAGGTGCTGTTGCCCTCATCTGTATTGCTGTTGCTGAGGCAAACCCACTGAAATCAATGAAG  
1► M E I K V L F A L I C I A V A E A K P T E I N E  
**AgeI (717)**  
1► M E I K V L F A L I C I A V A E A K P T E I N E  
**BglII (840)**  
801 ACCTCAATATAGCTGTGGCTCCAACCTTCCACCACAGATCTTGAGACTGACCTGTTCCACCACTGGGAGACCATGAATGTGATTAGCACTGACAC  
25► D L N I A A V A S N F A T T D L E T D L F T N W E T M N V I S T D T  
901 AGAGCAGGTGAACACAGATGCTGACAGGGCAAGCTGCCTGGCAAAAACCTCCCCAGATGCTCTGAGGGAGCTGGAGGCAATGCCAGAAGGGCTGGT  
58► E Q V N T D A D R G K L P G K K L P P D V L R E L E A N A R R A G  
1001 TGCACAAGAGGCTGCCTCATTGCTTCCACATTAAGTGCACCCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCCACACTTATGAAGGTGAAAAGG  
92► C T R G C L I C L S H I K C T P K M K K F I P G R C H T Y E G E K  
**EcoRV (1135)**  
1101 AGTCTGCTCAGGGAGGATTGGAGAGCAATTTGATATCCAGAGATTCTGGCTTCAAGGATAAGGAGCCACTGGACCAGTTTATTGCTCAAGTGGGA  
125► E S A Q G G I G E A I V D I P E I P G F K D K E P L D Q F I A Q V D  
1201 CCTCTGTGCTGATTGCACCACTGGCTGTCTGAAGGGCTTGCACATGTCAGTGTCTGACTCTCTGAAGAAGTGGCTTCCCGAGAGGTGACCACTTT  
158► L C A D C T T G C L K G L A N V Q C S D L L K K W L P Q R C T T F  
**NheI (1359)**  
1301 GCCAGCAAGATTGAGGTAGGGTGGACAAAATCAAGGTCTGGCTGGGACAGATGATAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTGG  
192► A S K I Q G R V D K I K G L A G D R •  
**HpaI (1497)**  
1401 ACAAACCACTAGAAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGAATAAACAGTT  
**EcoRI (1593)**  
1501 AACAAACAATTCGATTCATTTATGTTTCAGGTTGAGGGGAGGTGGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAATTC  
1601 AAAATACAGCATAGCAAATTTAACCCTCAAAATCAAGCCTTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAAT  
1701 GTGCATTAGCTGTTTGCAGCCTCACCTTCTTATGAGTGAATTAAGATAGTGTATTTCCCAAGGTTGAACTAGCTCTTCAATTTCTTATGTTTAAA  
1801 TGCACTGACCTCCACATTCCTTTTTAGTAAAATATTAGAAATAATTTAAATACATCATTGCAATGAAATAAATGTTTTTATTAGGAGAATCCAG  
1901 ATGCTCAAGGCCCTTATAATATCCCCAGTTTGTAGTGGACTTAGGAAACAAGGAACCTTAAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAG  
2001 CTTATCCTCAGTCTGCTCCTGTCACAAAGTGCACGAGTGGCGGCGGGTGCAGGCGGCAACTCCCGCCCCACGGCTGCTCGCGATCTCGGT  
125► D Q E E A V F H V C N G A P D R L A F E R G W P Q E G I E T  
2101 CATGGCCGGCCGGAGGCTCCCGAAGTTCGTGGACAGACCTCCGACCACTCGCGTACAGCTCGTCCAGGCGCGCACCCACACCCAGGCCAGGGTG  
94► M A P G S A D R F N T S V E S W E A Y L E D L G R V W V W A L T  
**SmaI (2288)**  
2201 TTGTCCGGCACCACTGGTCTGGACCGCTGATGAACAGGGTGCAGTCTGCTCCCGACACACCGCGAAGTCTGCTCCACGAAGTCCCGGAGAAC  
60► N D P V V Q D Q V A S I F L T V D D R V V G A F D D E V F D R S F G  
2301 CGAGCCGGTCCGAGAACTGACCGCTCCGCGACGTCGCGCGGGTGAACCGGAACGGCACTGGTCAACTTGGCCATGATGGCCCTCTATAGTG  
27► L R D T W F E V A G A V D R A T L V P V A S T L K A M  
**AseI (2436)**  
2401 AGTCGTATTATACTATGCCGATATCTATGCCGATGATTAATGTCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTCACTAAACGAGCTCT  
**SpeI (2591)**  
2501 GCTTATATAGACCTCCACCGTACACGCTACCGCCATTGGCGTCAATGGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCA  
2600 AAACAACTCCCATGACGTCATGGGGTGGAGACTTGGAAATCCCGTGAGTCAAACCGCTATCCAGCCATTGATGTACTGCCAAAACCGCATCATC  
**SnaBI (2719)**  
2700 ATGGTAATAGCGATGACTAATACGTAGATGACTGCCAAGTAGGAAAGTCCCATAAAGTTCATGACTGGGCATAATGCCAGGCGGGCCATTACCGTCAT  
**NdeI (2824)**  
2800 TGACGTCAATAGGGGGCTACTTGGCATATGATACACTTGTACTGCTGCAAGTGGGCGAGTTTACCGTAAATACTCCACCCATTGACGTCAATGAAAGT  
2900 CCCTATTGGCGTTACTATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGCTGTTGGCGGTCAGCCAGCGGGCCATTACCGTAAGTTATGTA  
**PacI (3010)**  
**PstI (3003)**  
**SdaI (3002)**  
**BspLU11I (3020)**  
3000 ACGCTGCGAGGTTAAATAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGCCGCTTGTGGCGTTTTTCCATAGGCTCCGC  
3098 CCCCTGACGAGCATCAAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGCGTTTCCCTGGAAAGCTCCCTCG  
3198 TGGCTCTCCTGTTCCGACCTGCGCTTACCGATACCTGTCCGCTTTCTCCCTTCGGAAGCGTGGCGTTTCTCATAGCTCACGCTGTAGGTATCT

3298 CAGTTCGGTGTAGGTCGTTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTTCAGCCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCC  
3398 AACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGT  
3498 GGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAA  
3598 ACAAAACACCGCTGGTAGCGGTGTTTTTTTTGTTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTCTACGGGG  
PacI (3750)  
3698 TCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATC A

# QUANTI-Luc™

A coelenterazine-based luminescence assay reagent

Catalog code: rep-qlc1, rep-qlc2

<https://www.invivogen.com/quantiluc>

For research use only

Version 19A04-MM

## PRODUCT INFORMATION

### Contents

QUANTI-Luc™ is provided as packs of individually sealed pouches.

- rep-qlc1: 2 pouches of QUANTI-Luc™
- rep-qlc2: 5 pouches of QUANTI-Luc™

Each pouch contains everything needed to prepare 25 ml of reagent allowing the preparation of 500 wells of a 96-well plate.

### Storage and Stability

- Store QUANTI-Luc™ pouches at -20°C for 12 months.
- Reconstituted QUANTI-Luc™ is stable for 1 week at 4°C and for 1 month at -20°C. Prepare aliquots to avoid repeated freeze-thaw cycles.

**Note:** This product is photosensitive and should be protected from light.

## DESCRIPTION

QUANTI-Luc™ is an assay reagent containing all the components required to quantitatively measure the activity of Lucia luciferase and other coelenterazine-utilizing luciferases. QUANTI-Luc™ contains the coelenterazine substrate and stabilizing agents for the luciferase reaction. The light signal produced is quantified using a luminometer and expressed as relative light units (RLU). The signal produced correlates to the amount of luciferase protein expressed, indicating promoter activity in the reporter assay.

QUANTI-Luc™ is optimized for use with Lucia luciferase reporter cell lines. Lucia luciferase is a secreted coelenterazine luciferase encoded by a synthetic gene. As Lucia luciferase is secreted, it can be directly measured in the cell culture medium using bioluminescent assays.

InvivoGen provides a recombinant Lucia luciferase protein (see Related Products) which is a positive control for QUANTI-Luc™. A dilution series of the recombinant Lucia luciferase protein can also be used to determine the linear range of the assay.

## METHODS

### Preparation of QUANTI-Luc™

1. Pour the pouch contents into a 50 ml screw cap tube.
2. Add 25 ml of sterile water.
3. Swirl product gently until powder is completely dissolved.
4. Use QUANTI-Luc™ assay solution immediately or store until required for use. Reconstituted QUANTI-Luc™ can be stored for 1 week at 4°C and for 1 month at -20°C. Prepare aliquots to avoid repeated freeze-thaw cycles.

**Note:** This product is photosensitive and should be protected from light.

### Detection of luciferase activity from cell culture medium

To obtain **end-point readings** using a luminometer **with an injector**.

1. Set the luminometer with the following parameters: 50 µl of injection, end-point measurement with a 4 second start time and 0.1 second reading time.
2. Pipet 10-20 µl of sample per well into a 96-well white (opaque) or black plate, or a luminometer tube.
3. Prime the injector with the QUANTI-Luc™ assay solution and proceed **immediately** with the measurement.

To obtain **end-point readings** using a luminometer **without injectors**.

1. Set the luminometer with a 0.1 second reading time.
2. Pipet 10-20 µl of sample per well into a 96-well white (opaque) or black plate, or a luminometer tube.
3. Add 50 µl of QUANTI-Luc™ assay solution to each well or tube.
4. Gently tap the plate several times to mix (do **not** vortex).
5. Proceed **immediately** with the measurement.

## RELATED PRODUCTS

Product	Catalog Code
QUANTI-Luc™ Gold (For standard and HTS assays)	rep-qlcg1
pSelect-zeo-Lucia™ (expression plasmid)	psetz-lucia
Recombinant Lucia luciferase protein	rec-lucia
<b>Reporter Cells</b>	
THP1-Dual™ (IRF-Lucia/NF-κB-SEAP) Cells	thpd-nfis
THP1-Lucia™ NF-κB Cells	thp1-nfkb

For a complete list of InvivoGen's Lucia luciferase Reporter Cell Lines visit <https://www.invivogen.com/lucia-reporter-cells>.

## TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873  
InvivoGen USA (International): +1 (858) 457-5873  
InvivoGen Europe: +33 (0) 5-62-71-69-39  
InvivoGen Hong Kong: +852 3622-3480  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)

