

# CRISPR/Cas9 genome-edited cells express nanoBRET-donor that monitors protein interaction and trafficking

Carl White<sup>1,2</sup>, Ethan See<sup>1,2</sup>, and Kevin Pfeleger<sup>1,2,3</sup>

<sup>1</sup>Molecular Endocrinology and Pharmacology, Harry Perkins Institute of Medical Research, Australia <sup>2</sup>Centre for Medical Research, The University of Western Australia, Australia <sup>3</sup>Dimerix Limited, Nedlands, Australia

- CRISPR/Cas9 genome editing enables endogenous expression and regulation of BRET donor-labelled proteins
- Two proteins fused to differing acceptor fluorophores monitor receptor trafficking
- CLARIOstar® exhibits high sensitivity in nanoBRET measurements using filters or LVMF monochromator™

## Introduction

G protein-coupled receptors (GPCRs) are membrane-spanning proteins that transmit extracellular stimuli to the inside of a cell. Activation by stimulating molecules (e.g. neurotransmitter, hormone or chemokine) starts a signaling cascade that results in a cellular response. An example is the chemokine CXCL12 that activates the receptor CXCR4, resulting in G protein activation and recruitment of  $\beta$ -arrestin2. Subsequently, the receptor is internalized for either recycling to the plasma membrane or lysosomal breakdown. GPCRs are important drug targets requiring receptor-protein interaction and trafficking studies to reveal how they function. Bioluminescence resonance energy transfer (BRET) is a versatile tool to study such interactions and trafficking. However it is limited by the ectopic expression of labelled interaction partners. CRISPR/Cas9 genome editing overcomes the limitation by enabling endogenous expression of luciferase-labelled proteins.

## Assay Principle

BRET uses a luciferase that in the presence of a substrate produces blue light. If a suitable fluorophore is in very close proximity and appropriate orientation, less light is emitted and resonance energy is transferred. Thus, interaction of fluorophore- and luciferase-labelled proteins can be monitored by relating fluorophore emission to luciferase signal.

The CRISPR/Cas9 genome editing method incises dsDNA at a precise target site. In the presence of a donor DNA template, homology-directed repair results in the insertion of the donor sequence.

CRISPR/Cas9 enabled insertion of the DNA coding for nanoluciferase (Nluc) into the endogenous genomic locus of CXCR4 of HEK293FT cells. The resulting CXCR4/Nluc fusion protein acts as a BRET donor and overcomes the need for exogenous donor expression.

## Materials & Methods

- White 96-well plate (Greiner)
- CLARIOstar® and LUMIstar® (BMG LABTECH)
- CXCL12 (Preprotech), AMD3100 (Sigma-Aldrich), Isoproterenol (Sigma-Aldrich)
- Furimazine (Promega)

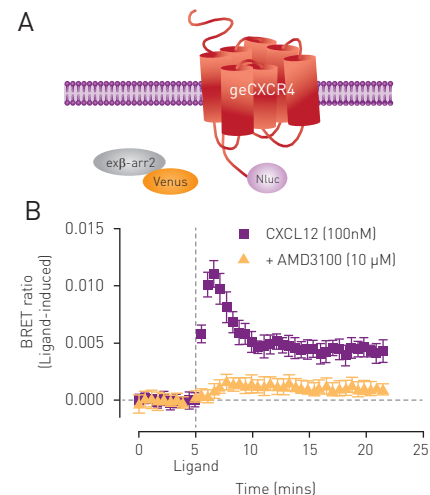
### Experimental procedure

For detailed descriptions please refer to White et al. (2017). Briefly, CRISPR/Cas9 genome-edited HEK293FT cells expressing CXCR4/Nluc were transfected with cDNA coding for the protein with acceptor-fluorophore

using Fugene®6 (Promega). 48 h after transfection cells were incubated with the luciferase substrate furimazine and filtered light emissions were analyzed on a CLARIOstar or LUMIstar Omega.

Settings stable through experiments		
Luminescence, top optic, plate mode		
Measurement interval time 1 s		
Incubation at 37 °C		
Optic settings (Filters or monochromator and [gains])		
Fig.1	Fig. 2	Fig. 3
CLARIOstar Filters	CLARIOstar Monochromator	LUMIstar Filters
Nluc 450-60 (3400)	Nluc 450-60 (3600)	Nluc 475-30 (3200)
Venus 570-100 (2800)	Venus 550-60 (3600)	Venus 535-30 (3600)
	HaloTag 660-100 (3200)	
Number of cycles / Cycle time		
40 / 33 s	30 / 49 s	60 / 31 s

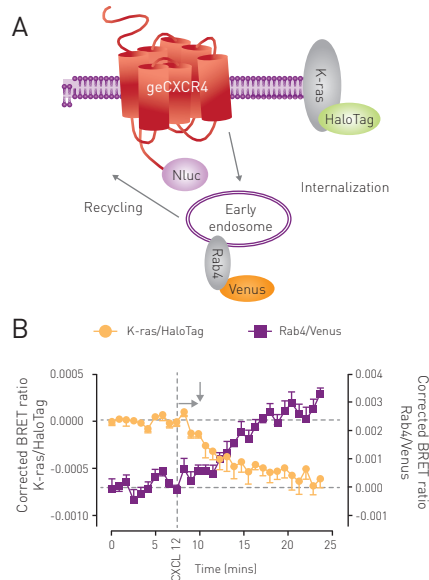
## Results & Discussion



**Fig. 1:** Monitoring  $\beta$ -arrestin2 recruitment to genome-edited CXCR4/Nluc using nanoBRET. HEK293FT cells expressing genome-edited CXCR4 fused to Nluc [geCXCR4/Nluc] transiently transfected with cDNA coding for  $\beta$ -arrestin2/Venus [ex $\beta$ -arr2/Venus] **(A)** were used to determine ligand-dependent CXCL12 (100 nM) recruitment of  $\beta$ -arrestin2 to CXCR4 in the absence or presence of the CXCR4 antagonist AMD3100 (10  $\mu$ M) **(B)**. Data previously published in White et al. (2017).

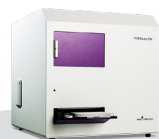


The suitability of genome-edited cells expressing CXCR4/Nluc to report on receptor interaction was studied by combining it with exogenously expressed  $\beta$ -arrestin-2/Venus [Fig. 1A]. Upon activation of CXCR4 with its ligand CXCL12,  $\beta$ -arrestin2 was recruited as reported by an increase in the BRET ratio [Fig. 1B, purple]. The CXCR4 antagonist AMD3100 inhibited CXCL12-induced recruitment of  $\beta$ -arrestin2 as seen by the reduction in energy transfer between CXCR4/Nluc and  $\beta$ -arrestin2/Venus [Fig. 1B, orange].



**Fig. 2:** HEK293FT cells expressing genome-edited CXCR4 fused to Nluc [geCXCR4/Nluc] transiently co-transfected with cDNA coding for HaloTag/K-ras and Rab4/Venus (A) were used to study genome-edited CXCR4/Nluc receptor internalization and trafficking induced by CXCL12 (30 nM) in the same cell using a BRET multiplex assay (B). Data previously published in White et al. [2017].

Genome-edited cells endogenously expressing CXCR4/Nluc were further used to study receptor internalization and trafficking. To this end, cells were transiently co-transfected with a K-ras fragment fused to HaloTag and a Rab4-Venus fusion protein serving as plasma membrane and early endosome marker, respectively [Fig. 2A]. As signals of Venus and HaloTag can be separated using the CLARIOstar's monochromator, BRET ratios of the respective proteins were determined in one experiment. Upon addition of the agonist, CXCR4 dissociates from the plasma membrane reported by a decrease in BRET ratio of the K-ras marker [Fig. 2B, orange]. At the same time, an increase in the BRET ratio and therefore proximity of the receptor and Rab4 is observed [Fig. 2B, purple]. The data indicate that the receptor is shuttled from the plasma membrane to the early endosome.



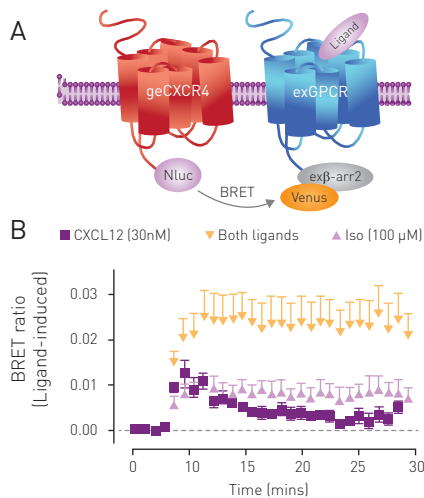
PHERAstar® FSX



CLARIOstar®



Omega Series



**Fig. 3:** GPCR-Heteromer Investigation Technology (HIT) BRET assay using genome-edited CXCR4. HEK293FT cells expressing genome-edited CXCR4/Nluc [geCXCR4/Nluc] were transiently transfected with cDNA coding for  $\beta$ -arrestin2/Venus and  $\beta$ -adrenoceptor (A) to carry out BRET assays using the GPCR-HIT configuration. Cells were stimulated with CXCL12 (30 nM, purple squares), isoprenaline (iso, 100  $\mu$ M, light purple triangles) or both ligands simultaneously (orange) (B).

The GPCR Heteromer Investigation Technology (GPCR-HIT, Dimerix) reports on GPCR heteromer formation by the recruitment of an interacting protein specifically to the heteromer complex. Cells expressing CXCR4/Nluc were transiently transfected with cDNA coding for  $\beta$ -adrenoceptor as well as the interacting protein  $\beta$ -arrestin2/Venus. Treating the cells with CXCL12 resulted in the expected recruitment of  $\beta$ -arrestin2/Venus to genome-edited CXCR4/Nluc, as did treatment with the  $\beta$ -adrenoceptor agonist isoprenaline, indicating the close proximity of  $\beta$ -adrenoceptor to CXCR4/Nluc. Applying both agonists resulted in a greater than additive BRET signal, again suggestive of heteromer formation.

## Conclusion

The novel CRISPR/Cas9 technique successfully fused the Nluc BRET donor to endogenously-expressed CXCR4. Luminescence generated by the resulting protein-luciferase fusion was sufficient to monitor receptor-protein interactions as well as trafficking. The multiplex internalization assay depends on two acceptor fluorophores whose selective detection was rendered possible by the CLARIOstar's monochromator.

## References

- White CW et al. (2017) *Sci Rep.* 7:3187. doi: 10.1038/s41598-017-03486-2