

**Human Hepatocyte Nuclear Factor 4 alpha
(NR2A1, HNF4 α)
Reporter Assay System**

384-well Format Assays
Product # IB36002

▪

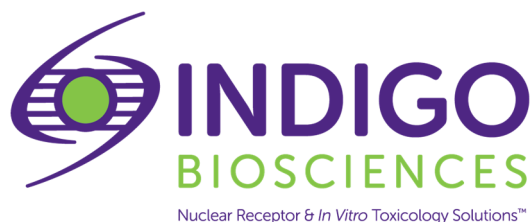
Technical Manual
(version 8.0i)

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

Technical Service:
814-234-1919
techserv@indigobiosciences.com



**Human HNF4 α Reporter Assay
 384-well Format Assays**

I. Description

- The Assay System.....3
- The Assay Chemistry.....3
- Considerations for the Preparation and Automated Dispensing of Test Compounds.....4
- Considerations for Automated Dispensing of Other Assay Reagents.....4
- Assay Scheme.....5
- Assay Performance.....6

II. Product Components & Storage Conditions7

III. Materials to be Supplied by the User.....7

IV. Assay Protocol

- *DAY 1 Assay Protocol*.....8
- *DAY 2 Assay Protocol*.....10

V. Literature Citations.....11

VI. Related Products.....11

VII. Limited Use Disclosures.....11

APPENDIX 1a: Example Scheme for Serial Dilution when using tip-based dispensing of test compounds.....12

APPENDIX 1b: Example Scheme for Serial Dilutions when using acoustic dispensing of test compounds.....13

I. Description

▪ The Assay System ▪

This nuclear receptor assay utilizes proprietary human cells engineered to provide constitutive, high-level expression of the **Human Hepatocyte Nuclear Factor 4 alpha** (NR2A1), a ligand-dependent transcription factor known as **HNF4 α** .

In INDIGO's Reporter Cells the HNF4 α receptor is expressed as a hybrid protein in which its native DNA binding domain (DBD) sequence has been exchanged with that of the DBD sequence from the yeast Gal4 gene. Additionally, the cells include the luciferase reporter gene functionally linked to the Gal4 upstream activation sequence (UAS) and a minimal promoter. HNF4 α 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 reporter gene. Thus, quantifying changes in luciferase expression in the treated reporter cells provides a specific and sensitive surrogate measure of the changes in HNF4 α activity. Unlike some other assay formats, 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*.

As is true *in vivo*, the expressed hybrid HNF4 α receptor displays a moderate level of constitutively activity. A true ligand interaction may provoke either *i.*) an *agonist* response that delivers an additional increase in receptor activity above its already high constitutive level of activity, or *ii.*) an *inverse-agonist* response in which a loss of receptor activity is observed. **Figures 2 and 3** demonstrates these responses. The principal application of this assay is in the screening of test compounds to quantify *inverse-agonist* or *agonist* activities that they may exert against the Human HNF4 α receptor.

These HNF4 α Reporter Cells are prepared using INDIGO's proprietary **CryoMite™** process. This cryo-preservation method yields high 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.

This kit is an all-inclusive assay system. As detailed on page 6, it includes cryopreserved HNF4 α Reporter Cells, two optimized media for use during initial cell culture and for preparing dilutions of test samples, a control agonist, luciferase detection reagents, a cell culture-ready assay plate, and a detailed assay protocol.

▪ The Assay Chemistry ▪

INDIGO's receptor reporter assays 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 O₂ and ATP as co-substrates, and yields as products oxyluciferin, AMP, PP_i, CO₂, 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 Assays feature a luciferase detection reagent specially formulated to provide stable light emission between 30 and 100+ minutes after initiating the luciferase reaction. Incorporating a 30-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.

▪ **Considerations for the Preparation and Automated Dispensing of Test compounds** ▪

Small molecule compounds are typically solvated at high concentration (ideally 1,000x-concentrated) in DMSO and stored frozen as master stocks. For **384-well format assays** these master stocks will be diluted by one of two alternative methods, the selection of which will be dictated by the type of dispensing instrument that is to be used. This Technical Manual provides detailed protocols for each of these two alternative methods:

- a.) Assay setups in which a conventional **tip-based** instrument is used to dispense test compounds into assay wells (in black text). Use **Compound Screening Medium (CSM)** to generate a series of **2x-concentration** test compound treatment media, as described in *Step 2a* of the **Assay Protocol**. The final concentration of DMSO carried over into assay reactions should not exceed 0.4%; strive to use 1,000x-concentrated stocks when they are prepared in any organic solvent.

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.

and,

- b.) Assay setups in which an **acoustic transfer** device is used to dispense test compounds into assay wells (text in blue text). Use DMSO to make a series of **1,000x-concentrated** test compound stocks that correspond to each desired final assay concentrations, as described in *Step 2b* of the **Assay Protocol**.

▪ **Considerations for Automated Dispensing of Other Assay Reagents** ▪

When dispensing into a small number of assay plates, first carefully consider the dead volume requirement of your tip-based 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 on a *per kit* basis. Always pool the individual reporter cell suspensions and all other respective assay kit reagents before processing multiple 384-well assay plates.

| Stock Reagent & Volume provided | Volume to be Dispensed (384-well plate) | Excess rgt. volume available for instrument dead volume |
|---|--|--|
| <i>when using tip dispensing of test cmpds</i> Reporter Cell Suspension 7.5 ml | 15 µl / well 5.8 ml / plate | ~ 1.7 ml |
| <i>when using acoustic dispensing of test cmpds</i> Reporter Cell Suspension 15 ml | 30 µl / well 11.5 ml / plate | ~ 3.4 ml |
| Detection Substrate 7.8 ml | 15 µl / well 5.8 ml / plate | ~ 2 ml |

▪ Assay Scheme ▪

The *Day 1* preparation, volumes, and chronology of dispensed cells and test compounds are different between assay setups using a *tip-based dispenser* (**1a**) and those using an *acoustic transfer device* (**1b**). Following 22 -24 hours incubation Detection Substrate is added. Light emission from each assay well is quantified using a plate-reading luminometer.

Figure 1a. Assay workflow if using conventional **tip-based** dispensing of test compounds.

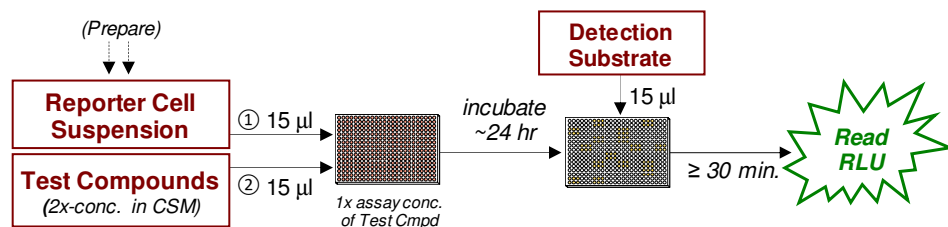
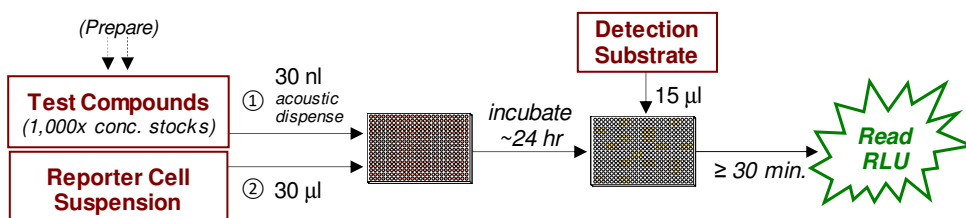


Figure 1b. Assay workflow if using **acoustic** dispensing of test compounds.



▪ Assay Performance ▪

Human HNF4 α , Agonist Assays
(Fold-Activation)

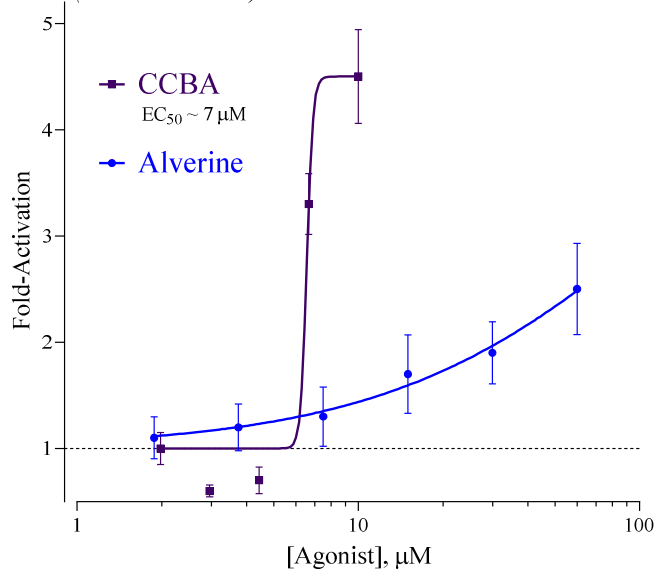


Figure 2. Agonist dose-response of Human HNF4 α .

Agonist dose-response of Human HNF4 α reporter cells using CCBA¹ (provided) and Alverine Citrate² (Millipore Sigma). Luminescence was quantified and average relative light units (RLU) and their corresponding values of standard deviation (SD) and Fold-Activation were determined for each treatment group (n = 3). GraphPad Prism software was used to curve-fit data using the least-squares method of non-linear regression, and EC₅₀ values were determined.

Human HNF4 α , Inhibition Assay
(Fold-Reduction)

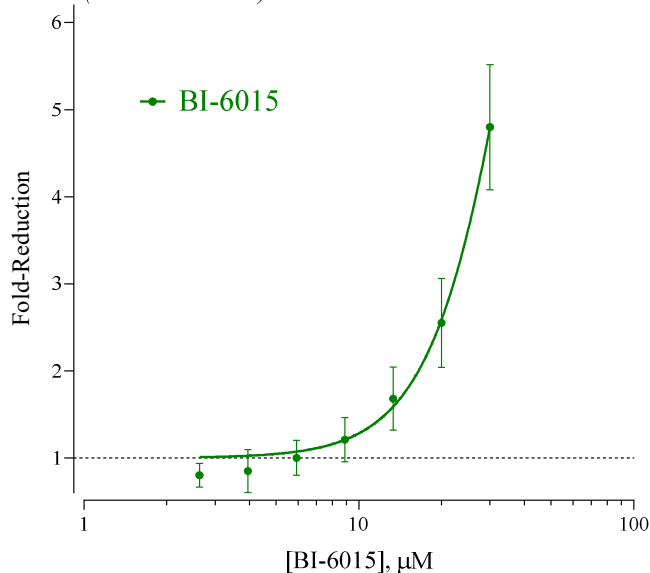


Figure 3. Inverse-Agonist dose-response of Human HNF4 α .

Inverse-agonist response of HNF4 α reporter cells using BI6015³ (Cayman Chem.). Luminescence was quantified and average relative light units (RLU) and their corresponding values of standard deviation (SD) and Fold-Reduction were determined for each treatment group (n = 3). INDIGO's Live Cell Multiplex assay was performed and confirmed the absence of compound-induced cytotoxicity (data not shown).

II. Product Components & Storage Conditions

This assay kit contains materials to perform assays in a single 384-well assay plate.

Cryopreserved mammalian 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 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 inventory the individual kit components be sure to first transfer and submerge the tube of cells in dry ice.

The aliquot of Reporter Cells is provided as a single-use reagent. Once thawed, the 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> |
|---|-----------------|----------------------|
| ▪ HNF4 α Reporter Cells | 1 x 1.0 mL | -80°C |
| ▪ Cell Recovery Medium (CRM) | 1 x 7 mL | -20°C |
| ▪ Compound Screening Medium (CSM) | 1 x 45 mL | -20°C |
| ▪ CCBA ¹ , 10 mM (in DMSO) (control agonist for HNF4 α) | 1 x 100 μ L | -20°C |
| ▪ Detection Substrate (Note: contains DTT) | 1 x 7.8 mL | -80°C |
| ▪ 384-well assay plate (white, sterile, cell-culture ready) | 1 | ambient |

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

- dry ice container
- cell culture-rated laminar flow hood.
- 37°C, humidified 5% CO₂ incubator for mammalian cell culture.
- 37°C water bath.
- 70% alcohol wipes
- 8-channel electronic, repeat-dispensing pipettes & tips suitable for dispensing 15 μ L.
- disposable media basins, sterile.
- sterile multi-channel media basins *or* deep-well plates, *or* appropriate similar vessel for generating dilution series of reference compound(s) and test compound(s).
- *Optional*: inverse-agonist/antagonist reference compound / antibody (*e.g.*, Fig. 3.)
- *Optional*: clear 384-well assay plate 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 1-8* are performed on **Day 1**, requiring less than 2 hours to complete. *Steps 9-13* are performed on **Day 2** and require less than 1 hour to complete.

DAY 1 Assay Protocol:

All steps should be performed using proper aseptic technique.

- 1.) Remove **Cell Recovery Medium (CRM)** and **Compound Screening Medium (CSM)** from freezer storage and thaw in a 37°C water bath.
- 2.) **Prepare dilutions of treatment compounds:** Prepare Test Compound treatment media for *Agonist-* or *Antagonist-mode* screens. NOTE that test and reference compounds will be prepared differently when using tip-dispensing vs. [acoustic dispensing](#). Regardless of the method, residual DMSO carried over into assay reactions should not exceed 0.4%.
 - a. *Tip dispensing method:* In *Step 6*, 15 µl / well of the prepared treatment media is added to the assay that has been pre-dispensed with 15 µl /well of Reporter Cells. Hence, to achieve the desired *final* assay concentrations one must prepare treatment media with a **2x**-concentration of the test and reference material(s). Use **CSM** to prepare the appropriate dilution series. Plan dilution volumes carefully; this assay kit provides 35 ml of CSM.
 - b. *Acoustic dispensing method:* In *Step 6*, 30 nl / well of **1,000x**-concentrated test compound solutions are added to the assay plate using an acoustic transfer device.

Preparing the positive control: This assay kit includes a 10 mM stock solution of CCBA, an agonist of HNF4α. The following 7-point treatment series, with concentrations presented in **1.5-fold** decrements, provides a complete dose-response: 10, 6.67, 4.44, 2.96, 1.98, 1.32 and 0.878 µM. Always include 'no treatment' (or 'vehicle') control wells.

APPENDIX 1a provides an example for generating such a dilution series to be used when *tip-dispensing* compound solutions prepared in CSM (15 µl / well).

APPENDIX 1b provides an example for generating such a series of 1,000x-concentrated solutions of compounds prepared in DMSO to be used when performing *acoustic dispensing* (transfer 30 nl / well).

When using *tip-based* instrumentation for dispensing test compounds ...

- 3.) *First*, retrieve the tube of **CRM** from the 37°C water bath, sanitize the outside with a 70% ethanol swab;
Second, retrieve **Reporter Cells** from -80°C storage and immerse in dry ice to transport the tube to a laminar flow hood. Perform a *rapid thaw* of the frozen cells by transferring a 6.5 ml volume of 37°C CRM into the tube of frozen cells. Recap the tube of Reporter Cells and place it in a 37°C water bath for 5 - 10 minutes. The resulting volume of cell suspension will be 7.5 ml.
- 4.) Retrieve the tube of Reporter Cell Suspension from the water bath. Sanitize the outside surface of the tube with a 70% alcohol swab, then transfer it into the cell culture hood.
- 5.) *Gently* invert the tube of cells several times to gain a homogenous suspension.
 - a. **for Agonist-mode assays:** Dispense **15 µl / well** of cell suspension into the assay plate.
- 6.) Dispense **15 µl / well** of 2x-concentrated treatment media (from *Step 2a*) into the assay plate.

(continued)

When using an *acoustic transfer* device for dispensing test compounds ...

3.) Dispense **30 nl / well** of the 1,000x-concentrated compounds (in DMSO solutions, from *Step 2b*) into the assay plate.

4.) *First*, retrieve the tube of **CRM** from the 37°C water bath, sanitize the outside with a 70% ethanol swab;

Second, retrieve **Reporter Cells** from -80°C storage and immerse in dry ice to transport the tube to a laminar flow hood. Perform a *rapid thaw* of the frozen cells by transferring a **6.5 ml** volume of 37°C CRM into the tube of frozen cells. Recap the tube of cells and place it in a 37°C water bath for 5 - 10 minutes. The resulting volume of cell suspension will be 7.5 ml.

5.) Retrieve the tube of cell suspension from the water bath. Sanitize the outside surface of the tube with a 70% alcohol swab. Add an additional **7.5 ml** of **CSM** to the tube. The resulting volume of cell suspension will be 15 ml.

6.) *Gently* invert the tube of cells several times to gain a homogenous cell suspension. Transfer into an appropriate media reservoir, then dispense **30 µl / well** of cell suspension into the Assay Plate that has been pre-dispensed with test compounds.

NOTE: Take special care to prevent cells from settling during the dispensing period. Allowing cells to settle during the transfer process, and/or lack of precision in dispensing uniform volumes across the assay plate *will* cause well-to-well variation (= increased Standard Deviation) in the assay.

NOTE: Following the dispensing of Reporter Cells and test compounds INDIGO recommends performing a *low-speed* spin of the assay plate (with lid) for ≤ 1 minute using a room temperature centrifuge fitted with counter-balanced plate carriers.

7.) Transfer the assay plate into a 37°C, humidified, 5% CO₂ 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.

8.) For greater convenience on *Day 2*, retrieve **Detection Substrate** from freezer storage and place 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 or in a **fume hood**.

9.) Approximately 30 minutes before intending to quantify receptor activity remove **Detection Substrate** from the refrigerator and place it in a low-light area so that it may equilibrate to room temperature. Gently invert the tube several times to ensure a homogenous solution.

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

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

11.) Following 22 - 24 hours of incubation dispense **15 µl / well** of **Detection Substrate** into all wells of the assay plate.

NOTE: 'Detection Substrate' contains a high concentration of DTT, which produces a strong odor that some users may find objectionable. It is advised to work in a **fume hood** when dispensing Detection Substrate into the assay plate and the following 'plate rest' period.

NOTE: Perform this reagent transfer carefully to avoid bubble formation! Scattered micro-bubbles will not pose a problem. However, bubbles covering the surface of the reaction mix, or large bubbles clinging to the side walls of the well, will cause lens-effects that will degrade the accuracy and precision of the assay data. INDIGO recommends performing a final *low-speed* spin of the assay plate (with lid) for ≤ 1 minute using a room temperature centrifuge fitted with counter-balanced plate carriers.

12.) Allow the plate(s) to rest at room temperature for 30 minutes. Do not shake the assay plate(s) during this period.

NOTE: the luminescent signal is unstable during the first 30 minutes of the luciferase reaction, however, after the initial 30-minute reaction period the luminescence signal achieves a stable emission output.

13.) Quantify luminescence.

14.) Data analyses.

V. Literature Citations

- ¹ 5-Chloro-n-(2-chlorophenyl)-2-hydroxybenzamide (CCBA; CAS #6626-92-2)
- ² Lee, S., *et al.*, (2013) Identification of Alverine and Benfluorex as HNF4 α Activators. *ACS Chem. Biol.*: **8**, 1730-1736.
- ³ Kiselyuk, A., *et al.* (2012) HNF4 α Antagonists Discovered by a High-Throughput Screen for Modulators of the Human Insulin Promoter. *Chemistry & Biology*:**19**, 806-818.
- ⁴ 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.

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

VI. Related Products

| Product No. | Product Descriptions |
|---|---|
| Human Hepatocyte Nuclear Factor 4 alpha | |
| IB36001 | Human Hepatocyte Nuclear Factor 4 alpha (HNF4 α) 1x 96-well format assay |
| IB36002 | Human Hepatocyte Nuclear Factor 4 alpha (HNF4 α) 1x 384-well format assays |
| LIVE Cell Multiplex (LCM) Assay | |
| LCM-01 | Reagent volumes sufficient to perform 96 Live Cell Assays in 1x96-well, or 2x48-well, or 3x32-well assay plate formats |
| LCM-05 | Reagent in 5x bulk volume to perform 480 Live Cell Assays performed in 5 x 96-well assay plates |
| LCM-10 | Reagent in 10x bulk volume to perform 960 Live Cell Assays performed in 10 x 96-well assay plates |
| INDIGlo Luciferase Detection Reagent (LDR) | |
| 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

VII. 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 TM 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 recently updated version available.

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

APPENDIX 1b for acoustic dispensing. Example scheme for the serial dilution of the CCBA using **DMSO** to generate **1,000x-concentrated** stocks. 30 nl / well of these prepared stocks are first dispensed into *empty* wells of the assay plate using an acoustic transfer device, followed by the dispensing of 30 μ l of HNF4 α reporter cells.

