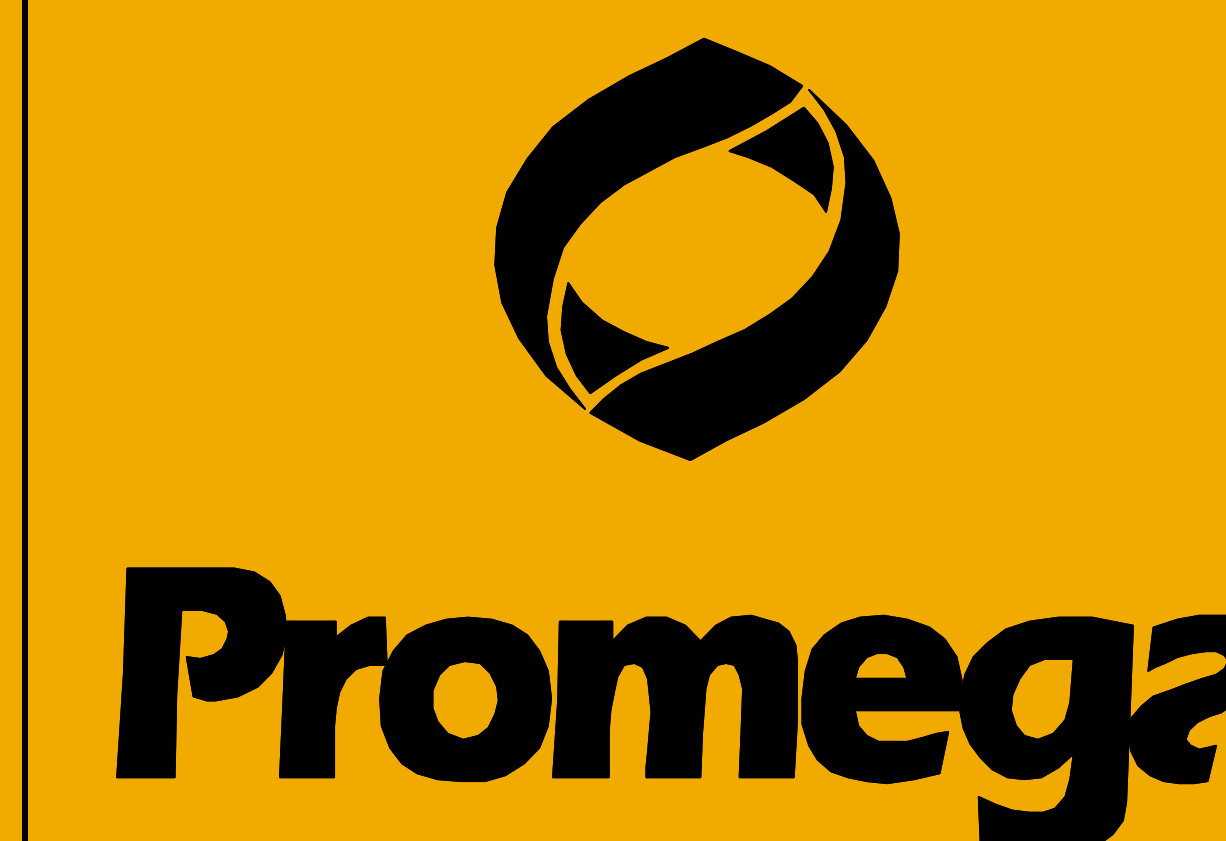


SRF-RE reporter assay for G₁₂-RhoA signaling pathway

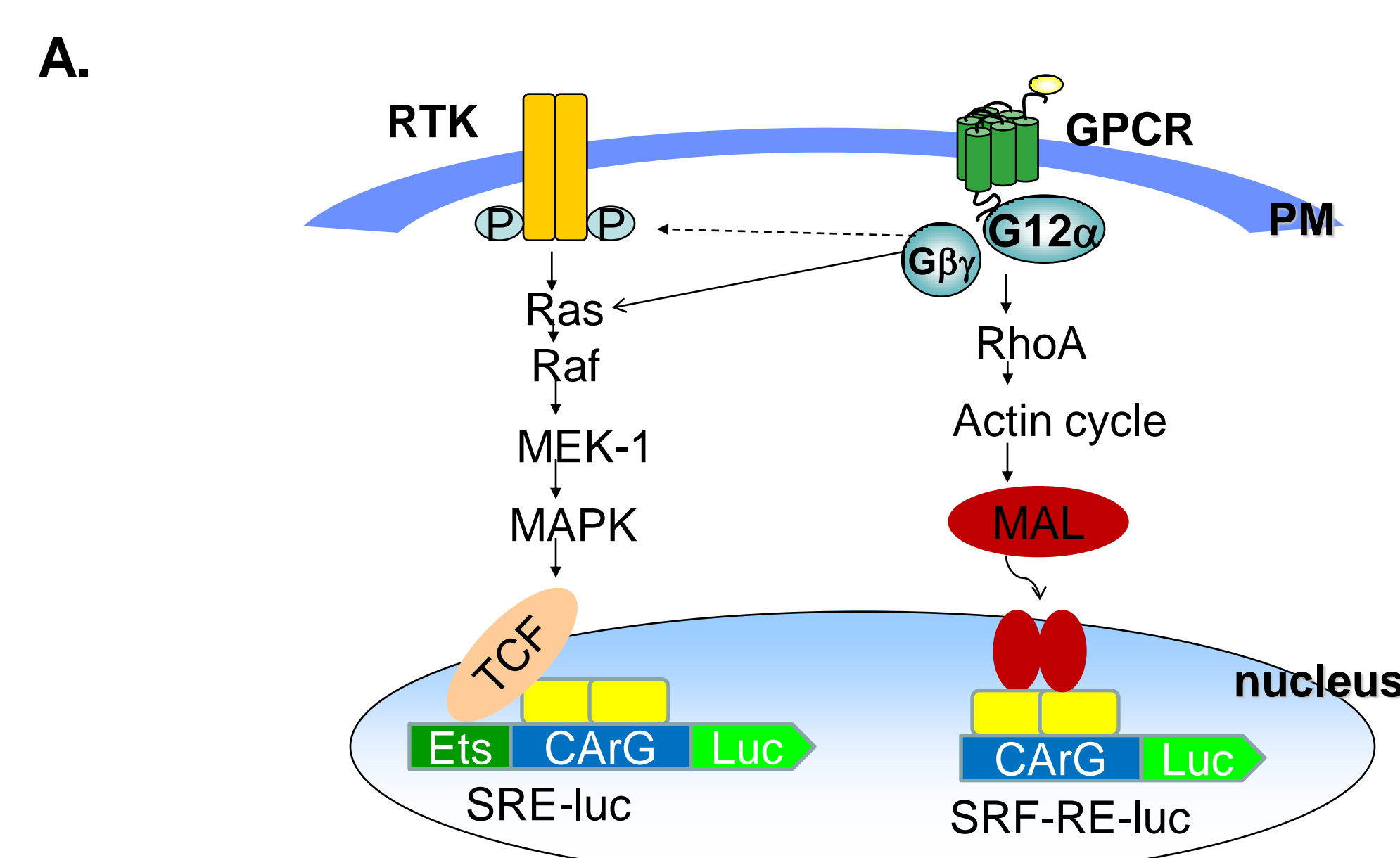
Jey Cheng, Denise Garvin, Pete Stecha, Keith Wood, and Frank Fan
Promega Corporation, 2800 Woods Hollow Road, Madison, WI 53711, USA



Abstract

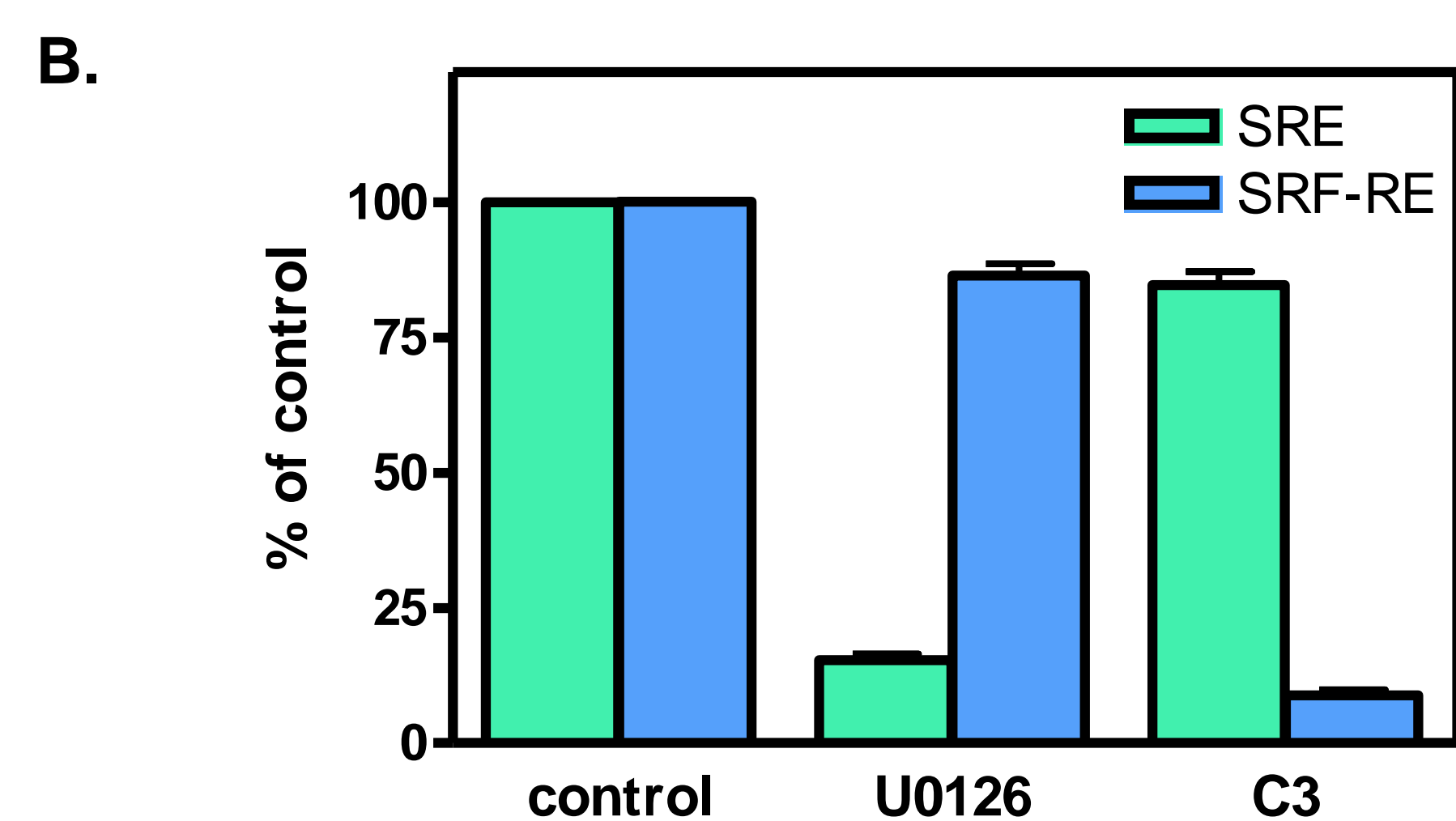
The G₁₂ subfamily of heterotrimeric G protein has been the subject of intense interest in both basic science and drug discovery research due to its involvement in regulating cell proliferation, cell migration and metastatic invasion, whereas not many drug-screening assays for G₁₂-linked pathway is available. Bioluminescent reporter assays are frequently used in high throughput screening due to their inherently high sensitivity, wide dynamic range and low susceptibility to compound interference. Here, We have incorporated the technical advancement of luciferase reporter genes and vectors to build an novel SRF-RE luciferase construct for RhoA signaling pathways. Large dynamic range of SRF-RE luciferase reporter assay allows efficient screening for GPCR modulators for G₁₂ pathways. The dual luciferase GPCR assay system, whereas the second plasmid expresses the target GPCR and Renilla luciferase as internal control, improves data quality by providing simultaneous monitoring of nonspecific effects such as cytotoxicity.

Dual Luciferase SRF-RE Reporter Assay for RhoA Signaling



SRF-RE-*luc2P* reporter responds to G_{12α}-linked RhoA signaling pathway.

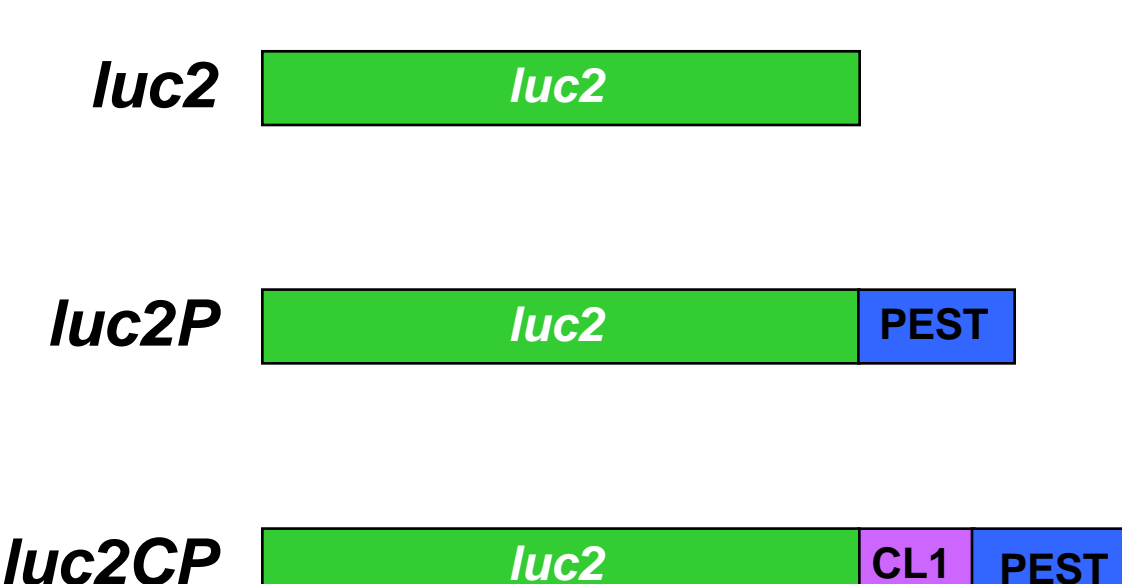
pf9A Vector expresses the GPCR of interest and Renilla-neo fusion control protein for normalization.



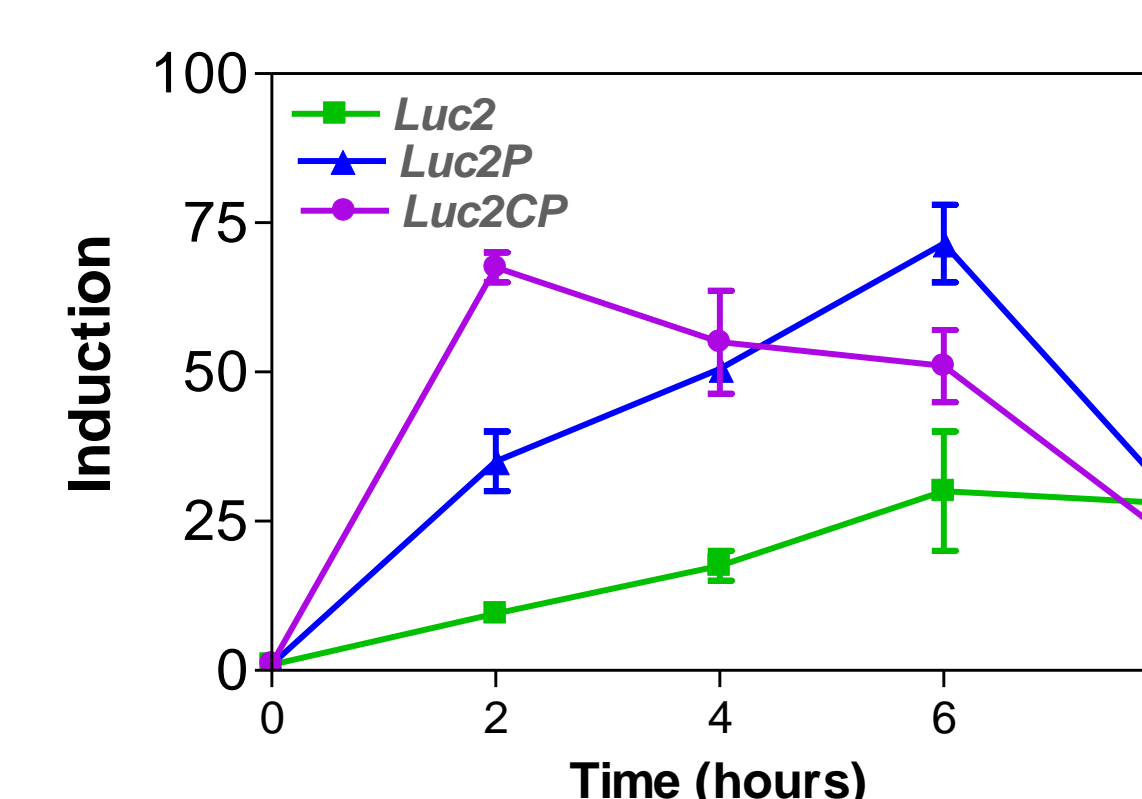
HEK293 cells were transiently transfected with SRE-*luc2P* or SRF-RE-*luc2P* together with Renilla luciferase in 96-well plates. Cells were pretreated with/without U0126 (MEK inhibitor) or C3 Transferase (RhoA inhibitor), then induced with 10ng/ml PMA for SRE or 20%FBS for SRF reporter assay. Firefly luciferase activity was quantified six hours after induction using the Dual-Glo™ Assay System.

Rapid Response™ Luciferase Enhances Reporter Dynamics

Rapid Response™ reporter genes

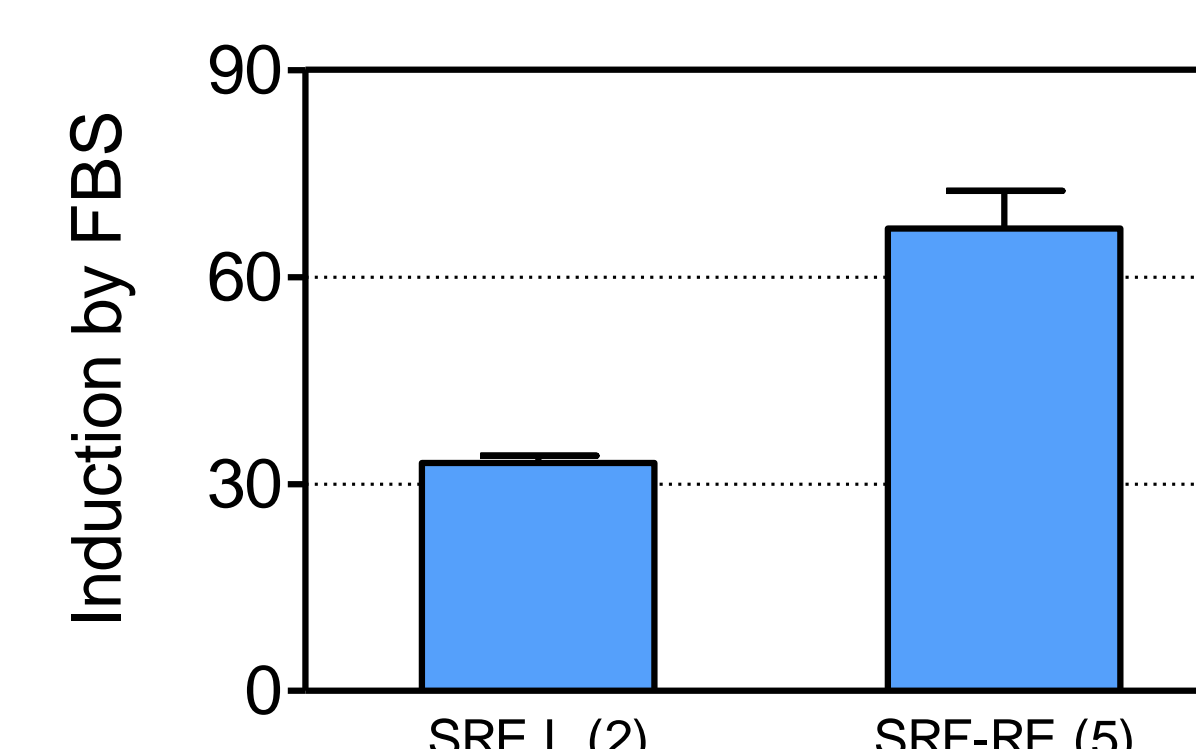
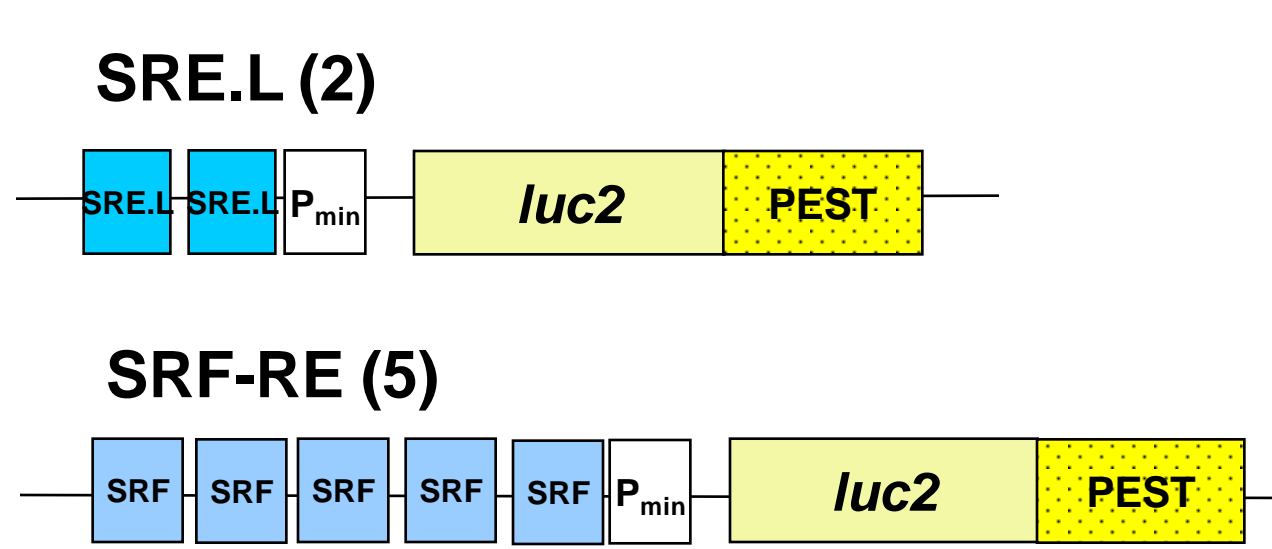


Induction of SRF-RE by FBS



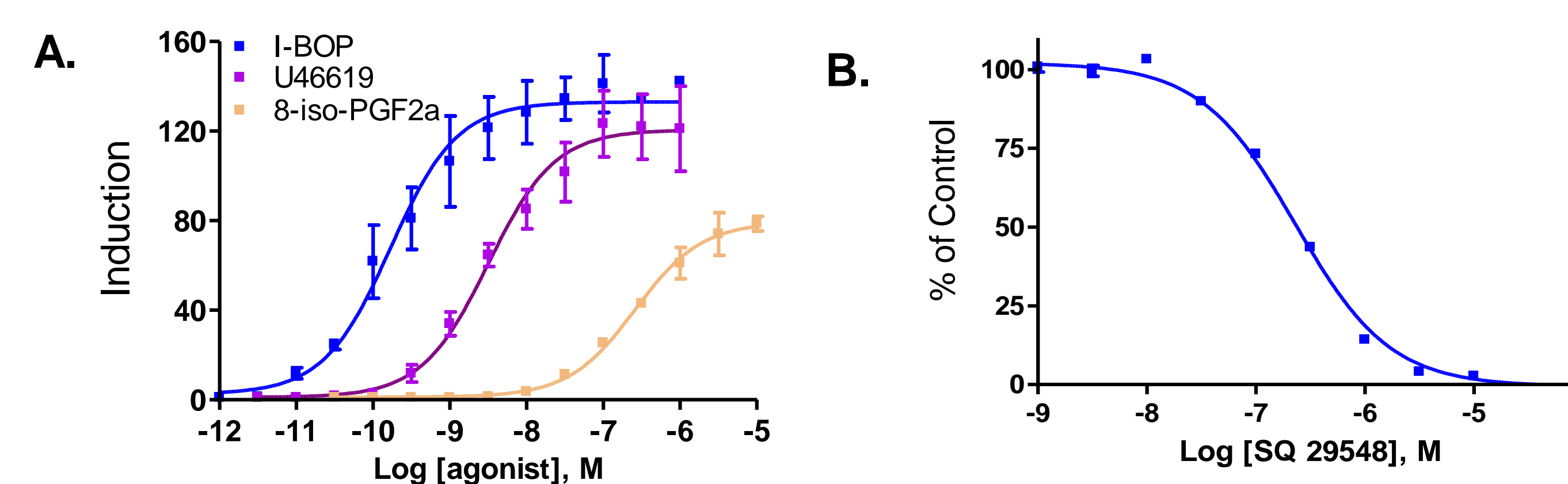
Versions of the *luc2* gene containing protein degradation sequences were used to enhance reporter dynamics. HEK293 cells transiently expressing SRF-RE-*luc2*, SRF-RE-*luc2P* or SRF-RE-*luc2CP* were induced with 20%FBS. Firefly luciferase activity was quantified every other hour for eight hours after induction using Luciferase Assay System.

Optimized Reporter Sequences/Repeats Further Improves Reporter Dynamics



HEK293 cells were transiently transfected with SRE.L(2)-*luc2P* or SRF(5)-RE-*luc2P* together with Renilla luciferase in 96-well plates. Cells were induced with 20%FBS. Firefly luciferase activity was quantified six hours after induction using the Dual-Glo™ Assay System.

Agonist and Antagonist Assay



HEK293 cells were transiently transfected with SRF-RE-*luc2P*, thromboxane A₂ receptor and Renilla luciferase, and serum starved overnight. Twenty four hours after transfection, firefly luciferase activity was measured after six hours induction with addition of 1:3 serial dilutions of agonists (A), or pretreatment with antagonists before induction with 0.1 μM U46619 (B) using the Dual-Glo™ Assay System in 96-well format. For agonist assay, fold Induction = induced Firefly RLU/uninduced Firefly RLU, normalized to *Renilla* RLU. For antagonist assay, % of Control = [Firefly RLU (antagonist+agonist)/Firefly RLU (agonist alone)] x100

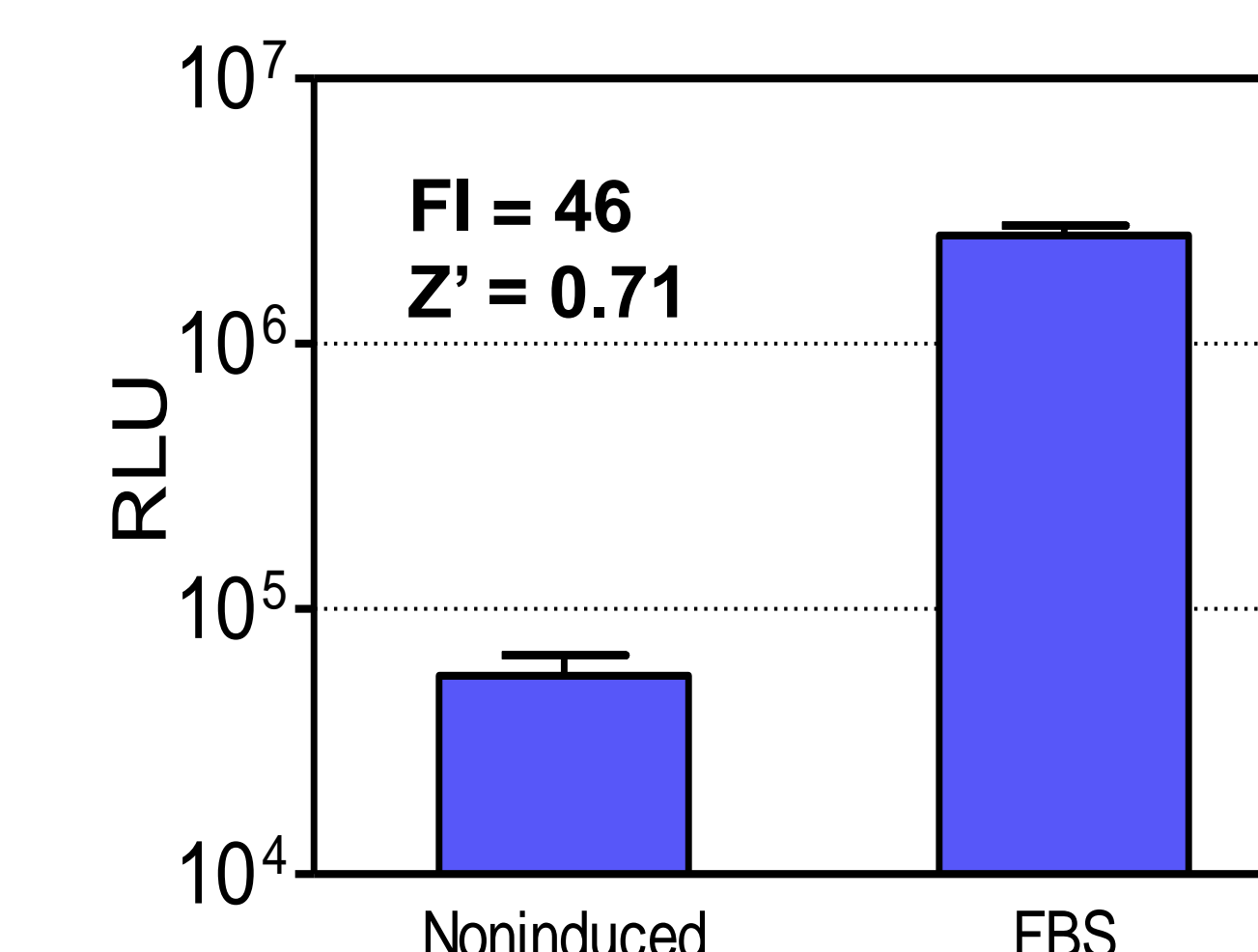
Summary of Fold change and EC₅₀ of Various GPCR/Agonist Pairs

GPCR	G protein subunits	agonist	Fold Induction	EC ₅₀ (M)
LPA receptor	G ₁₂ , G _q	LPA	27	3.0 × 10 ⁻⁷
Endothelin-B receptor	G ₁₂ , G _q	ET-1	11	8.0 × 10 ⁻¹⁰
Thromboxane A ₂ receptor	G ₁₂ , G _q	U46619	91	5.2 × 10 ⁻⁹

HEK293 cells transiently expressing SRF-RE-*luc2P*, Renilla luciferase and various endogenous (LPA receptor) or exogenous receptors (endothelin-B receptor and thromboxane A₂ receptor) were induced with agonists for six hours and analyzed in 96-well format. Fold Induction = induced Firefly RLU/uninduced Firefly RLU, normalized to *Renilla* RLU.

High Quality HTS assay

Day 1 Plate cells in T75 flask in complete medium
 Day 2 • Transfection in T75 flask
 • After 4 hrs, replate cells into 384 well plate, 20μl per well in serum-starving medium
 Day 3 • Add 5μl compounds (5) per well
 • 6hr induction
 • Add 25μl One-Glo reagent, wait 3-5 min
 • read



HEK293 cells transiently expressing SRF-RE-*luc2P* and Renilla luciferase were replated in serum-starving medium in 384-well plates. 24 hours after transfection, cells were induced with 20%FBS for six hours. Firefly luciferase activity was quantified using the Dual-Glo™ Assay System. Fold Induction = induced Firefly RLU/uninduced Firefly RLU, normalized to *Renilla* RLU. Z' was calculated as 1-[(3SD_{induced}+3SD_{uninduced})/(Ave_{induced}-Ave_{uninduced})]

Summary

Together the new SRF-RE luciferase reporter vector provide an excellent tool for G₁₂ pathway in research and drug discovery programs:

- ❖ Optimized SRF-RE sequence by deactivating the Ets binding domain of SRE allows specific detection of signaling via RhoA pathway.
- ❖ Use of Rapid Response™ *luc2P* gene improves reporter dynamics and reduces assay time.
- ❖ Optimized SRF-RE repeats further improves the dynamic response.
- ❖ Large dynamic range allows robust identification and ranking of GPCR modulators through G₁₂ pathways.
- ❖ Assay in HTS format shows good Z' value and dynamic range.