

Translating Genomes | Personalizing Medicine

X-MAN™ reporter cell lines: Tools to study endogenous promoter activity and protein dynamics

March 2014



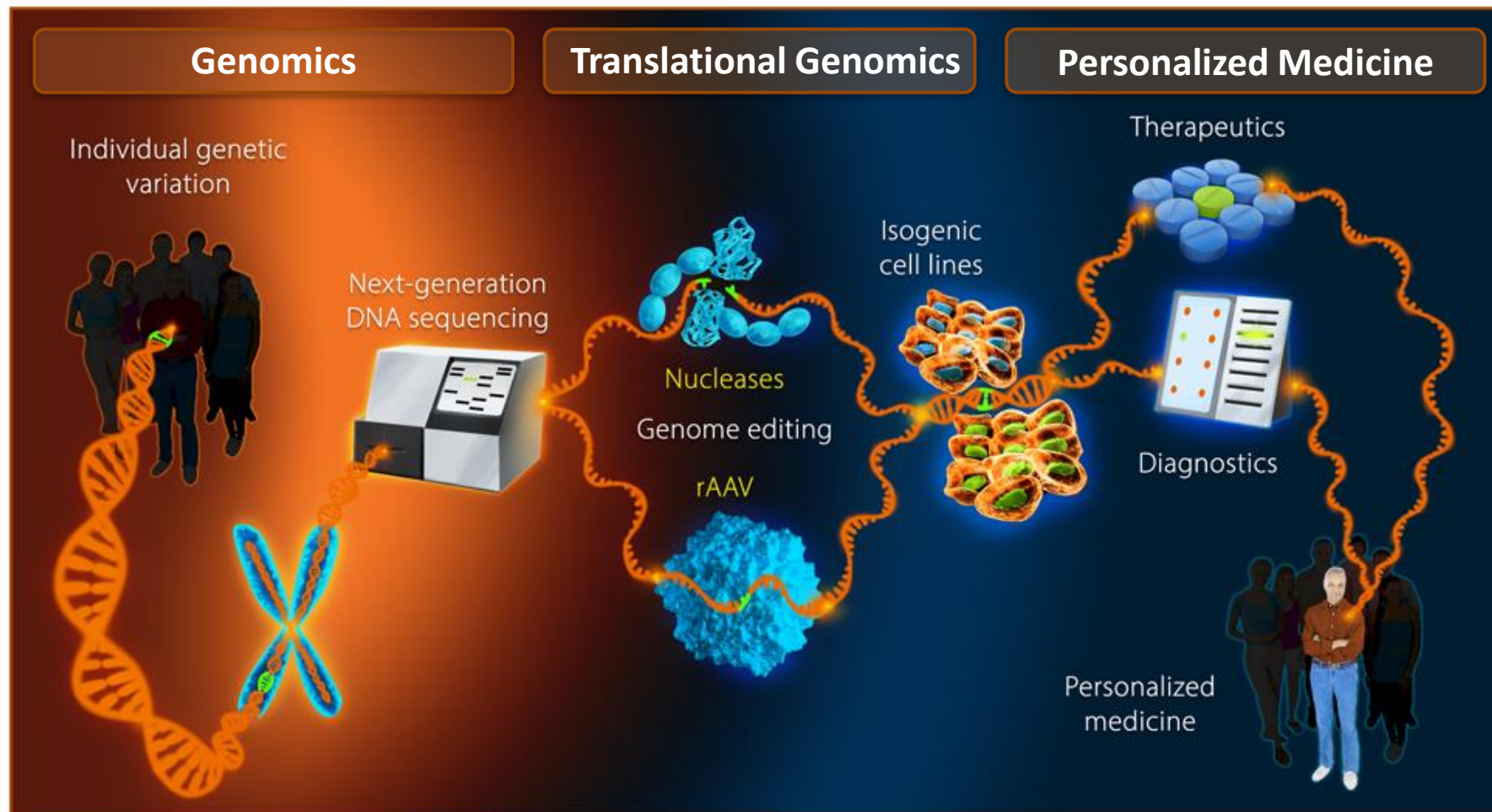
Agenda

➤ Introduction to Horizon's Technology

➤ X-MAN™ HaloTag® Reporter Kits

➤ X-MAN™ NanoLuc™ Reporter Kits

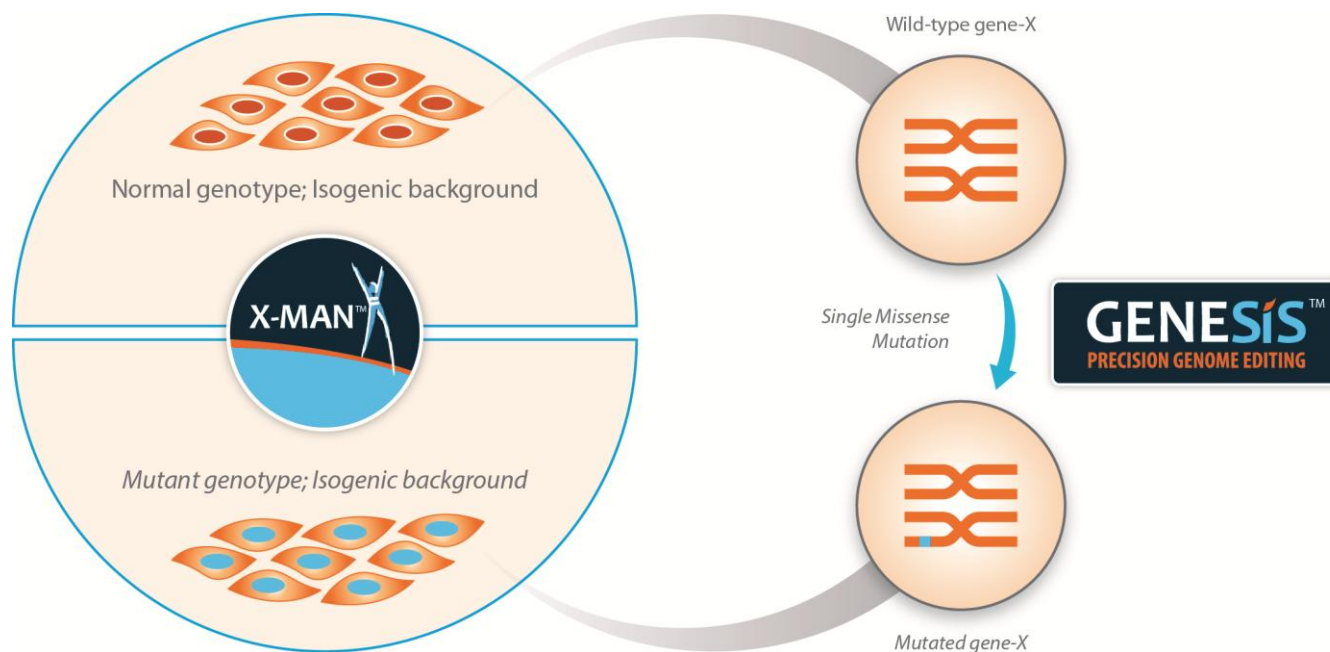
Entering the post-genomic era: Translating genetic information into personalised medicine



- Genome sequencing is now routine and emphasis is shifting to the ‘functional human genome’
- Horizon are experts in ‘Genome Editing’ in human/mammalian cells – **targeted gene KI/KOs**
- **An essential new tool to model patient genetics *in vitro* and discover novel targets, Tx’s and Dx’s**

GENESIS™: Horizon's gene editing platform & X-MAN™ cell lines

- GENESIS™ platform consists of **rAAV and nuclease-based** gene editing technologies
- Generated over 500 genetically-defined X-MAN™ isogenic human cell lines
 - **KOs** to study gene function or target validation
 - Accurate **KI-models** of target patient genetics “**surrogate patients**” + **perfect normal control**
- **Endogenous pathway reporter** lines

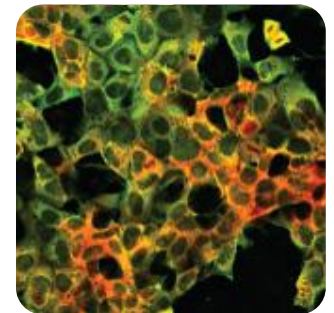
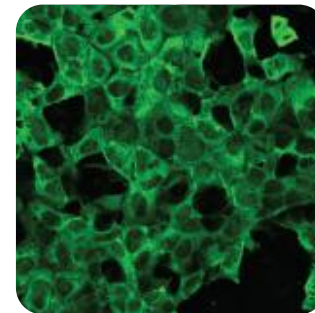
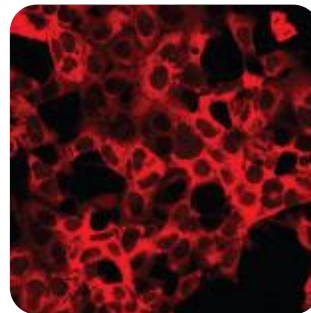


- **GENESIS™ Genome Editing:**
Make any mutation of interest
- Validate putative disease-causing mutations and SNPs
- Definitively study target biology and drug responses
- Create pathway reporter lines
- Build complex diseases from ground up, or sequentially correct them

X-MAN™ Reporter Kits: Live cell endogenous pathway reporters

- rAAV genome editing used to engineer endogenous tags and reporters
- Available in kit format from Horizon
 - **X-MAN™ NanoLuc™ Reporter Kits**
 - Knock-in of Luciferase reporter into endogenous locus for *in vitro* & *in vivo* pathway read-outs
 - **X-MAN™ HaloTag® Reporter Kits**
 - Knock-in of reporters/tag into endogenous locus for *in vitro* imaging / pull-downs / purification

Kits contain cell lines and all reagents necessary to perform a set of experiments



Agenda

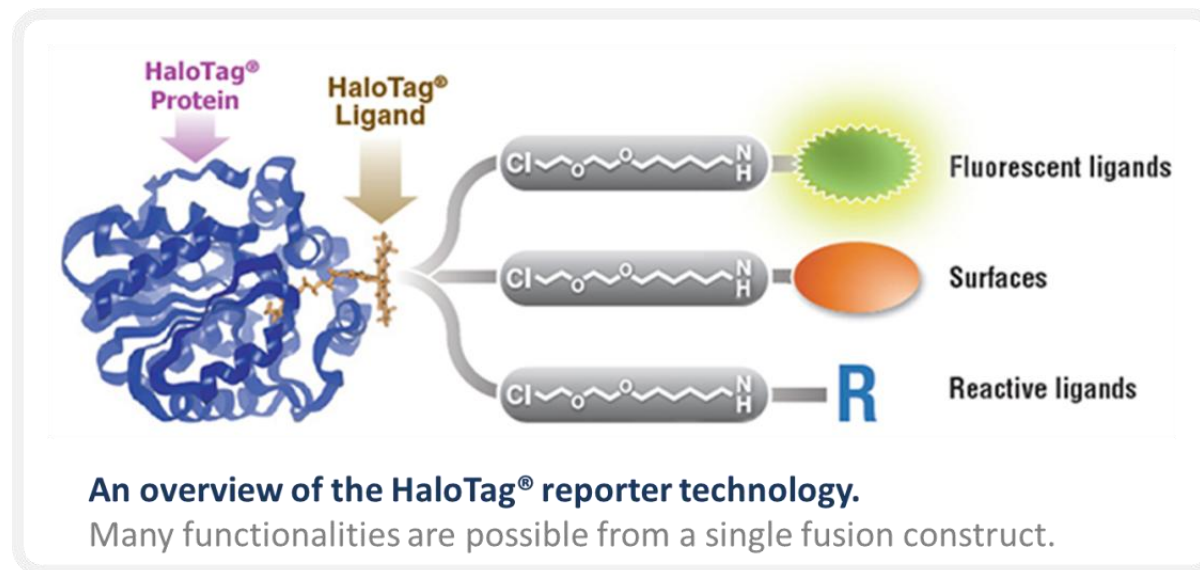
➤ Introduction to Horizon's Technology

➤ X-MAN™ HaloTag® Reporter Kits

➤ X-MAN™ NanoLuc™ Reporter Kits

X-MAN™ HaloTag® Reporter Lines: Technology Background

- X-MAN™ HaloTag® reporter cell lines have been generated using Horizon Discovery's GENESIS™ gene editing platform
- This technology enables introduction of HaloTag® as N- or C-terminal fusions to genes encoding proteins of interest at their native genetic loci, occurring at physiologically relevant levels
- HaloTag® reporter technology has many applications
 - No limit to the experiments that can be performed, only by the genes that you would target
 - Endogenous high-content screening

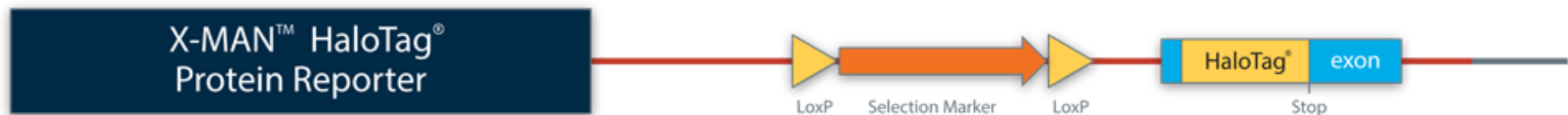


Flexible and efficient system Wider experimental range ... Reliable results

X-MAN™ HaloTag® Reporter Kit

X-MAN™ HaloTag® Protein Reporter Kit

- Engineered to generate N- or C-terminal fusion of HaloTag® to the endogenous protein of interest
- Ideal for studying a wide variety of properties of the protein of interest due to the broad range of HaloTag® ligands available



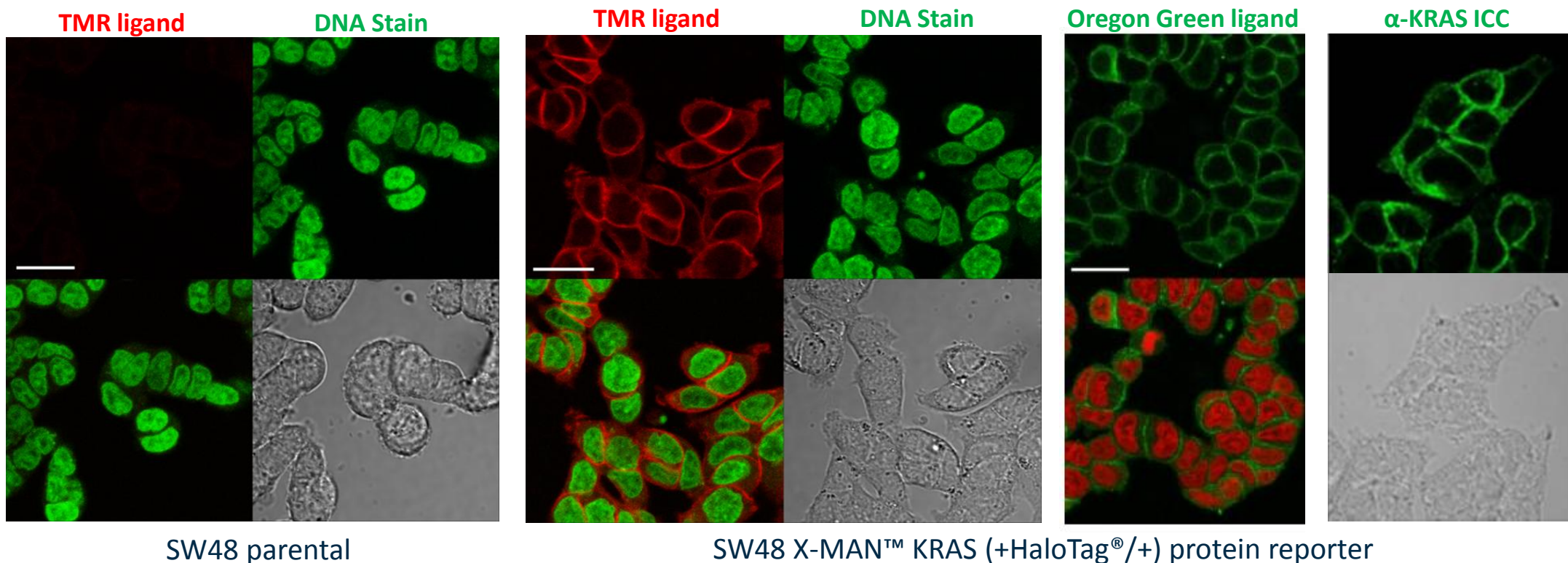
X-MAN™ HaloTag® Protein Reporter Control

- Can be used to demonstrate that the property is specific to the protein of interest, and not conferred by inherent properties of the HaloTag® protein
- Controls for protein reporter lines



HaloTag[®]: Live cell imaging of protein localisation (K-Ras)

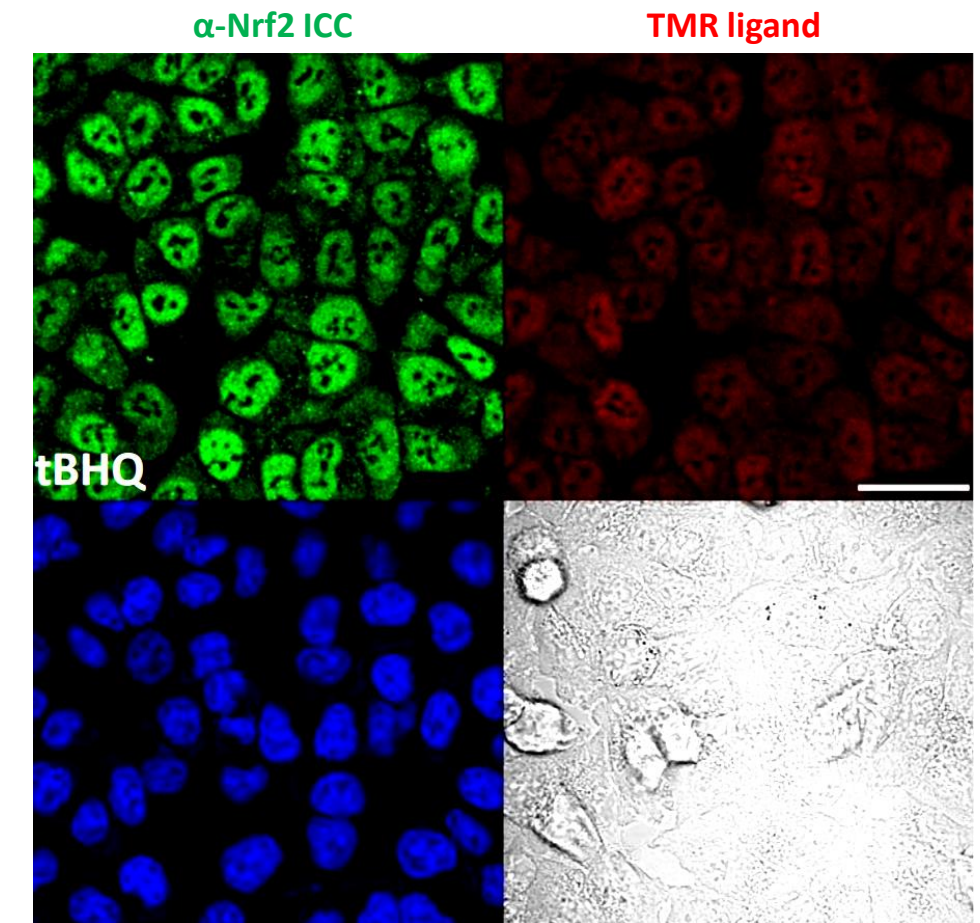
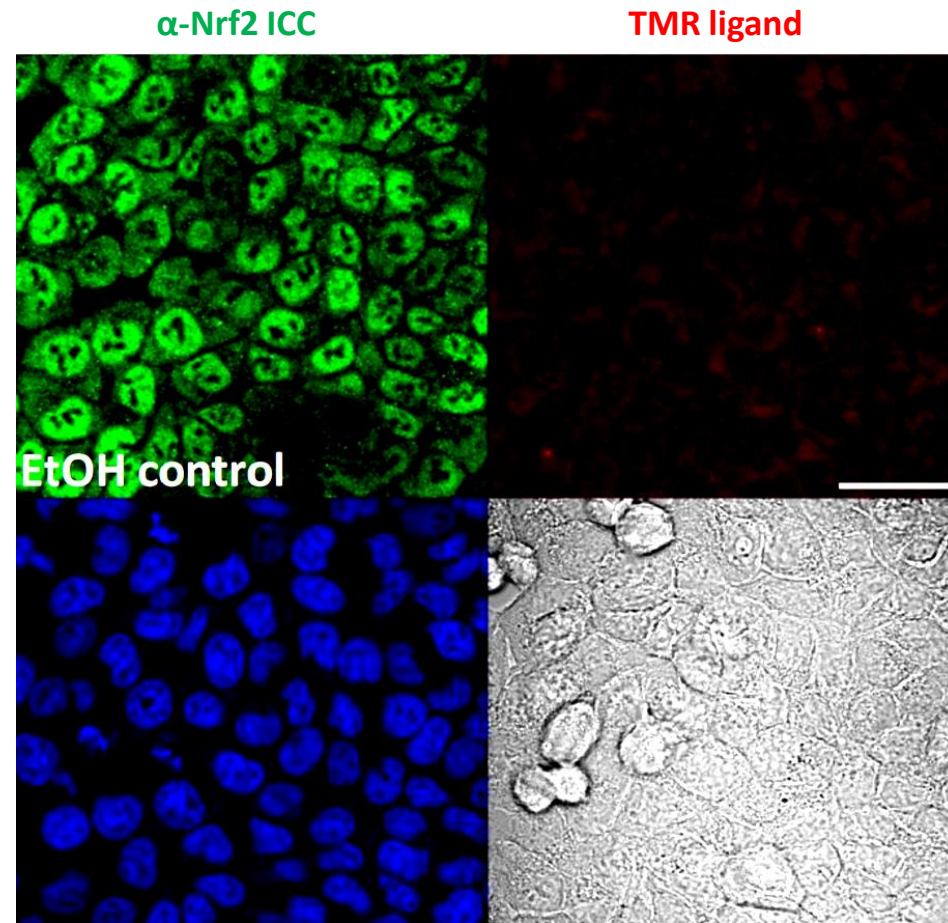
- Some targets/diseases you may want to screen directly for translocation events in live cells
- Easily assayed using HaloTag[®] and cell-permeable fluorescent HT-ligands, available in different colours
- Use of HaloTag[®] ligand shows selective staining for HaloTag[®] engineered lines over unmodified parental cells demonstrating specificity of signal
- HaloTag[®] ligand shows comparable results to ICC



Scale bars = 20 μ m

HaloTag®: Can provide superior alternative to ICC

- Well-characterised increase in stability and nuclear translocation of NFE2L2 (NRF2) in response to *tert*-butylhydroquinone (tBHQ) revealed using HaloTag® reporter cell line but not ICC



HCT116 X-MAN™ NFE2L2 (+HaloTag®/+) protein reporter

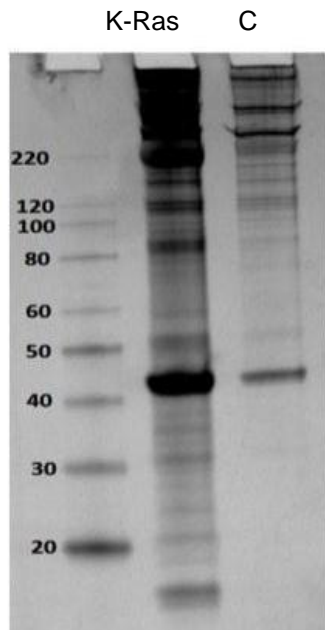
HCT116 X-MAN™ NFE2L2 (+HaloTag®/+) protein reporter

Scale bar = 20 μm

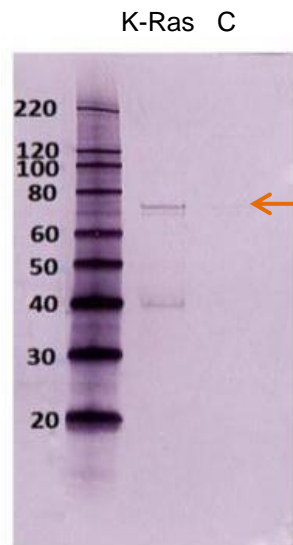
HaloTag®: Pull-down experiments with isogenic reporter lines demonstrate protein enrichment

Protein Enrichment Assays

K-Ras protein (Silver Stain)



Interacting protein (cRAF WB)



- HaloTag® ligand linked to resin enables stringent pull-downs
- High sensitivity via covalent interaction with matrix

Potential Interactors of KRAS Identified by LC/MS/MS from Halo-KRAS SW48 Horizon Cell line

Signaling Proteins And Exchange Factors

Rap1B
RhoA
RAC1
RAB8B
RhoAAct.
FARP1
RAB13
GNA11
RhoG
GNA13
ATP6V1G1
RheB
ATP6V1E1
HRAS
YWHAZ
ErbB2
RAB7L1
ARHGEF12

SpC
32
25
22
20
18
16
14
13
12
10
9
9
9
8
8
8
6
5

Cytochrome C Subunit II and III

COX4I1
UQCRC1
UQCRC2
COX6C
UQCRFS1
MT-CO2
COX2
CYC1
UQCRB
COX5A
COX5B
UQCRQ
UQCRH
COX6B1
COX7A2
COX7C
UQCR10
COX7A2L
CYB5B
COX11

SpC
94
86
83
46
41
37
29
26
25
25
23
13
9
6
4
3
3
3
2
2

NADH Dehydrogenase

NDUF51
NDUFA9
NDUF52
NDUFA13
NDUFV1
NDUFB10
NDUFA7
NDUFA6
NDUFA12
NDUF58
NDUF53
NDUFB5
NDUFA5
NDUFB9
NDUF57
NDUFV2
NDUFB6
NDUF55
NDUFA2
NDUFB4
NDUFAF3
NDUFA8
NDUF56
NDUFC2
NDUFB7
NDUFAF4
NDUFA10
NDUFA3
NDUF54
NDUFA4
NDUFB3
NDUFA11

SpC
77
41
38
34
33
28
26
26
24
21
19
17
17
16
15
14
12
11
10
10
10
9
8
8
7
6
6
5
5
4
3
2

X-MAN™ HaloTag® Kits Currently Available

NCBI Gene Name	Common name	X-MAN™ (+HaloTag®/+) protein reporter	X-MAN™ (+HaloTag®/+) protein reporter of mutated variants	X-MAN™ HaloTag® protein reporter control
HIF1A	HIF1α	x		x
NFE2L2	NRF2	x		x
KRAS	K-Ras	x	G12C	x
			G12D	
			G12V	
			G13D	

- Additional lines are constantly being developed by Horizon or can be developed as a custom project
- **X-MAN™ HaloTag® kits come with:**
 - A vial of the cell line of interest
 - Its parental cell line which can act as a control, and
 - All reagents and buffer necessary to perform a set of experiments.

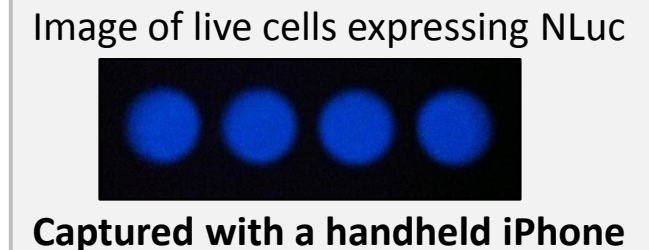
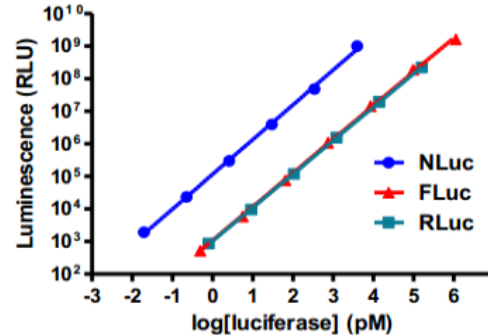
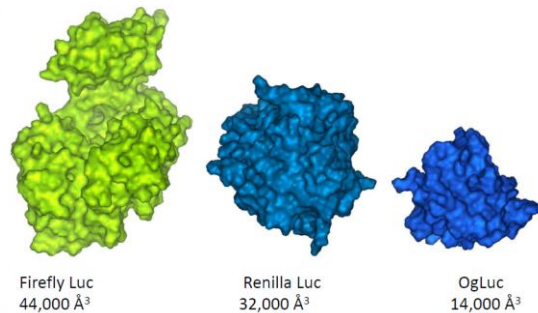
These products are governed by HaloTag® and Horizon Discovery Limited Use Label Licenses

Agenda

- Introduction to Horizon's Technology
- X-MAN™ HaloTag® Reporter Kits
- X-MAN™ NanoLuc™ Reporter Kits

X-MAN™ NanoLuc™ Reporter Lines: Technology Background

- NanoLuc™ derived from a new luciferase from Deep Ocean Shrimp (Oplophorus)
- Serially mutated to create a monomeric & 150-fold brighter luciferase vs FF and Renilla
- Secreted or intracellular; ATP independent light emission with novel substrates
- **Able to detect gene/protein levels at very low endogenous expression levels**
- Endogenous high-throughput screening now possible



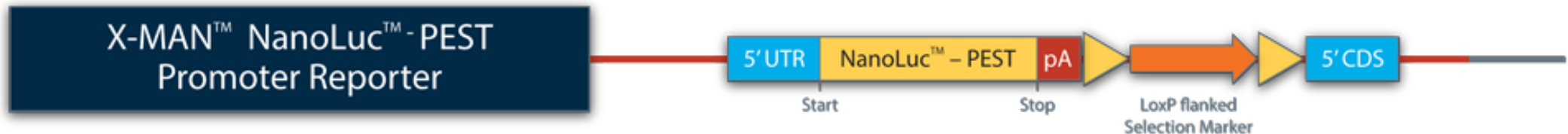
150 fold increased specific activity compared with firefly or Renilla luciferase

Better biology ... Endogenous HTS ... Biologically relevant results

X-MAN™ NanoLuc™ Kit Formats

X-MAN™ NanoLuc™-PEST promoter kit

- Promoter fusions with PEST-NanoLuc™ for rapid turnover and thus dynamic signal kinetics
- Short intracellular lifetime



X-MAN™ NanoLuc™ protein kit

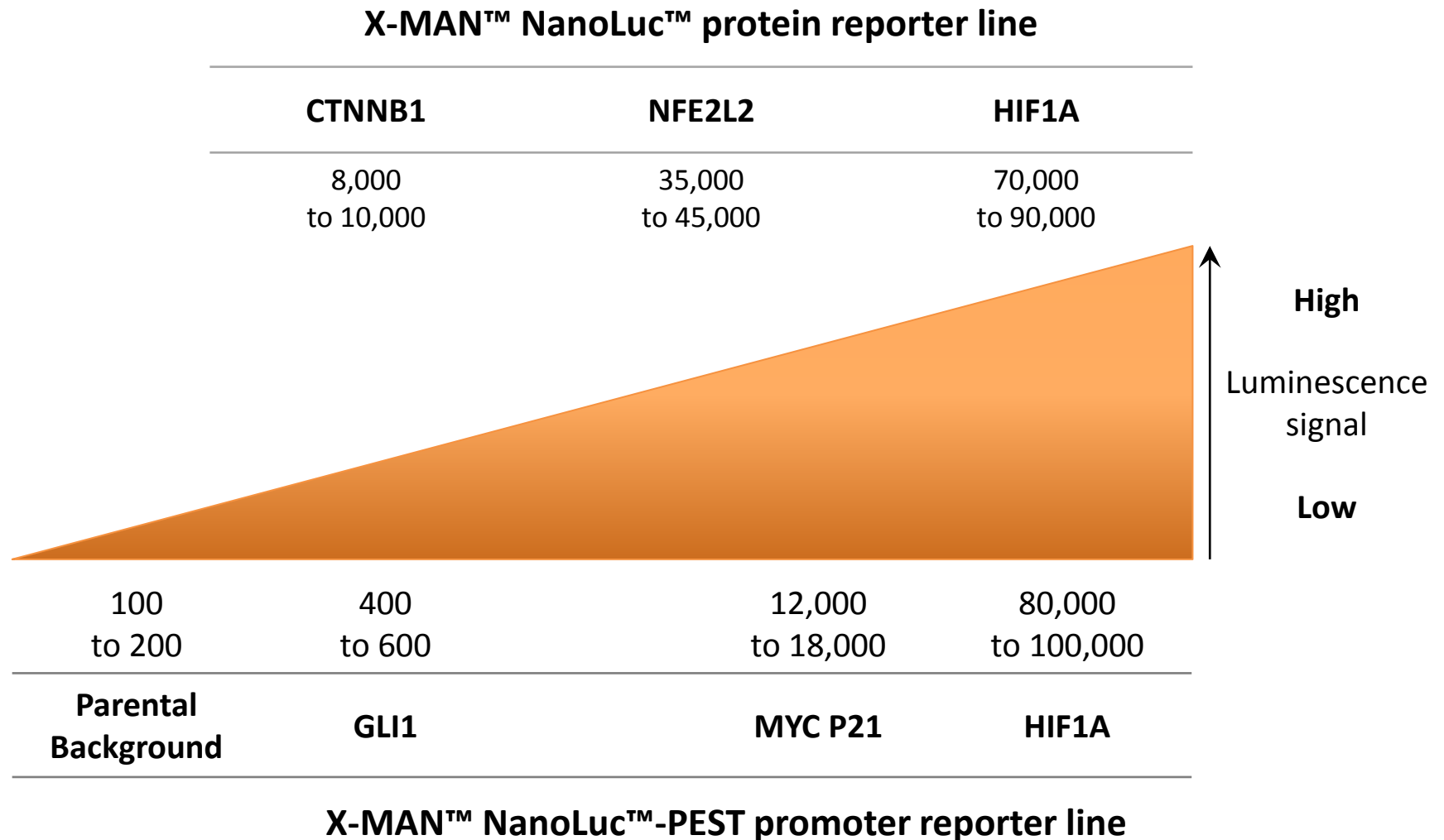
- Engineered to generate N- or C-terminal fusion of NanoLuc™ to the endogenous protein of interest
- Ideal for use in protein expression reporter assays



X-MAN™ NanoLuc™ protein reporter control cell lines are also available

NanoLuc™: Signal Range

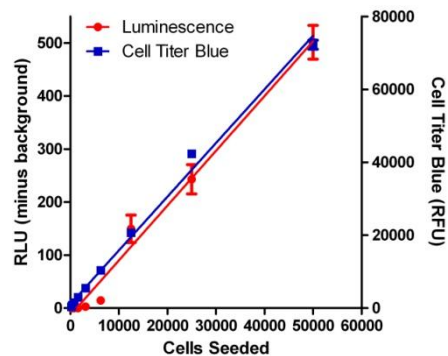
- Baseline luminescence signal is specific for each NanoLuc™ reporter line, dependent on gene or protein expression levels



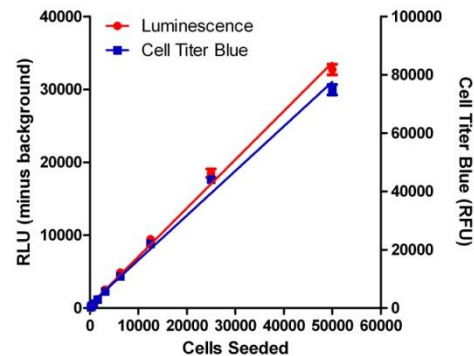
NanoLuc™: Signal Linearity

- Good signal linearity with increasing cell number
- Minimal background luciferase signal in parental cells, even at high cell densities

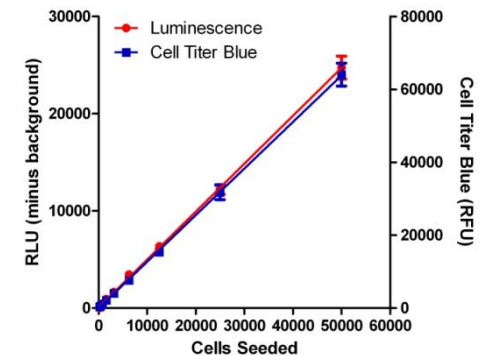
X-MAN™ GLI1 NanoLuc™-PEST Promoter Reporter



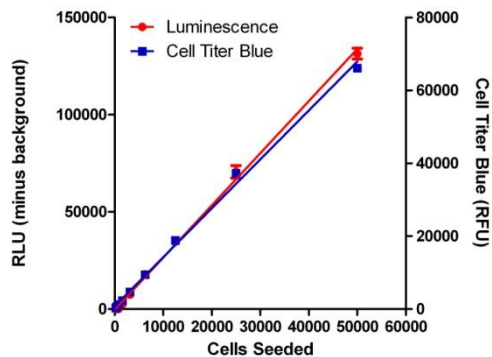
X-MAN™ MYC NanoLuc™-PEST Promoter Reporter



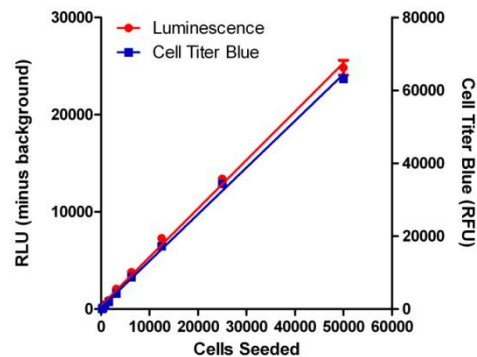
X-MAN™ P21 NanoLuc™-PEST Promoter Reporter



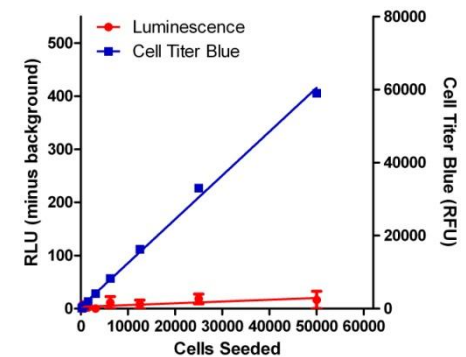
X-MAN™ HIF1A NanoLuc™-PEST Promoter Reporter



X-MAN™ HIF1A (+NanoLuc™/+) Protein Reporter



HCT116 Parental Cell Line

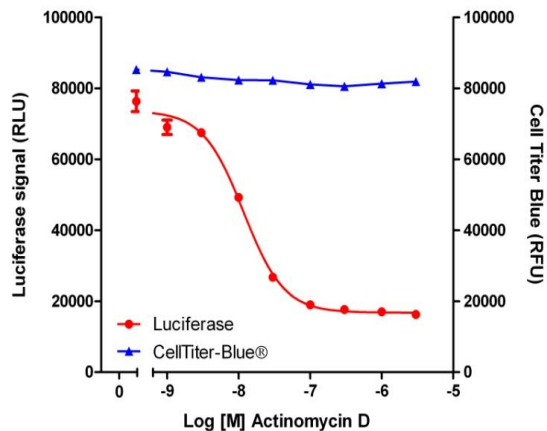


NanoLuc™: Excellent signal dynamics

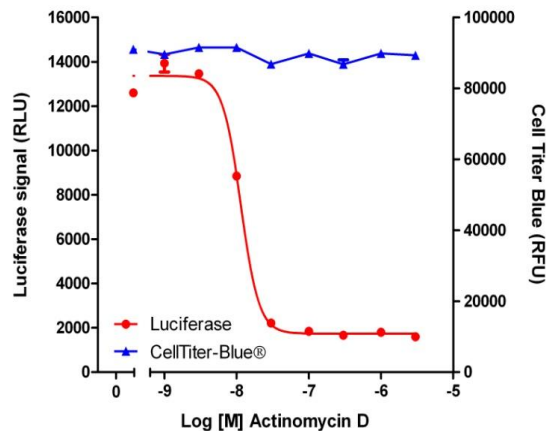
➤ Dynamics of promoter reporter lines:

- 6h treatment with a Actinomycin D
- Rapid and robust decreases in luciferase signal
- Good induction kinetics on inducible genes
- CellTiter-Blue® to control for effects on cell viability

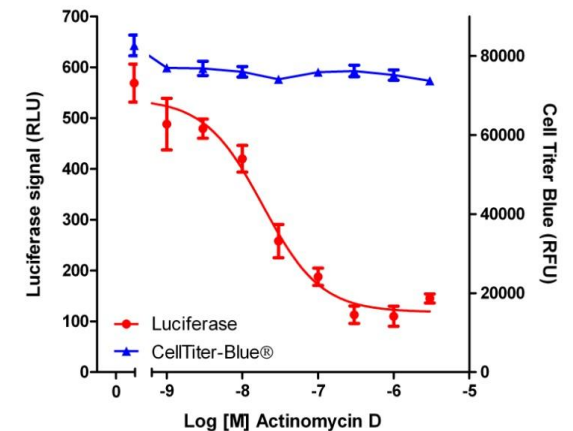
X-MAN™ HIF1A NanoLuc™-PEST Promoter Reporter



X-MAN™ MYC NanoLuc™-PEST Promoter Reporter



X-MAN™ GLI1 NanoLuc™-PEST Promoter Reporter



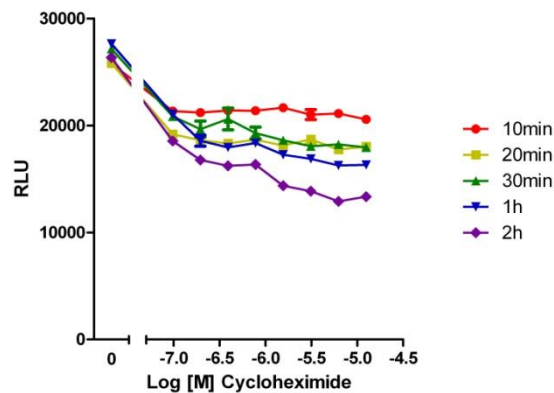
NanoLuc™: Excellent signal dynamics

➤ Dynamics of protein reporter lines:

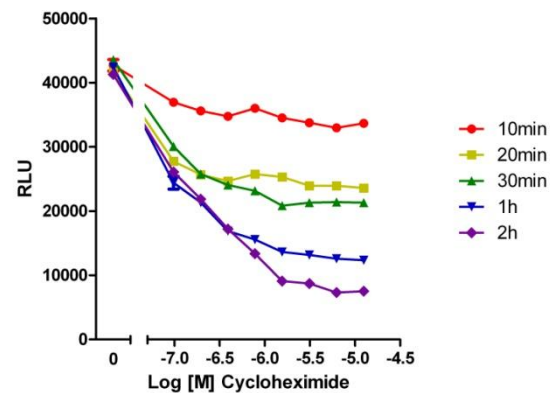
- Turnover: Treatment with Cycloheximide ($t_{1/2}$ dictated by target genes in this scenario)
- Accumulation: 4h treatment with bortezomib
- Good repression and accumulation kinetics of signals generally observed

Turnover

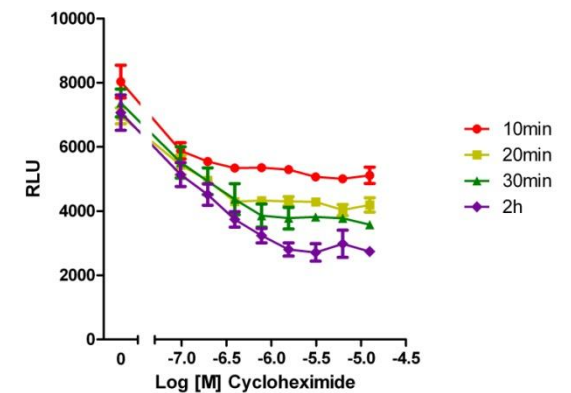
X-MAN™ HIF1A (+NanoLuc™/+) Protein Reporter



X-MAN™ NFE2L2 (+NanoLuc™/+) Protein Reporter

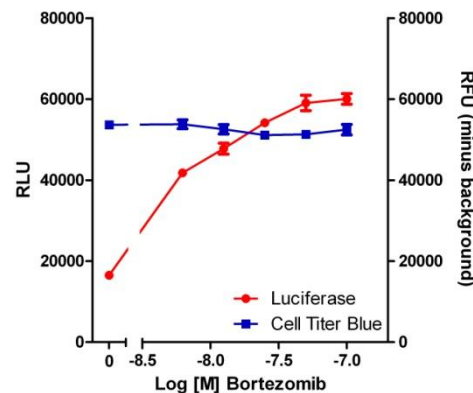


X-MAN™ CTNNB1 (+NanoLuc™/+) Protein Reporter

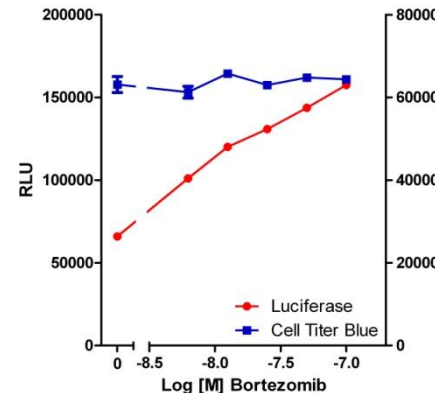


Accumulation

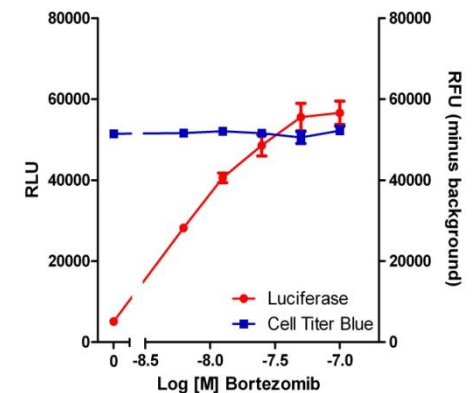
X-MAN™ HIF1A (+NanoLuc™/+) Protein Reporter



X-MAN™ NFE2L2 (+NanoLuc™/+) Protein Reporter



X-MAN™ CTNNB1 (+NanoLuc™/+) Protein Reporter

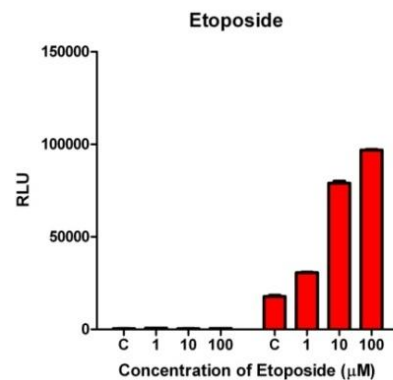
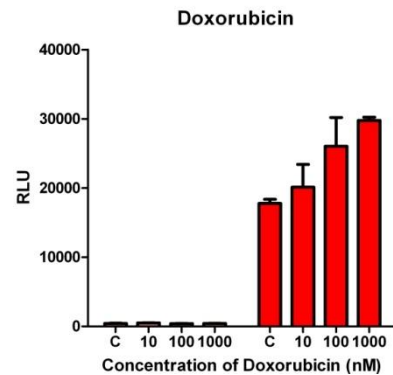


NanoLuc™ Functional Validation: p21 promoter reporter

➤ Monitor modulation of p21 transcription by DNA damaging agents

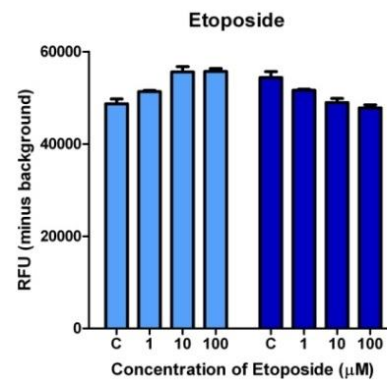
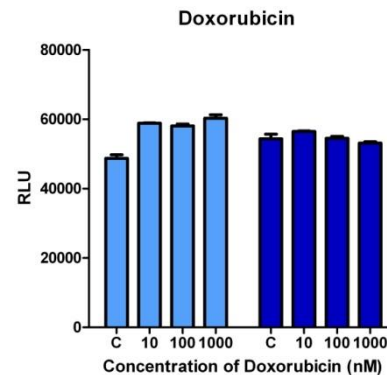
- 6h drug treatment, cell viability measured by multiplexing luciferase assay with CellTiter-Blue®

p21 Transcription
(NanoLuc™ activity measurement)



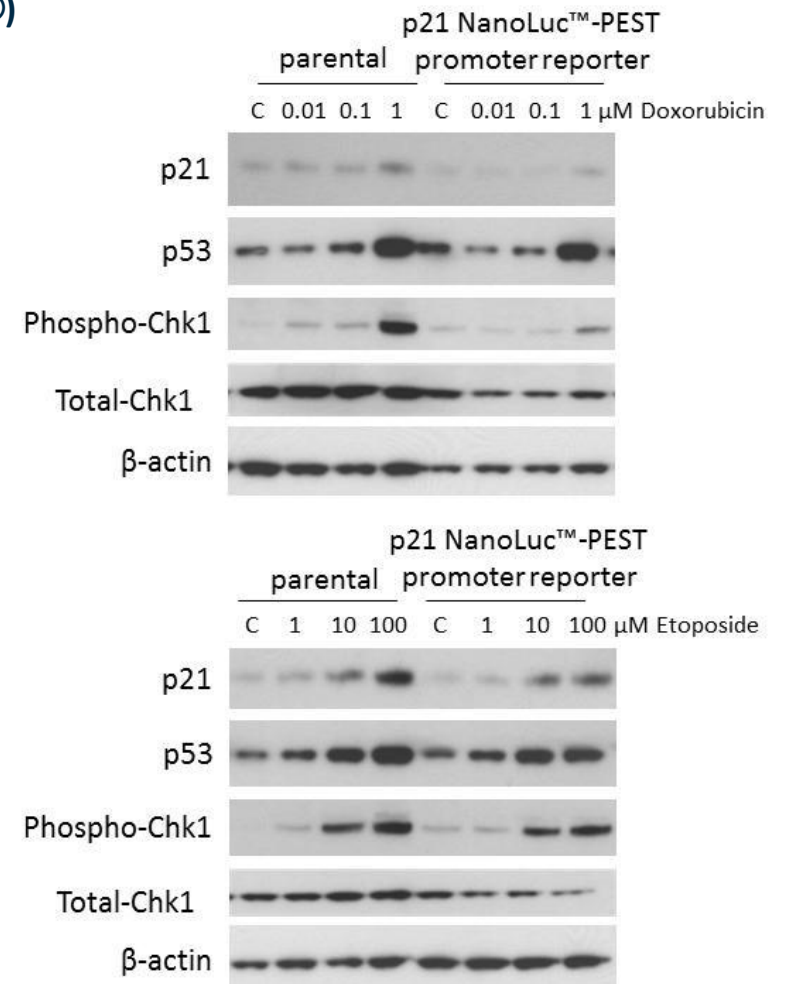
Parental
X-MAN™ p21 NanoLuc™-PEST promoter reporter

Cell Viability
(measured by CellTiter-Blue®)



Parental
X-MAN™ p21 NanoLuc™-PEST promoter reporter

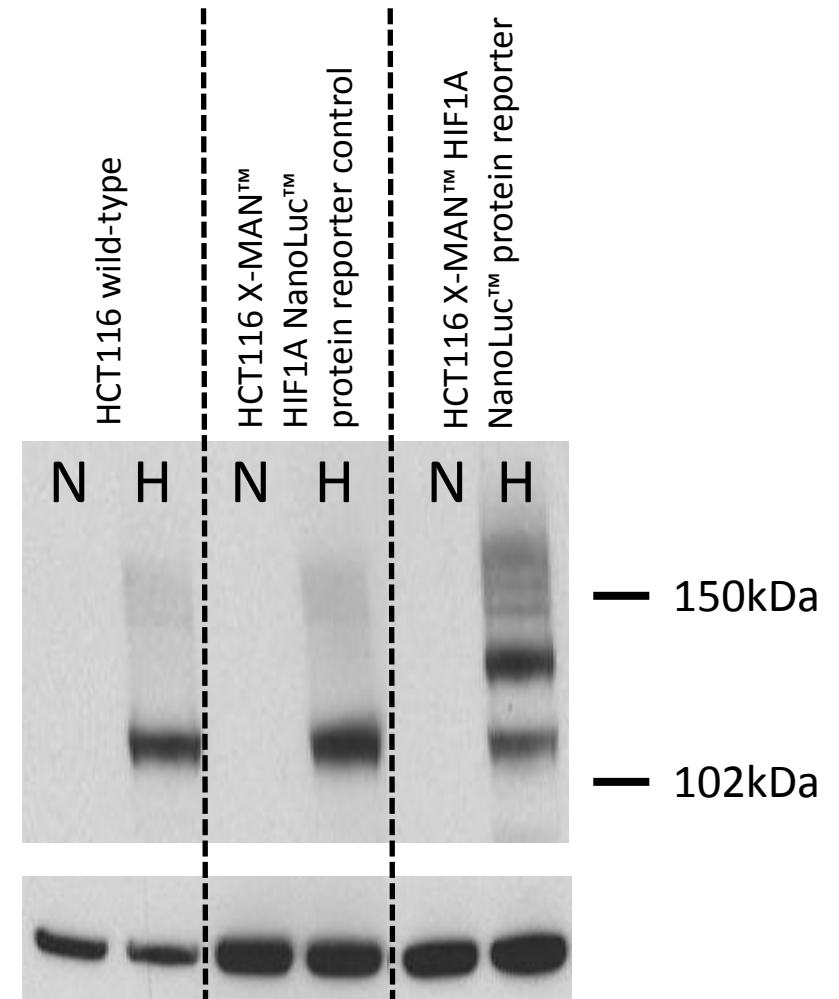
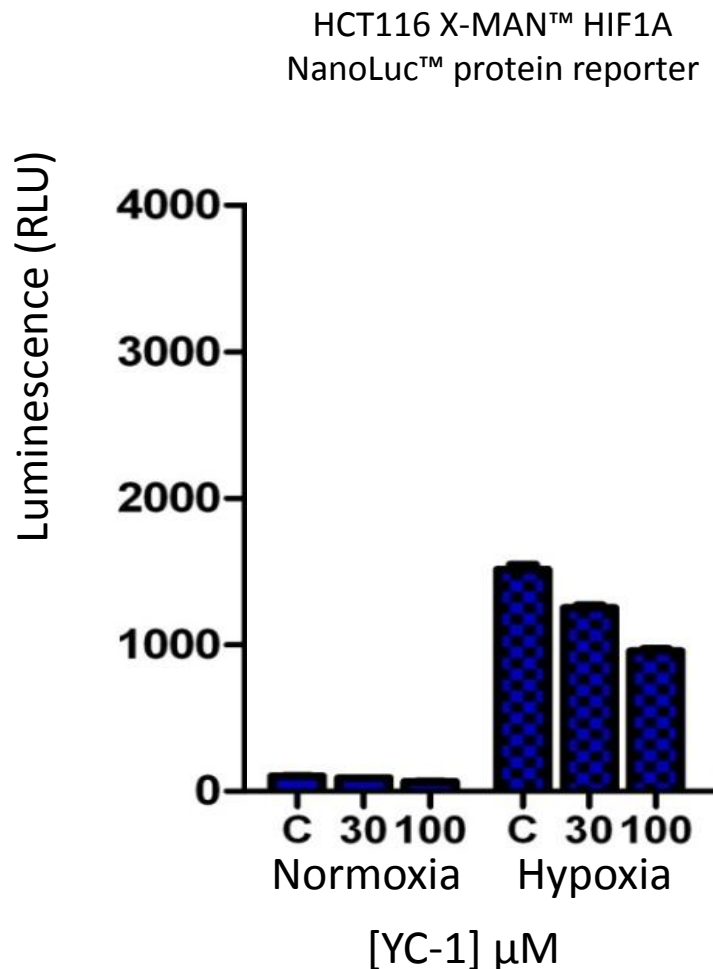
Protein Expression



NanoLuc™ Functional Validation: HIF1A protein reporter

➤ Monitor induction of HIF1A protein by hypoxia

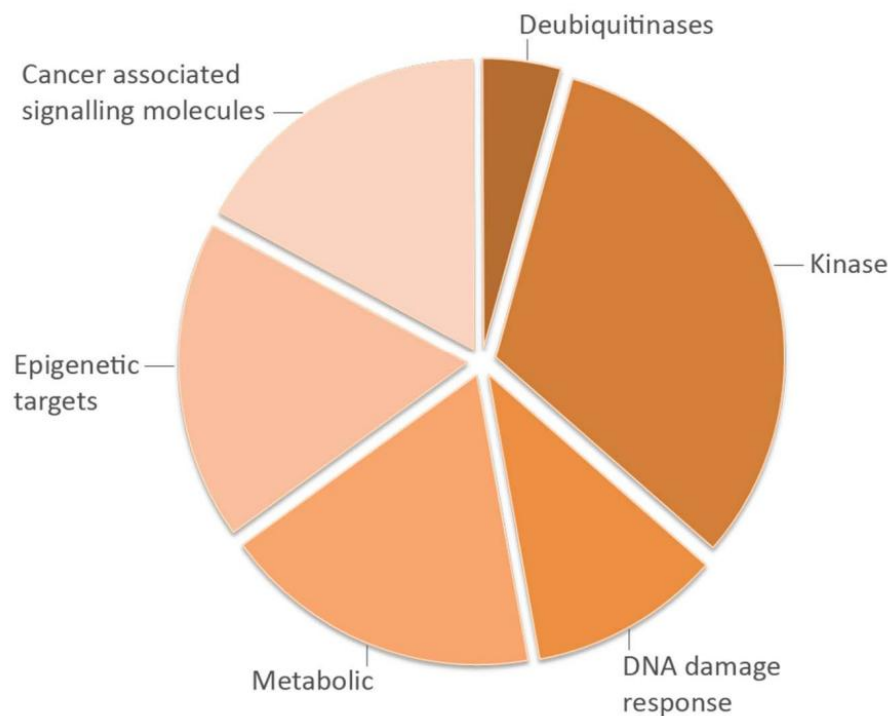
- Natural induction mechanisms preserved – 1% O₂ increases the HIF1A fusion protein
- Treatment with YC-1 HIF1A inhibitor decreases expression of HIF1A under hypoxia as determined by NanoLuc™ protein reporter



NanoLuc™ Screening Application: HIF1A protein reporter

➤ Used in 2 different screening formats

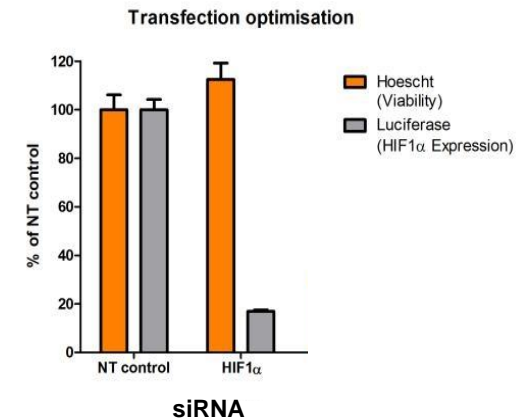
- siRNA screen at Horizon using our library of over 2000 druggable siRNAs
- Compound screen at NCATS, using the NCGC Pharmaceutical Collection of approved and investigational drugs



	siRNA Screen	Compound Screen
Plate format	384	1536
Number of Reagents	2235	2514
Z'	0.5	0.7
%CV	6.5%	6.1%

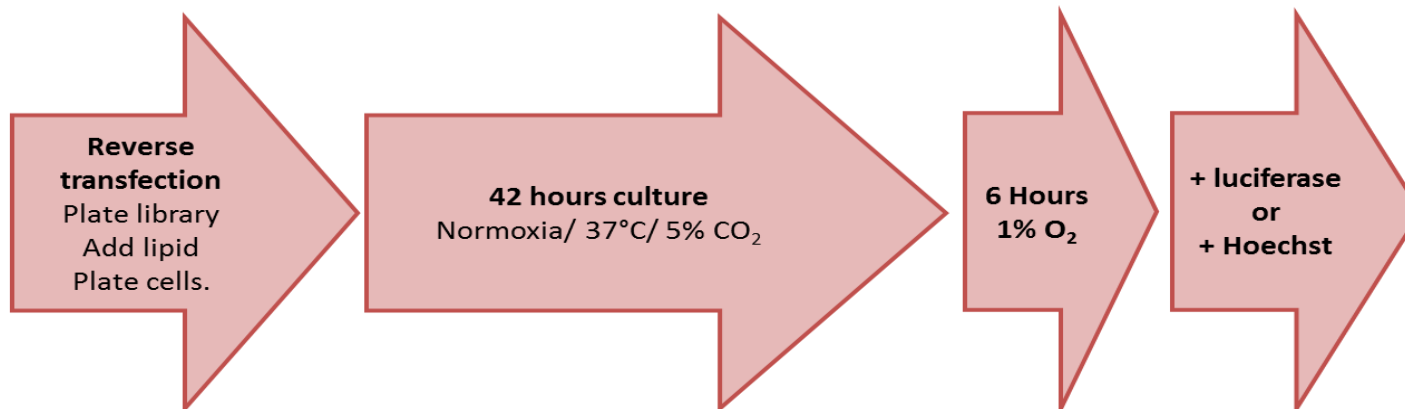
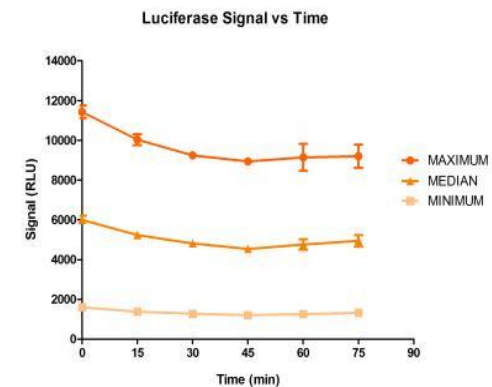
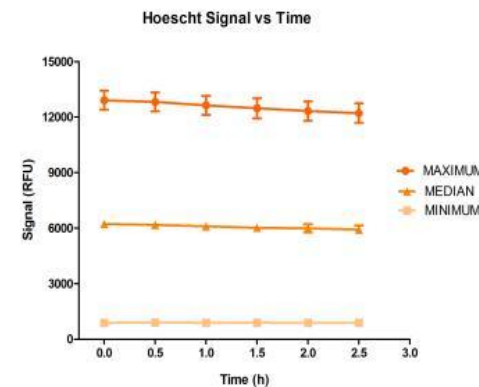
NanoLuc™ Screening Application: HIF1A siRNA Screen

➤ Transfection conditions were optimised to provide conditions that achieve optimal siRNA knockdown with minimal effects on viability.



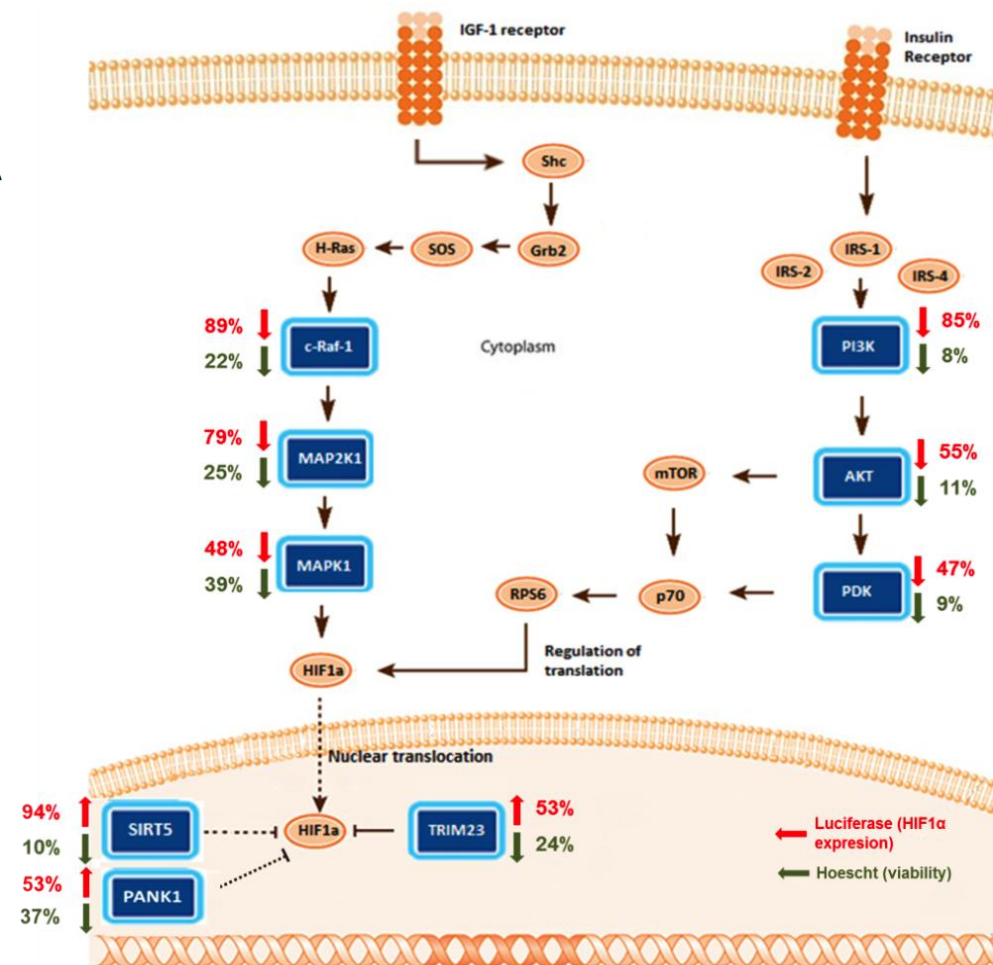
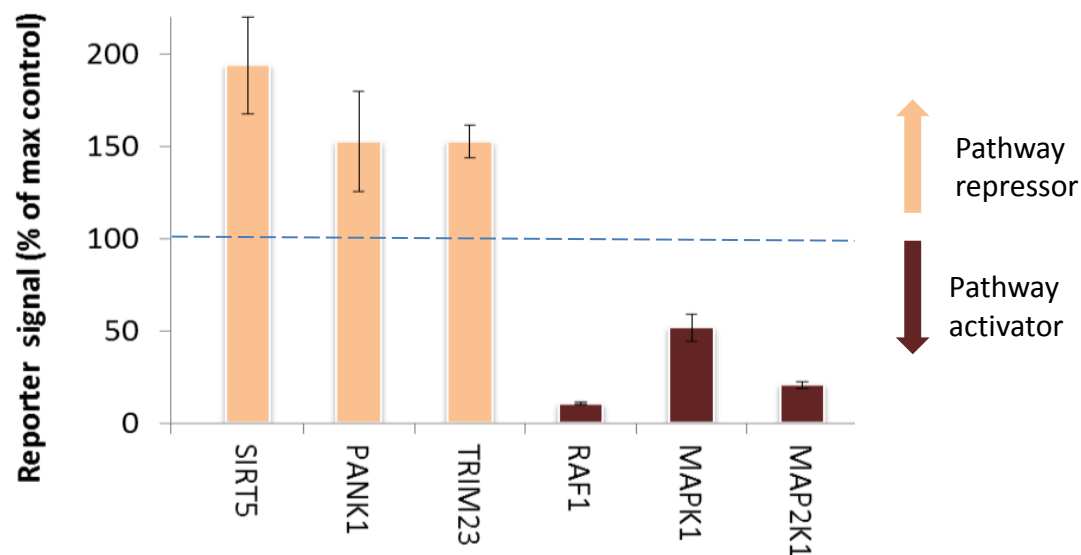
➤ We concurrently measured:

- Hoescht nuclear marker to assess viability
- NanoLuc™ luciferase signal to assess HIF1 α protein levels



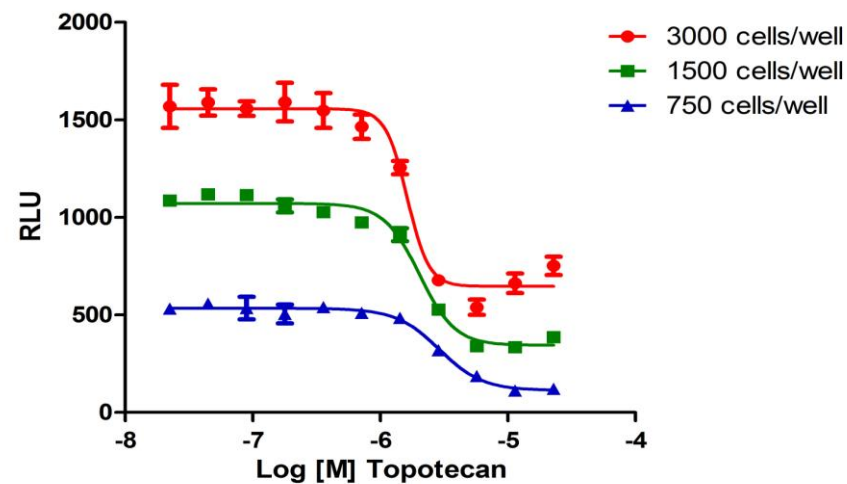
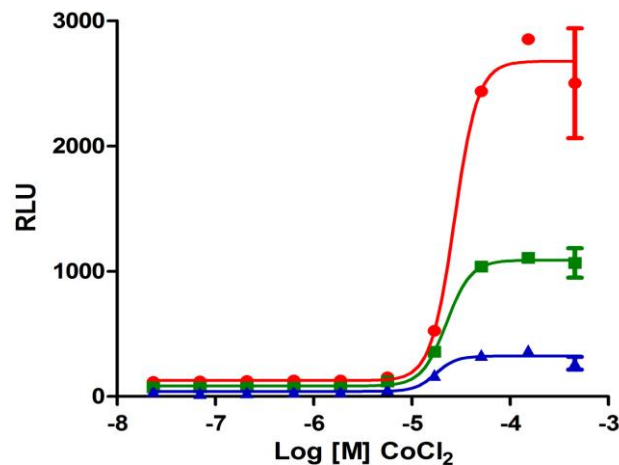
NanoLuc™ Screening Application: HIF1A siRNA Screen

- Subset of kinase, DUB & epigenome hits shown
- Known and novel transcription targets found
- Modulators both activating and repressing HIF1A



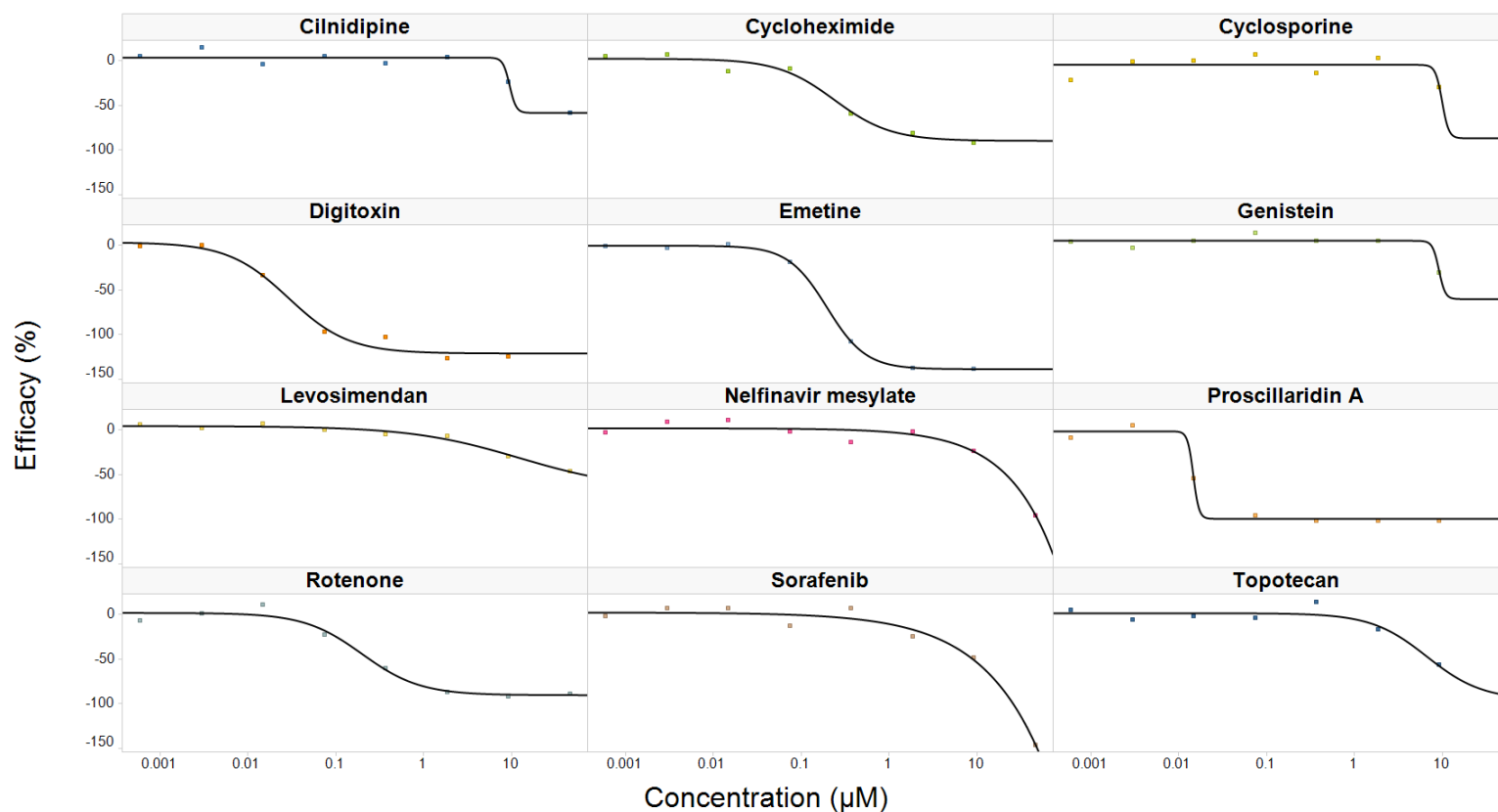
NanoLuc™ Screening Application: NCATS HIF1A Compound Screen

- Cells were seeded into 1536-well plates and induction of HIF1A performed using 1% oxygen or the hypoxic-mimetic CoCl_2
- Knockdown of HIF1A protein reporter signal was confirmed using Topotecan after a 16h drug exposure
 - 1500 cells/well selected for screen, giving a good assay window



NanoLuc™ Screening Application: NCATS HIF1A Compound Screen

- A high-throughput screen was run in 1536-well plate format using the NCGC Pharmaceutical Collection of approved and investigational drugs.
- 1500 cells per well, 18-hour compound incubation time, 1% O₂
 - 7 concentrations (0.6 nM-46 μM or 1.1 nM-92 μM), duplicate runs
 - The assay performed well returning strong hits for known regulators
 - QC parameters were good with Z' of 0.70 and CV of 6.1%.



X-MAN™ NanoLuc™ Kits Currently Available

NCBI Gene Name	Common name	X-MAN™ NanoLuc™- PEST promoter reporter	X-MAN™ (+NanoLuc™/+) protein reporter	X-MAN™ NanoLuc™ protein reporter control
HIF1A	HIF1α	x	x	x
NFE2L2	NRF2	---	x	x
CTNNB1	β-CATENIN	---	x	x
CDKN1A	p21, CIP1, P21 ^{CIP1}	x	---	---
MYC	c-MYC	x	---	---
GLI1	GLI	x	---	---

- Additional lines are constantly being developed by Horizon or can be developed as a custom project
- **X-MAN™ NanoLuc™ kits are supplied with:**
 - A vial of the cell line of interest
 - All reagents and buffer necessary to perform a set of experiments
- Kits are available in 10, 100, 500, 1,000 ml and custom reporter volume sizes
- These products are governed by NanoLuc™ and Horizon Discovery Limited Use Label Licenses

Selection of Customers / Collaborators (>650 unique partners)

INDUSTRY



ACADEMIA





Your Horizon Contact:

Holly Astley

Team Leader – Cell Line Production

h.astley@horizondiscovery.com

+44 (0) 1223 655 580 (Front desk)

Horizon Discovery Ltd, Building 7100, Cambridge Research Park, Waterbeach, Cambridge, CB25 9TL, United Kingdom

Tel: +44 (0) 1223 655 580 (Reception / Front desk) Fax: +44 (0) 1223 862 240 Email: info@horizondiscovery.com Web: www.horizondiscovery.com