



**Rabbit Progesterone Receptor  
(NR3C3, PGR, PR)  
Reporter Assay System**

**96-well Format Assays**  
Product # RB05001

■

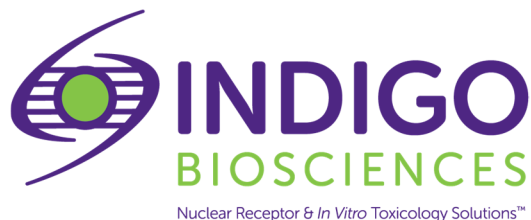
**Technical Manual**  
*(version 7.2i)*

**[www.indigobiosciences.com](http://www.indigobiosciences.com)**

3006 Research Drive, Suite A1, State College, PA 16801, USA

Customer Service:  
814-234-1919; FAX 814-272-0152  
[customerserv@indigobiosciences.com](mailto:customerserv@indigobiosciences.com)

Technical Service:  
814-234-1919  
[techserv@indigobiosciences.com](mailto:techserv@indigobiosciences.com)



## **Rabbit PGR Reporter Assay System 96-well Format Assays**

### **I. Description**

|                                                |   |
|------------------------------------------------|---|
| ▪ The Assay System.....                        | 3 |
| ▪ The Assay Chemistry.....                     | 3 |
| ▪ Preparation of Test Compounds.....           | 4 |
| ▪ Considerations for Automated Dispensing..... | 4 |
| ▪ Assay Scheme.....                            | 4 |
| ▪ Assay Performance.....                       | 5 |

### **II. Product Components & Storage Conditions .....7**

### **III. Materials to be Supplied by the User.....8**

### **IV. Assay Protocol**

|                                                         |   |
|---------------------------------------------------------|---|
| ▪ A word about <i>Antagonist</i> -mode assay setup..... | 7 |
| ▪ <i>DAY 1 Assay Protocol</i> .....                     | 7 |
| ▪ <i>DAY 2 Assay Protocol</i> .....                     | 9 |

### **V. Related Products.....10**

### **VI. Limited Use Disclosures.....11**

### **APPENDIX 1: Example Scheme for Serial Dilutions.....12**

## I. Description

### ▪ The Assay System ▪

This nuclear receptor assay utilizes proprietary human cells engineered to provide constitutive, high-level expression of the full-length **Rabbit** (*Oryctolagus cuniculus*) **Progesterone Receptor** (nr3c3), a ligand-dependent transcription factor referred to herein as **rbPGR**.

INDIGO's Reporter Cells include the luciferase reporter gene functionally linked to a PGR-responsive promoter. Thus, quantifying changes in luciferase expression in the treated reporter cells provides a sensitive surrogate measure of the changes in rbPGR activity. Luciferase gene expression occurs after ligand-bound PGR undergoes nuclear translocation, DNA binding, recruitment and assembly of the co-activators and accessory factors required to form a functional transcription complex, culminating in expression of the target gene. Unlike *in vitro* binding assays, and some other cell-based assay strategies, the readout from INDIGO's reporter cells demands the same orchestration of all intracellular molecular interactions and events that can be expected to occur *in vivo*.

rbPGR Reporter Cells are prepared using INDIGO's proprietary **CryoMite™** process. This cryo-preservation method yields exceptional cell viability post-thaw, and provides the convenience of immediately dispensing healthy, division-competent reporter cells into assay plates. There is no need for cumbersome intermediate treatment steps such as spin-and-rinse of cells, viability determinations, or cell titer adjustments prior to assay setup.

The principal application of this assay is in the screening of test samples to quantify functional activities, either agonist or antagonist, that they may exert against the rabbit progesterone receptor. It is an all-inclusive assay system that includes, in addition to rbPGR Reporter Cells, two optimized media for use during cell culture and for preparing dilutions of test samples, the reference agonist Progesterone, Luciferase Detection Reagent, a cell culture-ready assay plate, and a detailed protocol.

### ▪ The Assay Chemistry ▪

INDIGO's nuclear receptor assay kits capitalize on the extremely low background, high-sensitivity, and broad linear dynamic range of bio-luminescence reporter gene technology.

Reporter Cells incorporate the cDNA encoding beetle luciferase, a 62 kD protein originating from the North American firefly (*Photinus pyralis*). Luciferase catalyzes the mono-oxidation of D-luciferin in a  $Mg^{+2}$ -dependent reaction that consumes  $O_2$  and ATP as co-substrates, and yields as products oxyluciferin, AMP,  $PP_i$ ,  $CO_2$ , and photon emission. Luminescence intensity of the reaction is quantified using a luminometer and is reported in terms of Relative Light Units (RLU's).

INDIGO's Nuclear Receptor Assay kits feature a luciferase detection reagent specially formulated to provide stable light emission between 5 and 90+ minutes after initiating the luciferase reaction. Incorporating a 5-minute reaction-rest period ensures that light emission profiles attain maximal stability, thereby allowing assay plates to be processed in batch. By doing so, the signal output from all sample wells, from one plate to the next, may be directly compared within an experimental set.

### ▪ Preparation of Test Compounds ▪

Small molecule test compounds are typically solvated in DMSO at high concentrations; ideally 1,000x-concentrated stocks relative to the highest desired treatment concentration in the assay. Using high-concentration stocks minimizes DMSO carry-over into the assay plates. Immediately prior to setting up an assay, the master stocks are serially diluted using one of two alternative strategies:

1.) As described in *Step 7* and depicted in Appendix 1 for the reference agonist Progesterone, **Compound Screening Medium (CSM)** may be used as the diluent to make serial dilutions of test compounds to achieve the desired final assay concentration series.

Alternatively, if test compound solubility is expected to be problematic,

2.) DMSO may be used to make serial dilutions, thereby generating 1,000x-concentrated stocks for each independent test concentration. Treatment media are then prepared using CSM to make final 1,000-fold dilutions of the prepared DMSO dilution series.

Regardless of the dilution method used, the final concentration of total DMSO carried over into assay wells should not exceed 0.4%. Significant DMSO-induced cytotoxicity can be expected above 0.4%.

**NOTE:** CSM is formulated to help stabilize hydrophobic test compounds in the aqueous environment of the assay mixture. Nonetheless, high concentrations of extremely hydrophobic test compounds diluted in CSM may lack long-term stability and/or solubility, especially if further stored at low temperatures. Hence, it is recommended that test compound dilutions are prepared in CSM immediately prior to assay setup and are considered to be 'single-use' reagents.

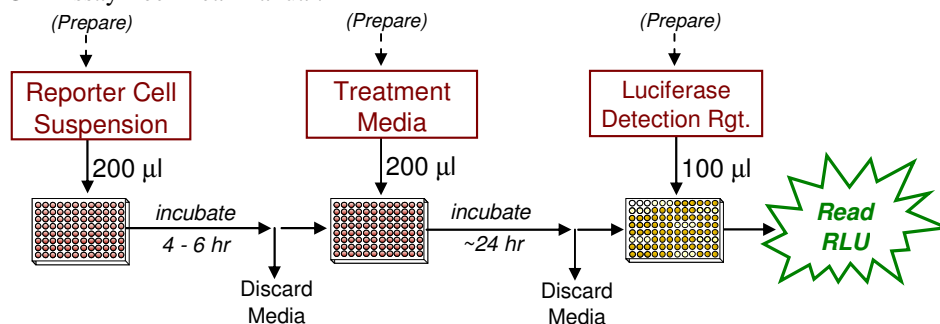
### ▪ Considerations for Automated Dispensing ▪

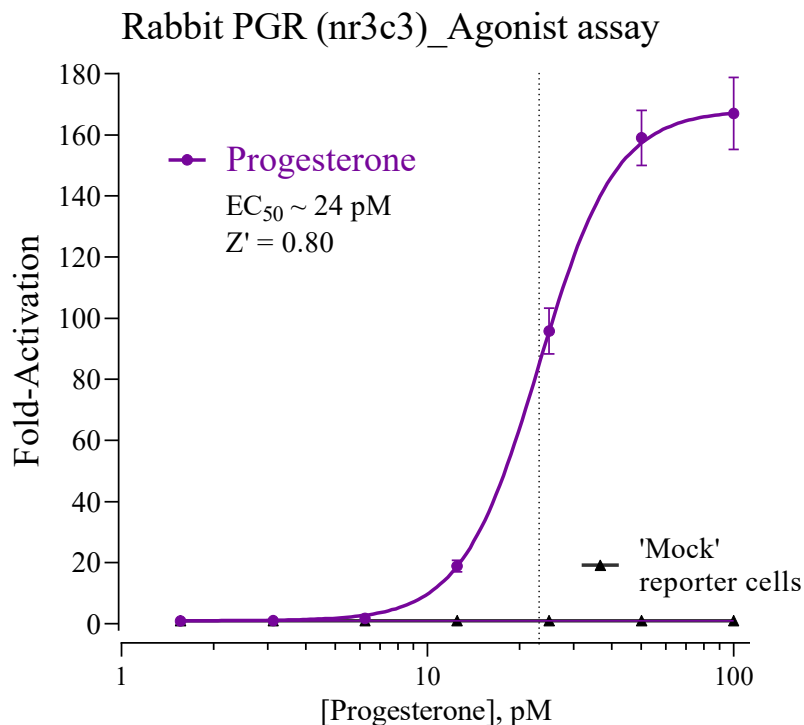
When processing a small number of assay plates, first carefully consider the dead volume requirement of your dispensing instrument before committing assay reagents to its setup. In essence, "dead volume" is the volume of reagent that is dedicated to the instrument; it will *not* be available for final dispensing into assay wells. The following Table provides information on reagent volume requirements, and available excesses.

| Stock Reagent<br>& Volume provided                                         | Volume to be<br>Dispensed<br>(96-well plate) | Excess rgt. volume<br>available for instrument<br>dead volume |
|----------------------------------------------------------------------------|----------------------------------------------|---------------------------------------------------------------|
| <b>Reporter Cell Suspension</b><br>21 ml<br>(prepared from kit components) | 200 µl / well<br>19.2 ml / plate             | ~ 1.8 ml                                                      |
| <b>LDR</b><br>12 ml<br>(prepared from kit components)                      | 100 µl / well<br>9.6 ml / plate              | ~ 2.4 ml                                                      |

### ▪ Assay Scheme ▪

**Figure 1.** Assay workflow. *In brief*, the Reporter Cell suspension is dispensed into assay wells and pre-incubated for 4-6 hours. Following the pre-incubation period, culture media are discarded, and the prepared treatment media are added. Following 22 - 24 hours incubation, treatment media are discarded, and Luciferase Detection Reagent is added. The intensity of light emission (in units of 'Relative Light Units'; RLU) from each assay well is quantified using a plate-reading luminometer. *Note:* If INDIGO's Live Cell Multiplex (LCM) Assay is to be incorporated, refer to the assay workflow schematic provided in the LCM Assay Technical manual.





**Figure 2. Agonist dose-response analyses of Rabbit PGR.**

Agonist dose-response analyses of PGR Reporter Cells using Progesterone (provided). In addition, to assess the level of background signal contributed by endogenously expressed PGR, or non-specific factors that may cause activation of the luciferase reporter gene, “mock” reporter cells were evaluated. (Mock reporter cells, which contain only the luciferase vector, are not provided with assay kits). The concentrated stock of progesterone was serially diluted in 2-fold decrements using CSM. Final assay concentrations for progesterone-treated cells ranged between 100 pM and 1.56 pM. Luminescence was quantified and average relative light units (RLU) and corresponding standard deviation (SD) and Fold-Activation values were determined for each treatment concentration (n = 3). Z' values were calculated as described by Zhang, *et al.* (1999)<sup>1</sup>. Non-linear regression and EC<sub>50</sub> analyses were performed using GraphPad Prism software.

<sup>1</sup> Zhang JH, Chung TD, Oldenburg KR. (1999) A Simple Statistical Parameter for Use in Evaluation and Validation of High Throughput Screening Assays. *J Biomol Screen.*:4(2), 67-73.

$$Z' = 1 - [3 * (SD^{Prog, EC100} + SD^{Bkg}) / (RLU^{Prog, EC100} - RLU^{Bkg})]$$

## II. Product Components & Storage Conditions

This rabbit PGR Assay kit contains materials to perform assays in a single collagen-coated 96-well assay plate.

**Reporter cells are temperature sensitive! To ensure maximal viability the tube of Reporter Cells must be maintained at -80°C until immediately prior to the rapid-thaw procedure described in Step 2 of this protocol.**

Assay kits are shipped on dry ice. Upon receipt of the kit transfer it to -80°C storage. If you wish to first inspect and inventory the individual kit components be sure to first transfer and submerge the tube of reporter cells in dry ice.

The aliquot of Reporter Cells is provided as a single-use reagent. Once thawed, reporter cells can NOT be refrozen, nor can they be maintained in extended culture with any hope of retaining downstream assay performance. Therefore, extra volumes of these reagents should be discarded after assay setup.

The date of product expiration is printed on the Product Qualification Insert (PQI) enclosed with each kit.

| <u>Kit Components</u>                                                                 | <u>Amount</u> | <u>Storage Temp.</u> |
|---------------------------------------------------------------------------------------|---------------|----------------------|
| ▪ Rabbit PGR Reporter Cells                                                           | 1 x 2.0 mL    | <b>-80°C</b>         |
| ▪ Cell Recovery Medium (CRM)                                                          | 2 x 10.5 mL   | -20°C                |
| ▪ Compound Screening Medium (CSM)                                                     | 1 x 45 mL     | -20°C                |
| ▪ Progesterone, 100 nM (in DMSO)<br>(reference agonist)                               | 1 x 30 µL     | -20°C                |
| ▪ Detection Substrate                                                                 | 1 x 6.0 mL    | <b>-80°C</b>         |
| ▪ Detection Buffer                                                                    | 1 x 6.0 mL    | -20°C                |
| ▪ 96-well, <i>collagen-coated</i> assay plate<br>(white, sterile, cell-culture ready) | 1             | <b>-20°C</b>         |

*NOTE:* This rbPGR Assay kit contains one 96-well assay plate in which the assay wells have been collagen-coated and dried; the assay plate should be stored frozen (-20°C or colder) until use.

## III. Materials to be Supplied by the User

The following materials must be provided by the user, and should be made ready prior to initiating the assay procedure:

### DAY 1

- container of dry ice (see Step 2)
- cell culture-rated laminar flow hood.
- 37°C, humidified 5% CO<sub>2</sub> incubator for mammalian cell culture.
- 37°C water bath.
- 70% alcohol wipes
- 8-channel electronic, repeat-dispensing pipettes & sterile tips
- disposable media basins, sterile.
- sterile multi-channel media basins (such as the Heathrow Scientific "Dual-Function Solution Basin"), or sterilized 96 deep-well blocks (e.g., Axygen Scientific, #P-2ML-SQ-C-S), or appropriate similar vessel for generating dilution series of reference and test compound(s).
- *Optional:* antagonist reference compound
- *Optional:* clear 96-well assay plate, collagen-coated, for viewing cells on Day 2.

**DAY 2** plate-reading luminometer.

## IV. Assay Protocol

Review the entire Assay Protocol before starting. Completing the assay requires an overnight incubation. *Steps 1-11* are performed on **Day 1**, requiring less than 2 hours of bench work to complete, but including a 4-hour incubation step. *Steps 12-17* are performed on **Day 2** and require less than 1 hour to complete.

### ▪ A word about Antagonist-mode assay setup ▪

Receptor inhibition assays expose the Reporter Cells to a constant, sub-maximal concentration (typically between EC<sub>50</sub> – EC<sub>85</sub>) of a known agonist AND the test compound(s) to be evaluated for antagonist activity. This rbPGR assay kit includes a 100 nM stock solution of Progesterone, a potent physiological agonist of PGR that may be used to setup antagonist-mode assays. 34 pM progesterone approximates EC<sub>70-80</sub> in this cell-based assay. Hence, it is an appropriate concentration of agonist to be used when screening test compounds for inhibitory activity.

Add the challenge agonist (progesterone) to a bulk volume of **CSM** at an EC<sub>50</sub> – EC<sub>85</sub> concentration. This medium is then used to prepare serial dilutions of test compounds to achieve the desired respective final assay concentrations. This is an efficient and precise method of setting up PGR antagonist assays, and it is the method presented in *Step 7b* of this protocol.

### DAY 1 Assay Protocol: All steps should be performed using aseptic technique.

**1.)** Remove the **2 tubes** of **Cell Recovery Medium (CRM)** from freezer storage, thaw and equilibrate to 37°C using a water bath.

**2.) Rapid Thaw of the Reporter Cells:** *First*, retrieve the two tubes of **CRM** from the 37°C water bath and sanitize their outside surfaces with a 70% ethanol swab.

*Second*, retrieve the tube of **Reporter Cells** from -80°C storage, place it directly into dry ice for transport to the laminar flow hood. When ready to begin, transfer the tube of reporter cells into a rack and, *without delay*, perform a rapid thaw of the cells by transferring 9.5 ml from **each of the 2 tubes** of 37°C CRM into the tube of frozen cells. Place the tube of Reporter Cells in a 37°C water bath for 5 minutes. The resulting volume of cell suspension will be **21 ml**.

**3.)** Retrieve the tube of Reporter Cell Suspension from the water bath and sanitize the outside surface with a 70% alcohol swab.

**4.)** Gently invert the tube of Reporter Cells several times to gain a homogenous cell suspension. Transfer the cell suspension into a reservoir. Using an 8-channel pipette, dispense **200 µl / well** of cell suspension into the assay plate.

*NOTE 4.1:* If INDIGO's Live Cell Multiplex Assay is to be incorporated, a minimum of 3 'cell blank' wells (meaning cell-free, but containing 'Compound Screening Media') must be included in the assay plate to allow quantification of plate-specific fluorescence background (refer to the LCMA Technical Manual).

*NOTE 4.2:* Increased well-to-well variation (= increased standard deviation!) will occur if care is not taken to prevent cells from settling in the reservoir during the dispensing period. Likewise, take care to ensure precision in dispensing exact volumes across the assay plate.

*NOTE 4.3:* Users sometimes wish to examine the reporter cells using a microscope. If so, the extra volume of cell suspension provided with each kit may be dispensed into a clear, preferably collagen-coated, 96-well assay plate. Continue to process this plate in identical manner to the white assay plate.

**5.) Pre-incubate reporter cells.** Place the assay plate into a cell culture incubator (37°C, ≥ 70% humidity, 5% CO<sub>2</sub>) for 4 - 6 hours.

6.) Near the end of the pre-incubation period remove Compound Screening Medium (CSM) from freezer storage and thaw in a 37°C water bath.

**7.) Prepare the Test Compound and Reference Compound treatment media at the desired final assay concentrations:** Use CSM to prepare an appropriate dilution series of the reference and test compound stocks. Prepare all treatment media at the desired final assay concentrations. In *Step 9*, the prepared treatment media will be dispensed at 200 µl / well into the assay plate. Manage dilution volumes carefully; this assay kit provides **45 ml** of CSM.

NOTE: Total DMSO carried over into assay reactions should not exceed 0.4%.

**a. Agonist-mode assays.** This rbPGR Assay kit includes a 100 nM stock solution of Progesterone, a potent reference agonist of rabbit PGR. The following 7-point treatment series, with concentrations presented in 2-fold decrements, provides a complete dose-response: 100, 50, 25, 12.5, 6.25, 3.13, and 1.56 pM. Always include 'no treatment' (or 'vehicle') controls. **APPENDIX 1** provides guidance for generating such a dilution series.

~ or ~

**b. Antagonist-mode assays.** When setting antagonist assays, first supplement a bulk volume of CSM with the challenge agonist Progesterone to achieve the desired final assay-concentration (refer to "*A word about antagonist-mode assay setup*", pg. 7). The agonist-supplemented CSM is then used to generate dilutions of test compound stocks to achieve their final assay concentrations.

8.) At the end of the pre-culture period, discard the media. The preferred method is to use a 'wrist flick' to eject media into an appropriate waste container. *Gently* tap the inverted plate onto a clean absorbent paper towel to remove residual droplets. Cells will remain tightly adhered to well bottoms.

9.) Dispense **200 µl / well** of each prepared treatment media into the assay plate.

NOTE: If well-to-well variation due to 'edge-effects' is a concern this problem may be mitigated by dispensing sterile liquid into the *inter-well* spaces of the assay plate. Simply remove 1 tip from the 8-channel dispenser and dispense 100 µl of sterile water into each of the seven inter-well spaces per column of wells.

10.) Transfer the assay plate into a cell culture incubator for 22 - 24 hours.

NOTE: Ensure a high-humidity (≥70%) environment within the cell culture incubator. This is critical to prevent the onset of deleterious "edge-effects" in the assay plate.

11.) For greater convenience on *Day 2*, retrieve **Detection Substrate and Detection Buffer** from freezer storage and place them in a dark refrigerator (4°C) to thaw overnight.



**DAY 2 Assay Protocol:** Subsequent manipulations do *not* require special regard for aseptic technique and may be performed on a bench top.

**12.)** Approximately 30 minutes before intending to quantify receptor activity, remove **Detection Substrate** and **Detection Buffer** from the refrigerator and place them in a low-light area so that they may equilibrate to room temperature.

*NOTE:* Do NOT actively warm Detection Substrate above room temperature. If these solutions were not allowed to thaw overnight at 4°C, a room temperature water bath may be used to expedite thawing.

**13.)** Set the plate-reader to "luminescence" mode. Set the instrument to perform a single 5 second "plate shake" prior to reading the first assay well. Read time may be set to 0.5 second (500 mSec) per well, *or less*.

**14.)** *Immediately before proceeding to Step 15*, transfer the entire volume of Detection Buffer into the vial of Detection Substrate, thereby generating a 12 ml volume of **Luciferase Detection Reagent (LDR)**. Mix gently to avoid foaming.

**15.)** Following 22 - 24 hours incubation in treatment media, discard the media contents by manually ejecting it into an appropriate waste container. *Gently* tap the inverted plate onto a clean absorbent paper towel to remove residual droplets.

**16.)** Add 100 µl of **LDR** to each well of the assay plate. Allow the assay plate to rest at room temperature for 5 - 10 minutes following the addition of LDR. Do not shake the assay plate during this period.

**17.)** Quantify luminescence.

## V. Related Products

| Product No.                                                                                                | Product Descriptions                                                                                                          |
|------------------------------------------------------------------------------------------------------------|-------------------------------------------------------------------------------------------------------------------------------|
| <b>Human PGR Assay Products</b>                                                                            |                                                                                                                               |
| IB05001-32                                                                                                 | Human PGR Reporter Assay System<br>3x 32 assays in 8-well strips (96-well plate format)                                       |
| IB05001                                                                                                    | Human PGR Reporter Assay System<br>1x 96-well format assay                                                                    |
| IB05002                                                                                                    | Human PGR Reporter Assay System<br>1x 384-well format assays                                                                  |
| <b>Rat PGR Assays</b>                                                                                      |                                                                                                                               |
| R05001-32                                                                                                  | Rat PGR Reporter Assay System<br>3x 32 assays in 8-well strips (96-well plate format)                                         |
| R05001                                                                                                     | Rat PGR Reporter Assay System<br>1x 96-well format assay                                                                      |
| <b>Rabbit PGR Assays</b>                                                                                   |                                                                                                                               |
| RB05001-32                                                                                                 | Rabbit PGR Reporter Assay System<br>3x 32 assays in 8-well strips (96-well plate format)                                      |
| RB05001                                                                                                    | Rabbit PGR Reporter Assay System<br>1x 96-well format assay                                                                   |
| <b>Monkey PGR Assays</b>                                                                                   |                                                                                                                               |
| C05001-32                                                                                                  | Cynomolgus Monkey PGR Reporter Assay System<br>3x 32 assays in 8-well strips (96-well plate format)                           |
| C05001                                                                                                     | Cynomolgus Monkey PGR Reporter Assay System<br>1x 96-well format assay                                                        |
| Bulk volumes of assay reagents may be custom manufactured to accommodate any scale of HTS. Please Inquire. |                                                                                                                               |
| <b>LIVE Cell Multiplex (LCM) Assay</b>                                                                     |                                                                                                                               |
| LCM-01                                                                                                     | Reagent volumes sufficient to perform <b>96</b> Live Cell Assays in 1x96-well, or 2x48-well, or 3x32-well assay plate formats |
| LCM-05                                                                                                     | Reagent in <b>5x bulk volume</b> to perform <b>480</b> Live Cell Assays performed in 5 x 96-well assay plates                 |
| LCM-10                                                                                                     | Reagent in <b>10x bulk volume</b> to perform <b>960</b> Live Cell Assays performed in 10 x 96-well assay plates               |
| <b>INDIGlo Luciferase Detection Reagent</b>                                                                |                                                                                                                               |
| LDR-10, -25, -50, -500                                                                                     | INDIGlo Luciferase Detection Reagents in 10 mL, 25 mL, 50 mL, and 500 mL volumes                                              |

Please refer to INDIGO Biosciences website for updated product offerings.

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

## ***VI. Limited Use Disclosures***

Products offered by INDIGO Biosciences, Inc. are for RESEARCH PURPOSES ONLY – not for therapeutic, diagnostic, or contact use in humans or animals.

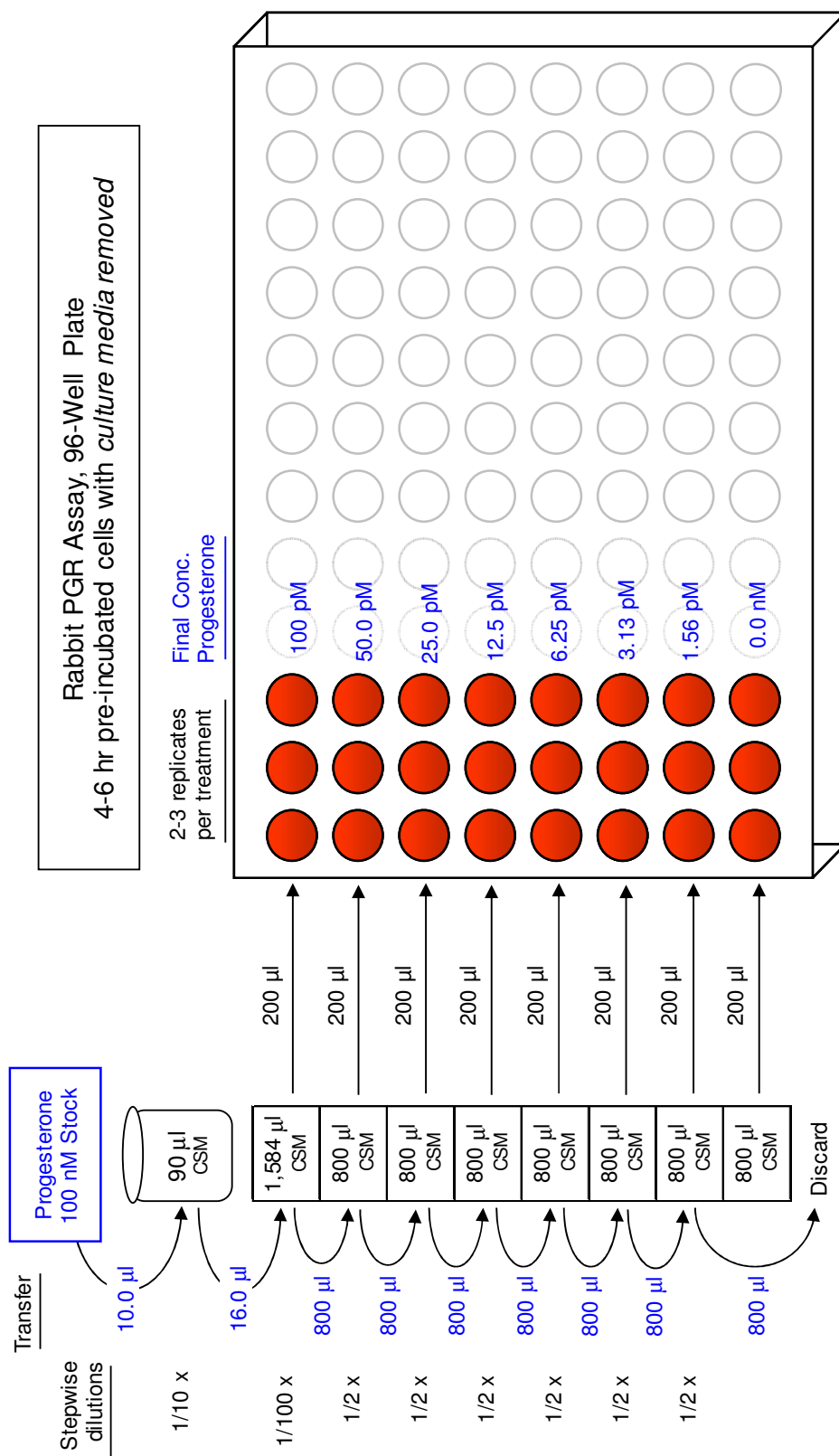
“CryoMite” is a Trademark <sup>TM</sup> of INDIGO Biosciences, Inc. (State College, PA, USA).

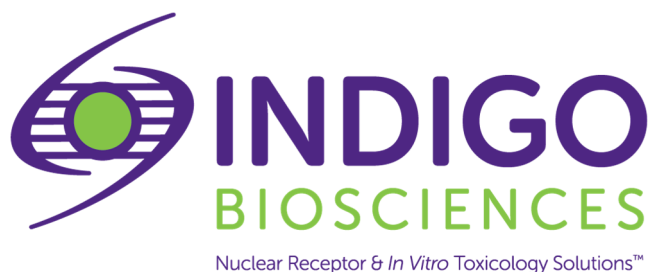
Product prices, availability, specifications, claims and technical protocols are subject to change without prior notice. The printed Technical Manual provided in the kit box will always be the most current version available.

Copyright © INDIGO Biosciences, Inc. (State College, PA, USA). All rights reserved.

## APPENDIX 1

Example scheme for the serial dilution of Progesterone reference agonist, and the setup of a rabbit PGR dose-response assay.





**Rabbit Progesterone Receptor  
(NR3C3, PGR, PR)  
Reporter Assay System**

**3x32 Assays in 96-well Plate Format**  
Product # RB05001-32

■

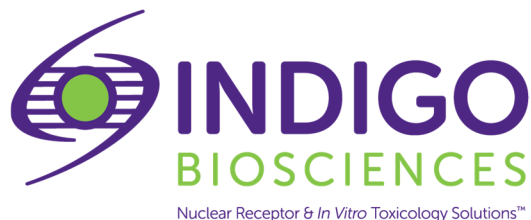
**Technical Manual**  
*(version 7.2i)*

**[www.indigobiosciences.com](http://www.indigobiosciences.com)**

3006 Research Drive, Suite A1, State College, PA 16801, USA

Customer Service:  
814-234-1919; FAX 814-272-0152  
[customerserv@indigobiosciences.com](mailto:customerserv@indigobiosciences.com)

Technical Service:  
814-234-1919  
[techserv@indigobiosciences.com](mailto:techserv@indigobiosciences.com)



## **Rabbit PGR Reporter Assay System** **3x 32 Assays in 96-well Plate Format**

### **I. Description**

- The Assay System.....3
- The Assay Chemistry.....3
- Preparation of Test Compounds.....4
- Assay Scheme.....4
- Assay Performance.....5

### **II. Product Components & Storage Conditions .....7**

### **III. Materials to be Supplied by the User.....8**

### **IV. Assay Protocol**

- A word about *Antagonist*-mode assay setup.....7
  - *DAY 1 Assay Protocol*.....7
  - *DAY 2 Assay Protocol*.....9

### **V. Related Products.....10**

### **VI. Limited Use Disclosures.....11**

### **APPENDIX 1: Example Scheme for Serial Dilutions.....12**

## I. Description

### ▪ The Assay System ▪

This nuclear receptor assay utilizes proprietary human cells engineered to provide constitutive, high-level expression of the full-length **Rabbit** (*Oryctolagus cuniculus*) **Progesterone Receptor** (nr3c3), a ligand-dependent transcription factor referred to herein as **rbPGR**.

INDIGO's Reporter Cells include the luciferase reporter gene functionally linked to a PGR-responsive promoter. Thus, quantifying changes in luciferase expression in the treated reporter cells provides a sensitive surrogate measure of the changes in rbPGR activity. Luciferase gene expression occurs after ligand-bound PGR undergoes nuclear translocation, DNA binding, recruitment and assembly of the co-activators and accessory factors required to form a functional transcription complex, culminating in expression of the target gene. Unlike *in vitro* binding assays, and some other cell-based assay strategies, the readout from INDIGO's reporter cells demands the same orchestration of all intracellular molecular interactions and events that can be expected to occur *in vivo*.

rbPGR Reporter Cells are prepared using INDIGO's proprietary **CryoMite™** process. This cryo-preservation method yields exceptional cell viability post-thaw, and provides the convenience of immediately dispensing healthy, division-competent reporter cells into assay plates. There is no need for cumbersome intermediate treatment steps such as spin-and-rinse of cells, viability determinations, or cell titer adjustments prior to assay setup.

The principal application of this assay is in the screening of test samples to quantify functional activities, either agonist or antagonist, that they may exert against the rabbit progesterone receptor. It is an all-inclusive assay system that includes, in addition to rbPGR Reporter Cells, two optimized media for use during cell culture and for preparing dilutions of test samples, the reference agonist Progesterone, Luciferase Detection Reagent, a cell culture-ready assay plate, and a detailed protocol.

### ▪ The Assay Chemistry ▪

INDIGO's nuclear receptor assay kits capitalize on the extremely low background, high-sensitivity, and broad linear dynamic range of bio-luminescence reporter gene technology.

Reporter Cells incorporate the cDNA encoding beetle luciferase, a 62 kD protein originating from the North American firefly (*Photinus pyralis*). Luciferase catalyzes the mono-oxidation of D-luciferin in a  $Mg^{+2}$ -dependent reaction that consumes  $O_2$  and ATP as co-substrates, and yields as products oxyluciferin, AMP,  $PP_i$ ,  $CO_2$ , and photon emission. Luminescence intensity of the reaction is quantified using a luminometer and is reported in terms of Relative Light Units (RLU's).

INDIGO's Nuclear Receptor Assay kits feature a luciferase detection reagent specially formulated to provide stable light emission between 5 and 90+ minutes after initiating the luciferase reaction. Incorporating a 5-minute reaction-rest period ensures that light emission profiles attain maximal stability, thereby allowing assay plates to be processed in batch. By doing so, the signal output from all sample wells, from one plate to the next, may be directly compared within an experimental set.

### ▪ Preparation of Test Compounds ▪

Small molecule test compounds are typically solvated in DMSO at high concentrations; ideally 1,000x-concentrated stocks relative to the highest desired treatment concentration in the assay. Using high-concentration stocks minimizes DMSO carry-over into the assay plates. Immediately prior to setting up an assay, the master stocks are serially diluted using one of two alternative strategies:

1.) As described in *Step 7* and depicted in Appendix 1 for the reference agonist Progesterone, **Compound Screening Medium (CSM)** may be used as the diluent to make serial dilutions of test compounds to achieve the desired final assay concentration series.

Alternatively, if test compound solubility is expected to be problematic,

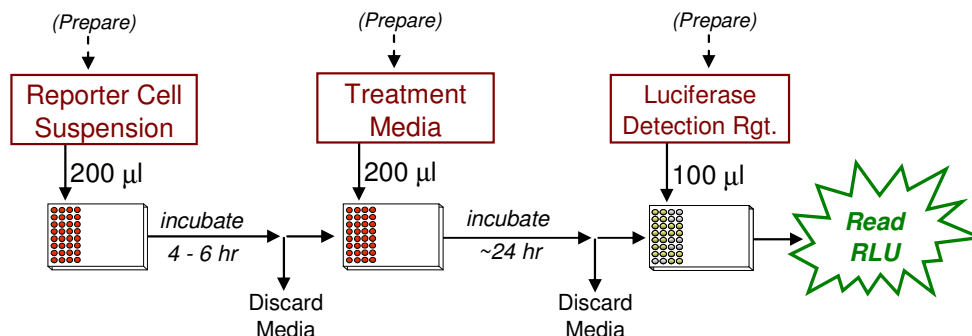
2.) DMSO may be used to make serial dilutions, thereby generating 1,000x-concentrated stocks for each independent test concentration. Treatment media are then prepared using CSM to make final 1,000-fold dilutions of the prepared DMSO dilution series.

Regardless of the dilution method used, the final concentration of total DMSO carried over into assay wells should not exceed 0.4%. Significant DMSO-induced cytotoxicity can be expected above 0.4%.

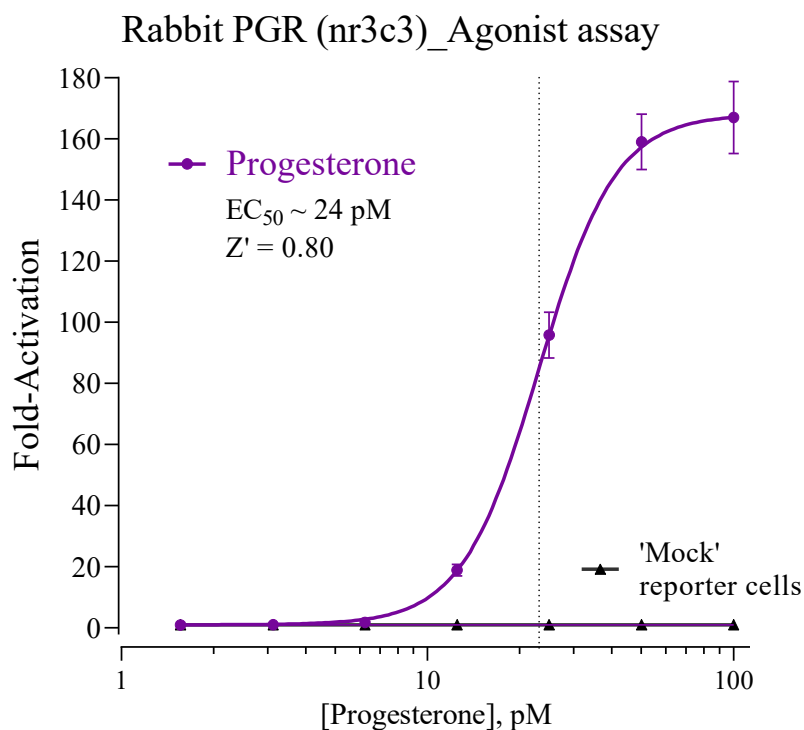
*NOTE:* CSM is formulated to help stabilize hydrophobic test compounds in the aqueous environment of the assay mixture. Nonetheless, high concentrations of extremely hydrophobic test compounds diluted in CSM may lack long-term stability and/or solubility, especially if further stored at low temperatures. Hence, it is recommended that test compound dilutions are prepared in CSM immediately prior to assay setup and are considered to be 'single-use' reagents.

### ▪ Assay Scheme ▪

**Figure 1.** Assay workflow. *In brief*, 200  $\mu$ l of Reporter Cells is dispensed into wells of the assay plate and pre-incubated for 4 - 6 hours. Following the pre-incubation period, culture media are discarded and 200  $\mu$ l/well of the prepared treatment media are added. Following 22 - 24 hours incubation, treatment media are discarded, and Luciferase Detection Reagent is added. The intensity of light emission (in terms of 'Relative Light Units'; RLU) from each assay well is quantified using a plate-reading luminometer.







**Figure 2. Agonist dose-response analyses of Rabbit PGR.**

Agonist dose-response analyses of PGR Reporter Cells using Progesterone (provided). In addition, to assess the level of background signal contributed by endogenously expressed PGR, or non-specific factors that may cause activation of the luciferase reporter gene, “mock” reporter cells were evaluated. (Mock reporter cells, which contain only the luciferase vector, are not provided with assay kits). The concentrated stock of progesterone was serially diluted in 2-fold decrements using CSM. Final assay concentrations for progesterone-treated cells ranged between 100 pM and 1.56 pM. Luminescence was quantified and average relative light units (RLU) and corresponding standard deviation (SD) and Fold-Activation values were determined for each treatment concentration (n = 3). Z' values were calculated as described by Zhang, *et al.* (1999)<sup>1</sup>. Non-linear regression and EC<sub>50</sub> analyses were performed using GraphPad Prism software.

<sup>1</sup> Zhang JH, Chung TD, Oldenburg KR. (1999) A Simple Statistical Parameter for Use in Evaluation and Validation of High Throughput Screening Assays. *J Biomol Screen.*:4(2), 67-73.

$$Z' = 1 - [3 * (SD^{Prog. EC100} + SD^{Bkg}) / (RLU^{Prog. EC100} - RLU^{Bkg})]$$

## II. Product Components & Storage Conditions

This rabbit PGR Assay kit contains materials to perform three distinct groups of assays in a 96-well plate format. Reagents are configured so that each group will comprise 32 assays. If desired, however, reagents may be combined to perform either 64 or 96 assays.

**Reporter cells are temperature sensitive! To ensure maximal viability the tube of Reporter Cells must be maintained at -80°C until immediately prior to the rapid-thaw procedure described in Step 2 of this protocol.**

Assay kits are shipped on dry ice. Upon receipt of the kit transfer it to -80°C storage. If you wish to first inspect and inventory the individual kit components be sure to first transfer and submerge the tube of reporter cells in dry ice.

The aliquot of Reporter Cells is provided as a single-use reagent. Once thawed, reporter cells can NOT be refrozen, nor can they be maintained in extended culture with any hope of retaining downstream assay performance. Therefore, extra volumes of these reagents should be discarded after assay setup.

The date of product expiration is printed on the Product Qualification Insert (PQI) enclosed with each kit.

| <u>Kit Components</u>                                               | <u>Amount</u> | <u>Storage Temp.</u> |
|---------------------------------------------------------------------|---------------|----------------------|
| ▪ Rabbit PGR Reporter Cells                                         | 3 x 0.6 mL    | <b>-80°C</b>         |
| ▪ Cell Recovery Medium (CRM)                                        | 2 x 10.5 mL   | -20°C                |
| ▪ Compound Screening Medium (CSM)                                   | 1 x 45 mL     | -20°C                |
| ▪ Progesterone, 100 nM (in DMSO)<br>(reference agonist)             | 1 x 30 µL     | -20°C                |
| ▪ Detection Substrate                                               | 3 x 2.0 mL    | <b>-80°C</b>         |
| ▪ Detection Buffer                                                  | 3 x 2.0 mL    | -20°C                |
| ▪ Plate frame                                                       | 1             | ambient              |
| ▪ Snap-in, 8-well strips<br>(white, sterile, collagen-coated wells) | 12            | <b>-20°C</b>         |

*NOTE:* This Assay kit contains 8-well strips that have been collagen-coated and dried; these strip wells should be stored frozen (-20°C or -80°C) until use.

## III. Materials to be Supplied by the User

The following materials must be provided by the user, and should be made ready prior to initiating the assay procedure:

### **DAY 1**

- container of dry ice (see *Step 2*)
- cell culture-rated laminar flow hood.
- 37°C, humidified 5% CO<sub>2</sub> incubator for mammalian cell culture.
- 37°C water bath.
- 70% alcohol wipes
- 8-channel electronic, repeat-dispensing pipettes & sterile tips
- disposable media basins, sterile.
- sterile multi-channel media basins (such as the Heathrow Scientific "Dual-Function Solution Basin"), *or* sterilized 96 deep-well blocks (*e.g.*, Axygen Scientific, #P-2ML-SQ-C-S), *or* appropriate similar vessel for generating dilution series of reference and test compound(s).
- *Optional:* antagonist reference compound
- *Optional:* clear 96-well assay plate, collagen-coated, for viewing cells on *Day 2*.

**DAY 2** plate-reading luminometer.

## IV. Assay Protocol

Review the entire Assay Protocol before starting. Completing the assay requires an overnight incubation. *Steps 1-11* are performed on **Day 1**, requiring less than 2 hours of bench work to complete, but including a 4-hour incubation step. *Steps 12-17* are performed on **Day 2** and require less than 1 hour to complete.

### ▪ A word about Antagonist-mode assay setup ▪

Receptor inhibition assays expose the Reporter Cells to a constant, sub-maximal concentration (typically between EC<sub>50</sub> – EC<sub>85</sub>) of a known agonist AND the test compound(s) to be evaluated for antagonist activity. This rbPGR assay kit includes a 100 nM stock solution of Progesterone, a potent physiological agonist of PGR that may be used to setup antagonist-mode assays. 34 pM progesterone approximates EC<sub>70-80</sub> in this cell-based assay. Hence, it is an appropriate concentration of agonist to be used when screening test compounds for inhibitory activity.

Add the challenge agonist (progesterone) to a bulk volume of **CSM** at an EC<sub>50</sub> – EC<sub>85</sub> concentration. This medium is then used to prepare serial dilutions of test compounds to achieve the desired respective final assay concentrations. This is an efficient and precise method of setting up PGR antagonist assays, and it is the method presented in *Step 7b* of this protocol.

### DAY 1 Assay Protocol: All steps should be performed using aseptic technique.

**1.)** Remove the **2 tubes** of **Cell Recovery Medium (CRM)** from freezer storage, thaw and equilibrate to 37°C using a water bath.

**2.) Rapid Thaw of the Reporter Cells:** *First*, retrieve the two tubes of **CRM** from the 37°C water bath and sanitize their outside surfaces with a 70% ethanol swab.

*Second*, retrieve tubes of **Reporter Cells** from -80°C storage and place them directly into dry ice for transport to the laminar flow hood: 1 tube for 32 assay wells, 2 tubes for 64 assay wells, or 3 tubes for 96 assay wells. When ready to begin, transfer the tube(s) of reporter cells into a rack and, *without delay*, perform a rapid thaw of the frozen cells by transferring **6.4 ml** of pre-warmed CRM into each tube of frozen cells. Recap the tube of Reporter Cells and immediately place it in a 37°C water bath for 5 - 10 minutes. The resulting volume of cell suspension will be **7.0 ml** per tube.

*Third*, during the 5 - 10 minutes incubation period, work in the cell culture hood to *carefully* mount four sterile 8-well strips into the blank assay plate frame. Strip-wells are fragile. Note that they have keyed ends (square and round), hence, they will fit into the plate frame in only one orientation.

**3.)** Retrieve the tube of Reporter Cell Suspension from the water bath and sanitize the outside surface with a 70% alcohol swab.

**4.)** If more than one tube of Reporter cells was thawed, combine them and gently invert several times to gain a homogenous cell suspension. Dispense **200 µl / well** of cell suspension into the assay plate.

*NOTE 4.1:* If INDIGO's Live Cell Multiplex Assay is to be incorporated, a minimum of 3 'cell blank' wells (meaning cell-free, but containing 'Compound Screening Media') must be included in the assay plate to allow quantification of plate-specific fluorescence background (refer to the LCMA Technical Manual).

*NOTE 4.2:* Increased well-to-well variation (= increased standard deviation!) will occur if care is not taken to prevent cells from settling in the reservoir during the dispensing period. Likewise, take care to ensure precision in dispensing exact volumes across the assay plate.

*NOTE 4.3:* Users sometimes wish to examine the reporter cells using a microscope. If so, the extra volume of cell suspension provided with each kit may be dispensed into a clear, preferably collagen-coated, 96-well assay plate. Continue to process this plate in identical manner to the white assay plate.

**5.) Pre-incubate reporter cells.** Place the assay plate into a cell culture incubator (37°C,  $\geq 70\%$  humidity, 5% CO<sub>2</sub>) for 4 - 6 hours.

**6.)** Near the end of the pre-incubation period remove Compound Screening Medium (CSM) from freezer storage and thaw in a 37°C water bath.

**7.) Prepare the Test Compound and Reference Compound treatment media at the desired final assay concentrations:** Use CSM to prepare an appropriate dilution series of the reference and test compound stocks. Prepare all treatment media at the desired final assay concentrations. In *Step 9*, **200  $\mu$ l / well** of the prepared treatment media are dispensed into the strip-wells of the assay plate. Manage dilution volumes carefully; this assay kit provides **45 ml** of CSM.

NOTE: Total DMSO carried over into assay reactions should not exceed 0.4%.

**a. Agonist-mode assays.** This rbPGR Assay kit includes a 100 nM stock solution of Progesterone, a potent reference agonist of rabbit PGR. The following 7-point treatment series, with concentrations presented in 2-fold decrements, provides a complete dose-response: 100, 50, 25, 12.5, 6.25, 3.13, and 1.56 pM. Always include 'no treatment' (or 'vehicle') controls. **APPENDIX 1** provides guidance for generating such a dilution series.

~ or ~

**b. Antagonist-mode assays.** When setting antagonist assays, first supplement a bulk volume of CSM with the challenge agonist Progesterone to achieve the desired final assay-concentration (refer to "*A word about antagonist-mode assay setup*", pg. 7). The agonist-supplemented CSM is then used to generate dilutions of test compound stocks to achieve their final assay concentrations.

**8.)** At the end of the cell pre-incubation period, discard the culture media. Because the assay plate is composed of a frame with snap-in strip-wells, the practice of physically ejecting media is NOT advised. Complete removal of the media is efficiently performed by tilting the plate on edge and aspirating media using either a single-tip, or an 8-pin manifold (*e.g.*, Wheaton Science Microtest Syringe Manifold, # 851381) affixed to a vacuum-trap apparatus. Do *not* touch the well bottom or run the tip of the aspiration device around the bottom circumference of the assay well. Such practices will result in destruction of the cells and greatly increased well-to-well variability.

**9.)** Dispense **200  $\mu$ l** of each treatment media into appropriate wells of the assay plate.

**10.)** Transfer the assay plate into a cell culture incubator for 22 - 24 hours.

NOTE: Ensure a high-humidity ( $\geq 70\%$ ) environment within the cell culture incubator. This is critical to prevent the onset of deleterious "edge-effects" in the assay plate.

**11.)** For greater convenience on *Day 2*, retrieve the appropriate number of vials of **Detection Substrate and Detection Buffer** from freezer storage and place them in a dark refrigerator (4°C) to thaw overnight.

**DAY 2 Assay Protocol:** Subsequent manipulations do *not* require special regard for aseptic technique and may be performed on a bench top.

**12.)** Approximately 30 minutes before intending to quantify receptor activity, remove **Detection Substrate** and **Detection Buffer** from the refrigerator and place them in a low-light area so that they may equilibrate to room temperature.

*NOTE:* Do NOT actively warm Detection Substrate above room temperature. If these solutions were not allowed to thaw overnight at 4°C, a room temperature water bath may be used to expedite thawing.

**13.)** Set the plate-reader to "luminescence" mode. Set the instrument to perform a single 5 second "plate shake" prior to reading the first assay well. Read time may be set to 0.5 second (500 mSec) per well, *or less*.

**14.)** *Immediately before proceeding to Step 15:* To read 32 assay wells, transfer the entire volume of 1 vial of Detection Buffer into 1 vial of Detection Substrate, thereby generating a 4 ml volume of **Luciferase Detection Reagent (LDR)**. Mix gently to avoid foaming.

**15.)** Following 22 - 24 hours incubation in treatment media, remove media contents from each well of the assay plate (as before in *Step 8*).

**16.)** Use an 8-channel pipette to dispense 100  $\mu$ l of **LDR** to each well of the assay plate. Allow the plate to rest at room temperature for 5 - 10 minutes following the addition of LDR. Do not shake the plate during this period.

**17.)** Quantify luminescence.

## V. Related Products

| Product No.                                                                                                | Product Descriptions                                                                                                          |
|------------------------------------------------------------------------------------------------------------|-------------------------------------------------------------------------------------------------------------------------------|
| <b>Human PGR Assay Products</b>                                                                            |                                                                                                                               |
| IB05001-32                                                                                                 | Human PGR Reporter Assay System<br>3x 32 assays in 8-well strips (96-well plate format)                                       |
| IB05001                                                                                                    | Human PGR Reporter Assay System<br>1x 96-well format assay                                                                    |
| IB05002                                                                                                    | Human PGR Reporter Assay System<br>1x 384-well format assays                                                                  |
| <b>Rat PGR Assays</b>                                                                                      |                                                                                                                               |
| R05001-32                                                                                                  | Rat PGR Reporter Assay System<br>3x 32 assays in 8-well strips (96-well plate format)                                         |
| R05001                                                                                                     | Rat PGR Reporter Assay System<br>1x 96-well format assay                                                                      |
| <b>Rabbit PGR Assays</b>                                                                                   |                                                                                                                               |
| RB05001-32                                                                                                 | Rabbit PGR Reporter Assay System<br>3x 32 assays in 8-well strips (96-well plate format)                                      |
| RB05001                                                                                                    | Rabbit PGR Reporter Assay System<br>1x 96-well format assay                                                                   |
| <b>Monkey PGR Assays</b>                                                                                   |                                                                                                                               |
| C05001-32                                                                                                  | Cynomolgus Monkey PGR Reporter Assay System<br>3x 32 assays in 8-well strips (96-well plate format)                           |
| C05001                                                                                                     | Cynomolgus Monkey PGR Reporter Assay System<br>1x 96-well format assay                                                        |
| Bulk volumes of assay reagents may be custom manufactured to accommodate any scale of HTS. Please Inquire. |                                                                                                                               |
| <b>LIVE Cell Multiplex (LCM) Assay</b>                                                                     |                                                                                                                               |
| LCM-01                                                                                                     | Reagent volumes sufficient to perform <b>96</b> Live Cell Assays in 1x96-well, or 2x48-well, or 3x32-well assay plate formats |
| LCM-05                                                                                                     | Reagent in <b>5x bulk volume</b> to perform <b>480</b> Live Cell Assays performed in 5 x 96-well assay plates                 |
| LCM-10                                                                                                     | Reagent in <b>10x bulk volume</b> to perform <b>960</b> Live Cell Assays performed in 10 x 96-well assay plates               |
| <b>INDIGlo Luciferase Detection Reagent</b>                                                                |                                                                                                                               |
| LDR-10, -25, -50, -500                                                                                     | INDIGlo Luciferase Detection Reagents in 10 mL, 25 mL, 50 mL, and 500 mL volumes                                              |

Please refer to INDIGO Biosciences website for updated product offerings.

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

## ***VI. Limited Use Disclosures***

Products offered by INDIGO Biosciences, Inc. are for RESEARCH PURPOSES ONLY – not for therapeutic, diagnostic, or contact use in humans or animals.

“CryoMite” is a Trademark <sup>TM</sup> of INDIGO Biosciences, Inc. (State College, PA, USA).

Product prices, availability, specifications, claims and technical protocols are subject to change without prior notice. The printed Technical Manual provided in the kit box will always be the most current version available.

Copyright © INDIGO Biosciences, Inc. (State College, PA, USA). All rights reserved.

Example scheme for the serial dilution of Progesterone reference agonist, and the setup of a rabbit PGR dose-response assay.

