

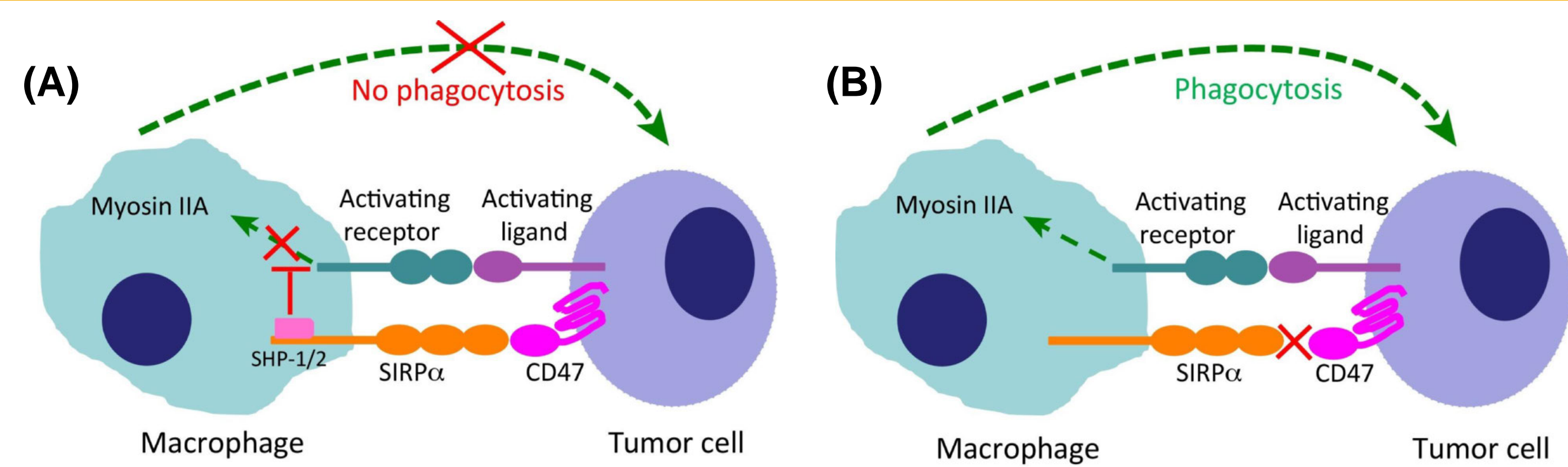
Cell-based Reporter Bioassays for Development of Fc-functional and Fc-silent SIRP α /CD47 Checkpoint Inhibitors

Jonathan Mitchell, Jamison Grailer, Jim Hartnett, Frank Fan, Mei Cong, Zhi-jie Jey Cheng
 Promega Corporation, 2800 Woods Hollow Rd, Madison, WI 53711

Abstract #3163



1. Introduction



Agent	Structure	Binds to		
		CD47	SIRP α	FcR
Anti-CD47 antibody - intact		+	-	+
Anti-CD47 antibody - F(ab) ₂		+	-	-
Anti-CD47 antibody - F(ab)		+	-	-
Anti-SIRP α antibody - intact		-	+	+
Anti-SIRP α antibody - F(ab) ₂		-	+	-
Anti-SIRP α antibody - F(ab)		-	+	-
Soluble SIRP α -Fc		+	-	+
Monomeric SIRP α		+	-	-
Anti-CD47 nanobody		+	-	-
Anti-CD47 single chain variable fragment		+	-	-

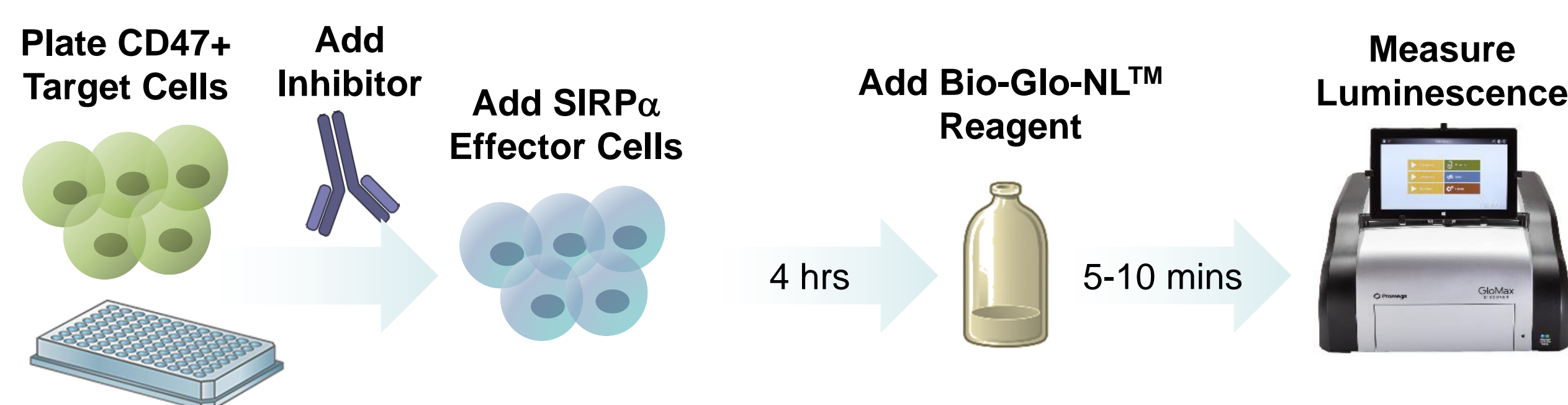
SIRP α /CD47 Immune Checkpoint.

(A) CD47 overexpressed on tumor cells engages the myeloid receptor SIRP α and delivers a “don’t eat me” signal that inhibits phagocytosis. (B) SIRP α /CD47 blockade promotes phagocytosis of tumor cells driven by pro-phagocytic ligands and/or Fc γ R engagement by Fc-functional antibodies (ADCP). (C) SIRP α /CD47 inhibitors vary in molecular structure and mechanism-of-action but can be broadly categorized as Fc-silent vs. Fc-functional.

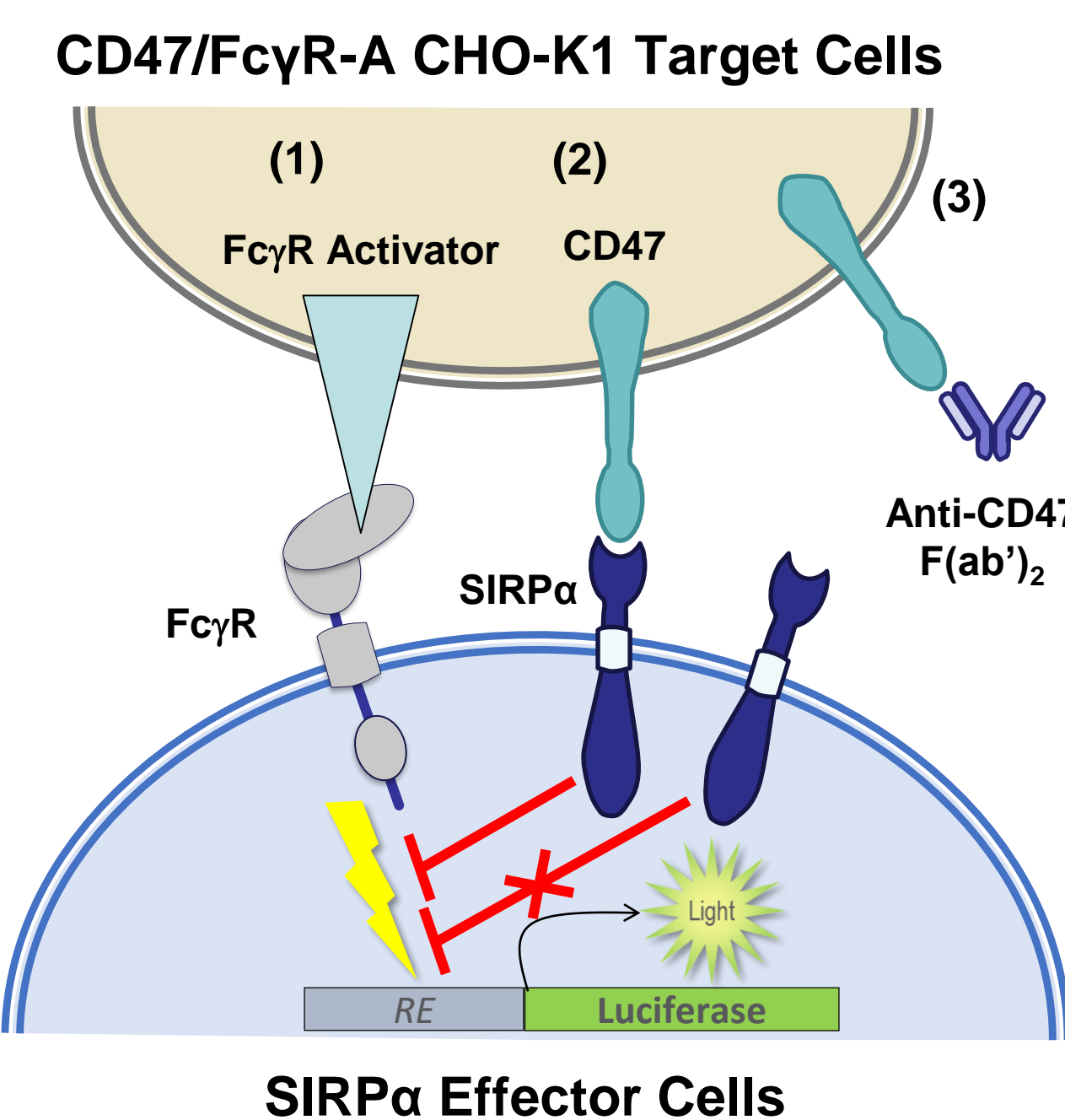
Figures from Veillette & Chen. 2018. Trends Immunol.

2. SIRP α /CD47 Blockade Bioassays: Design & Workflow

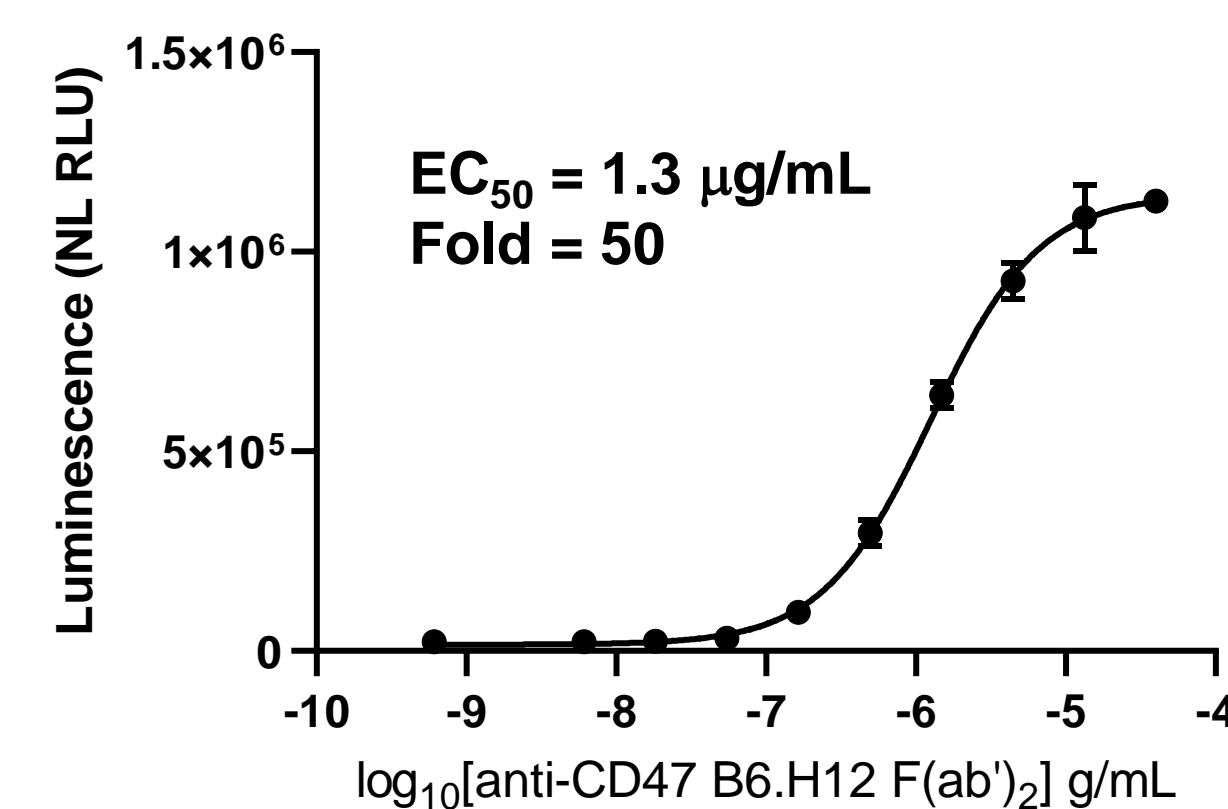
- We have developed a pair of reporter-based bioassays for measuring the activity of Fc-silent or Fc-functional SIRP α /CD47 inhibitors
- These SIRP α /CD47 Blockade Bioassays follow a simple, add-mix-read format



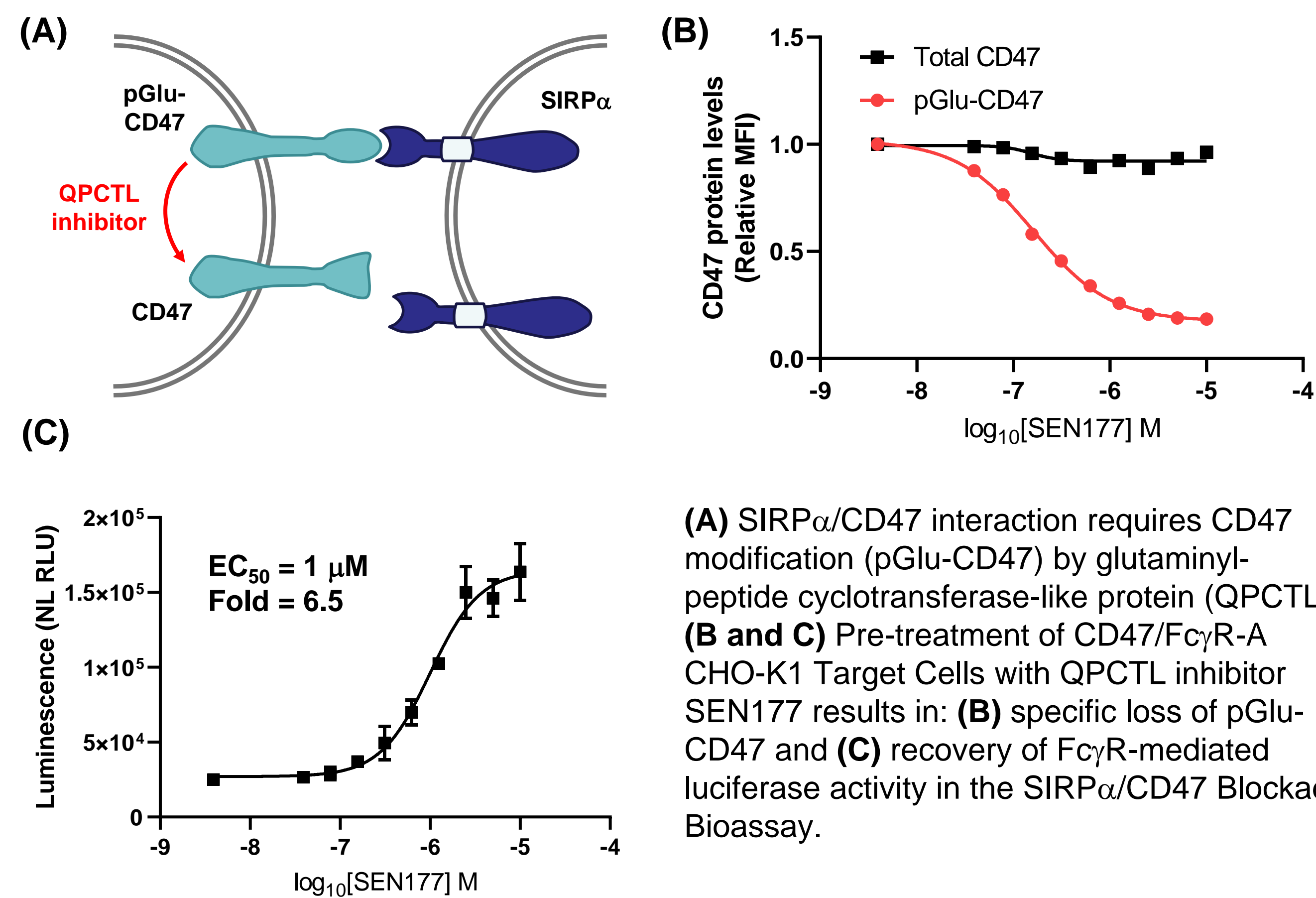
3. SIRP α /CD47 Blockade Bioassay Measures the Potency of Fc-silent SIRP α /CD47 Blocking Antibody



Luciferase activity indicating monocyte activation is:
 (1) induced by Fc γ R-A protein that engages Fc γ Rs
 (2) inhibited by co-engagement of SIRP α with CD47
 (3) restored by CD47 blockade using anti-CD47 F(ab)₂ fragment of blocking clone B6.H12



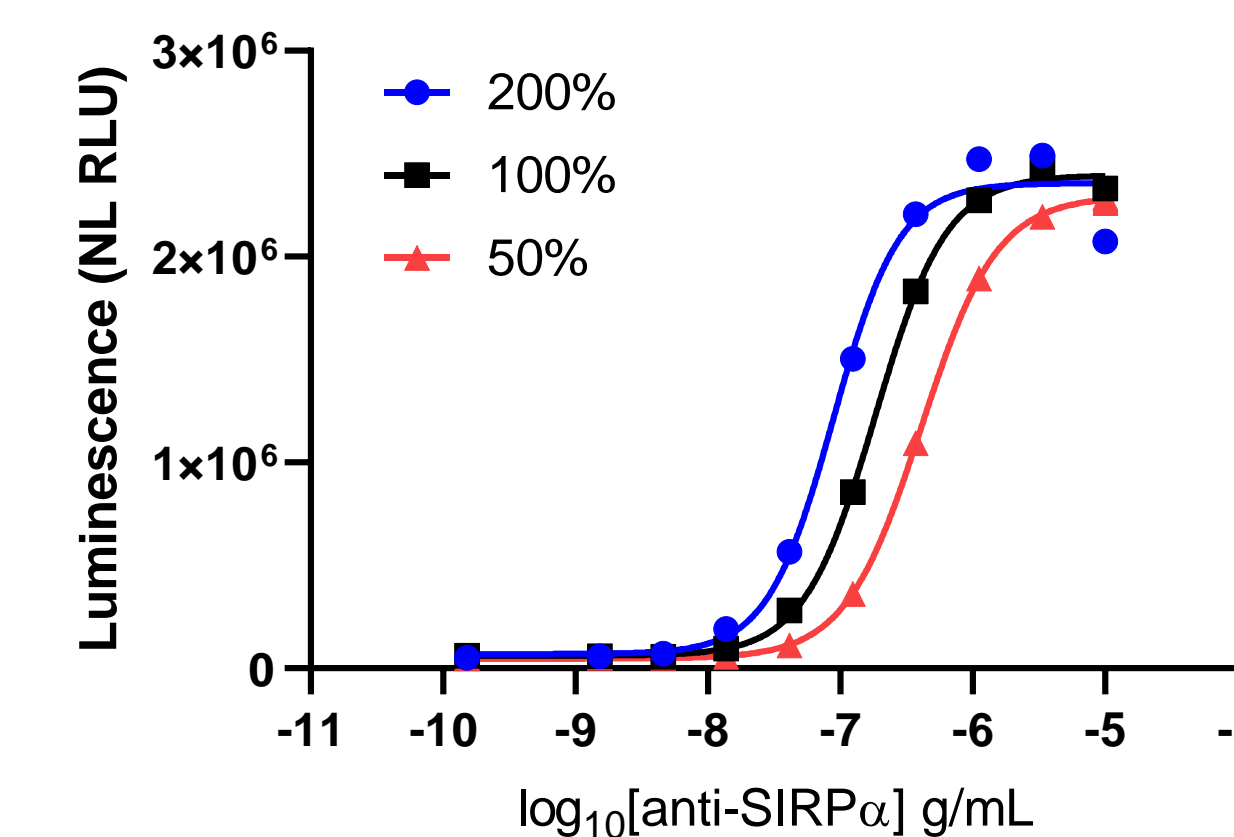
4. SIRP α /CD47 Blockade Bioassay Measures the Activity of Small Molecule SIRP α /CD47 Inhibitors



(A) SIRP α /CD47 interaction requires CD47 modification (pGlu-CD47) by glutaminyl-peptide cyclotransferase-like protein (QPCTL). (B and C) Pre-treatment of CD47/Fc γ R-A CHO-K1 Target Cells with QPCTL inhibitor SEN177 results in: (B) specific loss of pGlu-CD47 and (C) recovery of Fc γ R-mediated luciferase activity in the SIRP α /CD47 Blockade Bioassay.

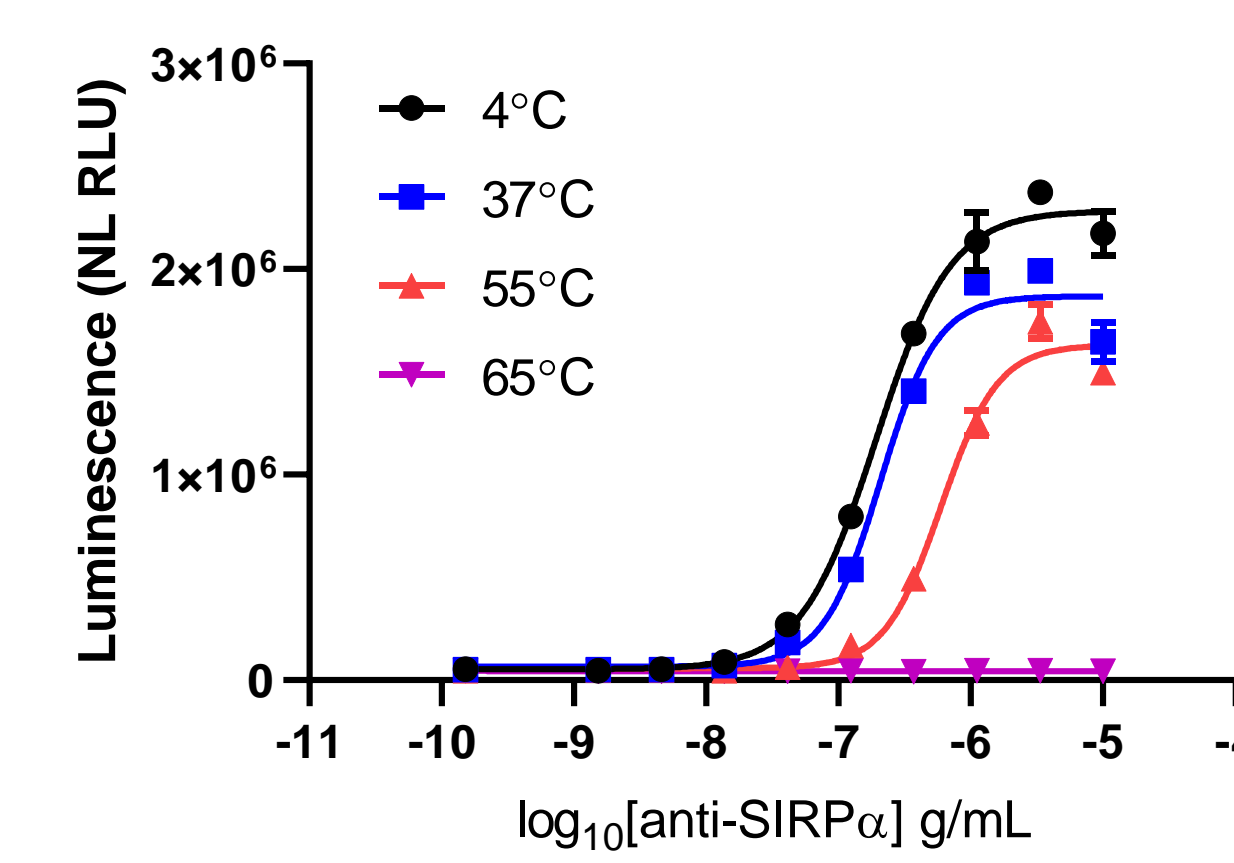
5. SIRP α /CD47 Blockade Bioassay can Measure Relative Potency and is Stability Indicating

(A) Relative potency of anti-SIRP α blocking Ab



Expected Relative Potency	Measured Relative Potency
50%	45.1%
200%	208%

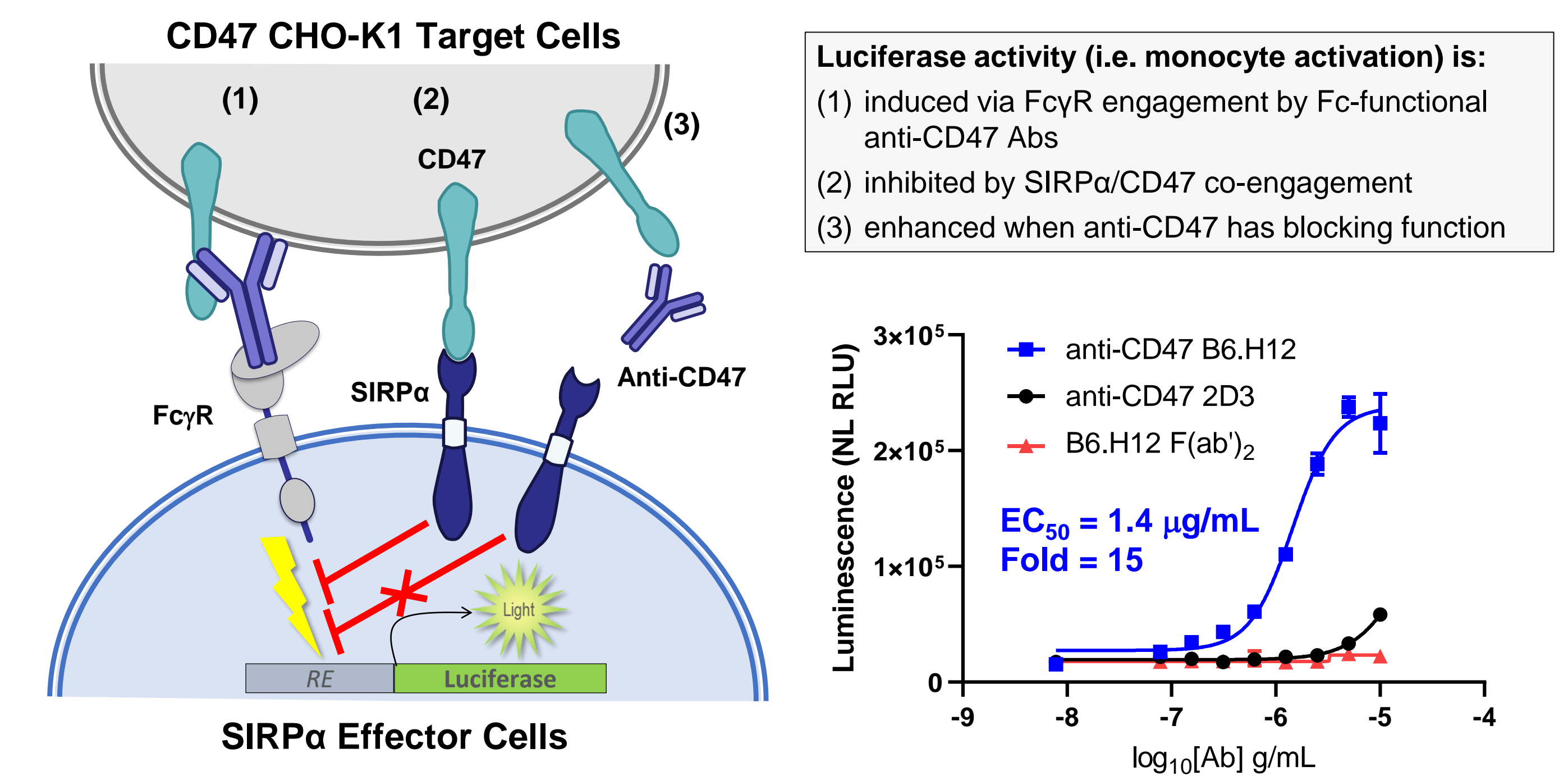
(B) Stability of anti-SIRP α blocking Ab



Temperature	EC ₅₀ (μg/mL)
4°C	0.19
37°C	0.21
55°C	0.59
65°C	No assay response

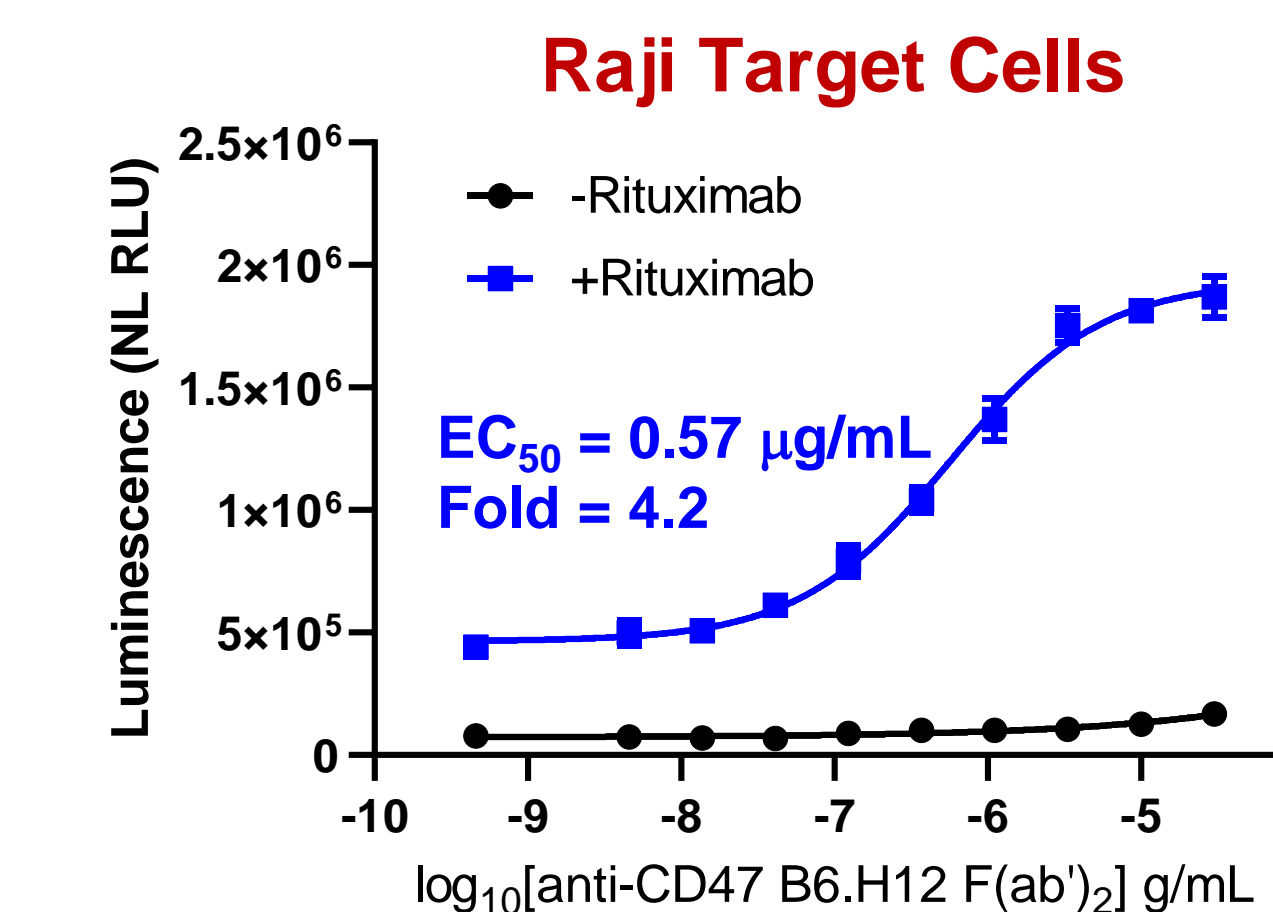
SIRP α /CD47 Blockade Bioassays were performed using anti-SIRP α blocking Ab (clone SE5A5): (A) Simulated potency series of anti-SIRP α Ab and measured relative potency compared to 100% reference. (B) Anti-SIRP α Ab was incubated at the indicated temperatures for 24 h prior to analysis in the SIRP α /CD47 Blockade Bioassay.

6. SIRP α /CD47 Blockade Bioassay, Fc-dependent Measures Potency of Fc-functional CD47 Blocking Abs



Fc γ R-mediated luciferase activity is observed in the SIRP α /CD47 Blockade Bioassay, Fc-dependent with full-length anti-CD47 blocking Ab (clone B6.H12, mouse IgG1 isotype), but not with B6.H12 F(ab)₂ fragment or a non-blocking anti-CD47 Ab (clone 2D3, mouse IgG1 isotype).

7. SIRP α /CD47 Blockade Bioassays Enable Testing of Drug Combinations using CD47+ Cancer Cells



SIRP α /CD47 Blockade Bioassay, Fc-dependent was performed using SIRP α Effector Cells and Raji target cells (human B-cell lymphoma, CD47+/CD20+). Anti-CD47 F(ab)₂ fragment was added at increasing concentrations in the presence or absence of anti-CD20 Ab (rituximab, EC₁₀₀). As expected, the anti-CD47 F(ab)₂ fragment enhanced rituximab-mediated luciferase activity. No response was observed with anti-CD47 F(ab)₂ fragment alone.

8. Conclusions

We have developed a pair of cell-based reporter gene assays for measuring biological activity of diverse SIRP α /CD47 inhibitors:

- SIRP α /CD47 Blockade Bioassay**
 - Suitable for Fc-silent CD47 blocking Abs, SIRP α blocking Abs, and small molecule inhibitors
- SIRP α /CD47 Blockade Bioassay, Fc-dependent**
 - Suitable for Fc-functional CD47 blocking Abs, drug combinations, and bispecific Abs

SIRP α /CD47 Blockade Bioassays offer a simple, high-throughput platform for drug development, lot release, and stability studies.

SIRP α /CD47 Blockade Bioassays can be performed using fresh cell cultures or thaw-and-use cells that eliminate the need for cell propagation.