



BCL2 inhibition by ABT-199 in T-cell acute lymphoblastic leukemia (T-ALL)

Pieter Van Vlierberghe

Bioluminescent Cell-based Assay Seminar Day
Monday 31 march 2014
UCL De Duve institute
Brussels, Belgium

Short Bio sketch



PhD 2003-2008
Pediatric Oncology
Erasmus Medical Center
Rotterdam, The Netherlands



Postdoc 2008-2013
Institute for Cancer Genetics
Columbia University Medical Center
New York, USA

Short Biosketch



**COLUMBIA UNIVERSITY
MEDICAL CENTER**

Discover. Educate. Care. Lead.



Turner BioSystems Modulus® II
Microplate Multimode Reader

Short Bio sketch



Odysseus type II grant
2013-2017
Center for Medical Genetics
Ghent University



GloMax®-Multi
Microplate Reader



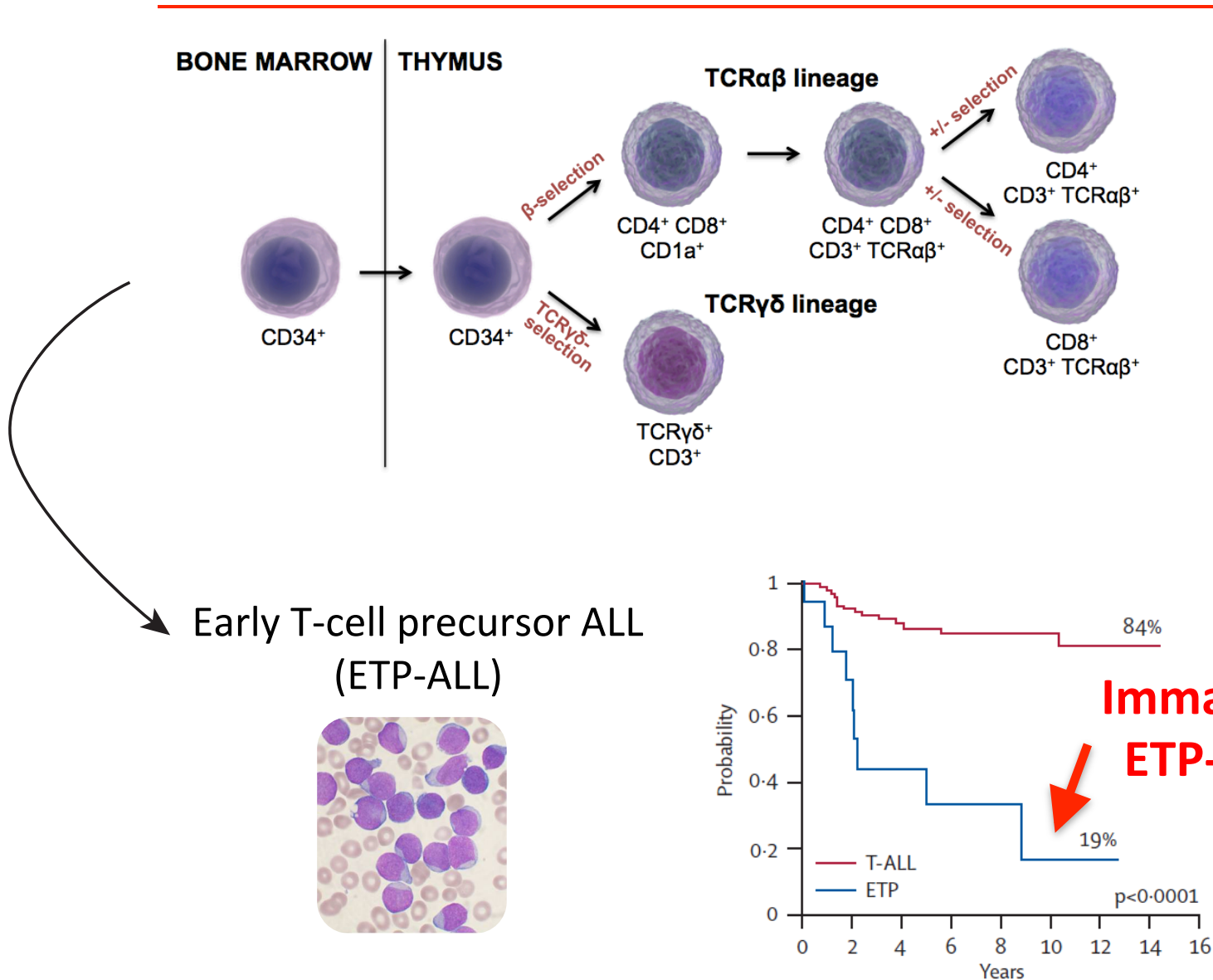
CellTiter-Glo®
Luminescent Cell Viability

Caspase-Glo® 3/7
Apoptosis Assay

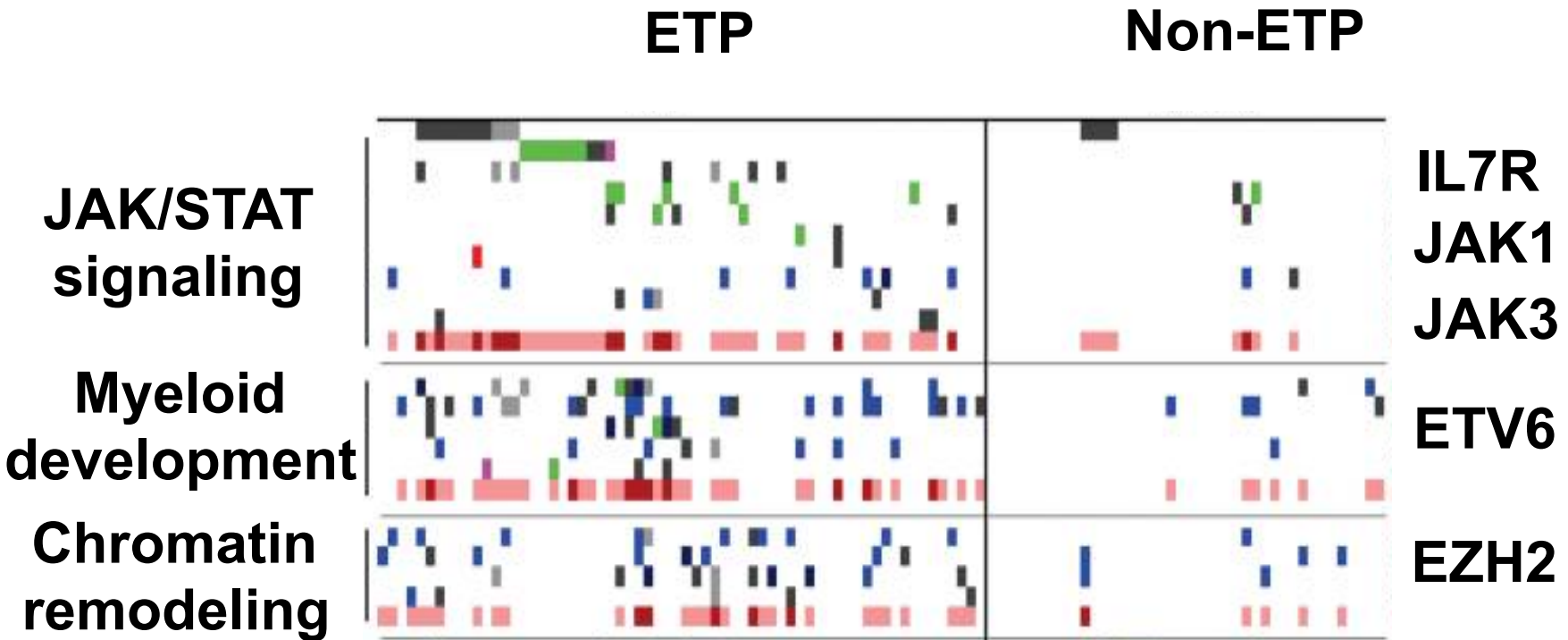
BCA protein assays

Luciferase assays
miRNA-mRNA regulation

From normal development to malignant transformation



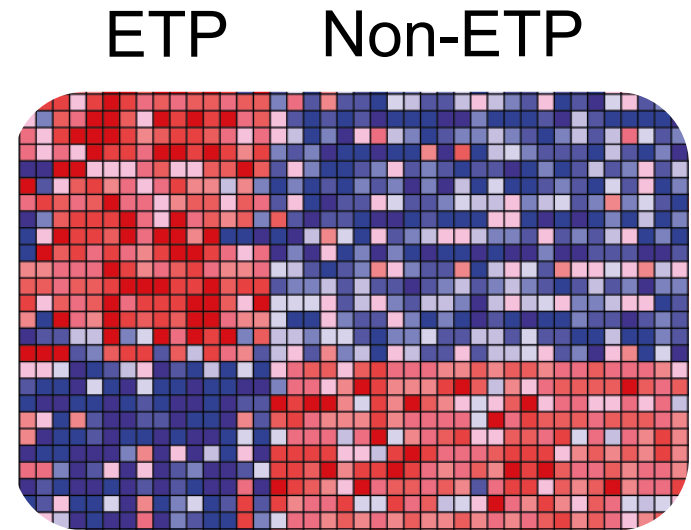
Genetic heterogeneity in human ETP-ALL



Genetic heterogeneity provides challenges for the implementation of a uniform targeted therapy in ETP-ALL

Unique features of human ETP-ALL

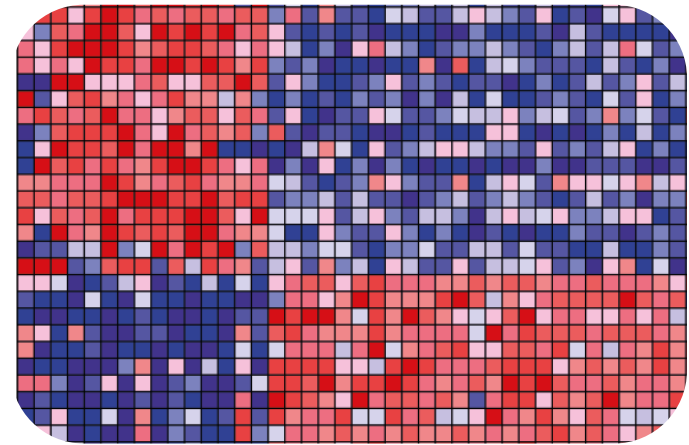
Human ETP-ALLs
have a unique
gene expression signature



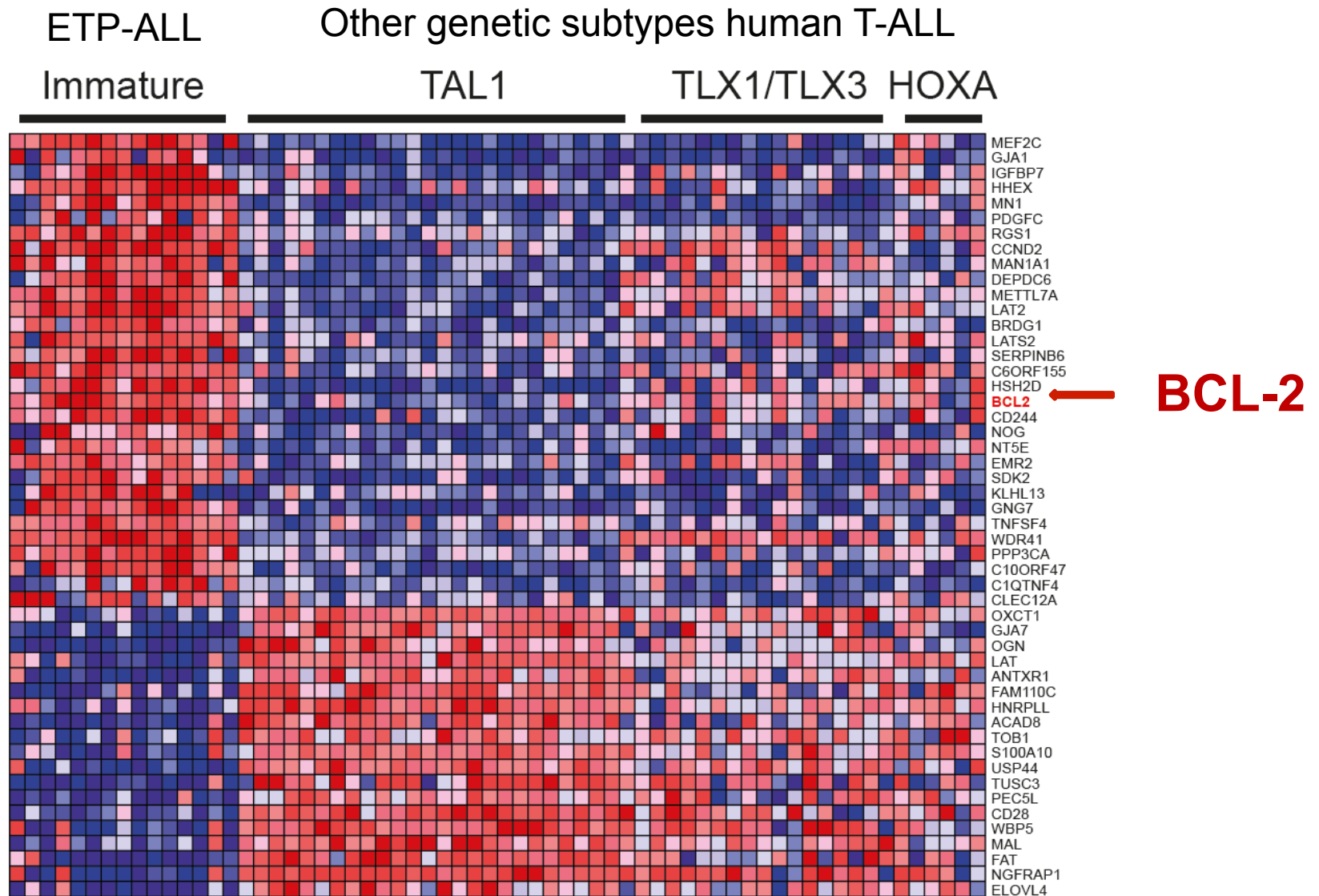
Therapeutic targets in human ETP-ALL gene expression profile



ETP Non-ETP

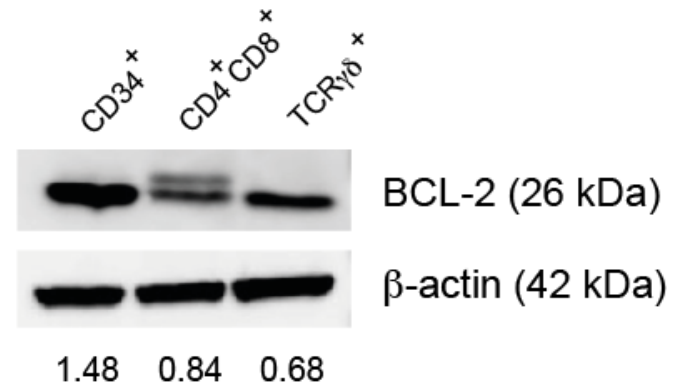
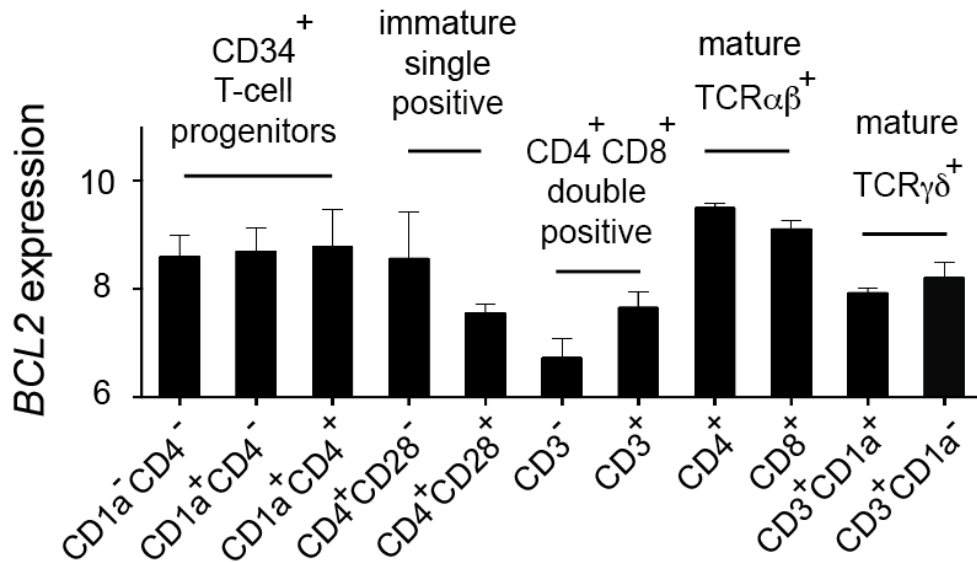
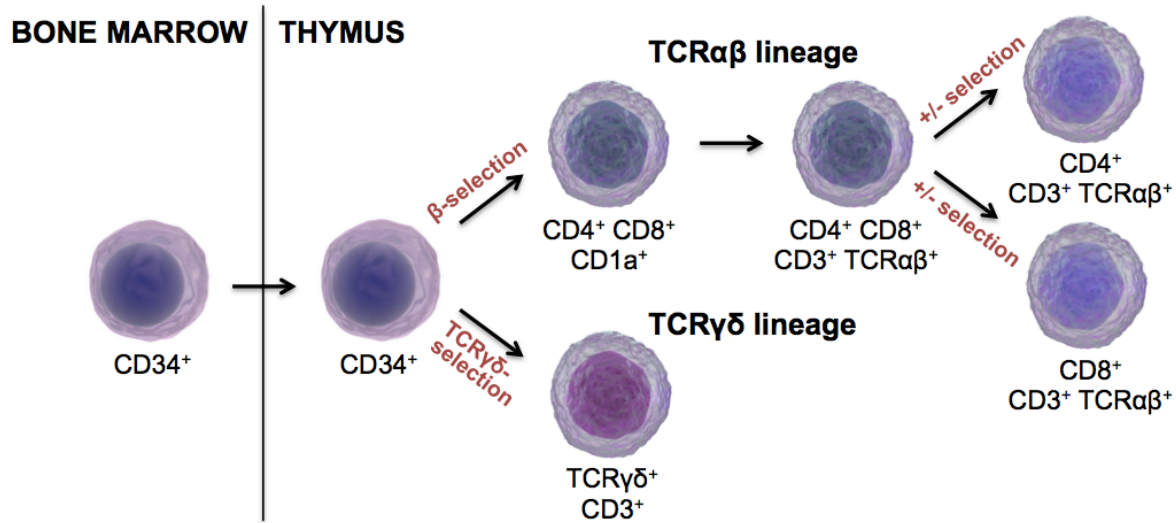


High *BCL2* expression in human ETP-ALL

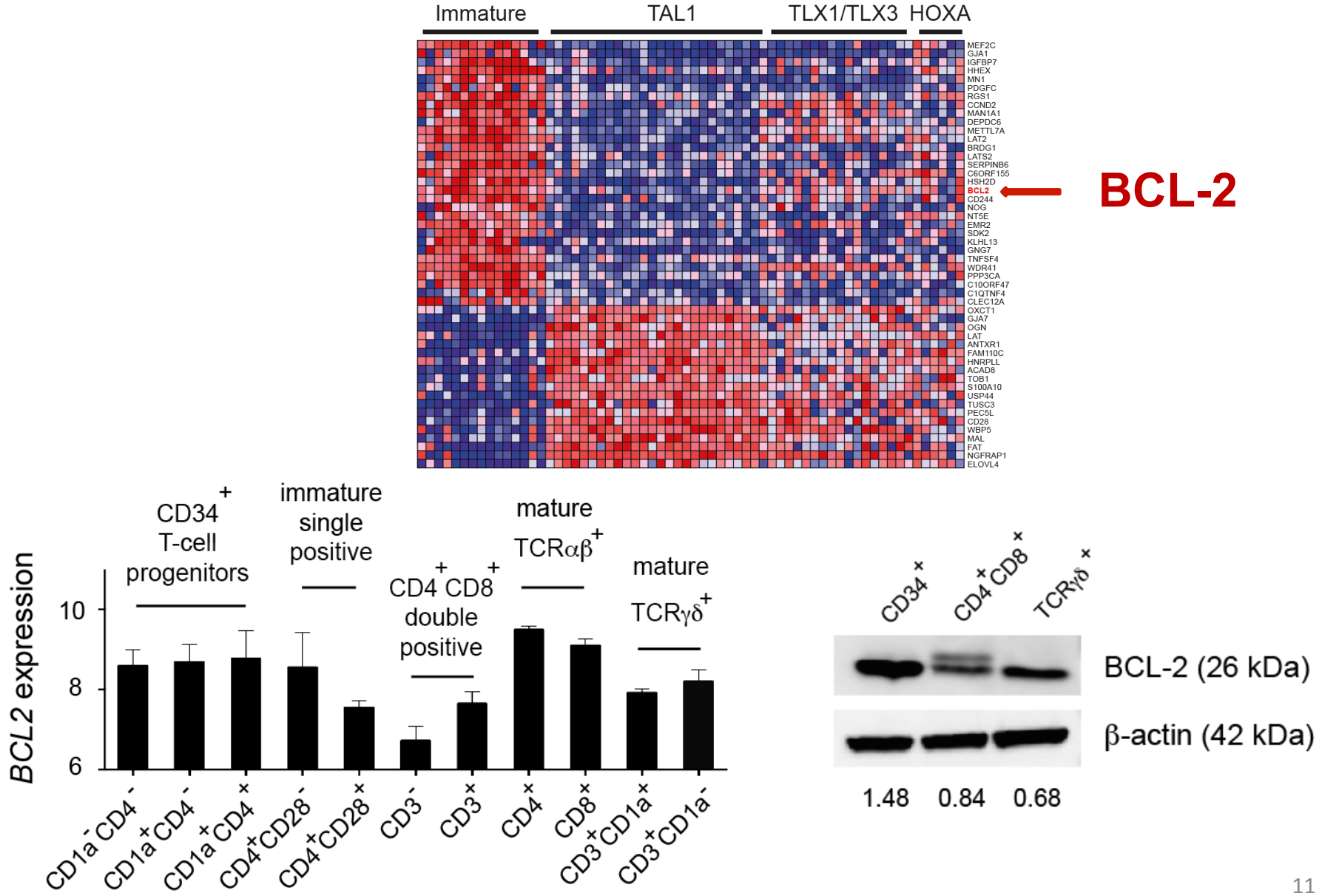


Fold change > 2; p<0.05

High *BCL2* expression a reflection of T cell differentiation?



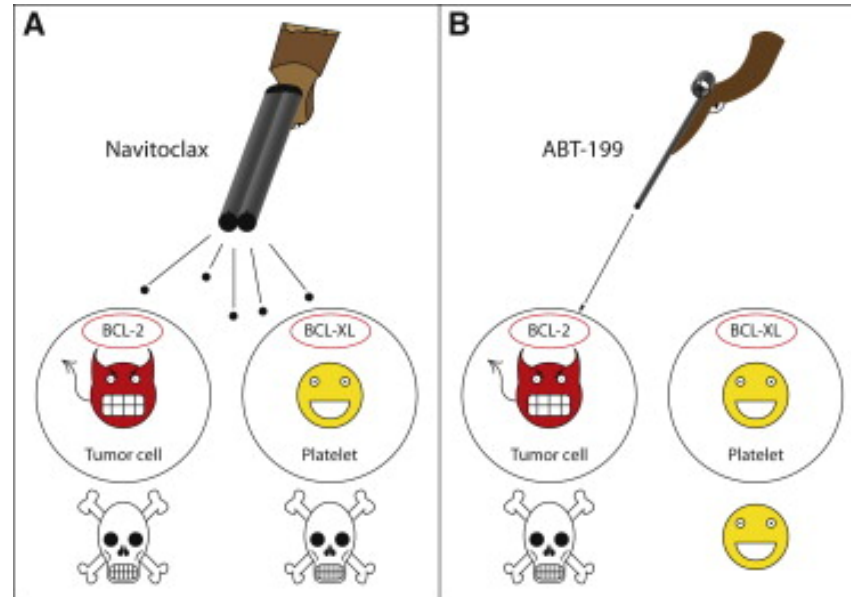
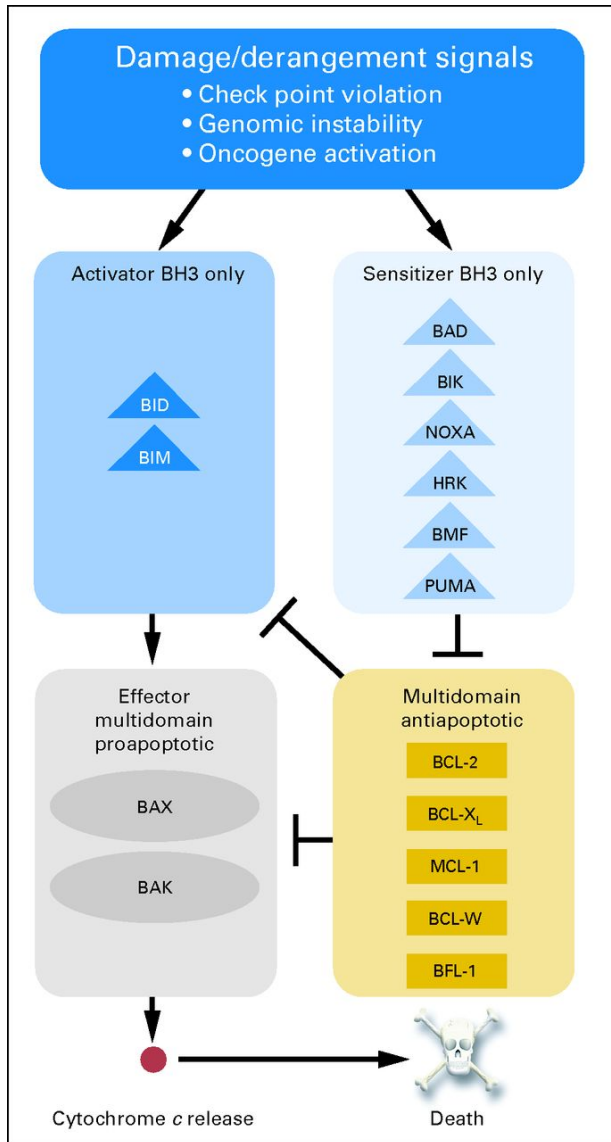
High *BCL2* expression a reflection of T cell differentiation?



BCL2 inhibition as a therapeutic strategy

abbvie

ABT-263 (Navitoclax) : BCL-2, BCL-xL
 ABT-199 : BCL-2 only



Dauids & Letai, Cancer Cell 2013

ABT-199 as a new therapeutic strategy in human cancer

**nature
medicine**

ABT-199, a potent and selective BCL-2 inhibitor, achieves antitumor activity while sparing platelets

January 2013

Andrew J Souers¹, Joel D Levenson¹, Erwin R Boghaert¹, Scott L Ackler¹, Nathaniel D Catron¹, Jun Chen¹, Brian D Dayton¹, Hong Ding¹, Sari H Enschede¹, Wayne J Fairbrother², David C S Huang^{3,4}, Sarah G Hymowitz², Sha Jin¹, Seong Lin Khaw^{3,4}, Peter J Kovar¹, Lloyd T Lam¹, Jackie Lee², Heather L Maecker², Kennan C Marsh¹, Kylie D Mason³⁻⁵, Michael J Mitten¹, Paul M Nimmer¹, Anatol Oleksijew¹, Chang H Park¹, Cheol-Min Park^{1,7}, Darren C Phillips¹, Andrew W Roberts³⁻⁵, Deepak Sampath², John F Seymour^{4,6}, Morey L Smith¹, Gerard M Sullivan¹, Stephen K Tahir¹, Chris Tse¹, Michael D Wendt¹, Yu Xiao¹, John C Xue¹, Haichao Zhang¹, Rod A Humerickhouse¹, Saul H Rosenberg¹ & Steven W Elmore¹

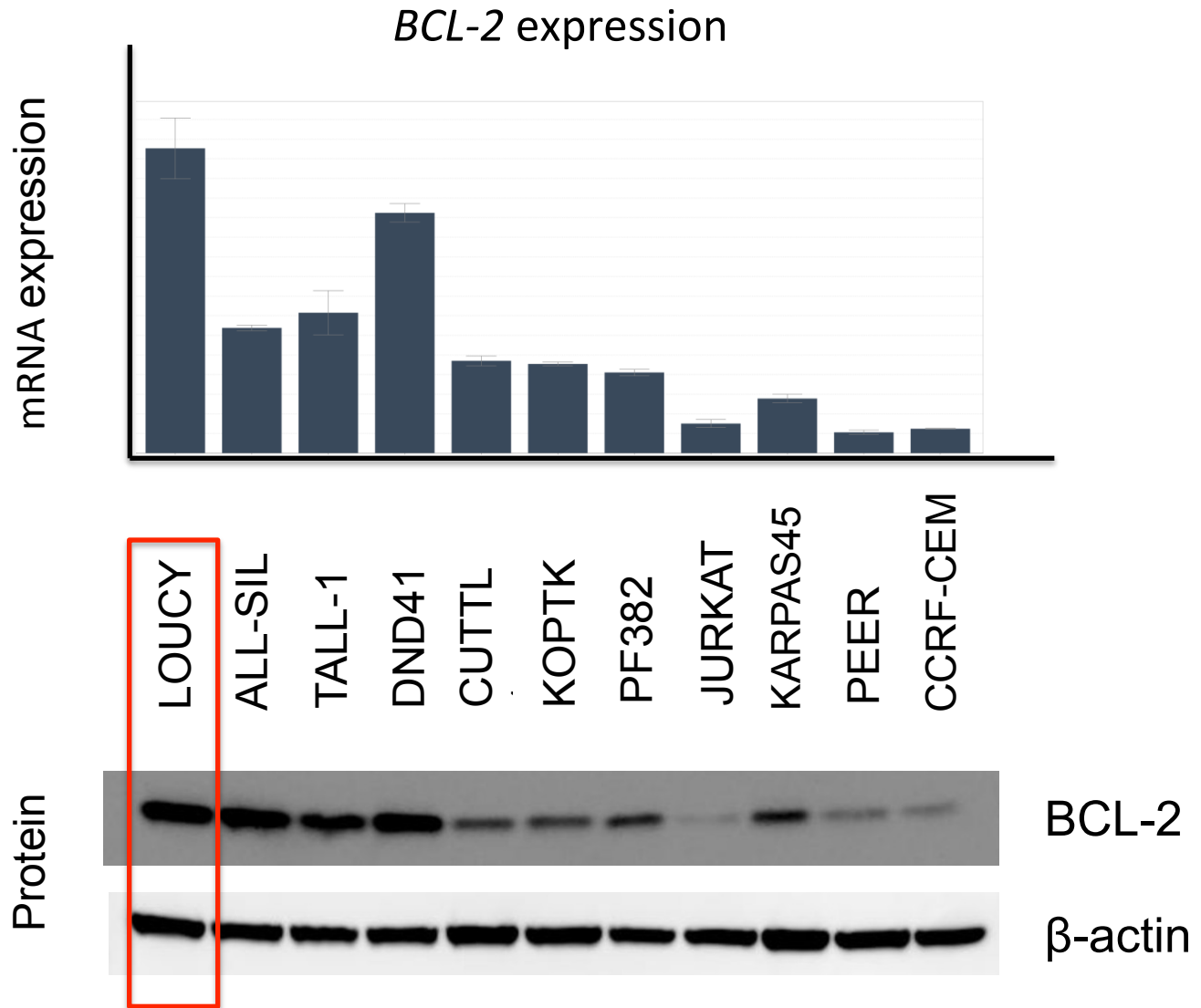
Clinical Phase I Trials in relapsed or treatment-resistant CLL

Tumor Lysis Syndrome

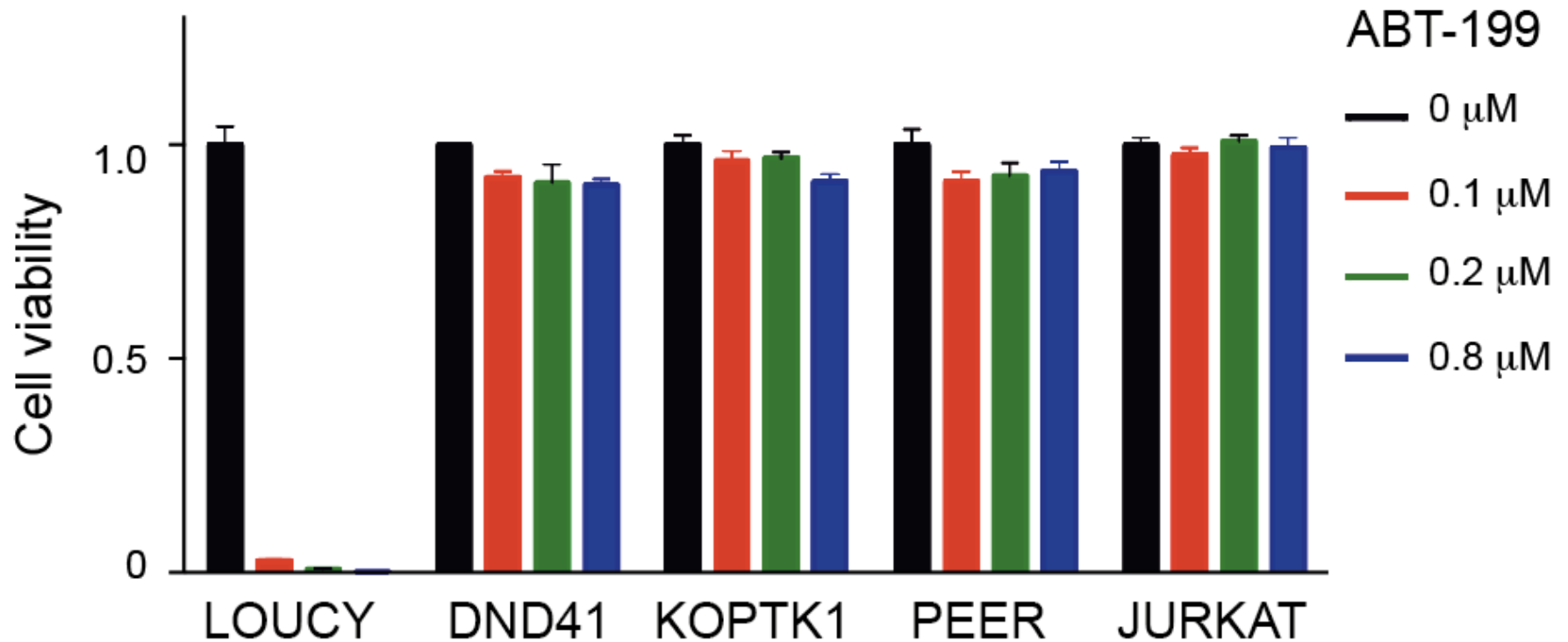
Reports on anti-tumoral effect in:

acute myeloid leukemia
multiple myeloma
estrogen receptor-positive breast cancer

ABT-199 expression in human T-ALL cell lines



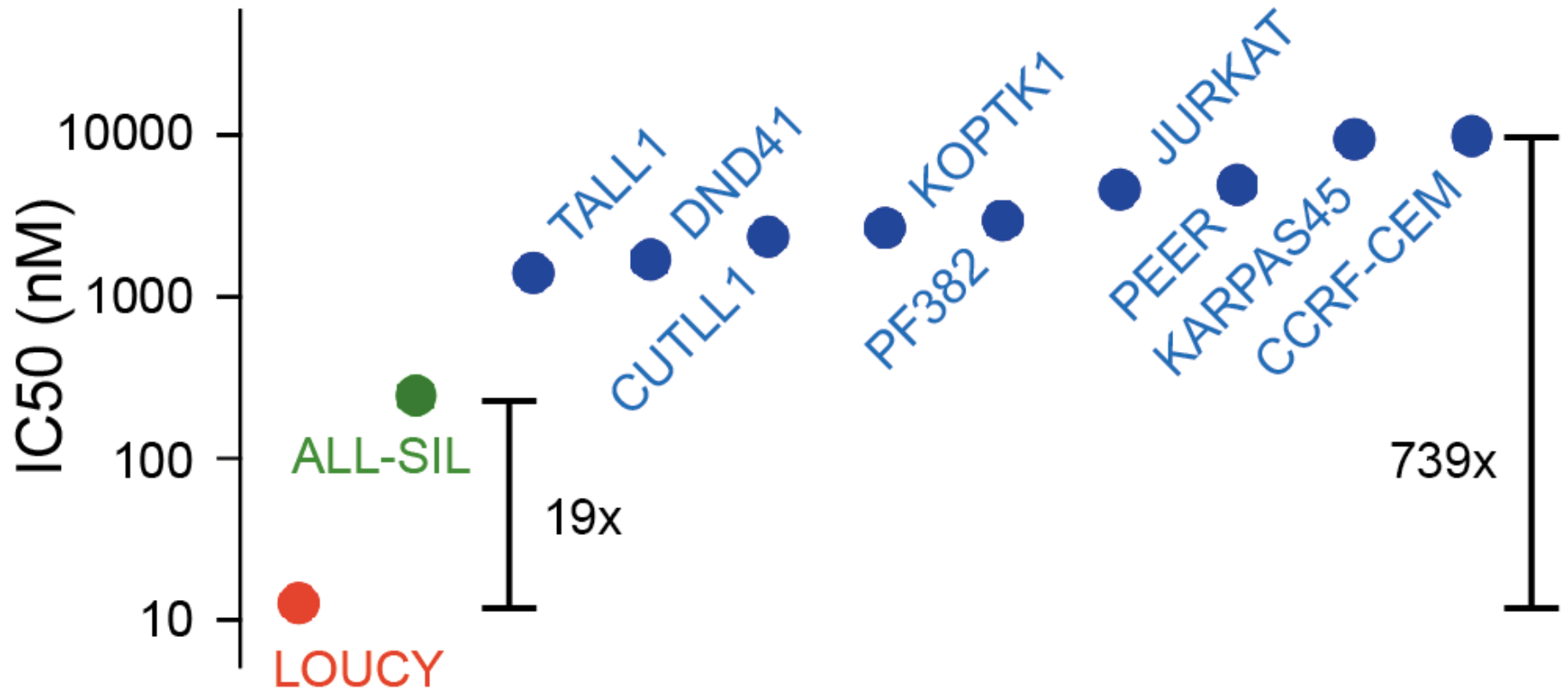
ABT-199 and cell viability in human T-ALL cell lines



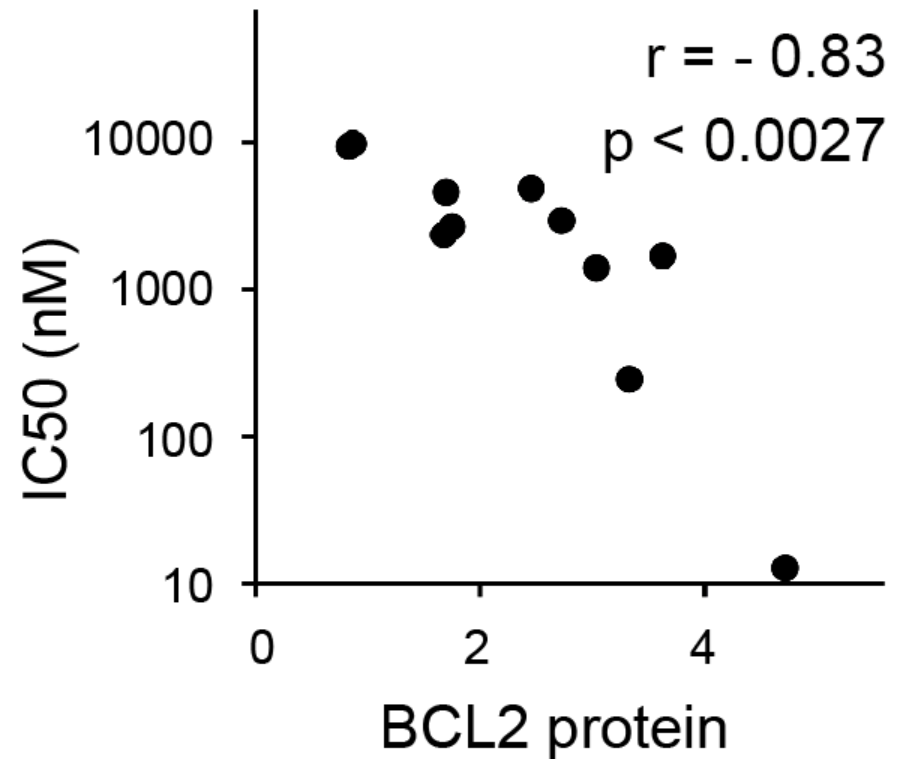
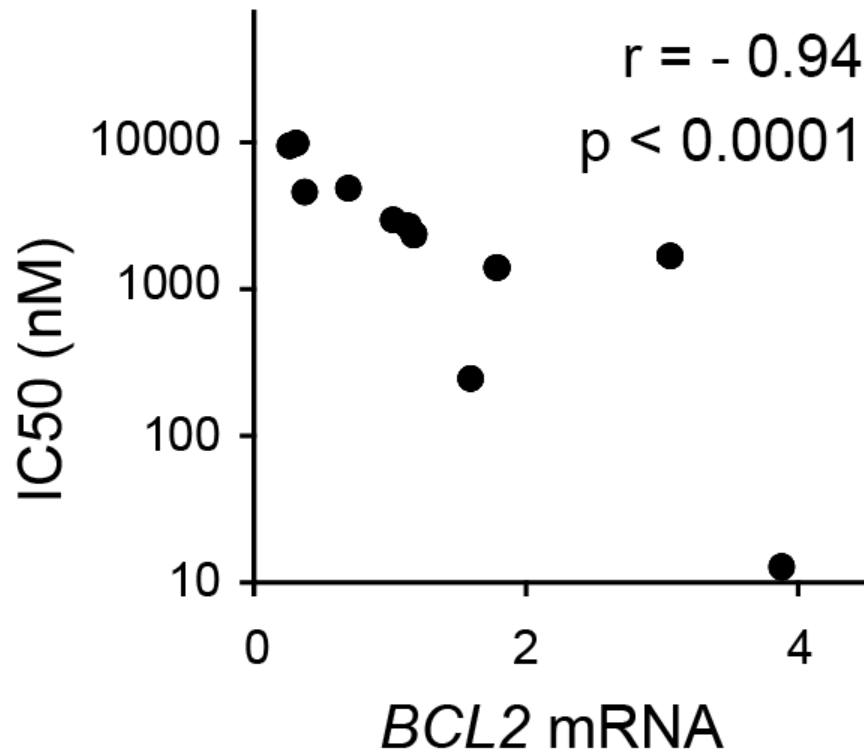
CellTiter-Glo® Luminescent Cell Viability

GloMax®-Multi Microplate Reader

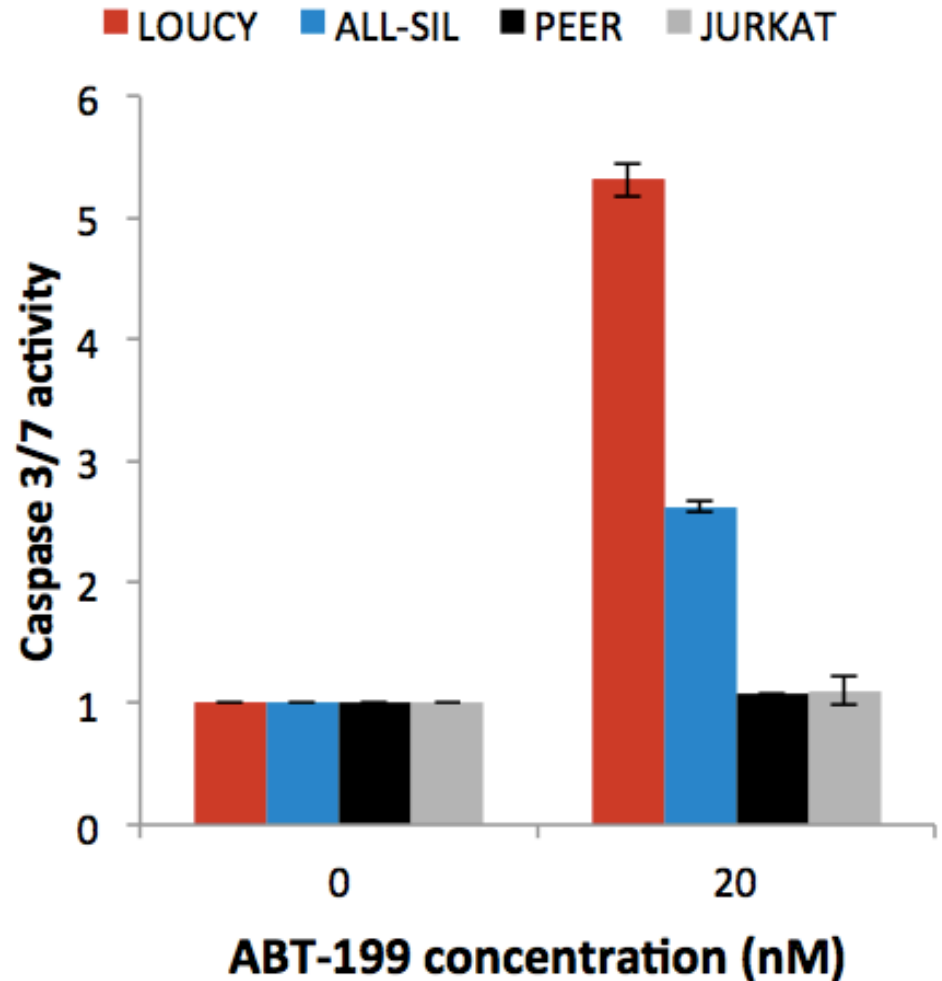
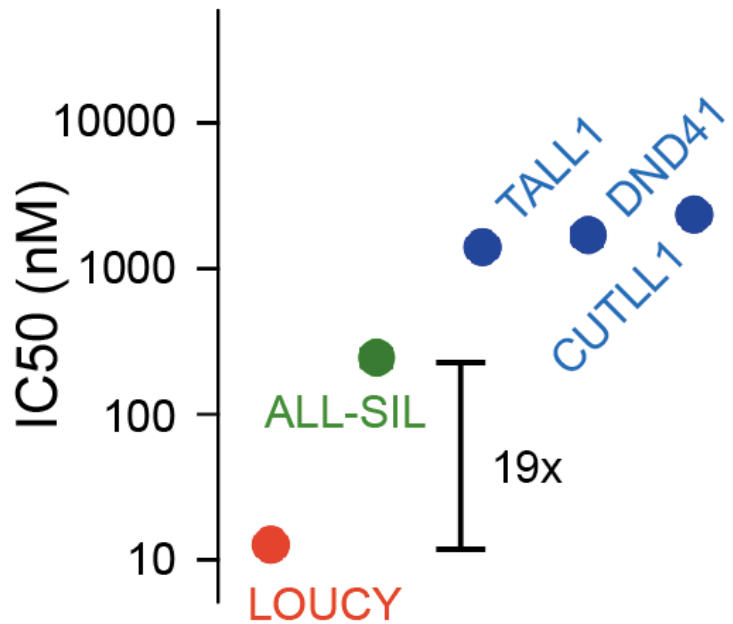
ABT-199 IC50 in human T-ALL cell lines



Correlation ABT-199 IC50 and BCL2 expression



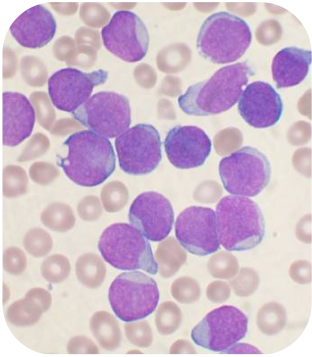
ABT-199 induces apoptosis



Caspase-Glo® 3/7 Apoptosis Assay
GloMax®-Multi Microplate Reader

In vivo xenograft model human T-ALL

Luciferase positive
LOUCY cells



Engraftment of cells
(6 weeks)

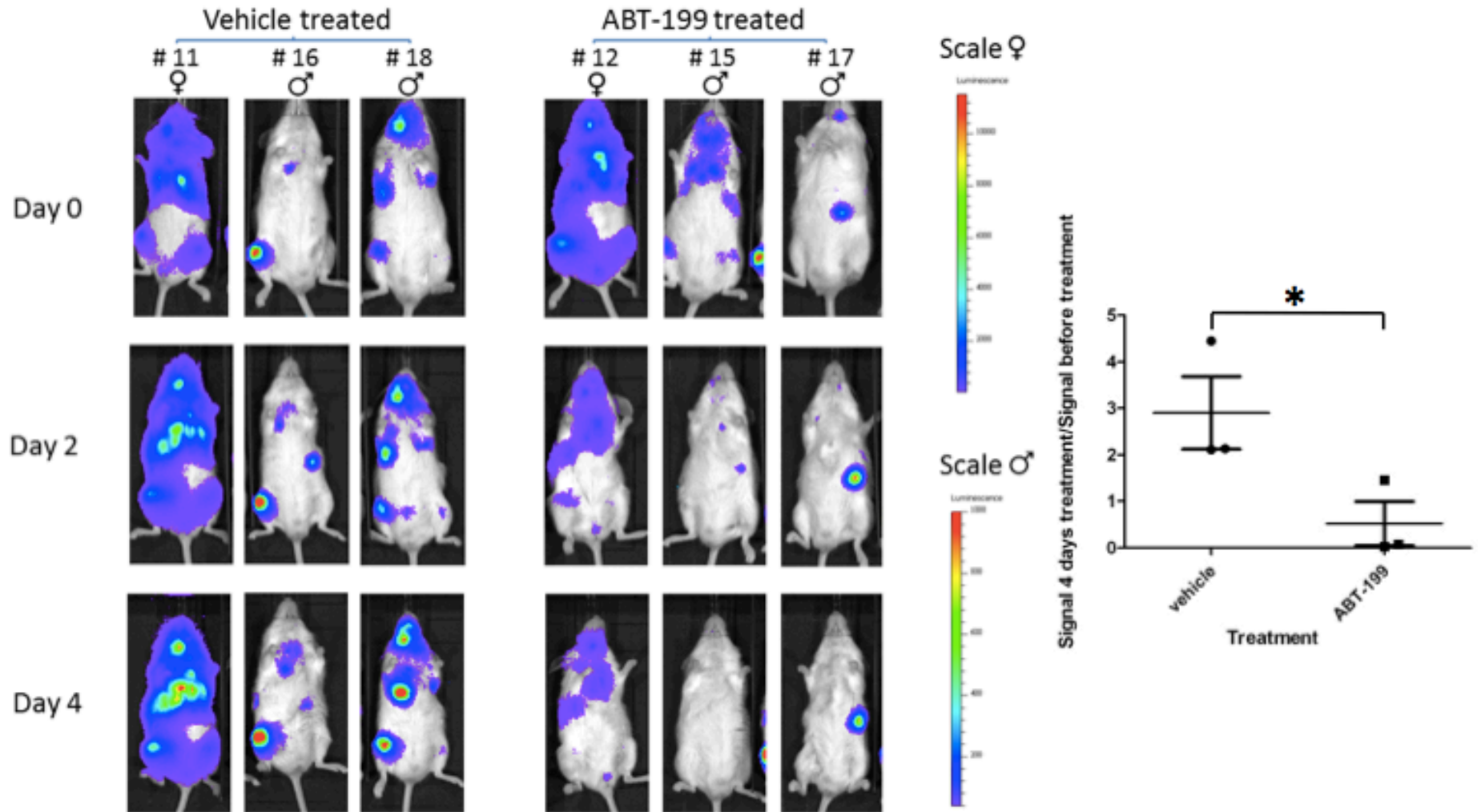


Measure luminescence on
different time points

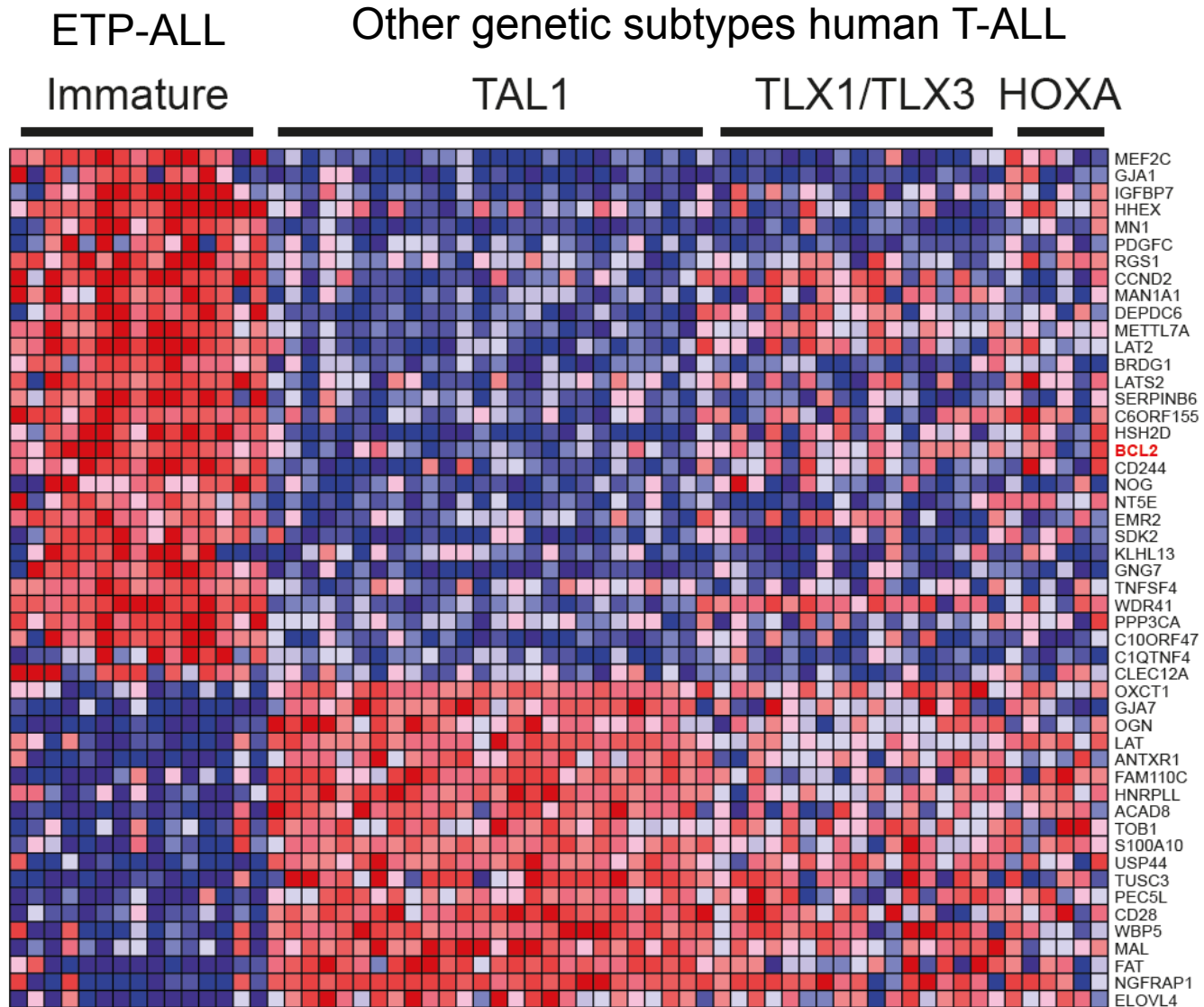


Oral gavage
100 mg ABT-199/kg
or vehicle

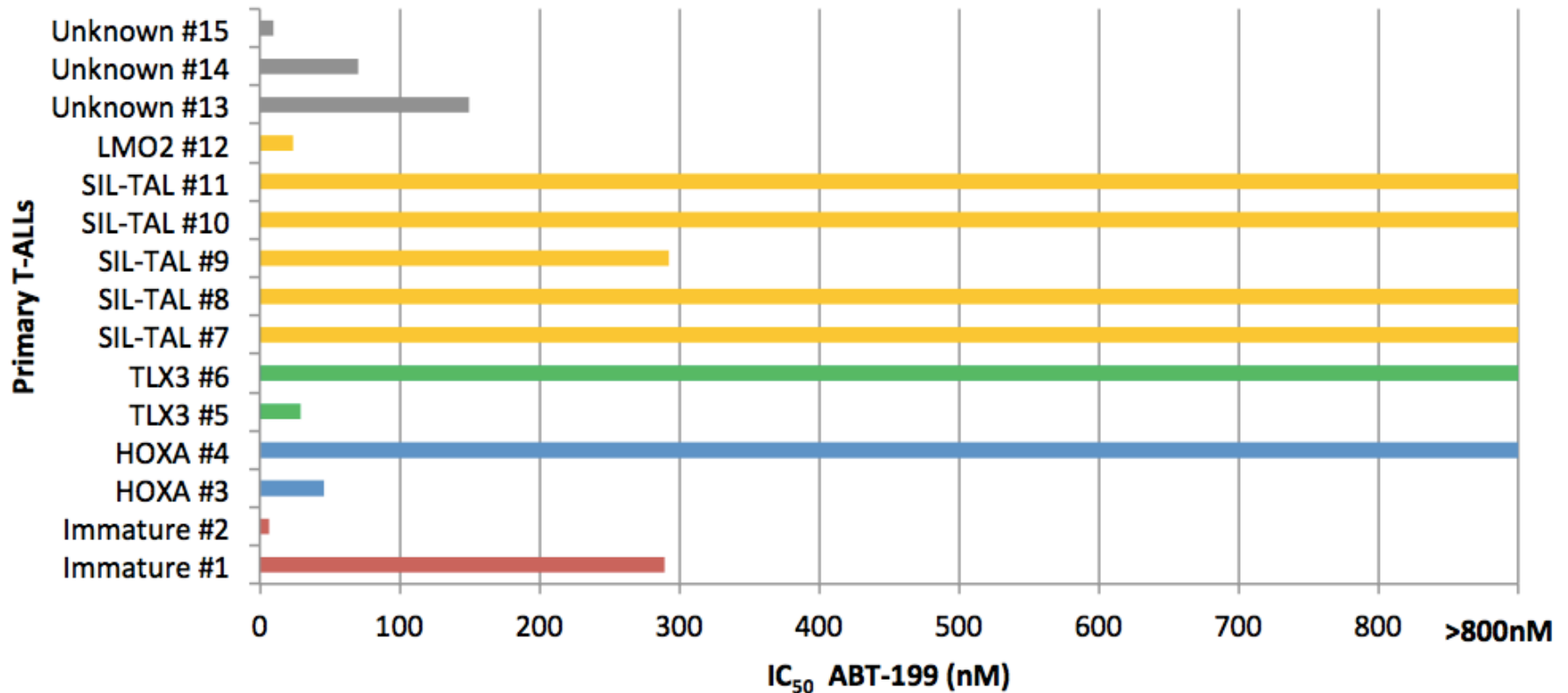
In vivo xenograft model human T-ALL



Genetic subtypes of human T-ALL

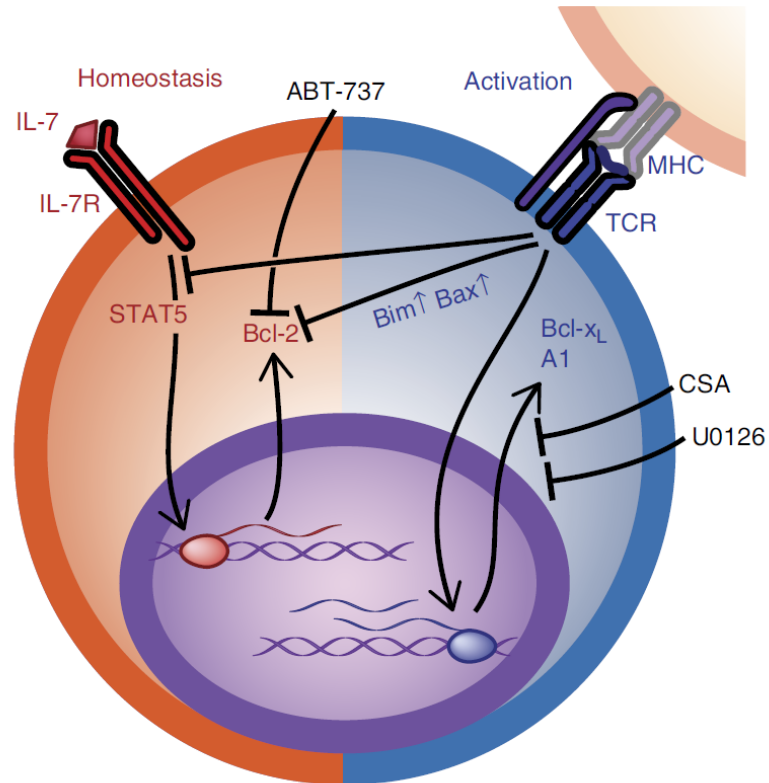
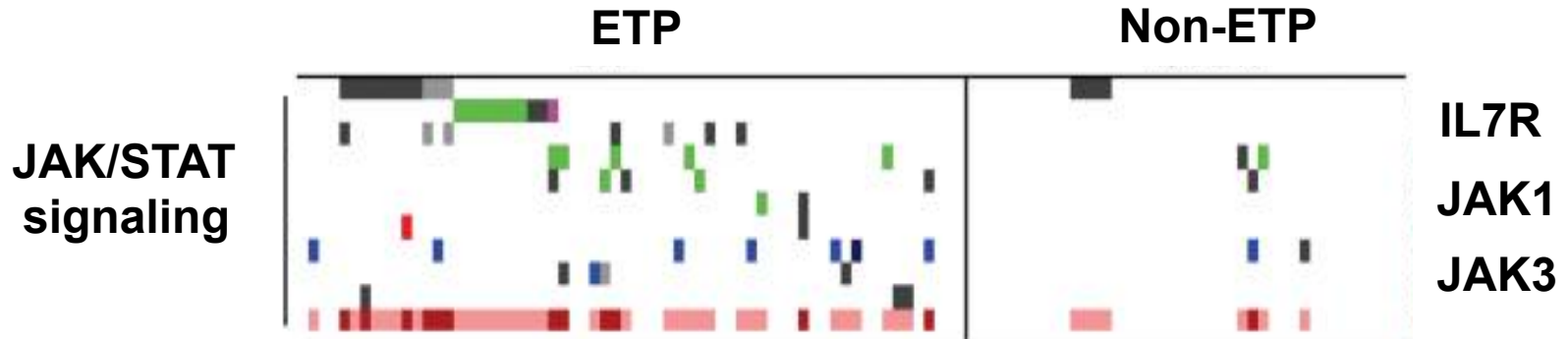


In vitro ABT-199 sensitivity primary T-ALL samples

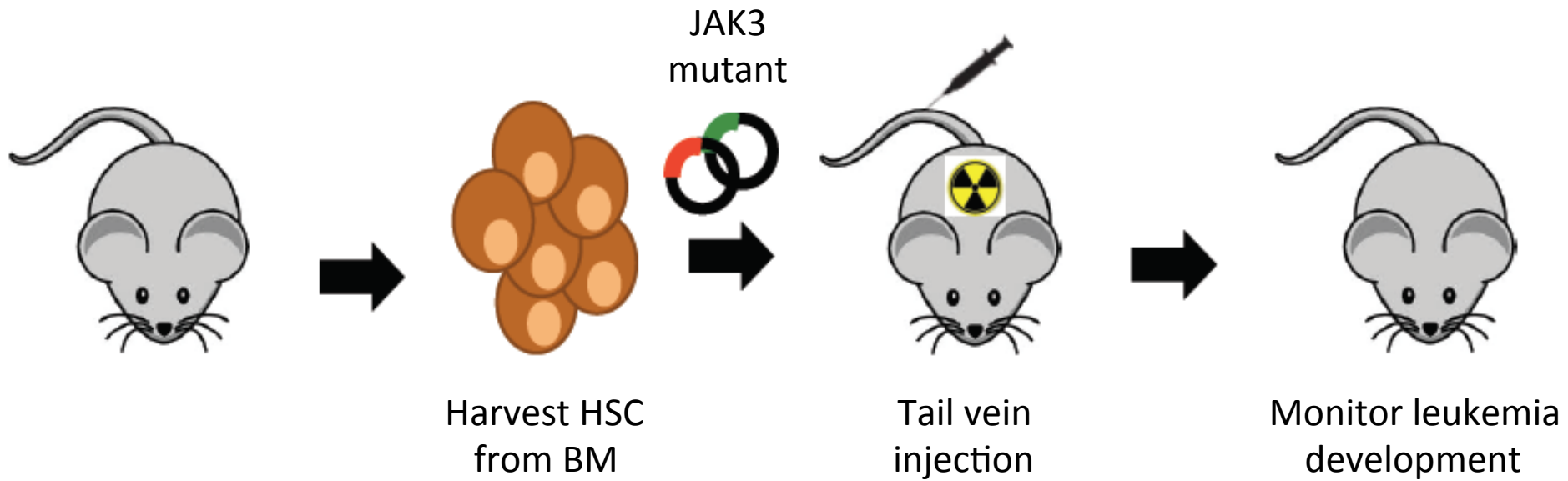


Hypothesis
The cell of origin and cooperative genetic defects define ABT-199 sensitivity in T-ALL

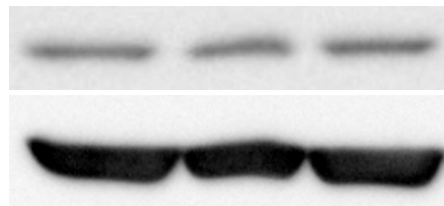
In vitro ABT-199 sensitivity human T-ALL



JAK3 mutant murine T-ALL



Control thymus

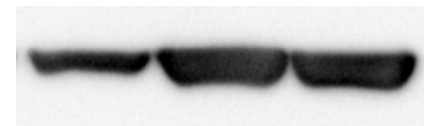


JAK3 mutant T-ALL tumors

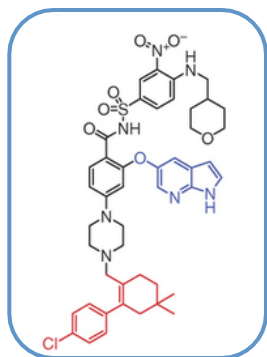
BCL-2



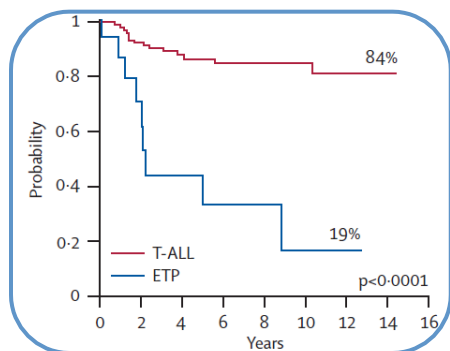
actin



Conclusions



ABT-199 is a promising new drug for human cancer



Our pre-clinical data suggest ABT-199 as a promising new drug for specific subtypes of pediatric T-ALL

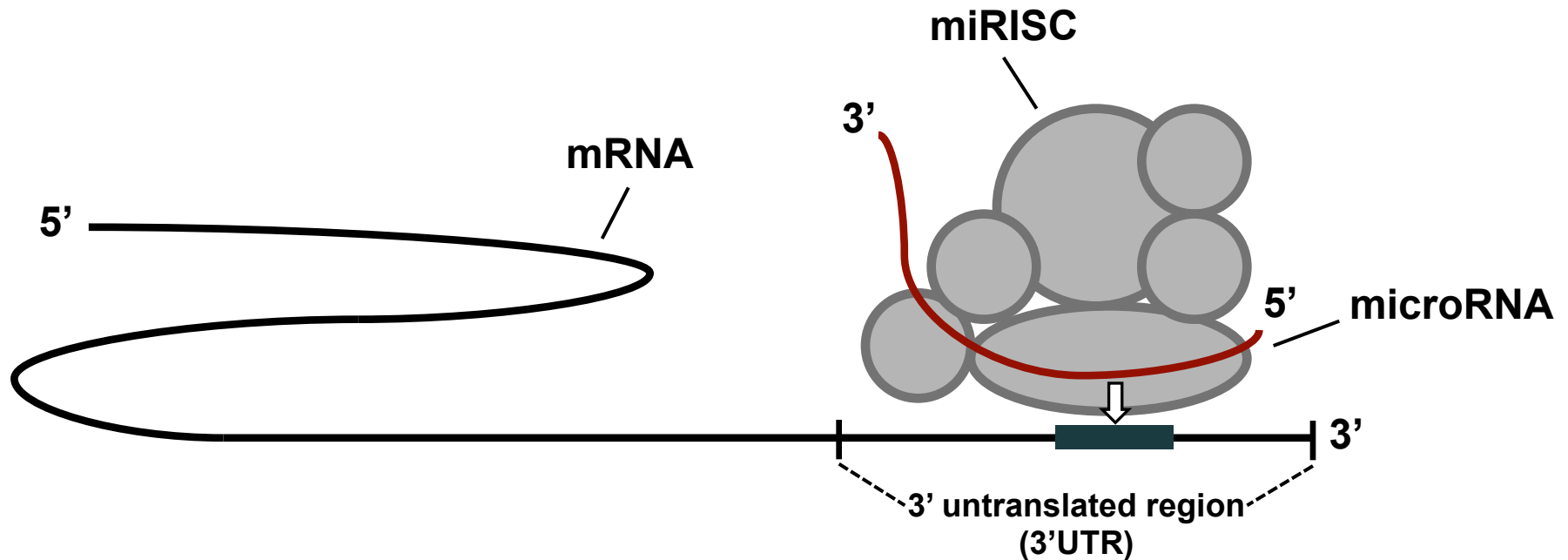
An unbiased high-throughput microRNA library screen identifies novel microRNA regulators of key oncogenes and tumor suppressor genes

Pieter Van Vlierberghe

Bioluminescent Cell-based Assay Seminar Day
Monday 31 march 2014
UCL De Duve institute
Brussels, Belgium

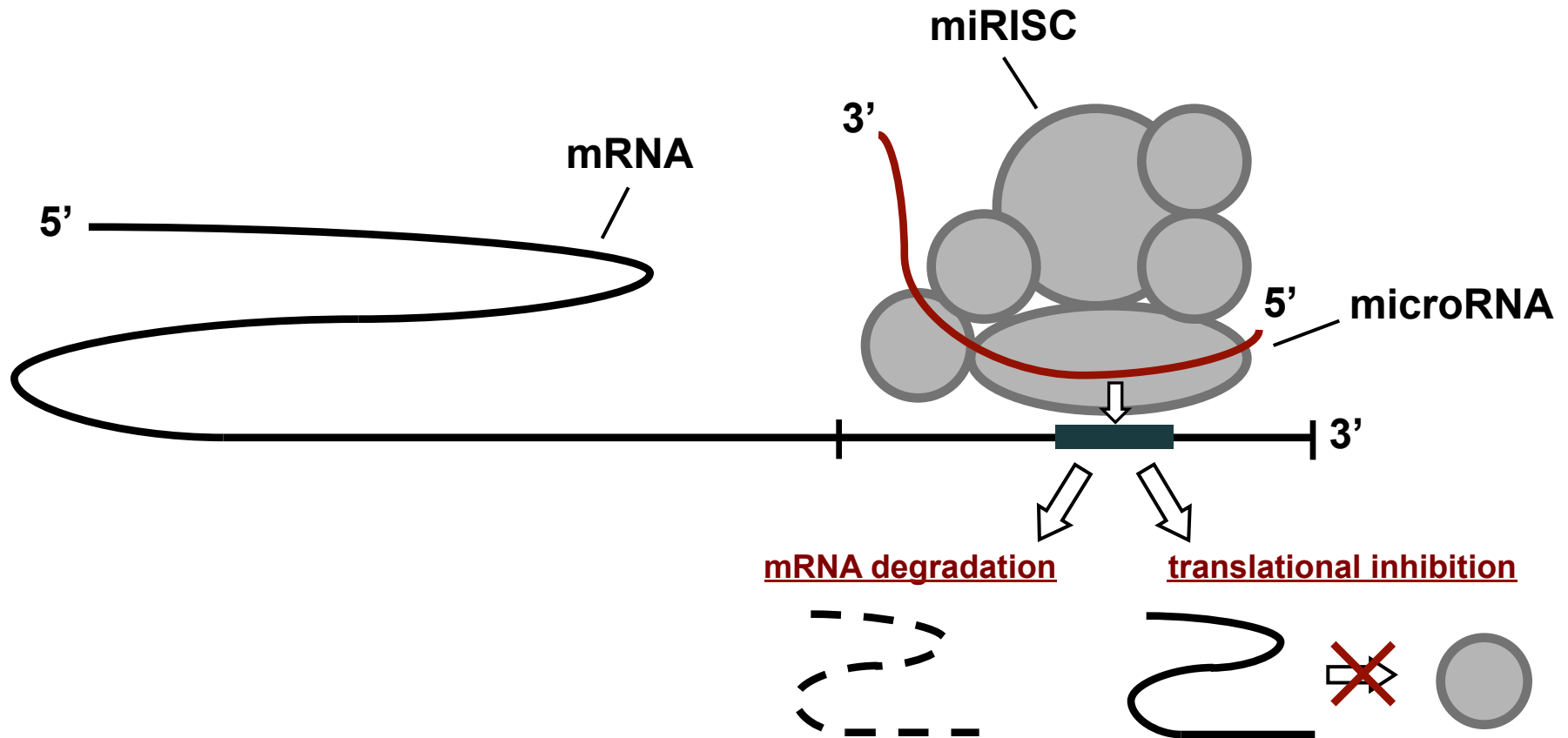
microRNAs

- small, non-coding RNA molecules
- post-transcriptional regulation of gene expression



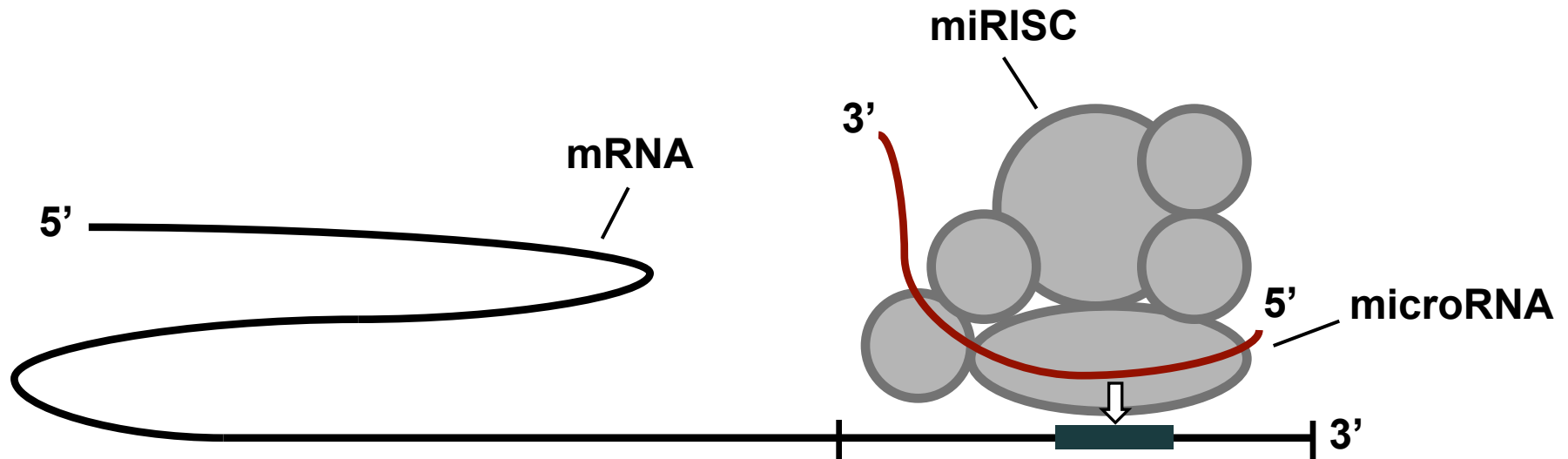
microRNAs

- small, non-coding RNA molecules
- post-transcriptional regulation of gene expression



microRNAs

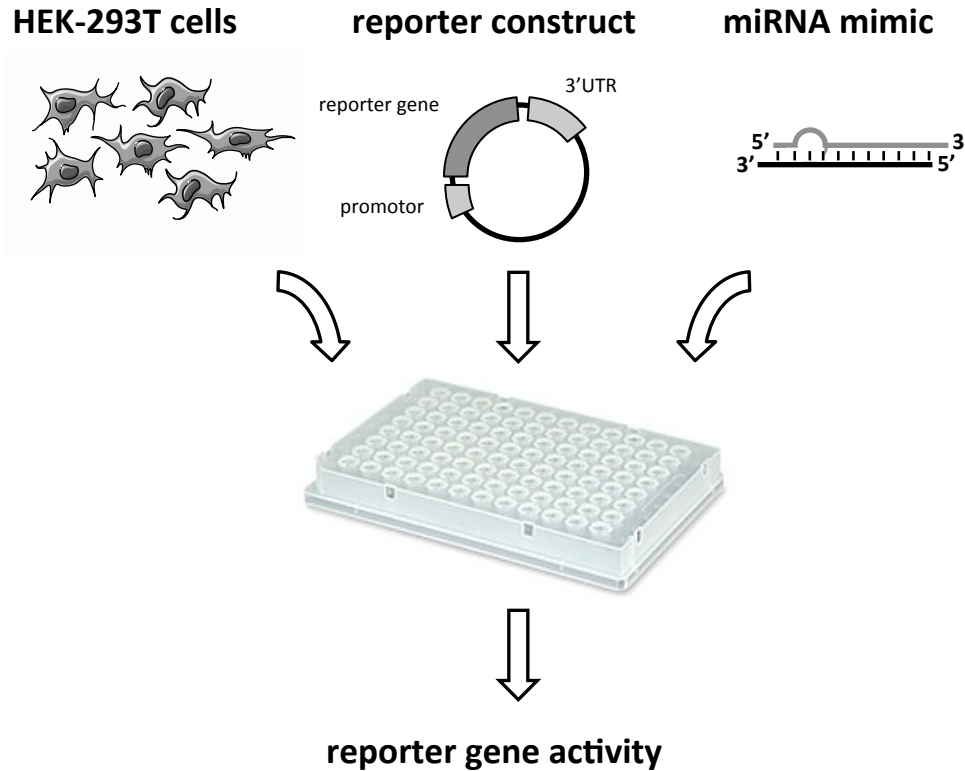
- small, non-coding RNA molecules
- post-transcriptional regulation of gene expression



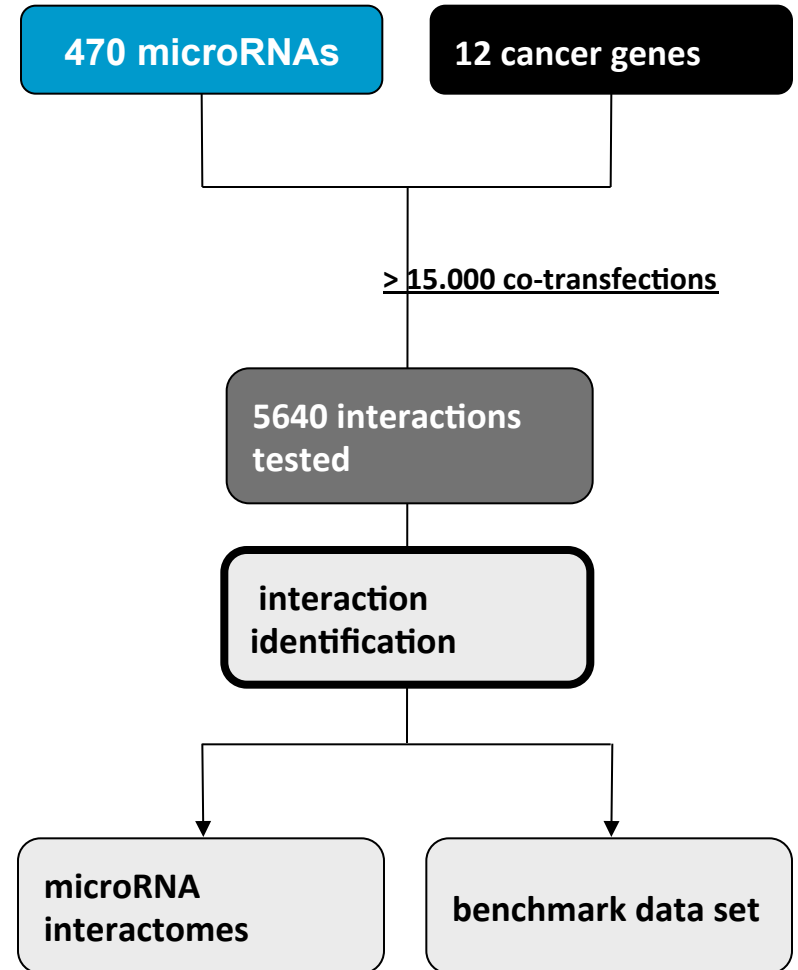
- microRNA function? microRNA targets!
- **microRNA target prediction algorithms!**

3'UTR microRNA library screen

3'UTR reporter assay



3'UTR microRNA library screen



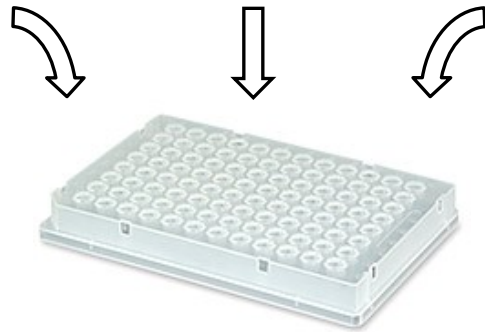
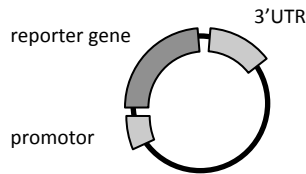
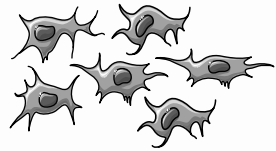
3'UTR microRNA library screen

3'UTR reporter assay

HEK-293T cells

reporter construct

miRNA mimic



reporter gene activity

microRNA interaction score

$$\Delta(\text{with } \frac{r_n - \text{median}_{r_n}}{MAD})$$

$$r_n = \log \frac{r_{\text{exp}}}{r_{\text{norm}}}$$

3'UTR microRNA library screen

470 microRNAs

12 cancer genes

> 15,000 co-transfections

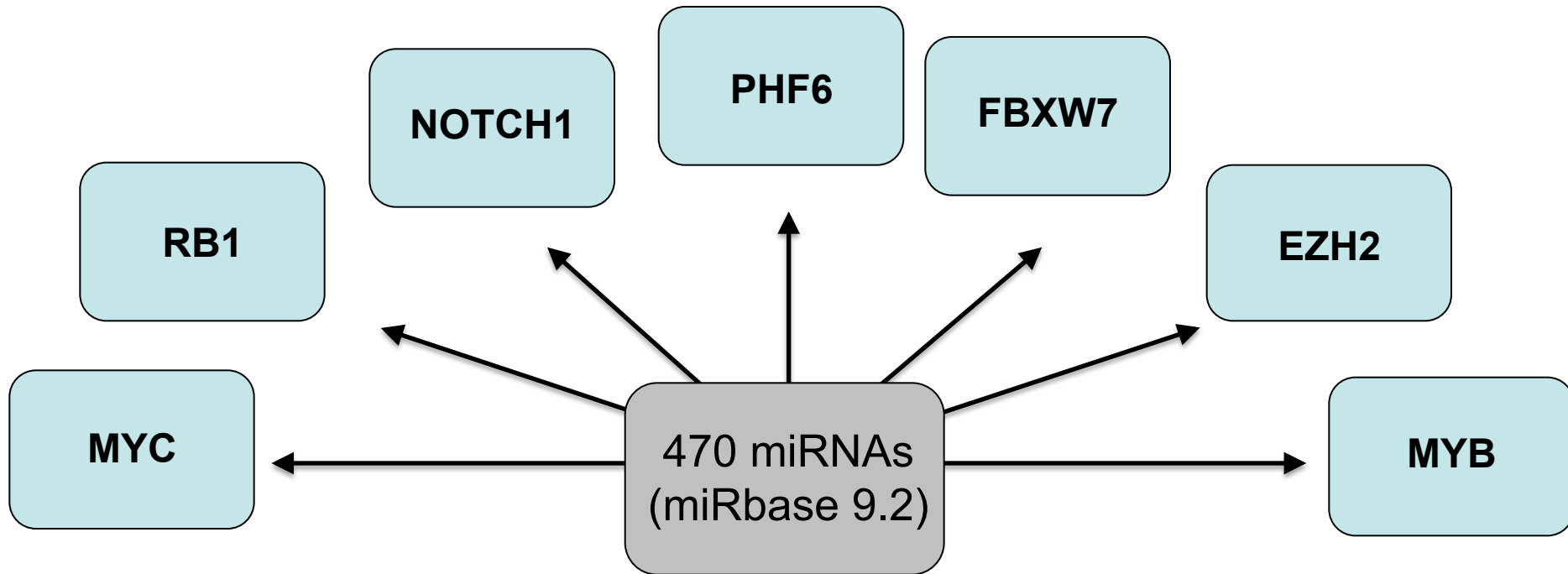
5640 interactions tested

interaction identification

microRNA interactomes

benchmark data set

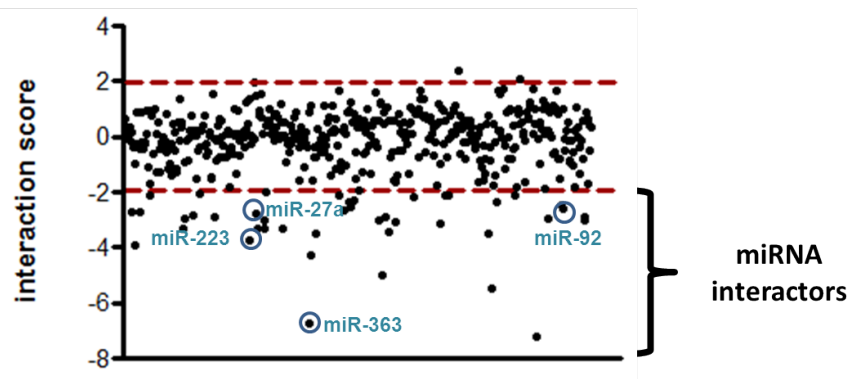
3'UTR library screen



Cancer gene microRNA interactomes

miRNA-FBXW7 interactome

← 470 miRNAs →



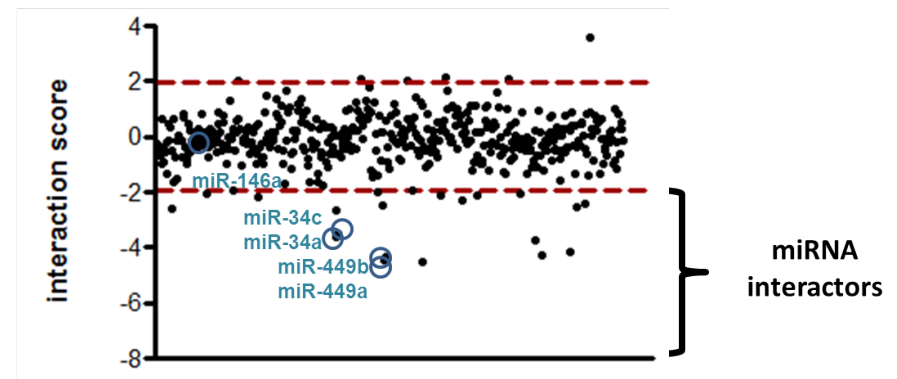
44 interactions identified

4 confirmed (already described in literature)

40 novel interactions (not yet described)

miRNA-NOTCH1 interactome

← 470 miRNAs →

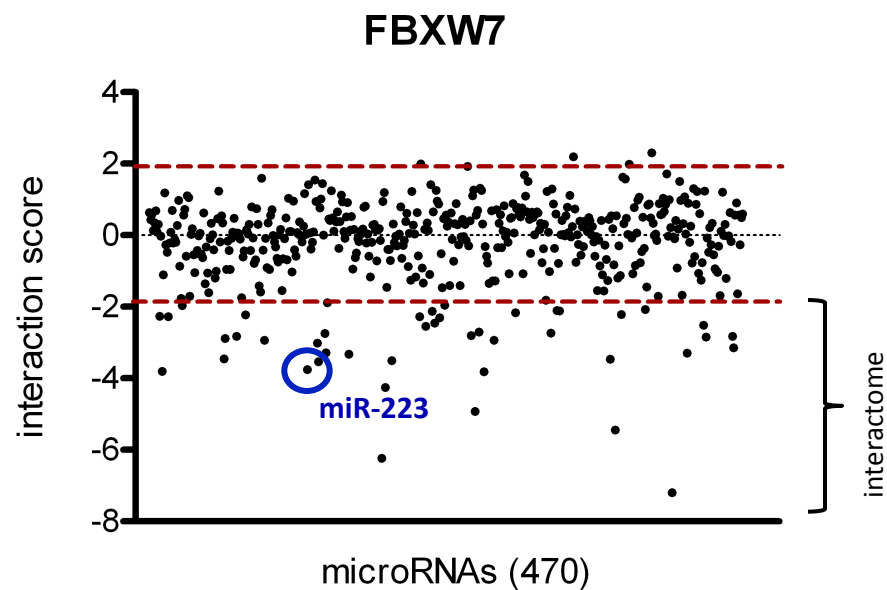
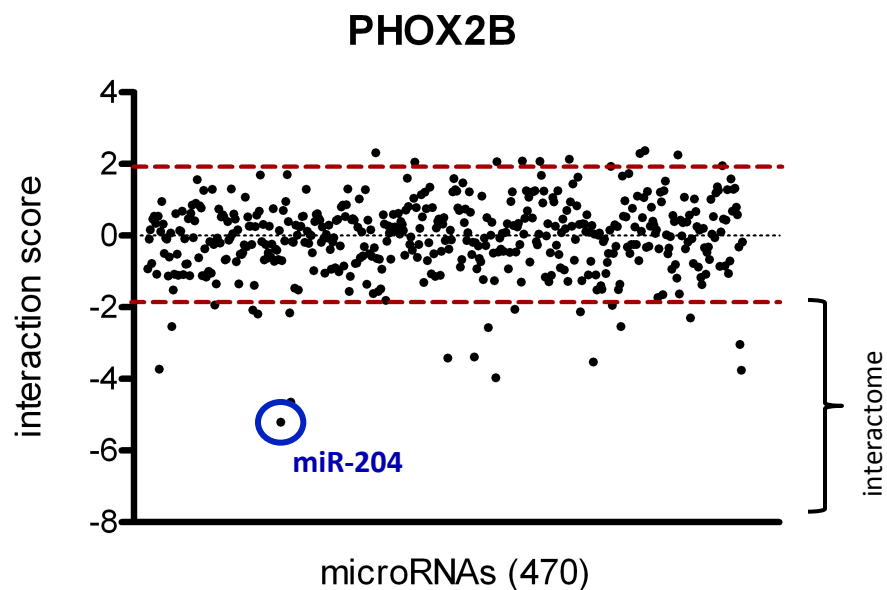


21 interactions identified

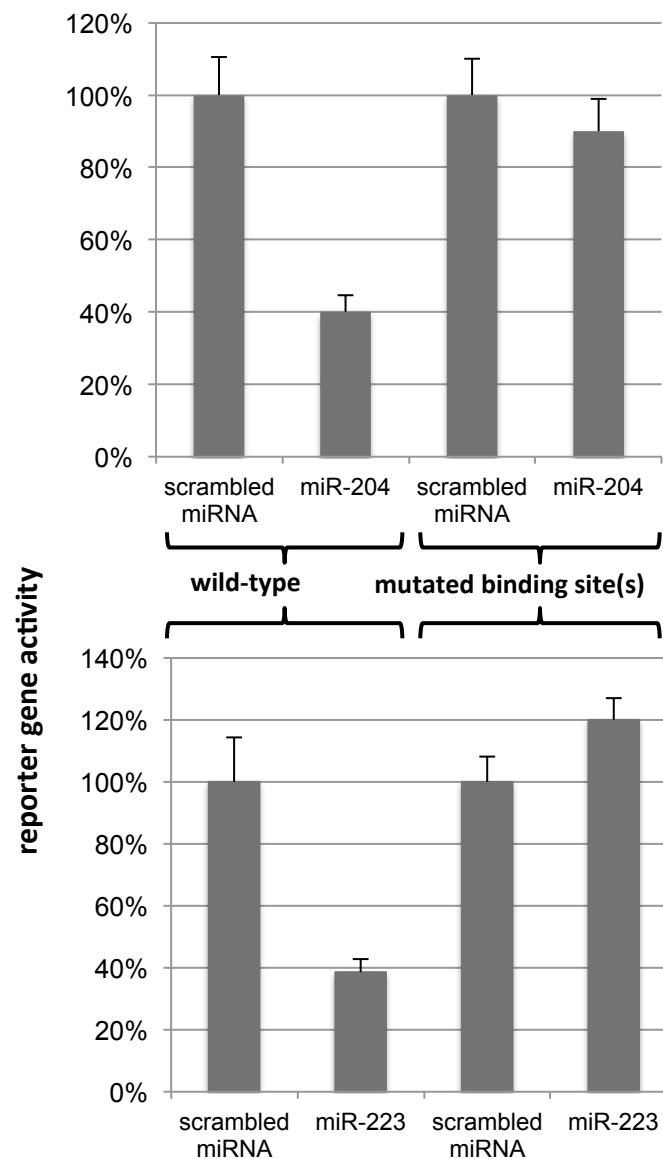
4 confirmed (already described in literature)

16 novel interactions (not yet described)

Cancer gene microRNA interactomes



rescue experiments!



Acknowledgements

Center for Medical Genetics (Ghent)

Sofie Peirs

Frank Speleman

Filip Matthijssens

Béatrice Lintermans

Evelien Van den Eeckhoudt

Department of Pediatric Hemato- Oncology (Ghent)

Tim Lammens

Barbara De Moerloose

Yves Benoit

Department of Clinical chemistry, microbiology and immunology (Ghent)

Tom Taghon

Inge Van de Walle

VIB Inflammation Research Center (Ghent)

Steven Goossens

Lab Jules Meijerink (Rotterdam)

Jules Meijerink

Kirsten Canté-Barrett

Lab Jean Soulier (Paris)

Jean Soulier

