

# Investigation of a Novel, High-throughput Method for Real-time Measurement of $\beta$ -arrestin Recruitment using a Live-cell Luminescence Based Assay System

P. Brescia<sup>1</sup>, P. Banks<sup>1</sup>, A. Paguio<sup>2</sup>, B. Binkowski<sup>2</sup> and A. Landreman<sup>2</sup>

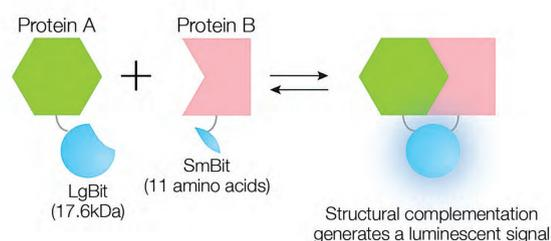
<sup>1</sup>BioTek Instruments, Inc. | Winooski, VT | USA • <sup>2</sup>Promega Corporation | Madison, WI | USA



## Abstract

GPCRs and their involvement in cellular signaling remain a focal point of concerted efforts into the investigation of druggable targets. Investigation of protein:protein interactions (PPIs) involved in GPCR signaling, such as the GPCR: $\beta$ -arrestin2 interaction, provides a means to better understand pathways involved with a variety of diseases such as cancer. The ability to perform live-cell measurements of these interactions remove many of the limitations of other approaches by providing real-time, kinetic results. The development of the NanoLuc<sup>®</sup> Binary Technology (NanoBiT<sup>®</sup>) in combination with Nano-Glo<sup>®</sup> Live Cell Reagent, allows real-time measurement of PPI dynamics in living cells using a simple bioluminescent readout. Here we show the ability to perform a GPCR:  $\beta$ -arrestin2 recruitment assay in a high-throughput, 1536-well microplate format using automated liquid handling. The NanoBiT CX3CR1/ARRB2 stable cell line was used to demonstrate this PPI when activated by addition of CX3CL1 ligand (fractalkine). A dose response titration of CX3CL1, as well as Z'-factor determination to assess assay performance, are demonstrated here.

## NanoBiT Overview



**Figure 1. The NanoBiT<sup>®</sup> assay.** LgBit and SmBit are fused to proteins A & B. A:B interaction facilitates LgBit:SmBit interaction, generating a bright luminescent enzyme. Non-lytic assay format allows real-time measurements of protein interaction dynamics for 1-2 hours.

## BioTek Instrumentation

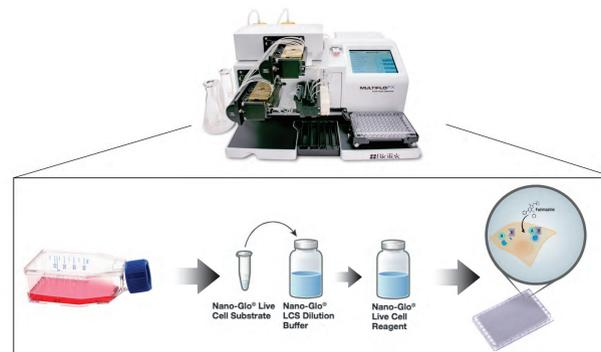


**Figure 2. Synergy<sup>™</sup> Neo2 Multi-Mode Microplate Reader.** The Synergy Neo2 Multi-Mode Microplate Reader is designed for speed and ultra high performance, incorporating BioTek's patented Hybrid Technology<sup>™</sup>, with its independent optical paths that ensure uncompromised performance. Continuously variable bandwidth quadruple monochromators, sensitive high transmission filter-based optics and up to 4 PMTs provide ultra-fast measurements with excellent results. Advanced environment controls, including available CO<sub>2</sub>/O<sub>2</sub> control, incubation to 65 °C and variable shaking support live cell assays; cell-based detection is optimized with direct bottom illumination.



**Figure 3. MultiFlo FX<sup>™</sup> Multimode Dispenser.** MultiFlo FX is an automated multi-mode reagent dispenser for 6- to 1536-well microplates. MultiFlo FX incorporates several unique technologies in its modular design, such as Parallel Dispense RAD<sup>™</sup> Random Access Dispense and the new, patent-pending AMX<sup>™</sup> Automated Media Exchange modules to facilitate a variety of liquid handling applications from 2D and 3D cell culture to concentration normalization assays, ELISA, bead-based assays and more. A fully configured MultiFlo FX replaces up to five liquid handlers, saving space, time and instrumentation budgets. MultiFlo FX integrates to the BioSpa<sup>™</sup> 8 Automated Incubator and a BioTek imager or multi-mode reader, for complete workflow automation for many cell imaging and biochemical applications.

## NanoBiT Procedure



Record luminescence at a single time point or continuously for up to 2 hours

**Figure 4. NanoBiT<sup>®</sup> Assay workflow.**

## Materials and Methods

NanoBiT<sup>®</sup> CX3CR1/ARRB2 cell line, HEK293 (PN:CS208104A, Promega Corp.) were cultured using standard tissue culture methods as per the manufacturers' recommendations. Blastocidin (PN:A11139-03, ThermoFisher) at 5 $\mu$ g/mL was used to maintain selective pressure. The cells were harvested at 80-90% confluence using TrypLE dissociation reagent (PN:12605036, Fisher Scientific) with gentle handling. The cells were collected by centrifugation and resuspended at the desired cell density in OptiMEM 1 (OM1) (PN:11058-021, ThermoFisher) for serum starvation during incubation at 37 °C, 5% CO<sub>2</sub> for 4-6 hours. Nano-Glo Live Cell Reagent (PN: N2012, Promega Corp.) was prepared as a 5X stock as per the manufacturers' recommendations and added to the cell suspension with mixing. Cells were plated in a volume of 5  $\mu$ L at a concentration of 600,000 cells/mL resulting in 3,000 cells/well using the MultiFlo FX. Baseline luminescence measurements were taken for ~10 minutes at ambient temperature using a Neo2. Fractalkine was prepared as an 8-point 1:3 serial dilution series in OM1 at 2x concentration and added as 8 replicates in equal 5  $\mu$ L volume. Luminescence measurements were taken using the minimal read time interval (~ every 90 seconds) for 60 minutes.

## Synergy Neo2 Reader Settings

Synergy Neo2 Reader Parameters	
Mode	Lum
Light Path	Single PMT
Optic Position	Top
Gain (PMT1, PMT2)	135
Integration Time	0.2 sec
Read Height	8 mm
Delay	0 msec

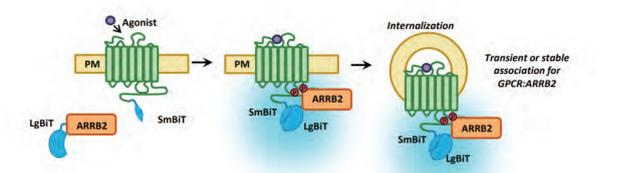
**Table 1. Synergy Neo2.** Synergy Neo2 equipped with a luminescence filter cube fitted with a 1536-well aperture was used, with the above settings, to rapidly capture kinetic measurements.

- The Synergy Neo2 was fitted with a 1536-well aperture (3.5 mm)
- The read mode selected was Luminescence in the Top position to take advantage of the high-sensitivity luminescence fiber-optic light-guide
- Photo-Multiplier Tube (PMT) gain setting was optimized and set to 135.
- The integration times was set to 0.2 seconds which showed sufficient signal-to-background.
- Read height was set to 8 mm for use with a 1536-well plate.

## NanoBiT<sup>®</sup> Calculations

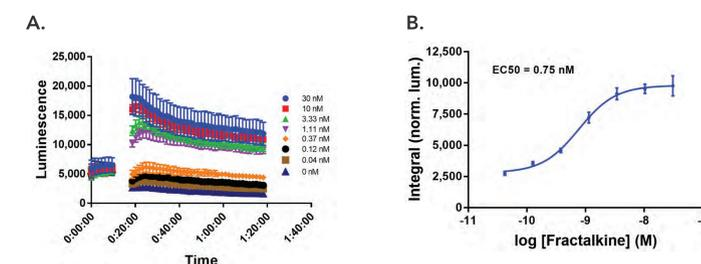
- Data normalization to account for cell plating differences was performed by dividing each data point, relative luminescence units (RLUs) post-fractalkine addition, by the final RLU measurement taken just prior to fractalkine addition.
- The integral under the curve was then calculated using kinetic measurement data post-fractalkine addition.
- Z'-factor was calculated using 36 replicate data points for negative and positive control wells using vehicle alone and the highest concentrating of fractalkine from the dose response titration, 0 and 30 nM, respectively.

## GPCR Activation Pathway



**Figure 5. NanoBiT<sup>®</sup> CX3CR1/ARRB2 cell line.** Stable expression of CX3CR1/ARRB2 in HEK293 cells allow the real-time investigation of GPCR: $\beta$ -arrestin2 recruitment.

## Fractalkine Titration and Z'-Factor

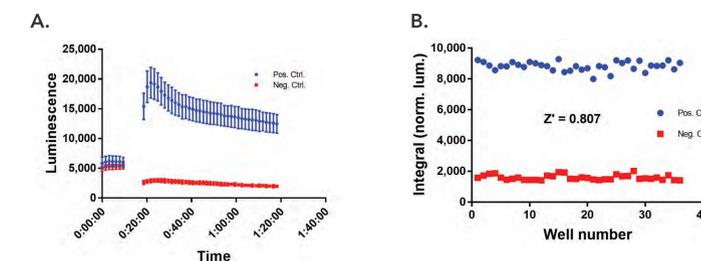


**Figure 6. Agonist titration.** (A) Raw kinetic luminescence data. (B) Fractalkine titration dose response curve based on integral of area under the kinetic curve post fractalkine addition.

Fractalkine EC <sub>50</sub> (nM)	
1536-well assay	0.75

**Table 3. EC<sub>50</sub> Value.** The EC<sub>50</sub> value for the agonist Fractalkine was determined in a high-throughput, 1536-well assay format.

- Cells were plated as 8 replicates for each agonist concentration in a 1536-well, solid white microplate.
- The luminescence baseline signal was measured on the Synergy Neo2 for ~10 mins. prior to agonist addition.
- The plate was removed and an 8-point, 1:3 serial dilution series of the CX3CR1 agonist fractalkine, including a zero compound point, was added in an equal volume using the MultiFlo FX.
- The luminescence signal was read kinetically for 60 minutes immediately following agonist addition.
- The EC<sub>50</sub> concentration was determined using a four-parameter dose-response curve fit in Prism (GraphPad Software, Inc., La Jolla, CA)(Figure 6).
- An EC<sub>50</sub> value of 0.75 nM correlate well with previously published data (Table 3).



**Figure 7. Z' Factor determination.** (A) Raw kinetic luminescence data. (B) Control measurement based on integral of area under the kinetic curve post fractalkine addition.

Z'-Factor	
1536-well assay	0.83

**Table 4. Z'- Factor.**

- The Z factor was calculated using 36 replicate measurement of +/- fractalkine (30 nM).
- The assay resulted in Z' factor of 0.83 indicative of very robust assay performance with low variability.

## Conclusions

- The live-cell NanoBiT<sup>®</sup> assay allowed for the investigation of dynamic receptor recruitment within a biologically relevant environment in a format amenable to high-throughput screening efforts.
- A dose response of fractalkine was performed using CX3CR1-LgBit:SmBit-AARB2 in a HEK293 background resulted in a EC<sub>50</sub> = 0.75 which correlates with previous reported values.
- The Synergy Neo2 provides rapid detection which is necessary for high-throughput assay formats. Read time was < 8 min for 1536-wells with the above reader parameters while still providing a highly robust assay (z' = 0.83).
- The combination of assay and instrumentation provide an ideal solution for high-throughput detection of protein:protein interactions.