

A Versatile Platform for Cell Therapy Cytotoxicity Assays Using mRNA Transfection

Brock Binkowski, Julia Gilden, Nicholas Hess, Aileen Paguio, Mei Cong & Marjeta Urh

Promega Corporation, 2800 Woods Hollow Rd, Madison WI, 53711

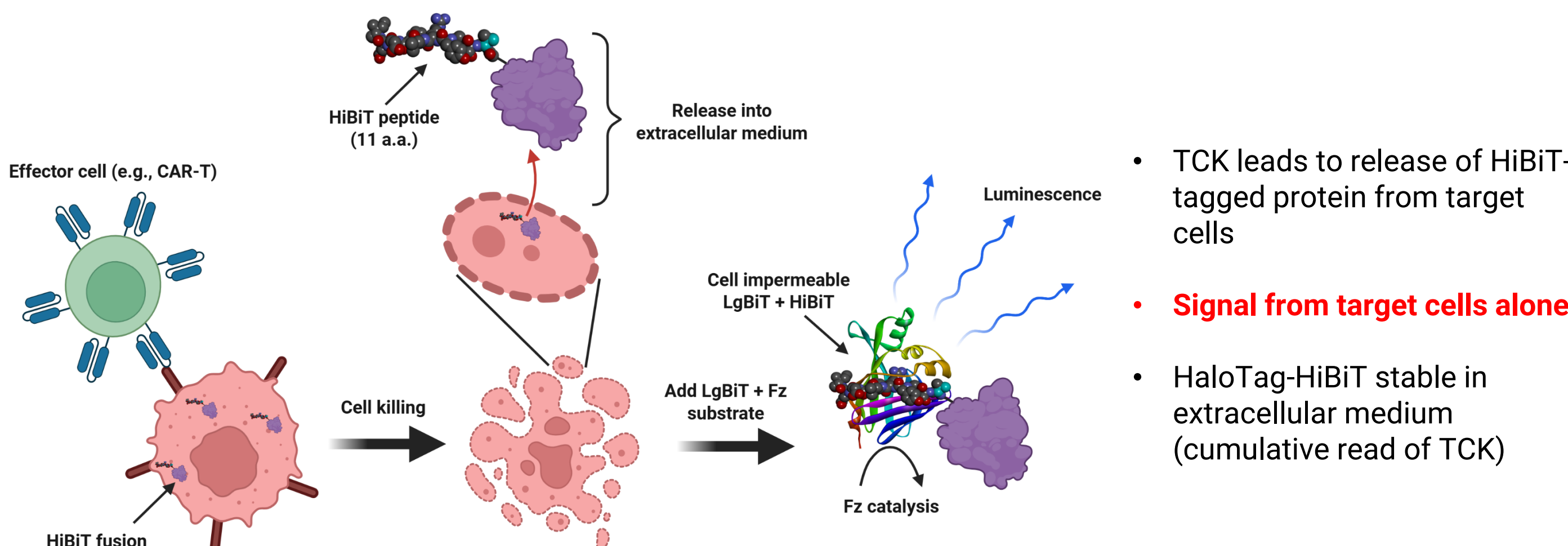


1. Introduction

The HiBiT Target Cell Killing Bioassay (HiBiT TCK Bioassay) is a sensitive and selective method for quantifying target cell lysis in coculture experiments with effector cells. In this assay, target cells engineered to express a HiBiT-tagged protein are cocultured with effector cells, where target cell death promotes release of the HiBiT-tagged protein into the extracellular medium. Once there, HiBiT can bind to cell-impermeable LgBiT to generate a bright, luminescent enzyme in the presence of furimazine substrate. In this manner, a non-lytic detection reagent containing LgBiT and furimazine substrate can be used to quantify target cell death in a gain-of-signal assay format with a cumulative readout. To date, this approach has been used exclusively with stable expression of the HiBiT-tagged protein, requiring time and effort to make stable clones.

To expand beyond the use of stable clones, we developed a novel mRNA transfection reagent that enables rapid and uniform expression of HaloTag-HiBiT in a wide range of target cell types. This approach dramatically reduces the workload for creating HiBiT target cells, allowing users to readily label target cell type(s) of interest. We demonstrate labeling of a variety of target cell types and their use in the HiBiT TCK Bioassay. Additionally, we demonstrate the use of mRNA transfection to achieve uniform and highly tunable expression of target receptors in cells stably expressing a HiBiT-tagged protein, allowing drug potency/efficacy testing across a broad range of receptor expression levels in a single cell type.

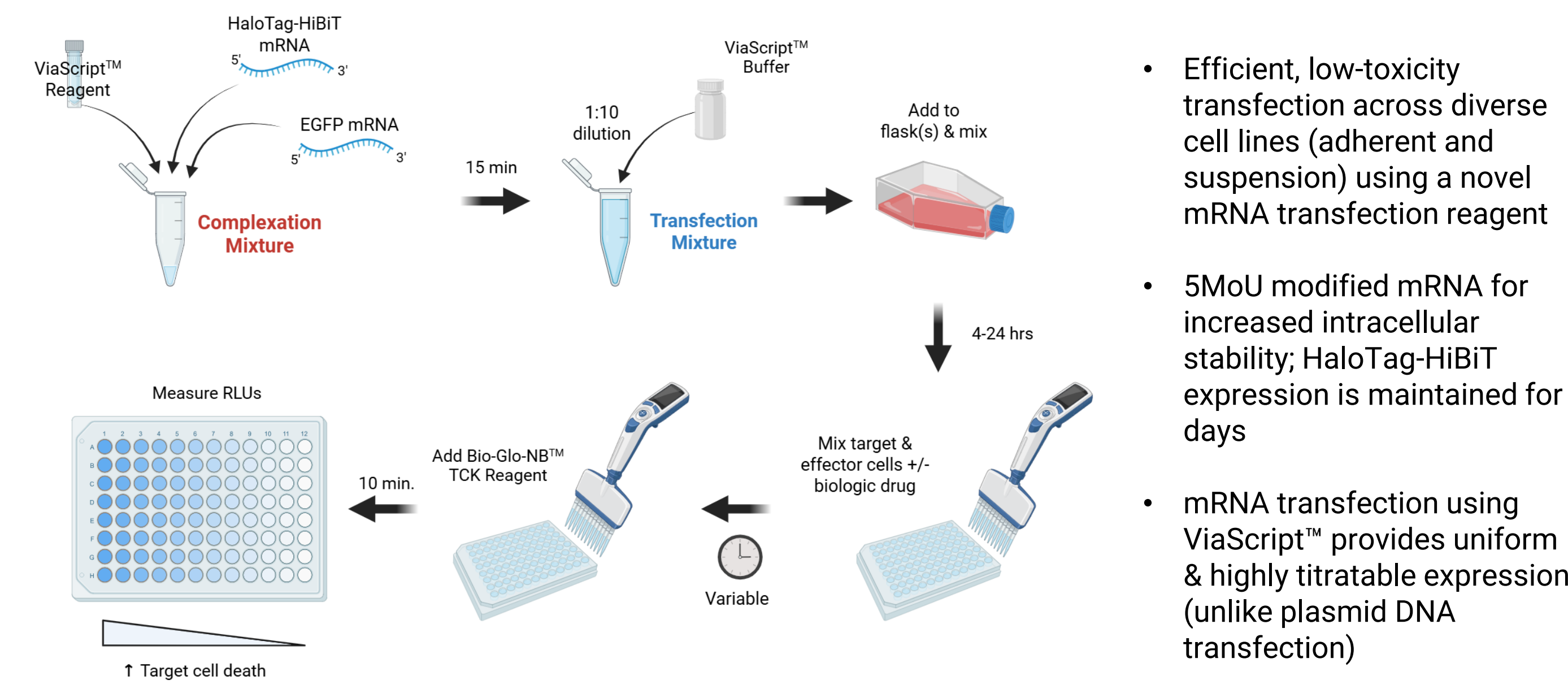
2. HiBiT Target Cell Killing (TCK) Bioassay



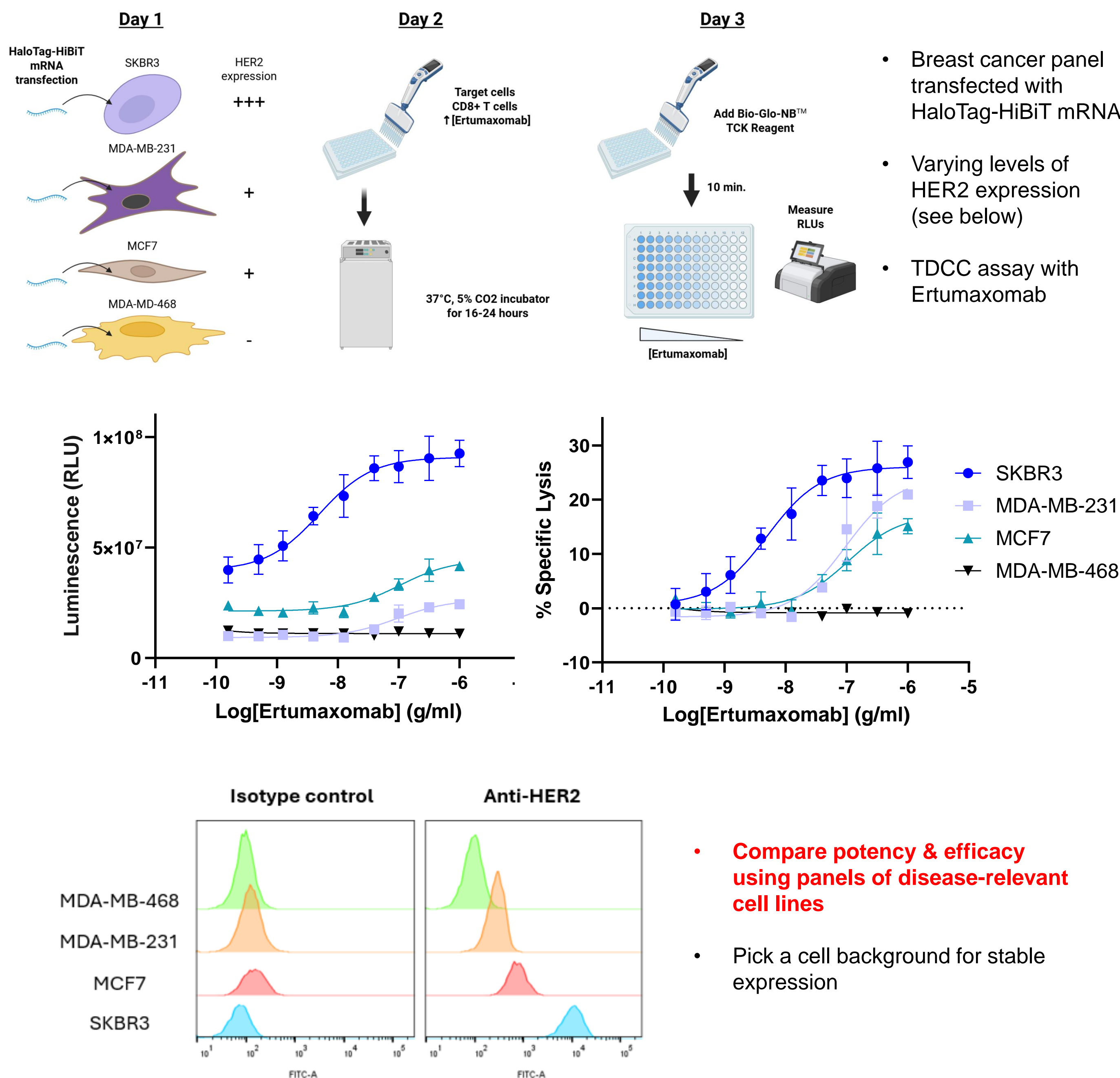
For a complete listing of available target cell lines, including target receptor KO's, visit:
<https://www.promega.com/products/reporter-bioassays/target-cell-killing-bioassays/>

3. ViaScript™ TCK Bioassay

Extending the HiBiT TCK Bioassay beyond use of stable clones

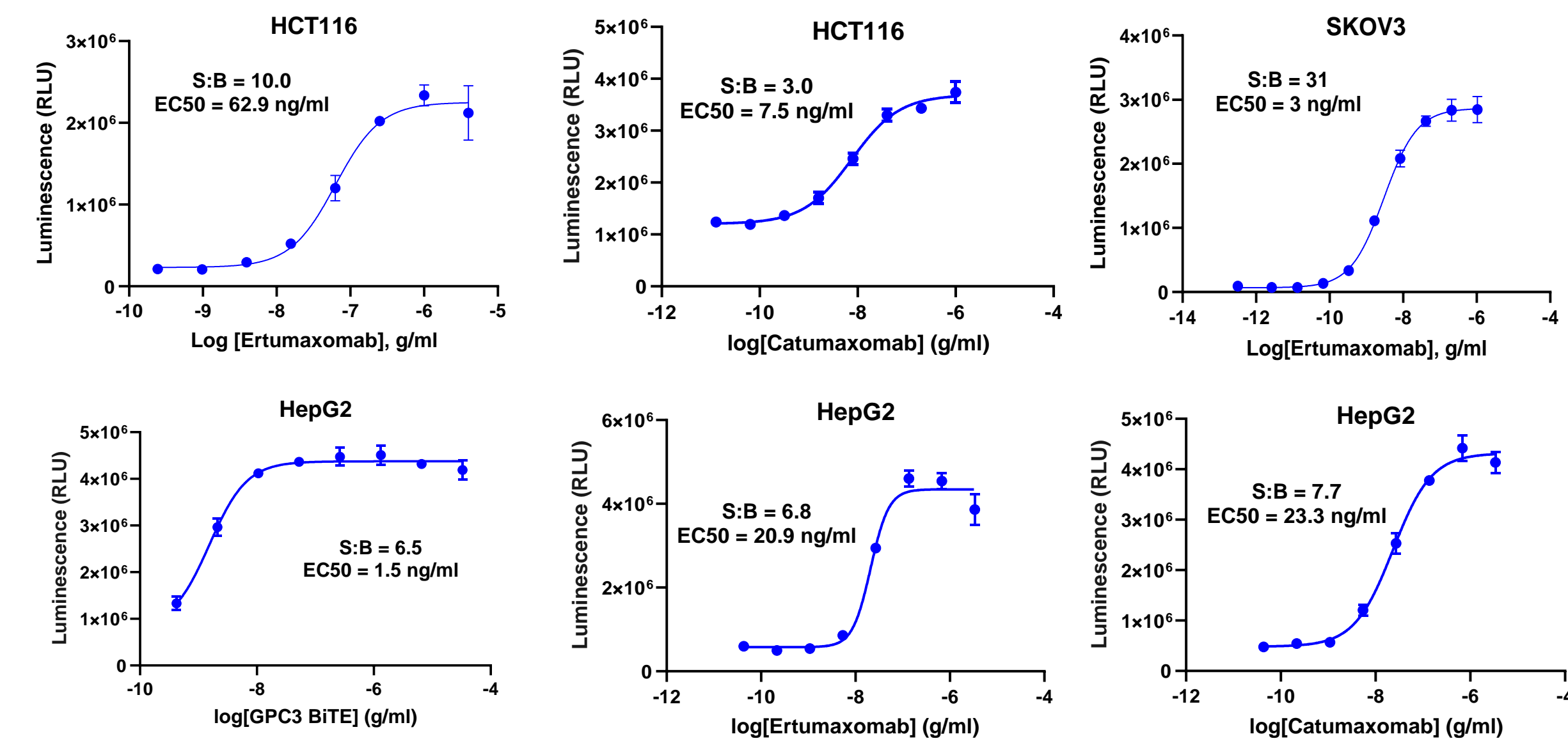


4. Easily Screen Panels of Target Cells

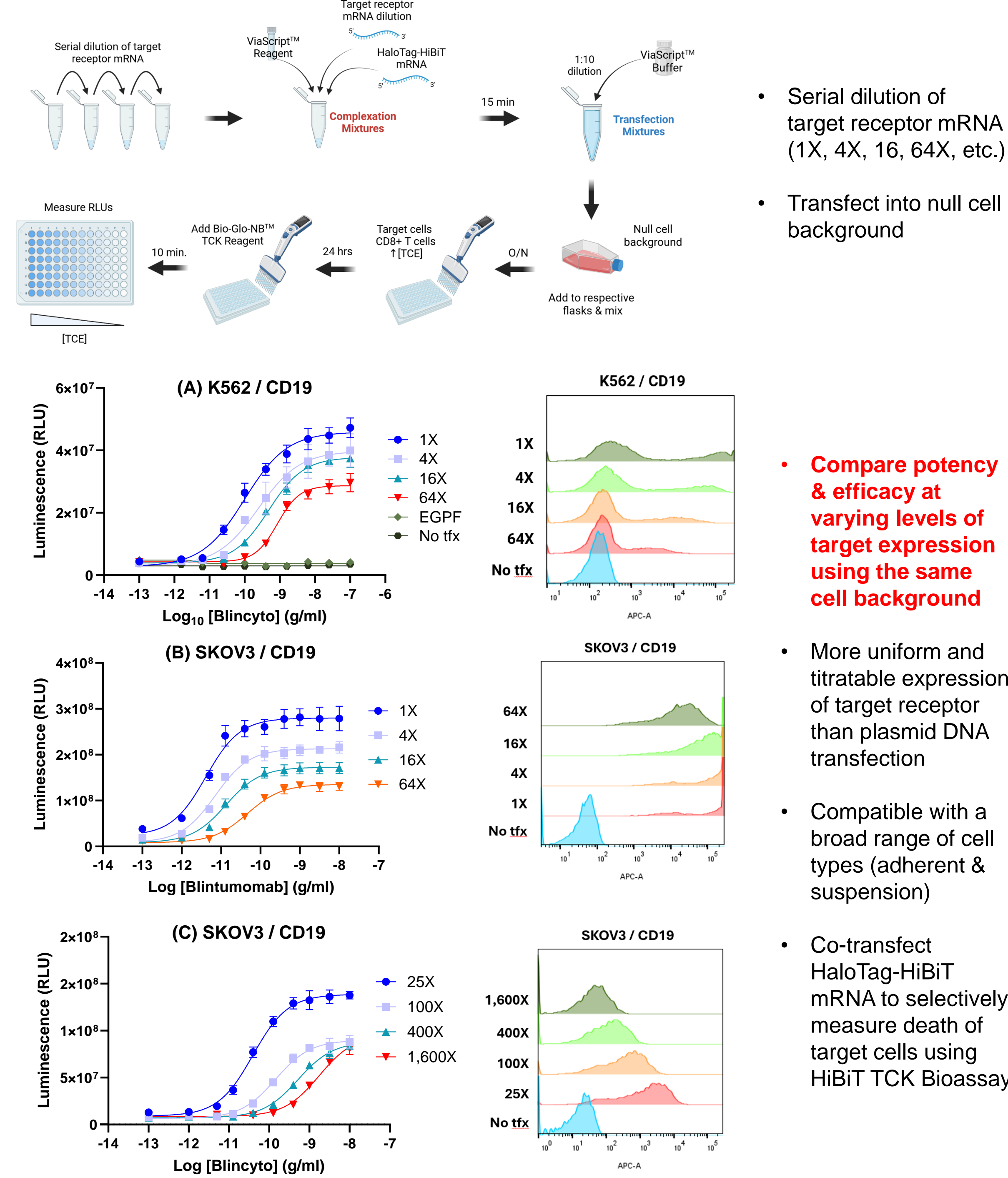


ViaScript TCK Approach Applied to a Breast Cancer Panel. Day 1, transfection of HaloTag-HiBiT mRNA using ViaScript™ Reagent. Day 2, mixing of CD8+ T cells with target cells (10:1 E:T ratio) in the presence of varying concentrations of Ertumaxomab (HER2 X CD3). Day 3, addition of Bio-Glo-NB™ TCK Reagent followed by luminescence measurement.

6. Further POC with Adherent Cells



7. Titrate Target Receptor Expression in the Same Cell Background



Modulating CD19 Expression in K562 & SKOV3. Day 1, CD19 mRNA transfection using ViaScript™ of a K562 (HaloTag-HiBiT) stable cell line or parental SKOV3 (also transfected with HaloTag-HiBiT). Day 2, mixing of CD8+ T cells with target cells (10:1 E:T ratio) in the presence of varying concentrations of Blinicyto (CD19 X CD3). Day 3, addition of Bio-Glo-NB™ TCK Reagent followed by luminescence measurement. Panels A, B & C are independent experiments.

8. Summary

- HiBiT TCK Bioassay:** Release of HiBiT-tagged protein from dead/dying cells leads to signal increase; no signal contribution from effector cells
- ViaScript™ TCK Bioassay:**
 - Transfection of HaloTag-HiBiT mRNA using a novel reagent
 - Easily apply to panels of disease-relevant cell lines
- Modulate target receptor expression in the same cell background:**
 - Highly titratable (far better than use of plasmid DNA)
 - Determine drug product potency & efficacy as a function of target receptor molecules per cell
 - HiBiT TCK Bioassay readout when co-transfecting HaloTag-HiBiT mRNA
- Available for purchase via Promega's Early Access Program. Contact nicholas.hess@promega.com.

Corresponding author: brock.binkowski@promega.com
For Research Use Only. Not for use in diagnostic procedures.