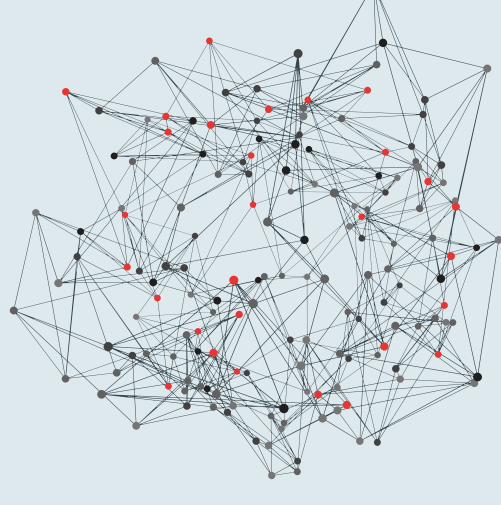


# Interrogating Protein Interactions with the NanoBRET™ Assay

The number of protein interactions in human cells has been estimated to be upwards of **650,000** (Stumpf *et al.*, 2008<sup>1</sup>). Some of the processes in which protein interactions are involved include:

- Normal cellular processes
- Cellular response to stress
- Pathological Processes & Cancer
- \_\_\_\_ (fill in the blank with the interaction you're studying)

## In your protein interactions studies, perhaps you've asked questions such as:

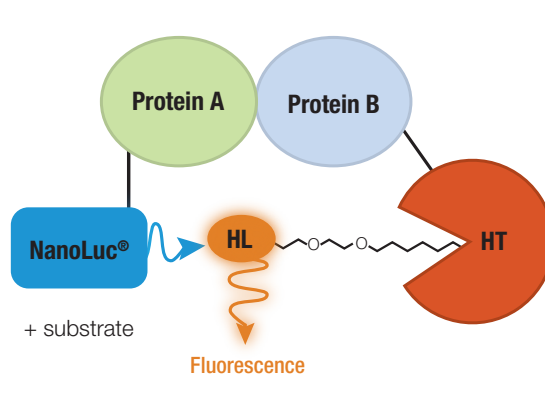


- Can I study my PPI within its natural biological context?
- Can I monitor PPI dynamics in real-time?
- Could inhibition of my PPI be important for disease treatments?
- Will my inhibitor compound be effective on full-length proteins in natural cellular context?
- How do mutations impact my PPI?

**NanoBRET™ PPI Assays can help you answer these questions.**

The new BRET-based protein:protein interaction (PPI) assay uses NanoLuc® Luciferase as the BRET energy donor and HaloTag® protein, labeled with the NanoBRET™ 618 fluorophore, as the energy acceptor to measure the interaction of specific protein pairs.

**The improvements to this BRET system make it easier and more reliable.**

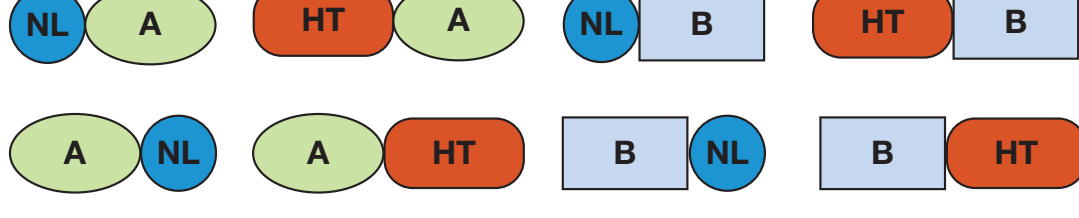


## How do you design a NanoBRET™ Assay to monitor your protein interactions?

1

### Generate Fusion Proteins

Prepare up to 8 possible clones containing combinations of protein A or protein B fused to NanoLuc (NL) and HaloTag (HT) at the amino (N) or carboxy (C) terminus of the protein.



*Did you know?*

### Find My Gene™ makes this easy:

- Search a library of nearly 10,000 constructs to find genes encoding your protein of interest
- Receive your ORFs cloned into a Flexi® Vector that already has the HaloTag® protein appended to the N-terminus of the fusion protein
- Transfer an ORF to the remaining HaloTag® (C-terminal) and NanoLuc® (N- and C-terminal) fusion vectors using Flexi® cloning

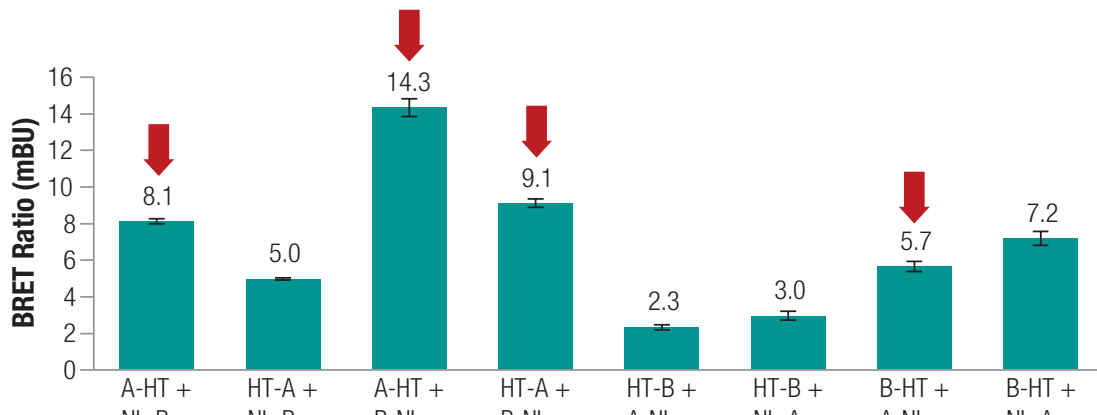
FindMyGene provides a rapid, efficient, high-fidelity method to create fusions, so that there is no need to resequence. For more information, [click here »](#)

Alternatively, NanoBRET™ Starter Systems are available, and provide the vectors required to create NanoLuc® Luciferase and HaloTag® protein fusions to target proteins of interest. For more information, [click here »](#)

2

### Test the Combinations

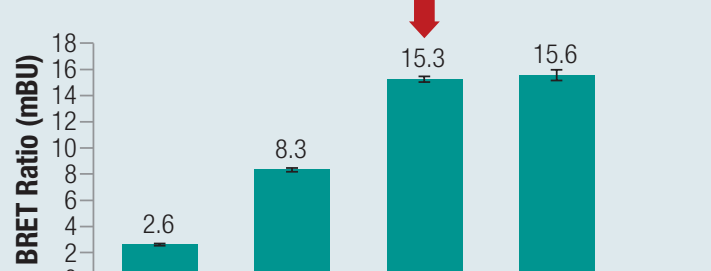
Using your clones, test these combinations to find the best energy transfer.



3

### Optimize Transfection

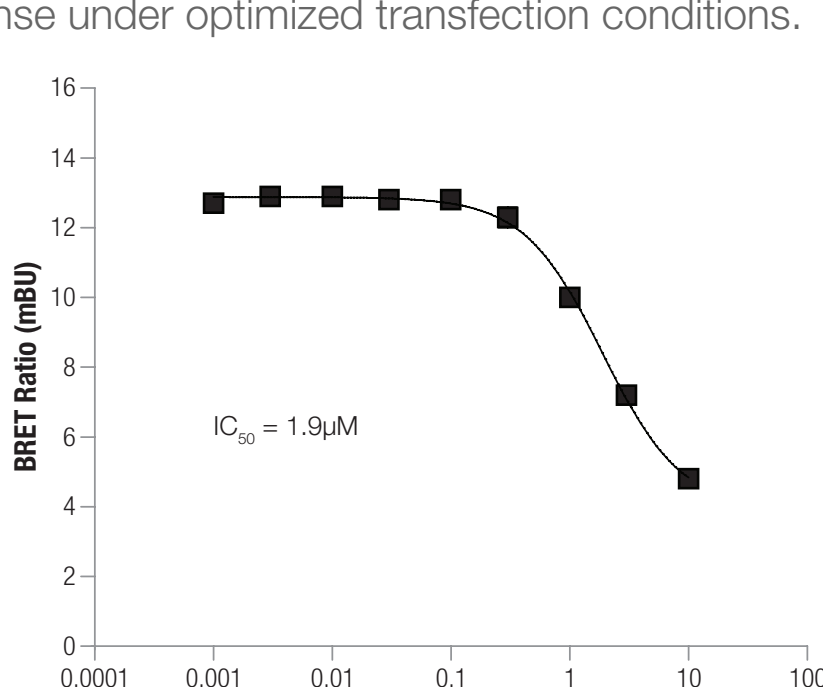
Find optimal donor to acceptor DNA ratio to minimize unbound donor and maximize dynamic range.



4

### Validate

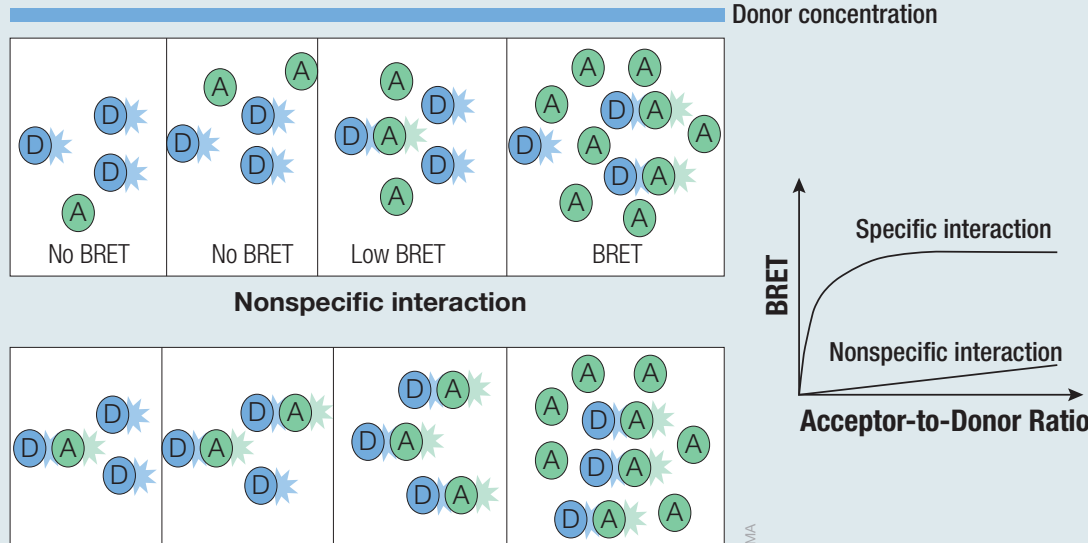
Use a known inhibitor or activator to test for a specific response under optimized transfection conditions.



*Alternatively*

### Test Specificity

using a donor saturation assay.



For more information and other tools for your PPI studies visit:

[www.promega.com/Try-NanoBRET](http://www.promega.com/Try-NanoBRET)

<sup>1</sup>Stumpf M.P.H. et al., (2008) Estimating the size of the human interactome. PNAS 105 (19): 6959-6964.