



LightSwitch™ Dual Assay System

DA010 (100 assays)

DESCRIPTION

The LightSwitch Dual Assay System is used to normalize for variation between transfection replicates in cell lines that are particularly difficult to transfect by utilizing a co-transfection control. Please read the [Do I Need a Co-transfection Control?](#) technical note to help determine whether your project needs a co-transfection control. Please note that the use of a co-transfection control is not sufficient for normalizing between conditions, treatments, or cell lines. For normalizing between experimental conditions, use LightSwitch promoter or 3'UTR control constructs.

All LightSwitch promoter and 3'UTR reporter constructs utilize RenSP, a novel luciferase developed by SwitchGear Genomics for reporter gene assays. The optimized luciferase protein has a half-life of ~1 hour enabling a detailed analysis of kinetic responses with a highly robust signal. RenSP reporter signal is detected using LightSwitch Luciferase Assay Reagents. The Cypridina TK control construct (pTK-CLuc Vector) contains the Cypridina luciferase reporter gene driven by an HSV-TK constitutive promoter. Its signal is detected using the BioLux® Cypridina Luciferase Assay Kit developed by New England Biolabs.

The Cypridina TK control construct and LightSwitch GoClone experimental constructs contain different luciferase genes that use different substrates, so there is no cross-reaction between the two reporter genes and their substrates. As an added advantage, the Cypridina reporter protein is secreted into the culture media while the RenSP reporter protein remains within the cell. Normalizing the LightSwitch signal using the Cypridina reporter signal may decrease variation between transfection replicates in difficult cell lines.

See the [LightSwitch Luciferase Assay System Manual](#) and [BioLux® Cypridina Luciferase Assay Kit Manual](#) for more details.

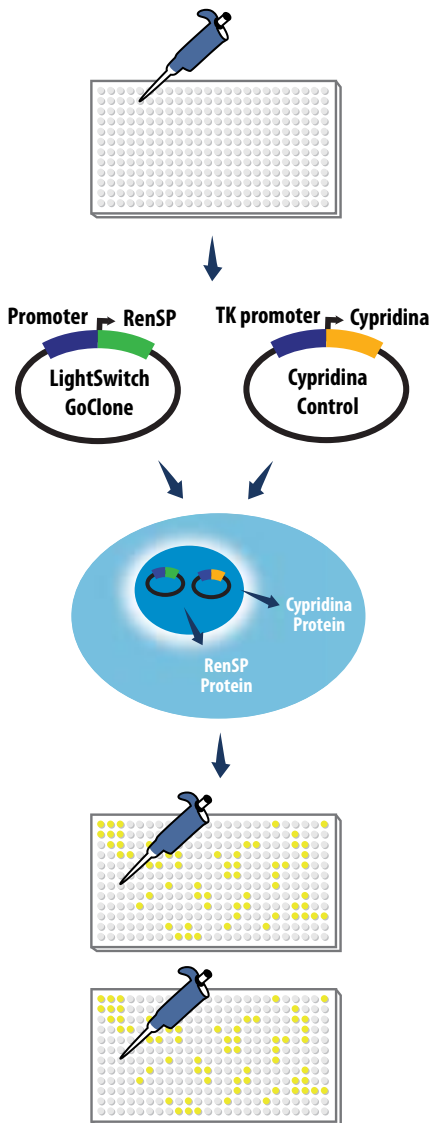
Components & Storage (100 assays)

1 tube	LightSwitch Assay Substrate (100X)	-20C for 6 months
150 uL	LightSwitch Substrate Solvent	room temp
10 mL	LightSwitch Assay Buffer	-20C for 6 months
1 tube	BioLux Cypridina Luciferase Substrate	-20C for 6 months
500 uL	BioLux Cypridina Luciferase Substrate Solvent	room temp
5 mL	BioLux Cypridina Luciferase Assay Buffer (1X)	-20C for 6 months

! Additional SwitchGear Product Needed:

Cypridina TK Control Construct (Cat # SN0322S)

(Cypridina luciferase gene regulated by the HSV-TK constitutive promoter)



LightSwitch Dual Assay Workflow

Step 1: Seed cells in plate format.

Step 2: Co-transfect a SwitchGear GoClone Construct with the Cypridina TK control construct into cells using FuGENE HD. Apply stimulus of interest.

Step 3: Transfer an aliquot of media to secondary assay plate.

Step 4: Add LightSwitch Luciferase Assay Reagent to primary assay plate and read on luminometer.

Step 5: Add BioLux Assay Reagent to secondary assay plate and read on luminometer.

GETTING STARTED

- ▶ Use low passage number cells (highly passaged cells generally yield noisy data).
- ▶ Optimize the transfection protocol for your specific cell line and experimental design using the SwitchGear LightSwitch Transfection Optimization Kit (TFXOPT).
 - Important variables to optimize for your transfection:
 1. FuGENE to plasmid ratio 3:1 or 6:1 (nL FuGENE : ng plasmid)
 2. Cells seeded per well 5,000 or 15,000 cells per well (96 well format)
 3. Duration of transfection 24 or 48 hrs
 4. Amount of supernatant removed for secondary assay

DAY 1

Goal: Seed cells to yield 40-80% confluence after 24 hours

1. Seed the appropriate number of cells for the assay in white 96 or 384-well TC plates (see Table 1).
2. Seed additional wells in a clear TC plate in parallel for assessing confluence.

Table 1

Number of wells in plate	Total volume per well	Number of cells per well (confluency of 40-80%)
96	100µl	5,000-15,000
384	50ul	2,500-10,000

DAY 2

Goal: Prepare GoClone reporter constructs and Cypridina TK control construct and transfect into the seeded cells

Prepare constructs and reagents

1. Thaw GoClone constructs (plasmid DNA) and Cypridina TK control construct DNA at room temperature and mix well.
2. Centrifuge the tubes or plates of DNA to remove condensation from lids.
3. Bring FuGENE HD transfection reagent to room temperature, mix well.
4. Pre-warm Opti-MEM media in a 37°C water bath.

Transfections

- Combine the reagents in Table 2 for each transfection. Conduct at least three replicate transfections per construct in each condition (e.g. untreated and treated). Volumes listed in the table are for a single replicate transfection in a single well. We recommend at least 50ng of GoClone plasmid DNA and 10ng of Cypridina TK control construct DNA per well for 96-well experiments.

Co-transfection Assay Component	Per well (96-well)	Per well (384-well)
FuGENE HD* Transfection Reagent	0.18µl	0.12µl
Opti-MEM (serum free media)	2.15µl	1.38µl
GoClone construct (30ng/µl)	1.67µl	1.00µl
Cypridina control construct (10ng/µl)	1.00µl	0.5µl
TOTAL	5.00µl	3.00µl

- * The FuGENE HD to plasmid ratio shown in this example is 3:1 (180nl FuGENE HD to 60ng total plasmid DNA)
- Mix DNA, FuGENE HD, OptiMEM combination well. Let sit at room temperature for 30 minutes.
 - Gently drip 5µL or 3µL (96- or 384-well formats) onto seeded cells.
 - Shake plate gently, cover with lid or breathable sealing tape and return to incubator.
 - Incubate for 24 to 48 hours before assaying.
 - Alternatively, incubate cells with transfection mix for 16 to 24 hours before changing conditions or applying a stimulus.

DAY 3

Goal: Measure LightSwitch GoClone luciferase activity with LightSwitch Assay Reagents

Note: Luciferase assays may be conducted immediately or the plates may be frozen at -80°C (freezing generally increases cell lysis and luciferase signal). If using frozen plates, thaw and bring to room temperature before assaying.

- Remove plate from incubator and bring to room temperature.
- Transfer 20µl supernatant to a second white 96-well plate and either set plate aside or freeze at -80C.
- Prepare LightSwitch Assay Reagents
 - Reconstitute 100X Substrate by adding 100µL of Substrate Solvent to tube of lyophilized Assay Substrate. *Protect from light and minimize time at room temperature. 100X Substrate may be stored at -20C for 2-3 weeks. For best results, use freshly reconstituted substrate.*

B. Prepare Assay Solution by thawing 10mL bottle of Assay Buffer in room temperature water bath and add 100 μ L of reconstituted 100X Substrate prior to use. *For best results, avoid additional freeze/thaw cycles. To thaw re-frozen buffer, incubate in a warm (37C) water bath for at least 1 hour and mix well to ensure that all components go back into solution.*

4. Use a multi-channel pipettor to add 100 μ L LightSwitch Assay Solution (buffer+substrate) directly to each sample well in a white 96-well plate. *If cells were grown in another plate or flask format, transfer samples to a white 96-well plate in 100 μ L total volume (media or PBS).*
5. Cover plate, protect from light, and incubate for 30 minutes at room temperature. *If assaying more than one plate, stagger addition of assay solution so that each plate incubates for 30 minutes before reading.*
6. Read each well for 2 seconds in a plate luminometer (SpectraMax L or equivalent).

Goal: Measure Cypridina control construct luciferase activity with BioLux Assay Reagents

1. Thaw BioLux Assay Buffer (1X) completely to room temperature (protect from light) and mix well before use.
2. If using frozen supernatant, allow plate to thaw completely and bring to room temperature.
3. Prepare BioLux Assay Solution
 - A. Add 0.5mL of absolute ethanol (not included) to the BioLux Substrate Solvent and mix well.
 - B. Carefully lift the rubber cap of the lyophilized BioLux Substrate vial and add 60 μ L of the ethanol-solvent mixture to the lyophilized BioLux Substrate.
 - C. Gently mix to dissolve the substrate (do not vortex, do not create bubbles by pipeting) and incubate at room temperature for 30 minutes (protect from light). Note: Be sure to dissolve any residual lyophilized powder on the rubber cap.
 - D. Aliquot the reconstituted substrate (100x solution) and store at -80°C if desired.
 - E. Prepare BioLux Assay Solution by adding 50 μ L of reconstituted substrate to 5mL of BioLux Assay Buffer.
4. Add 50 μ L of the BioLux assay solution to each well of the assay plate that contains a Cypridina supernatant aliquot.
5. Cover plate, protect from light, and incubate for 30 minutes at room temperature. *If assaying more than one plate, stagger addition of assay solution so that each plate incubates for 30 minutes before reading.*
6. Read each well for 2 seconds in a plate luminometer (SpectraMax L or equivalent).

Goal: Conduct normalization calculations

Compare well-to-well variation between LightSwitch Reagent only data and the data created using the Cypridina co-transfection control normalization to determine whether the co-transfection control will be needed in subsequent experiments. To determine whether the co-transfection decreases well-to-well variation, divide the raw signal from the LightSwitch reagent data by the raw signal from the corresponding BioLux data. Repeat the data analysis on these new values that account for differences in the Cypridina signal and establish whether the correction decreases variation in your experimental setup.

RELATED PRODUCTS

Item	Catalog No.
LightSwitch Assay Reagent™ (1000 assays)	LS100
GoClones™ Promoters	S7- - - - -
GoClones™ Promoter Controls	S79- - - -
GoClones™ 3'UTRs	S8- - - - -
GoClones™ 3'UTR Controls	S89- - - -
LightSwitch Transfection Optimization Kit	TFXOPT
FuGENE HD® Transfection Reagent	F500

REFERENCES

- Nakajima, Y. et al. (2004) *Biosci. Biotechnol. Biochem.* 63:565-570.
- Yamagishi, K., Enomoto, T. and Ohmiya, Y. (2006) *Anal. Biochemistry*, 354:15-21.
- Wu, C., Suzuki-Ogoh, C. and Ohmiya, Y. (2007) *Biotechniques*, 42:290-292.
- Lorenz, W., Cormier, M., O'Kane, D., Hua, D., Escher, A., and Szalay, A. (1996) *J. Biolumin. Chemilumin.* 11:31-37.
- Haddock, S., Moline, M., and Case, J. (2010) *Ann. Rev. Mar. Sci.* 2:293-343.
- Matthews, J., Hori, K., and Cormier, M. (1977) *Biochemistry*. 16:85-91.
- Lorenz, W., McCann, R., Longiaru, M., and Cormier, M. (1991) *Proc. Nat. Acad. Sci.* 88:4438-4442.
- Hori, K., Wampler, J., Matthews, J., and Cormier, M. (1973) *Biochemistry*. 12:4463-4468.

LIMITED USE LICENSE

For research use only. This product and/or its use is subject to one or more of the following SwitchGear Genomics patents: U.S. Pat. Pub. Nos. US2007/0161031, US2009/0018031, US2008/0220983, US61/397340 and various corresponding patent applications and issued patents. BY USE OF THIS PRODUCT, RESEARCHER AGREES TO BE BOUND BY THE TERMS OF THIS LIMITED USE STATEMENT. If the researcher is not willing to accept the conditions of this limited use statement, and the product is unused, SwitchGear Genomics will accept return of the unused product and provide the researcher with a full refund. Researchers may use this information and/or product in their research, but express written consent from SwitchGear Genomics is required for the transfer of any information including, but not limited to, the DNA sequences or genomic coordinates of these fragments or the products or their derivatives for use by others. Researchers may not reverse engineer, amplify from or transform to prepare more DNA from SwitchGear Genomics constructs without express written consent from SwitchGear Genomics. Researchers shall have no right to modify or otherwise create variations of the nucleotide sequence of the reporter gene without prior expressed written consent of SwitchGear Genomics. In addition, researchers must either: (1) use luminescent assay reagents purchased from SwitchGear Genomics for all determinations of luminescence activity of this product and its derivatives; or (2) contact SwitchGear Genomics to obtain a license for use of the product and its derivatives.

Unless otherwise agreed, no reach-through payments shall be owed to SwitchGear Genomics relating to an organization's commercialization of products that are the discoveries resulting from the research use of this product or its derivatives, provided that such products do not fall within the scope of the valid claims of any issued patents assigned or licensed to SwitchGear Genomics, or that such commercialization would not be a violation of the terms of this license. No other uses including, but not limited to commercial use, of this information, products or their derivatives is authorized without the express written consent of SwitchGear Genomics. Commercial Use means any and all uses of this product and derivatives by a party for monetary or other consideration and may include, but is not limited to use in: (1) product development or manufacture; (2) to provide a service utilizing SwitchGear information or products; and (3) the resale of the information, products or their derivatives, regardless of their intended use. With respect to such Commercial Use, or any diagnostic, therapeutic or prophylactic uses, please contact SwitchGear Genomics for supply and licensing information.

SWITCHGEAR GENOMICS MAKES NO REPRESENTATIONS OR WARRANTIES OF ANY KIND, EITHER EXPRESSED OR IMPLIED, INCLUDING FOR MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE WITH REGARDS TO THE PRODUCT. The terms of this agreement shall be governed under the laws of the State of California, USA.

Licensed under certain patents and patent applications from the National Institute of Advanced Industrial Science and Technology ("AIST") for Research and Development Purposes. U.S. Patent Nos. 7,718,389 and 7,989,621; U.S. Appln. Serial Nos. 12/588,671 and 13/067,565; Japanese Patent Nos. 4,761,150 and 4,484,429; Japanese Appln. Serial No.: 2006-280827; 2007-536587; 2009-257631; EPO Appln. Serial No.: 06 810 525.3; Chinese Appln. Serial No.: 200680035410.3 For use of the Biolux® Cypridina Luciferase Assay Kit, or associated assay reagents, in human diagnosis and measurement in relation to human health, contact busdev@neb.com.