

What can you do with a luciferase Reporter Assay?

Signaling Pathway Analysis Applications

Presented Fall 2009



Click the icon in the upper left hand corner to view speaker notes for slides.

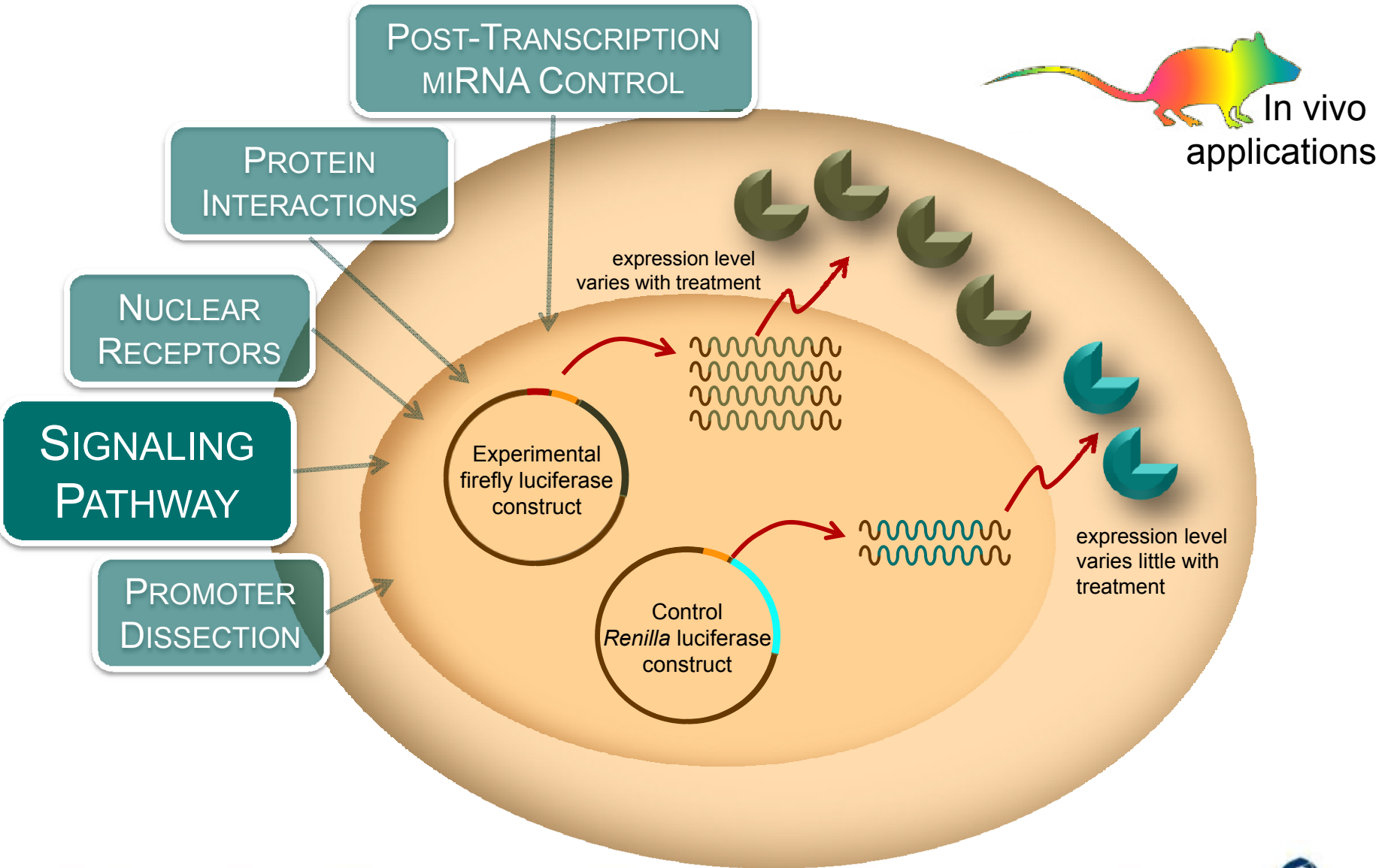


Have a question?
Ask a Scientist



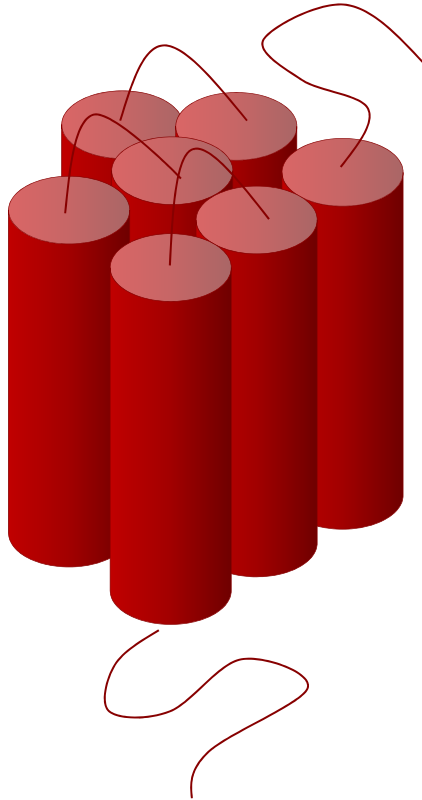
Apo-ONE, CellTiter-Fluor, and One-Glo are trademarks and GloMax is a registered trademark of Promega Corporation. HighWire Press is a registered trademark of the Board of Trustees of the Leland Stanford Junior University

Application Overview





What signaling pathway does my protein use?

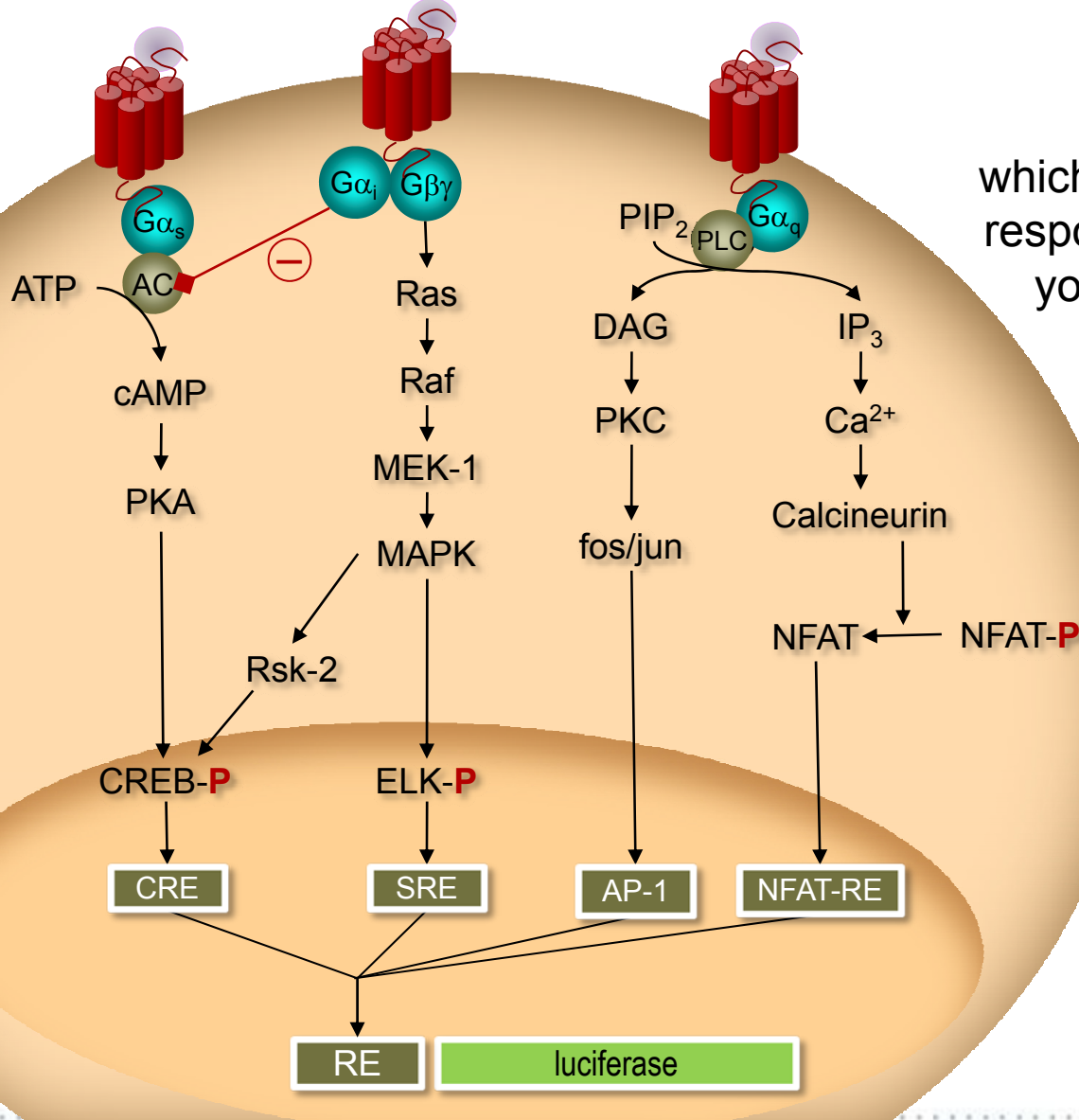


You've identified a novel GPCR receptor.

- Which pathway does it use for signal transduction?
- Which G protein does it utilize?

Synthetic promoter constructs can help you identify which pathways are involved

Investigate signal transduction pathways

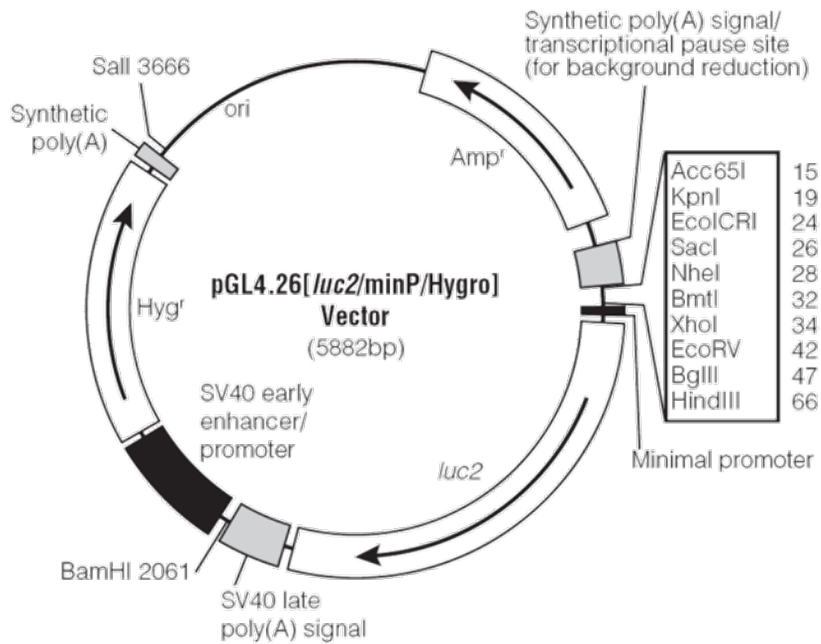


The reporter assay activation can tell you which signaling pathway is responding to activation of your receptor of interest

Now you can use inhibitors, siRNA, etc., to define which factors are involved



Easily create response element reporters

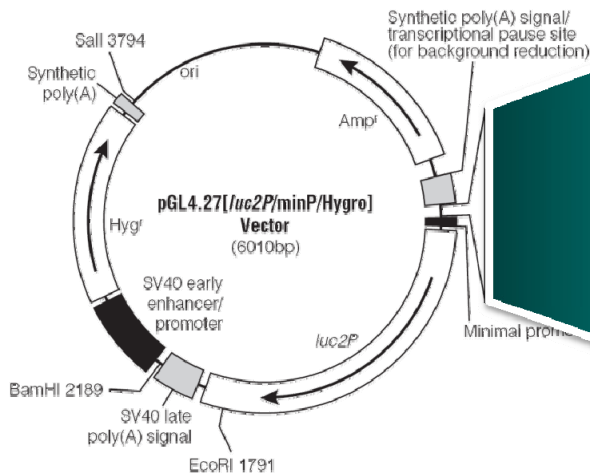


5951MA

- Response elements for many transcription factors are known
- Response elements can be placed upstream of a minimal promoter to create synthetic reporter constructs for monitoring signal transduction

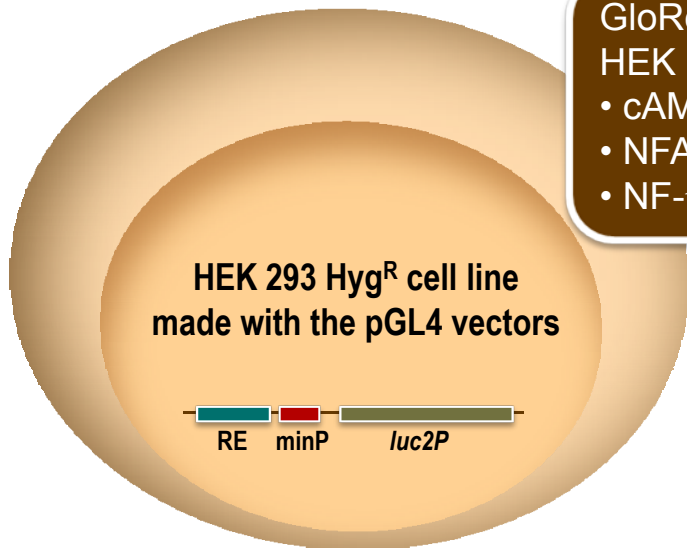
RapidResponse™ versions available

Tools to speed your research



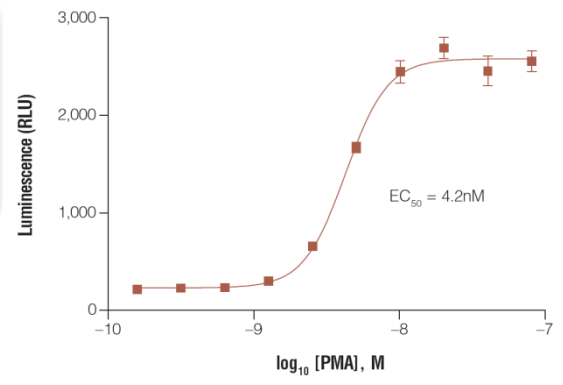
Pre-designed, ready-to-use response element pGL4 Vectors

Response Element	Signaling Pathway	pGL4 Vector
cAMP	cAMP/PKA	pGL4.29
NFAT Nuclear Factor of Activated T-cells	Ca ²⁺ /Calcineurin	pGL4.30
NF-κB	NF-κB	pGL4.32
Serum Response	MAP/ERK	pGL4.33
Serum Response Factor	RhoA	pGL4.34



GloResponse™ HEK 293 cell lines:

- cAMP RE
- NFAT RE
- NF-κB RE



Dose response curve
GloResponse NFAT-RE-luc2P
HEK293 Cell Line

Case Study: A20 addition controls lymphoma

nature

Vol 459 | 4 June 2009 | doi:10.1038/nature07169

LETTERS

Frequent inactivation of A20 in B-cell lymphomas

Motohiro Kato^{1,2}, Masashi Sanada^{1,5}, Itaru Kato⁶, Yasuharu Sato⁷, Junko Takita^{1,2,3}, Kengo Takeuchi⁸, Akira Niwa⁴, Yuyan Chen^{1,2}, Kumi Nakazaki^{1,4,5}, Junko Nomoto⁹, Yoshitaka Asakura⁹, Satsuki Muto¹, Azusa Tamura¹, Mitsuru Iio¹, Yoshiki Akatsuka¹, Yasuhide Hayashi^{1,2}, Hiraku Mori^{1,3}, Takashi Igarashi¹⁰, Mineo Kurokawa¹, Shigeru Chiba¹, Shigeo Mori¹⁴, Yuichi Ishikawa⁹, Koji Okamoto¹⁰, Kensei Tobinai¹¹, Hitoshi Nakagawa¹⁰, Tatsutoshi Nakahata⁹, Tadashi Yoshino⁷, Yukio Kobayashi⁷ & Seishi Ogawa^{1,2}

A20 is a negative regulator of the NF- κ B pathway and was initially identified as being rapidly induced after tumour-necrosis factor- α stimulation¹. It has a pivotal role in regulation of the immune response and prevents excessive activation of NF- κ B in response to a variety of external stimuli^{2,3}. Recent genetic studies have disclosed putative associations of polymorphic A20 (also called TNFAIP3) alleles with autoimmune disease risk^{4,5}. However, the involvement of A20 in the development of human cancers is unknown. Here we show, using a genome-wide analysis of genetic lesions in 238 B-cell lymphomas, that A20 is a common genetic target in B-lineage lymphomas. A20 is frequently inactivated by somatic mutations and/or deletions in mucosa-associated tissue lymphoma (18 out of 87; 21.8%) and Hodgkin's lymphoma of nodular sclerosing histology (5 out of 15; 33.3%), and, to a lesser extent, in other B-lineage lymphomas. When re-expressed in a lymphoma-derived cell line with no functional A20 alleles, wild-type A20, but not mutant A20, resulted in suppression of cell growth and in duration of apoptosis, accompanied by down regulation of NF- κ B activation. The A20-deficient cells stably generated tumours in immunodeficient mice, whereas the tumorigenicity was effectively suppressed by re-expression of A20. In A20-deficient cells, suppression of both cell growth and NF- κ B activity due to re-expression of A20 depended, at least partly, on cell-surface-receptor signalling, including the tumour-necrosis factor receptor. Considering the physiological function of A20 in the negative modulation of NF- κ B activation induced by multiple upstream stimuli, our findings indicate that uncontrolled signalling of NF- κ B caused by loss of A20 function is involved in the pathogenesis of subsets of B-lineage lymphomas.

Malignant lymphomas of B-cell lineage are mature lymphoid neoplasms that arise from various lymphoid tissues^{6,7}. To obtain a comprehensive registry of genetic lesions in B-lineage lymphomas, we performed a single nucleotide polymorphism (SNP) array analysis of 238 primary B-cell lymphoma specimens of different histologies, including 64 samples of diffuse large B-cell lymphomas (DLBCLs), 52 follicular lymphomas, 35 mantle cell lymphomas (MCLs), and 87 mucosa-associated tissue (MALT) lymphomas (Supplementary Table 1). Three Hodgkin's lymphoma-derived cell lines were also analysed. Interrogating more than 250,000 SNP sites, this platform permitted the identification of copy number changes at an average resolution of less than 12 kilobases (kb). The use of large numbers of

SNP-specific probes is a unique feature of this platform, and combined with the CNAG/ACNAR software, enabled accurate determination of 'allele-specific' copy numbers, and thus allowed for sensitive detection of loss of heterozygosity (LOH) even without apparent copy-number reduction, in the presence of up to 70–80% normal cell contamination^{8,9}.

Lymphoma genomes underwent a wide range of genetic changes, including numerical chromosomal abnormalities and segmental gains and losses of chromosomal material (Supplementary Fig. 1), as well as copy-number-neutral LOH, or uniparental disomy (Supplementary Fig. 2). Each histology type had a unique genomic signature, including a distinctive underlying molecular pathogenesis for different histology types (Fig. 1a and Supplementary Fig. 3). On the basis of the genomic signatures, the initial pathological diagnosis of MCL was re-evaluated and corrected to DLBCL in two cases. Although most copy number changes involved large chromosomal segments, a number of regions showed focal gain and deletion, accelerating identification of their candidate gene targets. After excluding known copy number variations, we identified 46 loci showing focal gain (19 loci) or deletion (27 loci) (Supplementary Tables 2 and 3 and Supplementary Fig. 4).

Genetic lesions on the NF- κ B pathway were common in B-cell lymphomas and found in approximately 40% of the cases (Supplementary Table 1), underpinning the importance of aberrant NF- κ B activation in lymphomagenesis^{10,11} in a genome-wide fashion. They included focal gain/amplification at the *REL* locus (16.4%) (Fig. 1b) and *TRAF6* locus (5.9%), as well as focal deletions at the *PTEN* locus (5.5%) (Supplementary Figs 1 and 4). However, the most striking finding was the common deletion at 6q23.3 involving a 103-kb segment. It exclusively contained the A20 gene (also called TNFAIP3), a negative regulator of NF- κ B activation^{12,13} (Fig. 1b), which was previously reported as a candidate target of 6q23 deletions in ocular lymphoma¹⁴. LOH involving the A20 locus was found in 50 cases, of which 12 showed homozygous deletions as determined by the loss of both alleles in an allele-specific copy number analysis (Fig. 1b, Table 1 and Supplementary Table 4). On the basis of this finding, we searched for possible tumour-specific mutations of A20 by genomic DNA sequencing of entire coding exons of the gene in the same series of lymphoma samples (Supplementary Fig. 5). Because two out of the three Hodgkin's lymphoma-derived cell lines had biallelic A20 deletions/mutations (Supplementary Fig. 6), 24 primary samples from Hodgkin's lymphoma were also analysed for mutations, where

Cancer Genome Project, Department of Pathology, Cell Therapy and Transplantation Medicine, and Hematology and Oncology, Graduate School of Medicine, University of Tokyo, 7-3-1 Hongo, Bunkyo-ku, Tokyo 113-8655, Japan. ²Core Research for Evolutional Science and Technology, Japan Science and Technology Agency, 4-1-8, Honcho, Kawaguchi-shi, Saitama 332-0012, Japan. ³Department of Pediatrics, Graduate School of Medicine, Kyoto University, Sakyo-ku, Yoshino-cho, Shogoin, Sakyo-ku, Kyoto 606-8507, Japan. ⁴Department of Pathology, Okayama University Graduate School of Medicine, Dentistry and Pharmaceutical Sciences, 2-5-1 Shikata-cho, Kita-ku, Okayama 700-8558, Japan. ⁵Division of Pathology, The Cancer Institute of Japanese Foundation for Cancer Research, Japan, 3-10-6 Ariake, Koto-ku, Tokyo 135-8550, Japan. ⁶Hematology Division, Hospital, and ⁷Early Oncogenesis Research Project, Research Institute, National Cancer Center, 5-1-7 Tsukiji, Chuo-ku, Tokyo 104-0045, Japan. ⁸Division of Immunology, Aichi Cancer Center Research Institute, 1-1 Kanokosaka Chikusa-ku, Nagoya 464-8688, Japan. ⁹Quincy Children's Medical Center, 779 Shinobu-cho, Hokkaido-machi, Ohtakawa 207-8577, Japan. ¹⁰Division of Hematology, Internal Medicine, Shizuoka University Fujigaoka Hospital, 1-30, Fujigaoka, Aoba-ku, Yokohama-shi, Kanagawa 227-8501, Japan. ¹¹Department of Pathology, Teikyo University School of Medicine, 2-11-4 Kaga, Itabashi-ku, Tokyo 173-8505, Japan.

712

©2009 Macmillan Publishers Limited. All rights reserved

- Genetic analysis for SNPs and deletions in B-cell lymphomas
- Large number of hits on the NF- κ B regulator, A20
- Examined the ability of wild-type and cancer-associated A20 mutants to control NF- κ B activity.
- Utilized the pGL4.32 [Luc2P/NF- κ B-RE/Hygro] Vector to assay NF- κ B activity.

Kato, M., et al. (2009) *Nature* **459**, 712-716.

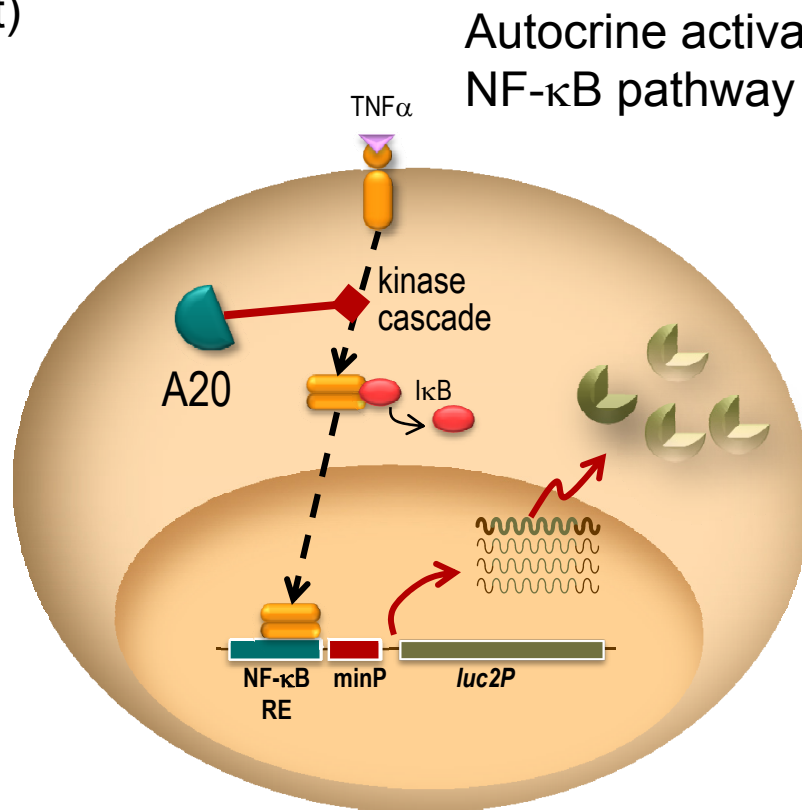
Case Study:
Kato, M., et al.

A20 expression stops NF- κ B pathway

KM-H2 cells do not
express A20
(double mutant)

A20 expressed
through on/off
system

Lymphoma
mutant A20's
expression
through on/off
system



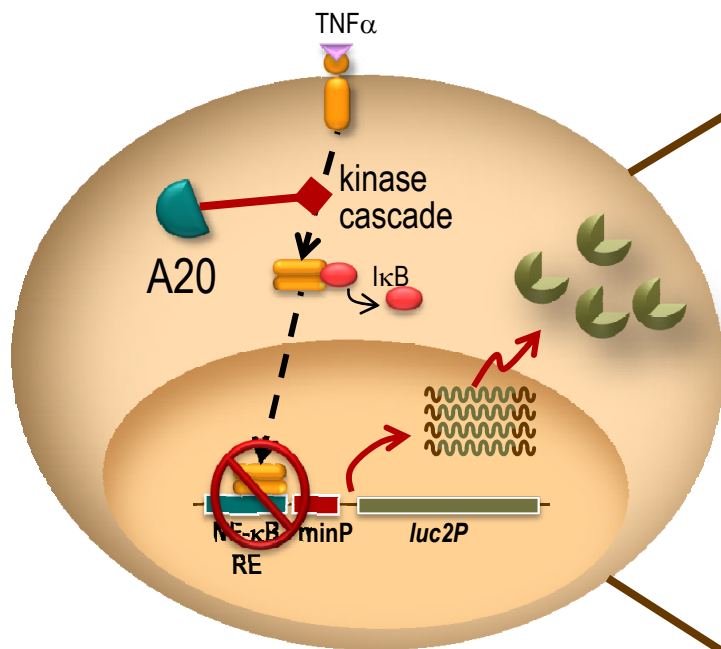
A20 on
↓ Luc Activity
↓ Viability
↑ Apoptosis

A20 Mutant on
= Luc Activity
↑ Viability
X Apoptosis

KM-H2 lymphoma cells stably transfected with
pGL4.32[luc2P/NF- κ B RE/Hygro] Vector

Case Study:
Kato, M., et al.

How the authors determined the cell state



3 Separate Assays
3 Separate sets of cells

Assay for apoptosis with an Annexin V Assay

10⁶ cells/assay
Flow cytometry

Assay for remaining viable cells by measuring reducing potential

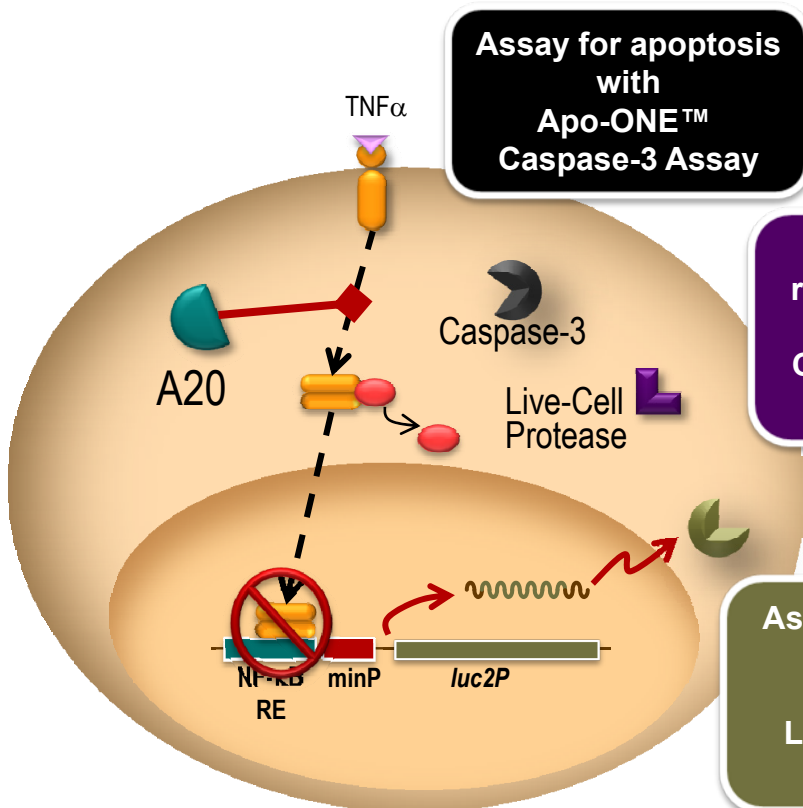
96-well assays
Spectrophotometer

Assay for luciferase activity with Luciferase assay requiring production of cell lysate

96-well assays
Luminometer

Case Study:
Kato, M., et al.

The assays could have been streamlined...



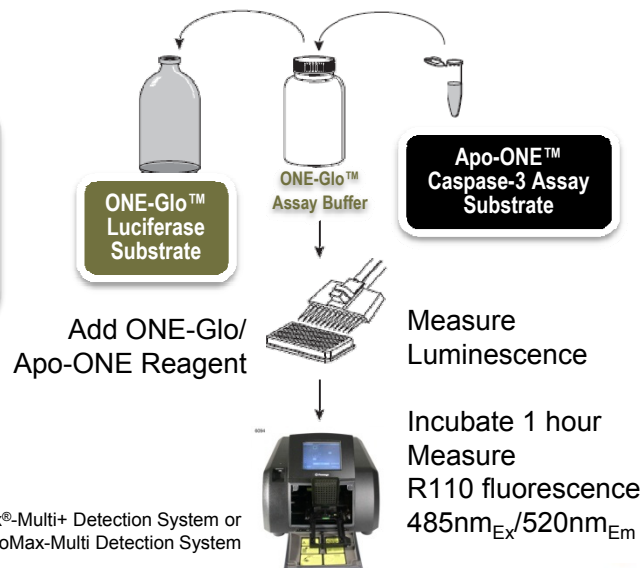
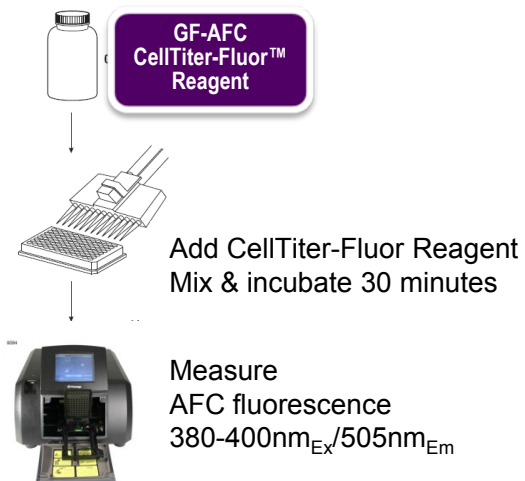
Assay for apoptosis with Apo-ONE™ Caspase-3 Assay

Assay for remaining viable cells with CellTiter-Fluor™ Assay

Assay for luciferase activity with ONE-Glo™ Luciferase Assay System

Measure viable cell number, apoptosis & luciferase activity

- on the same cells
- in the same well
- in the same experiment



GloMax®-Multi+ Detection System or GloMax-Multi Detection System

More Information



[pGL4.29 \[*luc2P*/CRE/Hygro\] Vector Product Protocol](#)



[GloResponse™ CRE-*luc2P* HEK293 Cell Line Technical Bulletin](#)



[pGL4.30 \[*luc2P*/NFAT-RE/Hygro\] Vector Product Protocol](#)



[GloResponse™ NFAT-RE-*luc2P* HEK293 Cell Line Technical Bulletin](#)



[pGL4.32 \[*luc2P*/NF- \$\kappa\$ B-RE/Hygro\] Vector Product Protocol](#)



[GloResponse™ NF- \$\kappa\$ B-RE-*luc2P* HEK293 Cell Line Technical Bulletin](#)



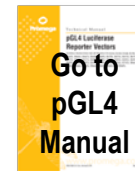
[pGL4.33 \[*luc2P*/SRE/Hygro\] Vector Product Protocol](#)



[pGL4.34 \[*luc2P*/SRF-RE/Hygro\] Vector Product Protocol](#)

pGL4.23-pGL4.28
citations from
HighWire Press®

pGL4.29-pGL4.34
citations from
HighWire Press®



Schagat, T., Paguio, A. and Kopish, K. (2007) Normalizing genetic reporter assays: Approaches and considerations for increasing consistency and statistical significance. *Cell Notes* **17**, 9-12.



Shultz, S., *et al.* (2008) Expanding applications of bioluminescence technology: Predictive bioassays for TNF α biologicals potency and dose-standardization studies. *Promega Notes* **100**, 19-21.



Paguio, A., *et al.* (2006) Using luciferase reporter assays to screen for GPCR modulators. *Cell Notes* **16**, 22-25.



Schumacher, M., Dicato, M. and Diederich, M. (2008) Improved response ratio for NF- κ B inhibition analysis using new luciferase reporter vector (pGL4.32 [*luc2P*/NF- κ B-RE/Hygro]). *Cell Notes* **21**, 18-19.