

# A Role for ABC Transporters and Porphyrin Metabolism Leukemia ?

John D. Schuetz

St. Jude Children's Research Hospital

Memphis, TN 38105

September 2019

Solvo Biotechnology Transporter Conference

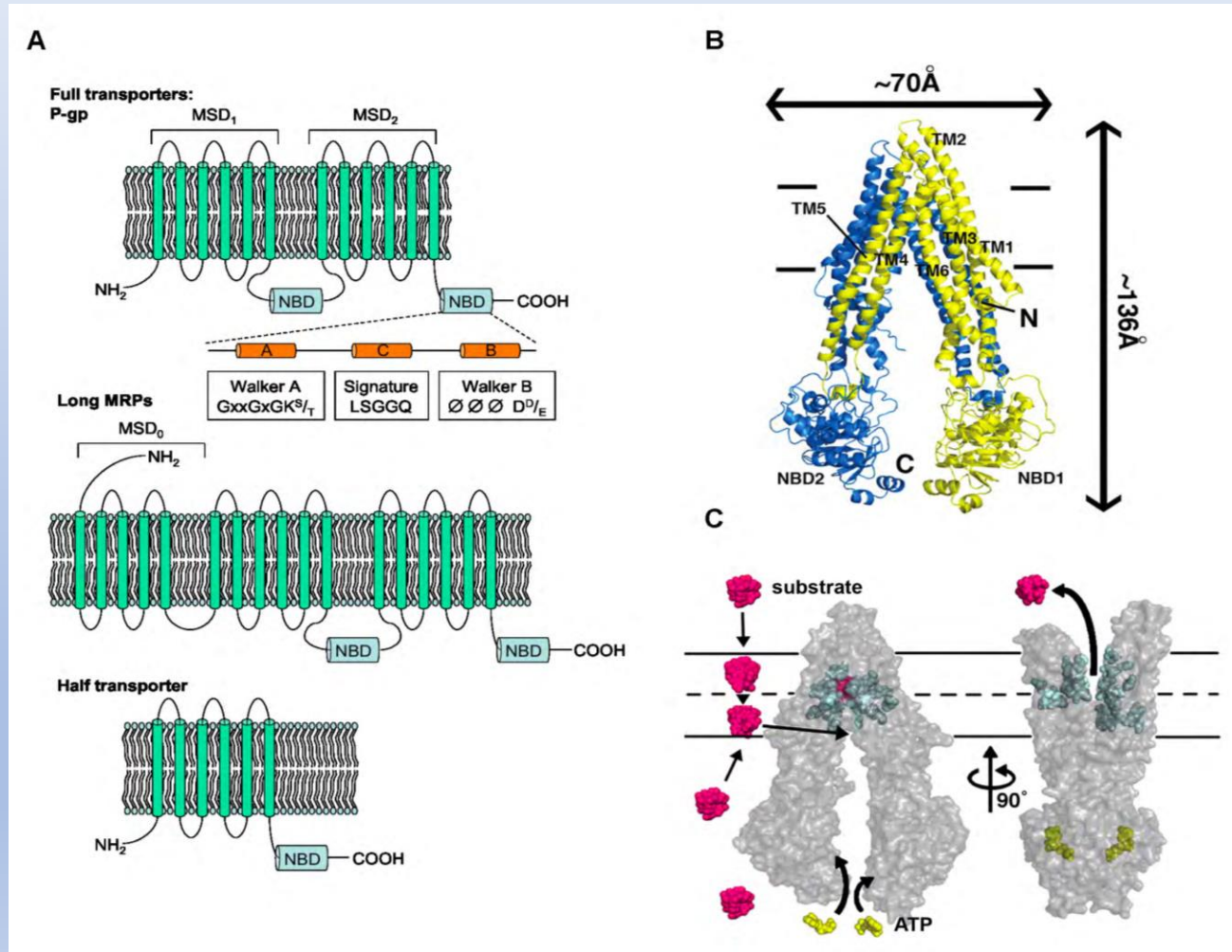
Cambridge, MA



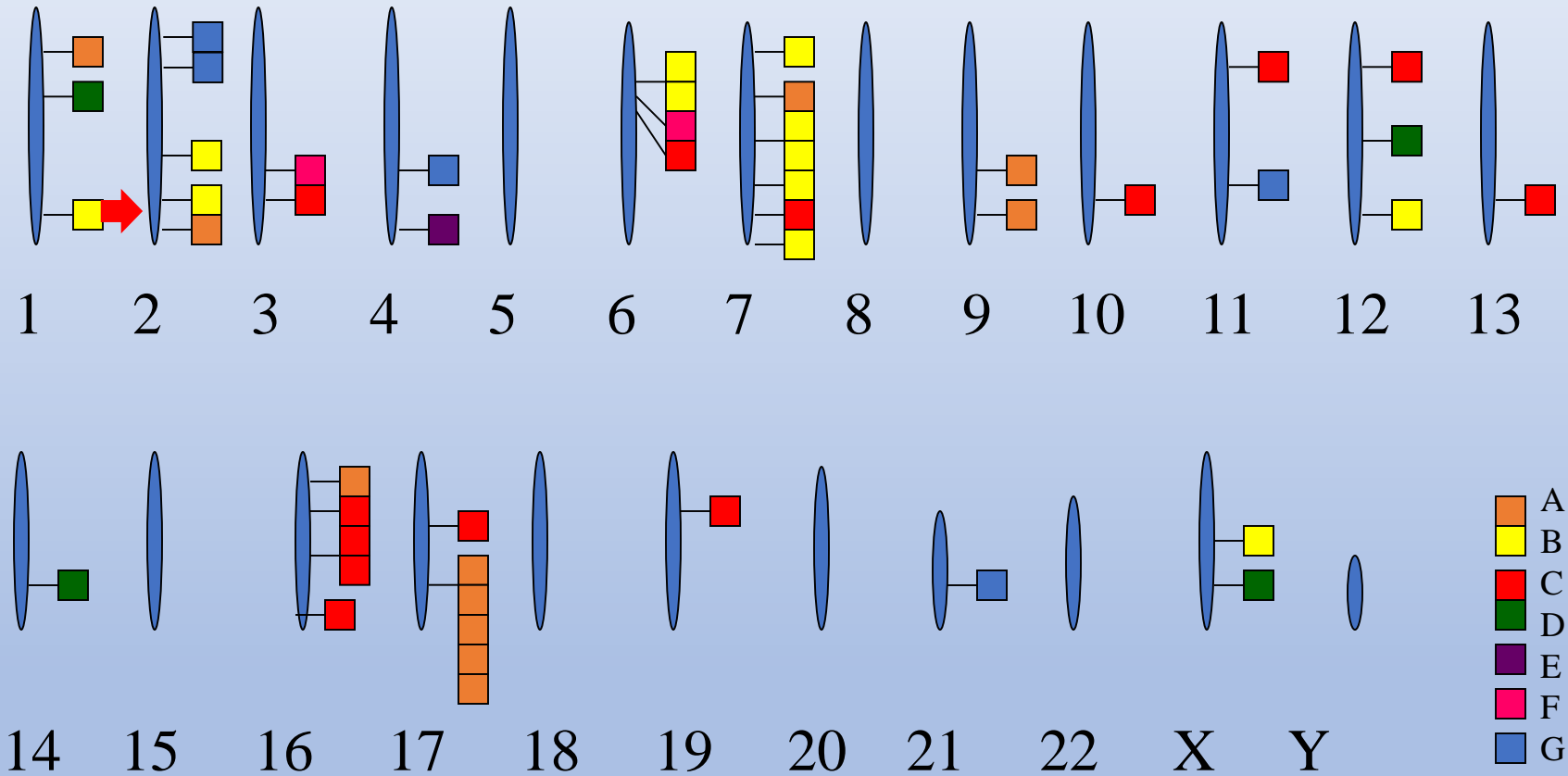
# Outline

- Intro-ABC transporters
- ABCG2 in acute myeloid leukemia (AML)
- Heme/porphyrin synthesis
- PPIX mechanism of death
- Identification of dual function ABCG2 inhibitors

# ABC transporters have two primary functional domains



# 48 Human ABC Genes



# Acute myeloid leukemia (AML) is a disease of altered hematopoiesis in myeloid lineage

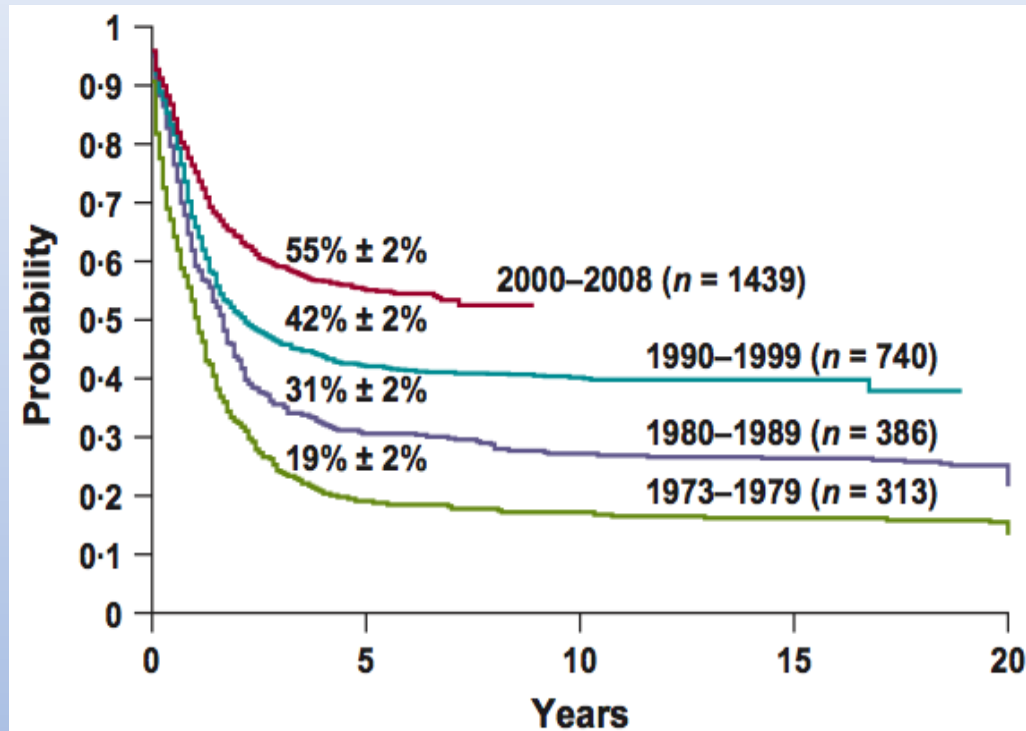
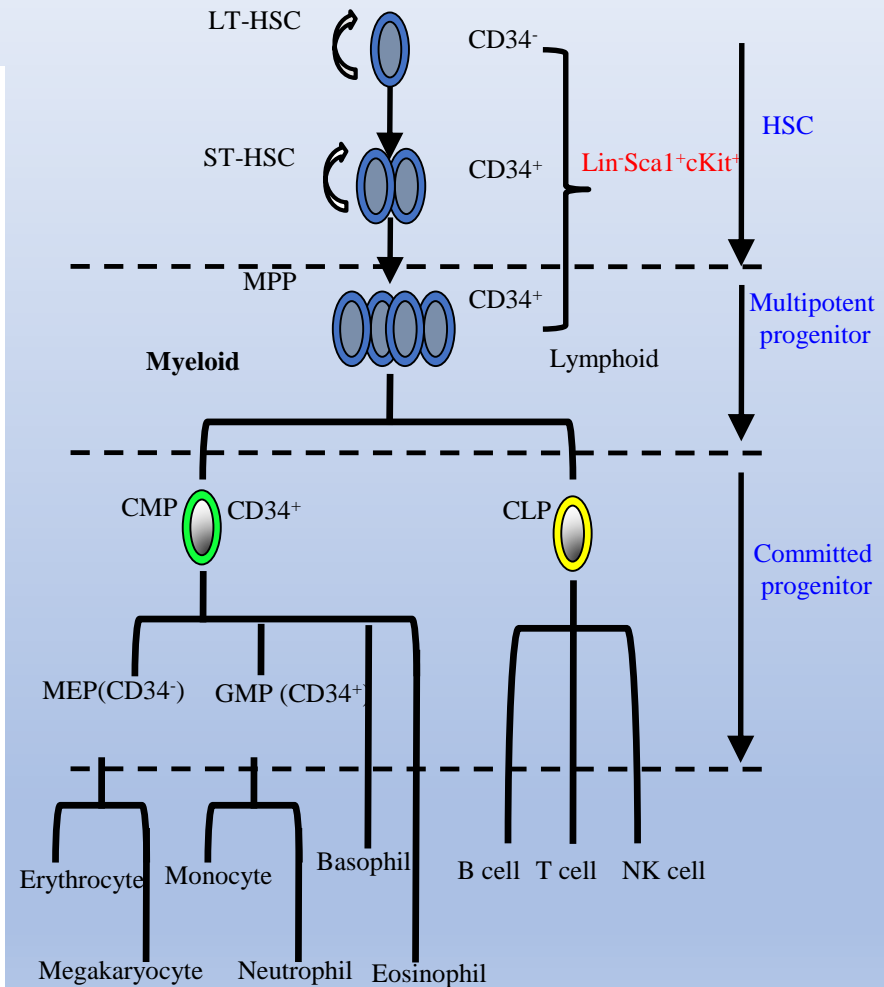
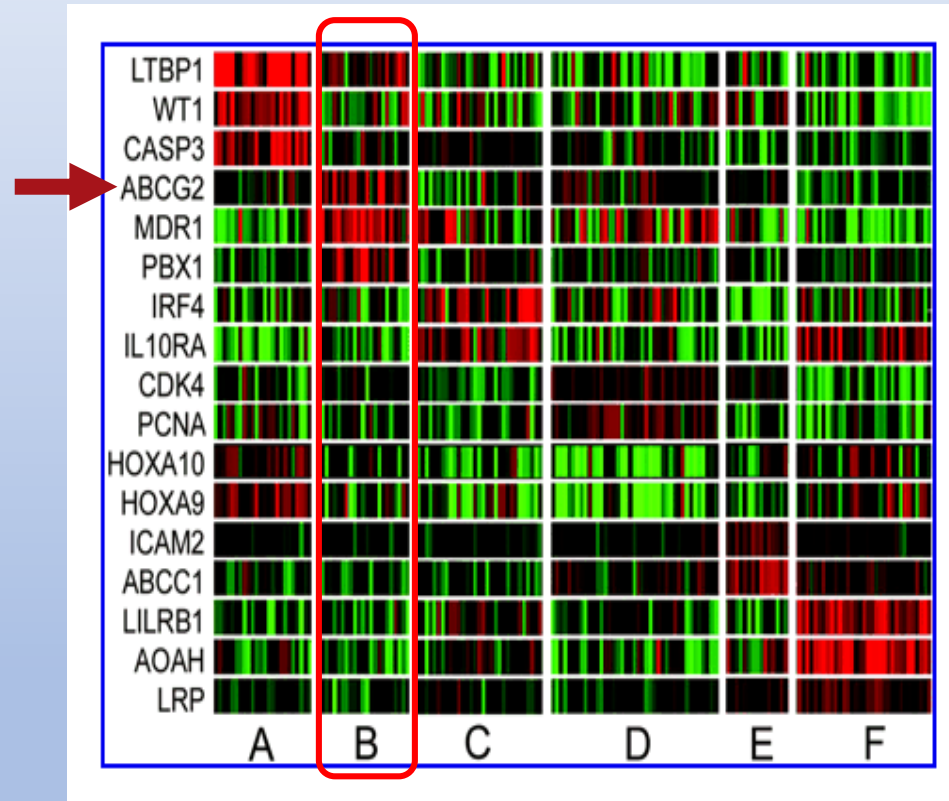


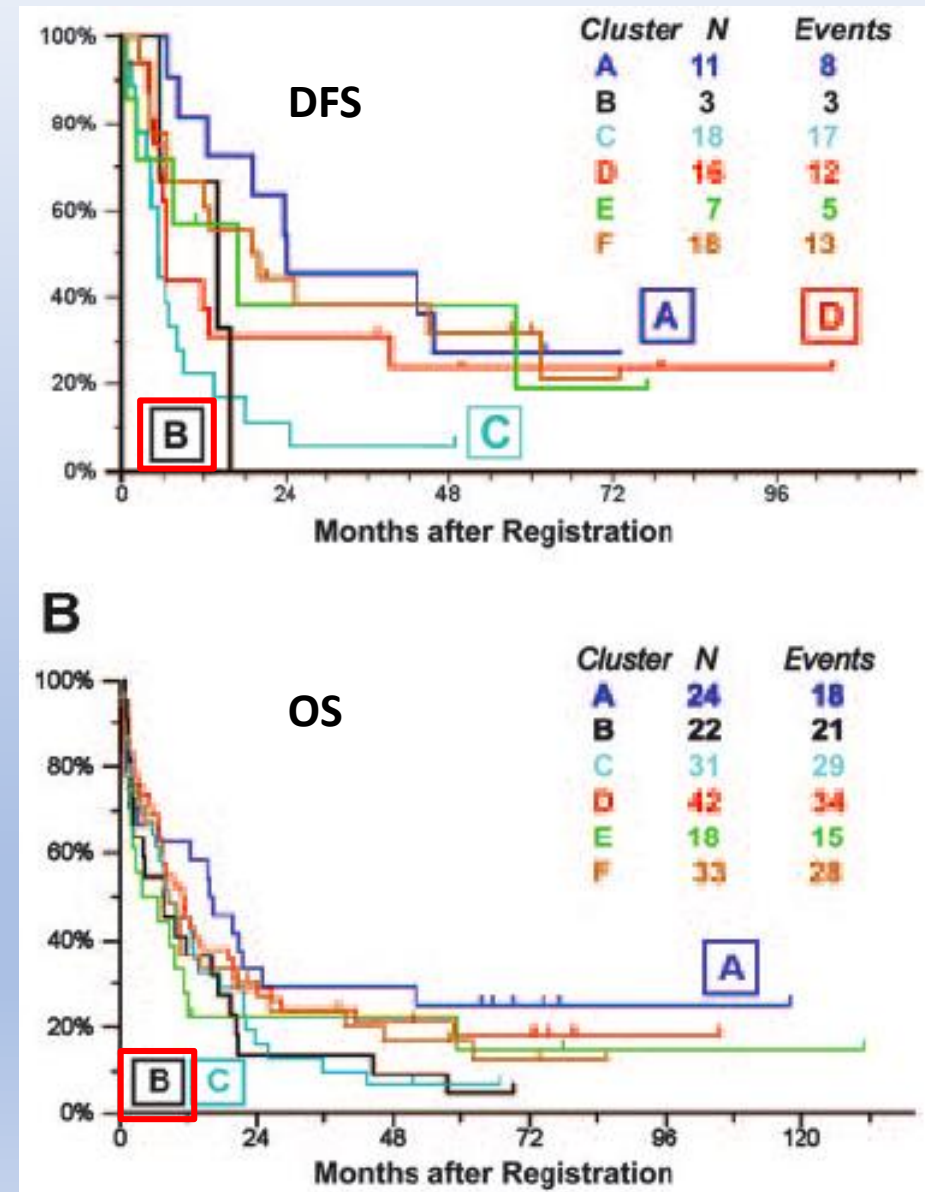
Fig 1. Overall survival of patients who were <20 years of age and diagnosed with AML during the time periods indicated. The data were obtained from [www.seer.cancer.gov/popdata](http://www.seer.cancer.gov/popdata).



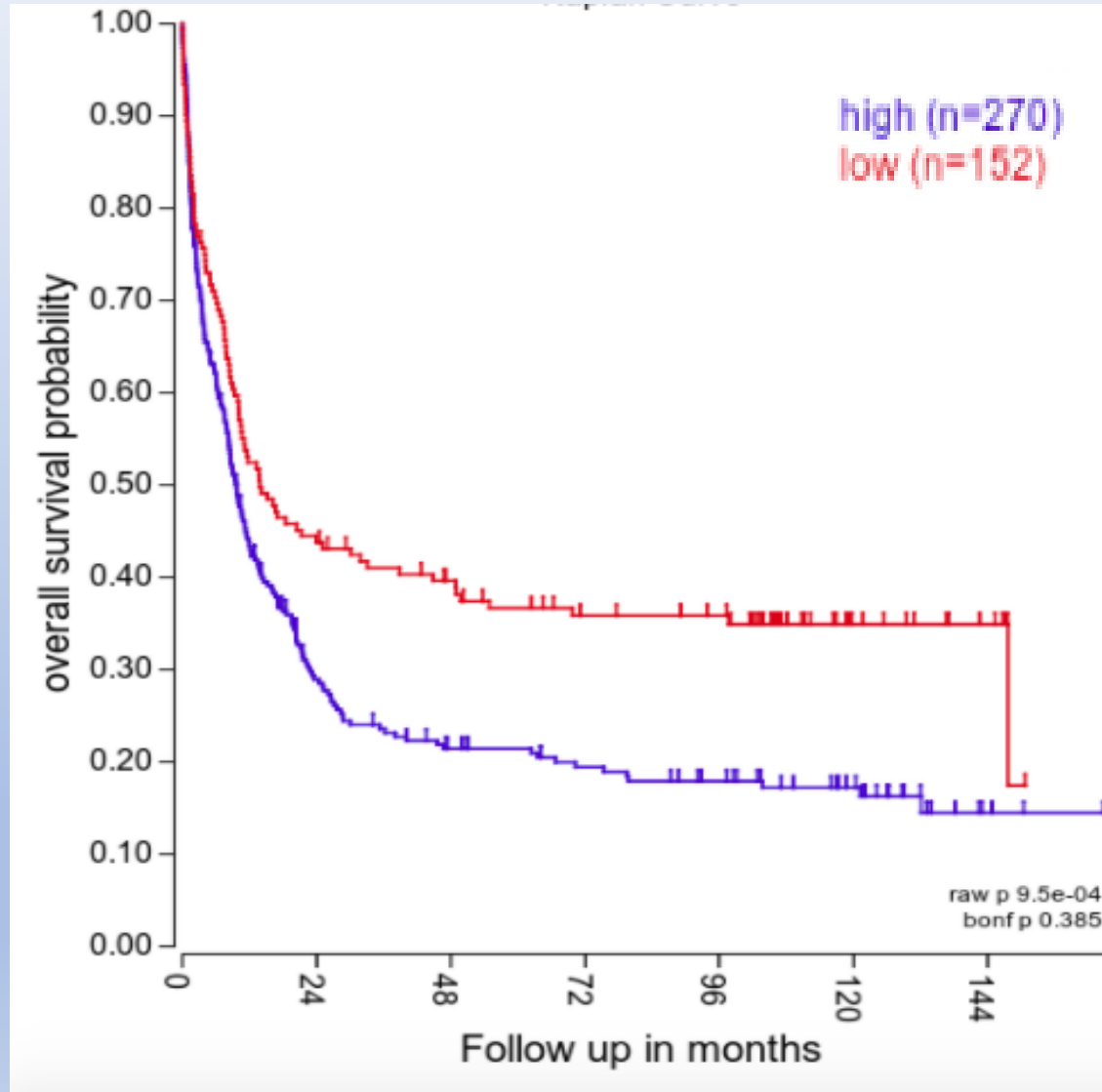
# ABCG2 is highly expressed among AML patients with poor disease-free and overall survival



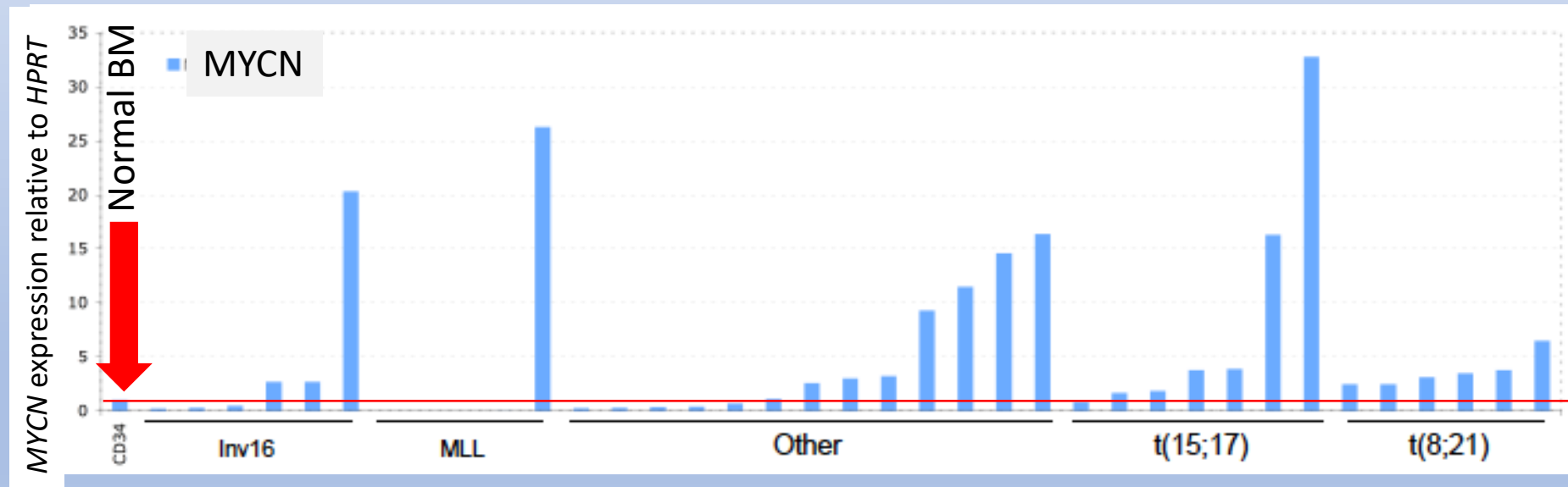
Blood 2006



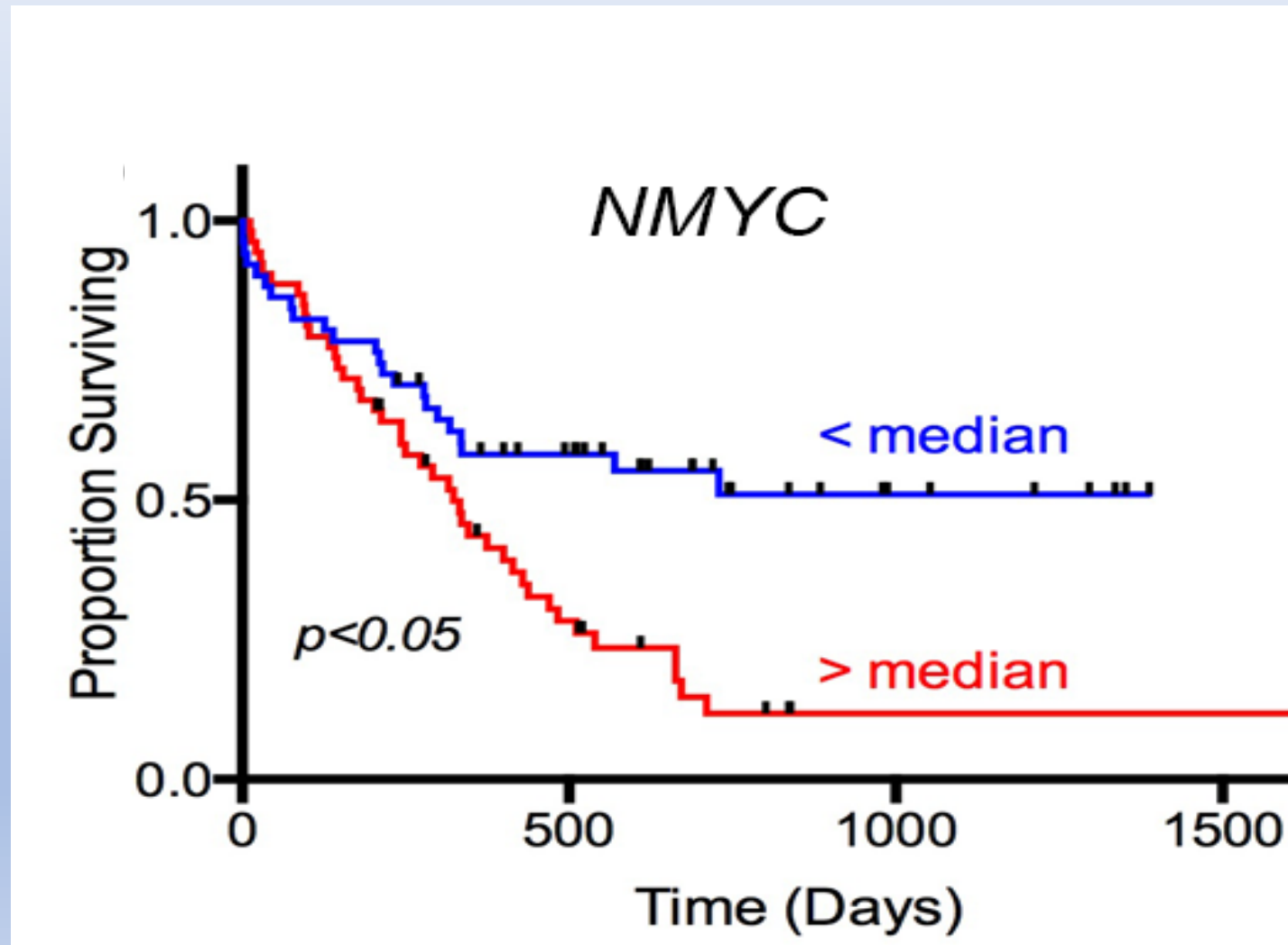
# ABCG2 alone is a poor prognostic factor in AML



# MYCN is a transcription factor highly expressed in pediatric AML

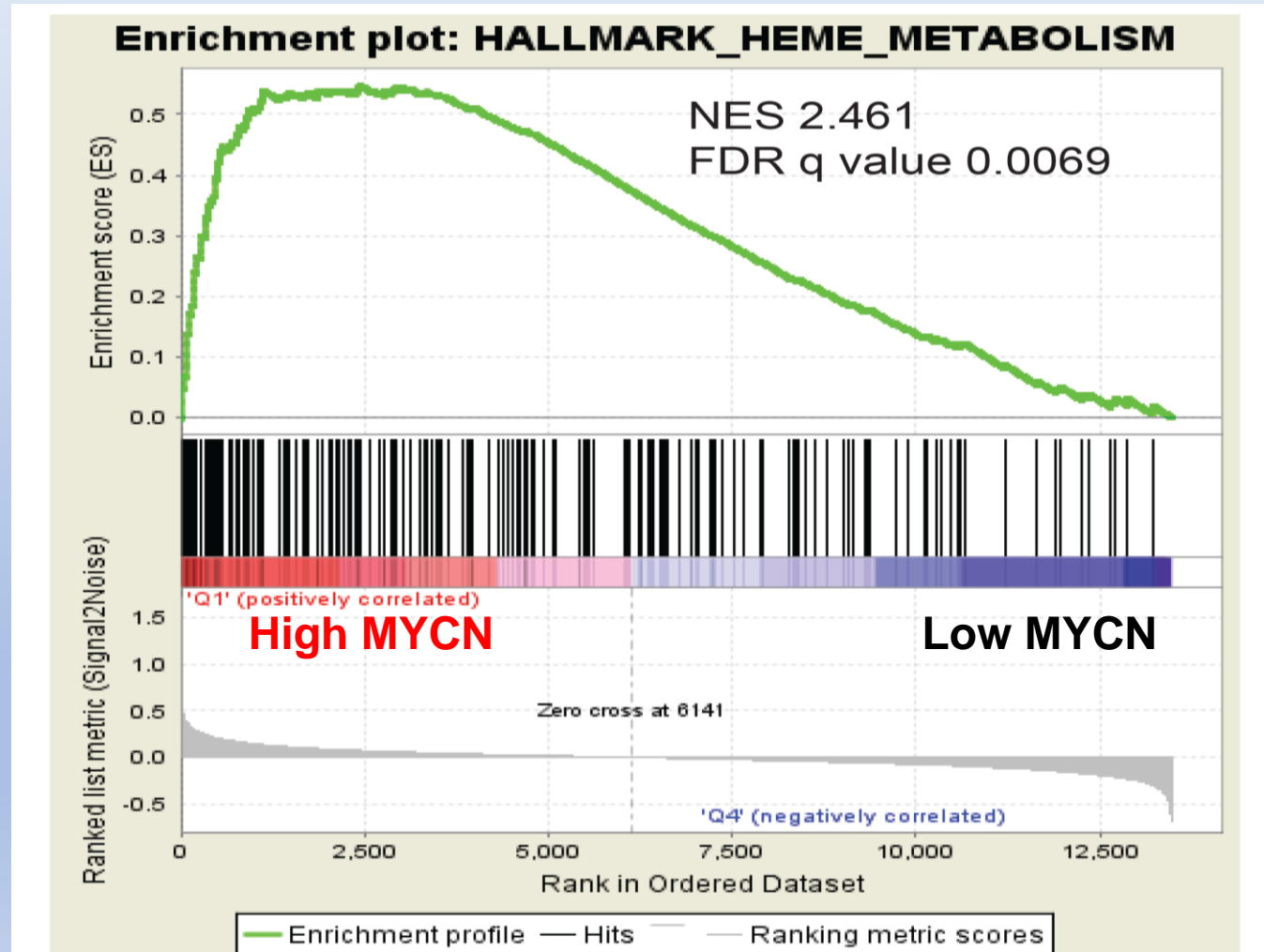


# MYCN is a poor prognostic factor in AML

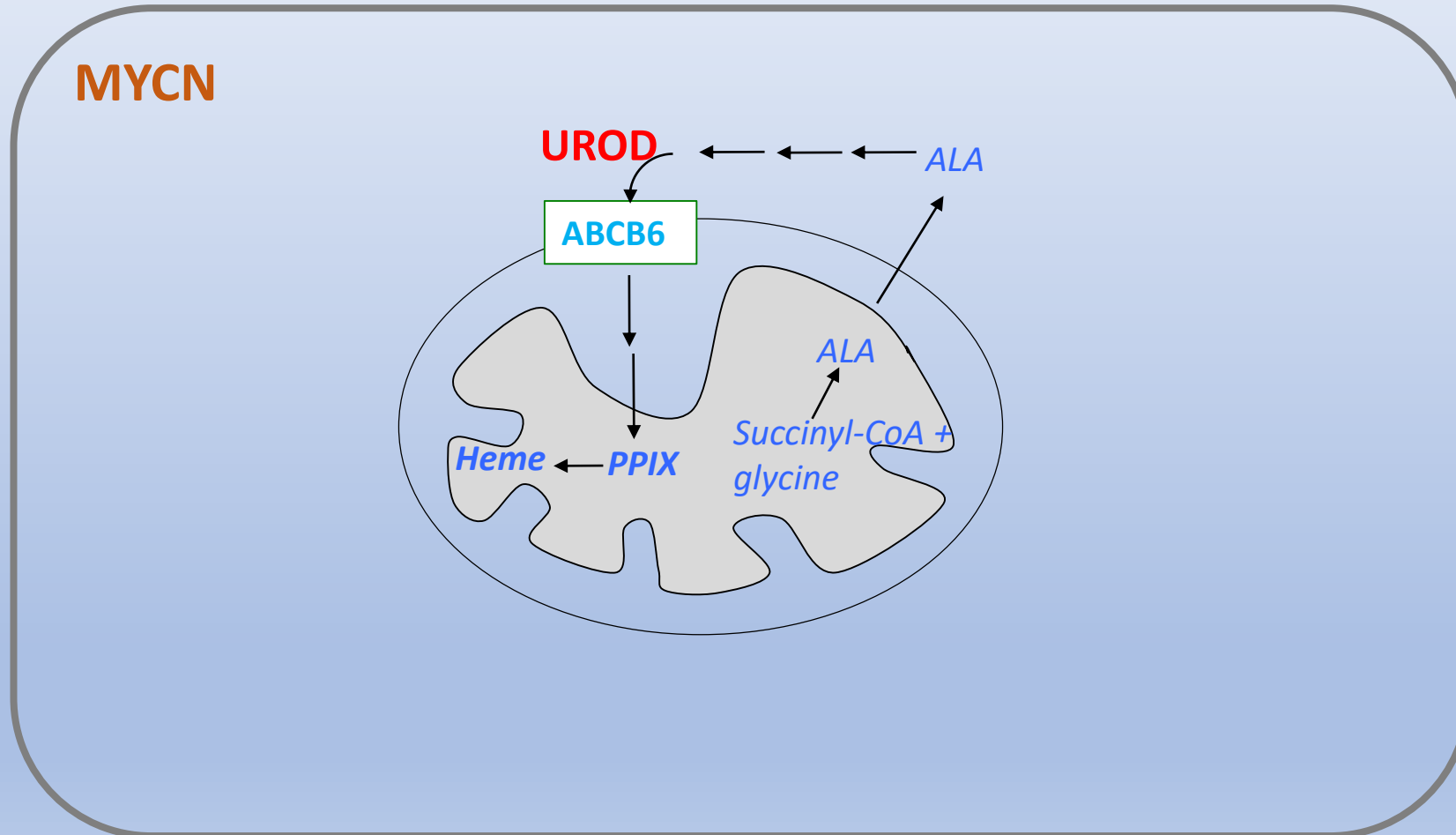


Does MYCN produce metabolic effects?

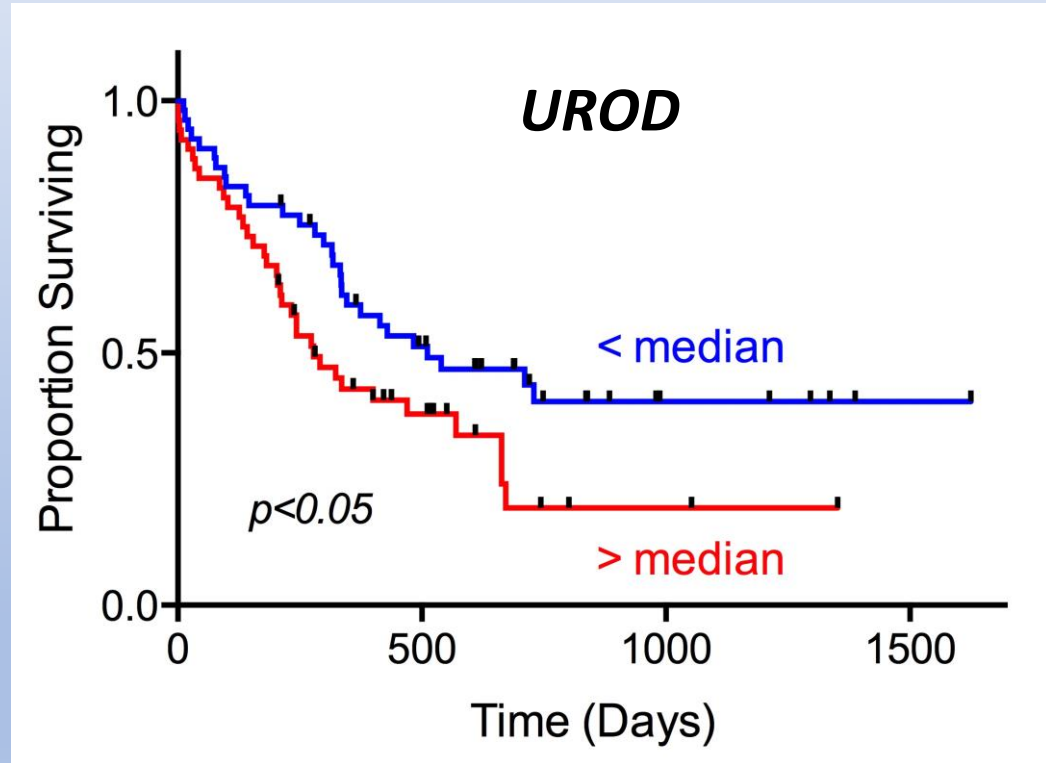
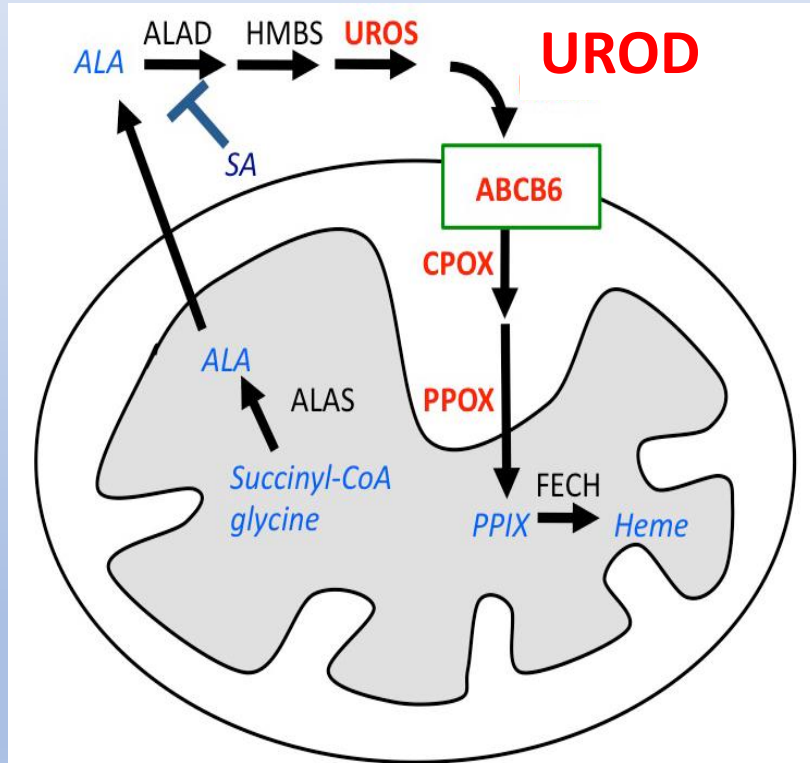
# GSEA reveals upregulated heme metabolism in pediatric AML with high MYCN expression



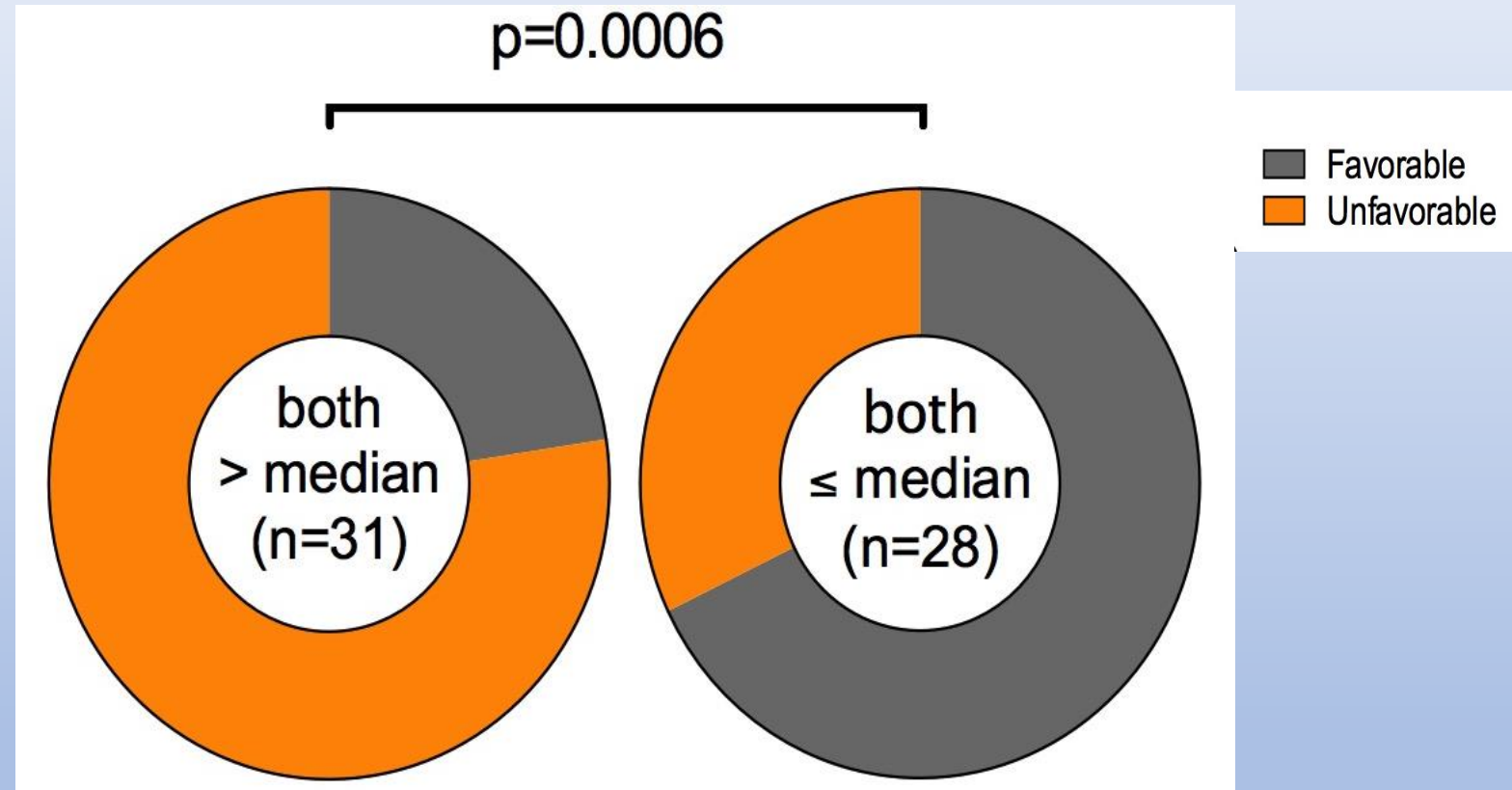
UROD was the most elevated heme pathway gene in High MYCN adult and pediatric AML



# Porphyria biosynthesis is a poor prognostic factor in adult AML



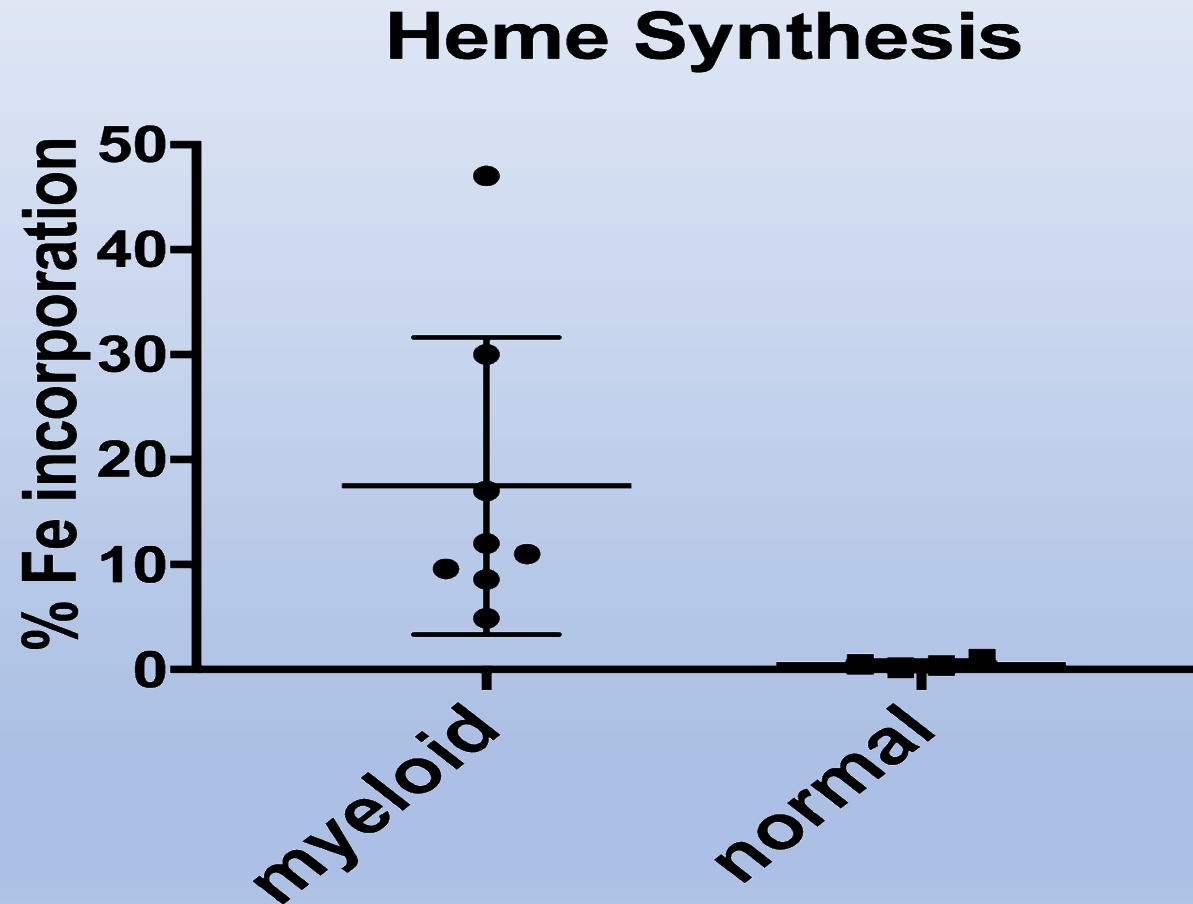
# High *MYCN* and *UROD* expression increases unfavorable outcome in adult AML



**Favorable:** complete remission

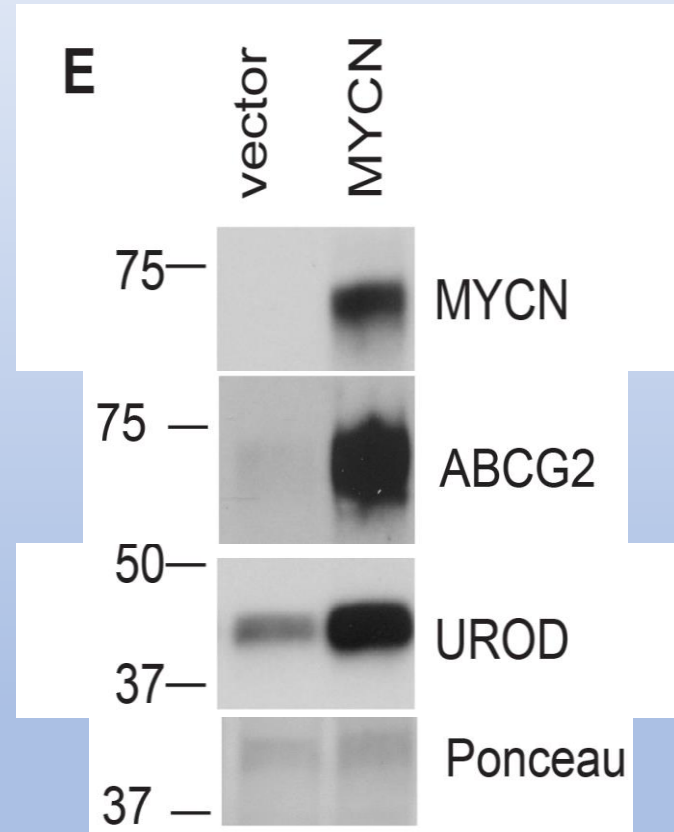
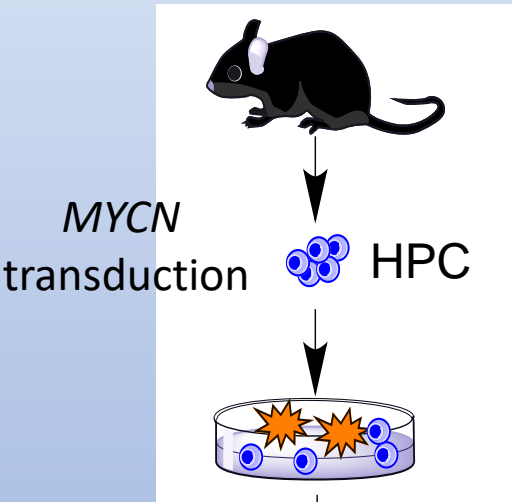
**Unfavorable:** relapse and refractory disease

# AML patients have increased heme synthesis

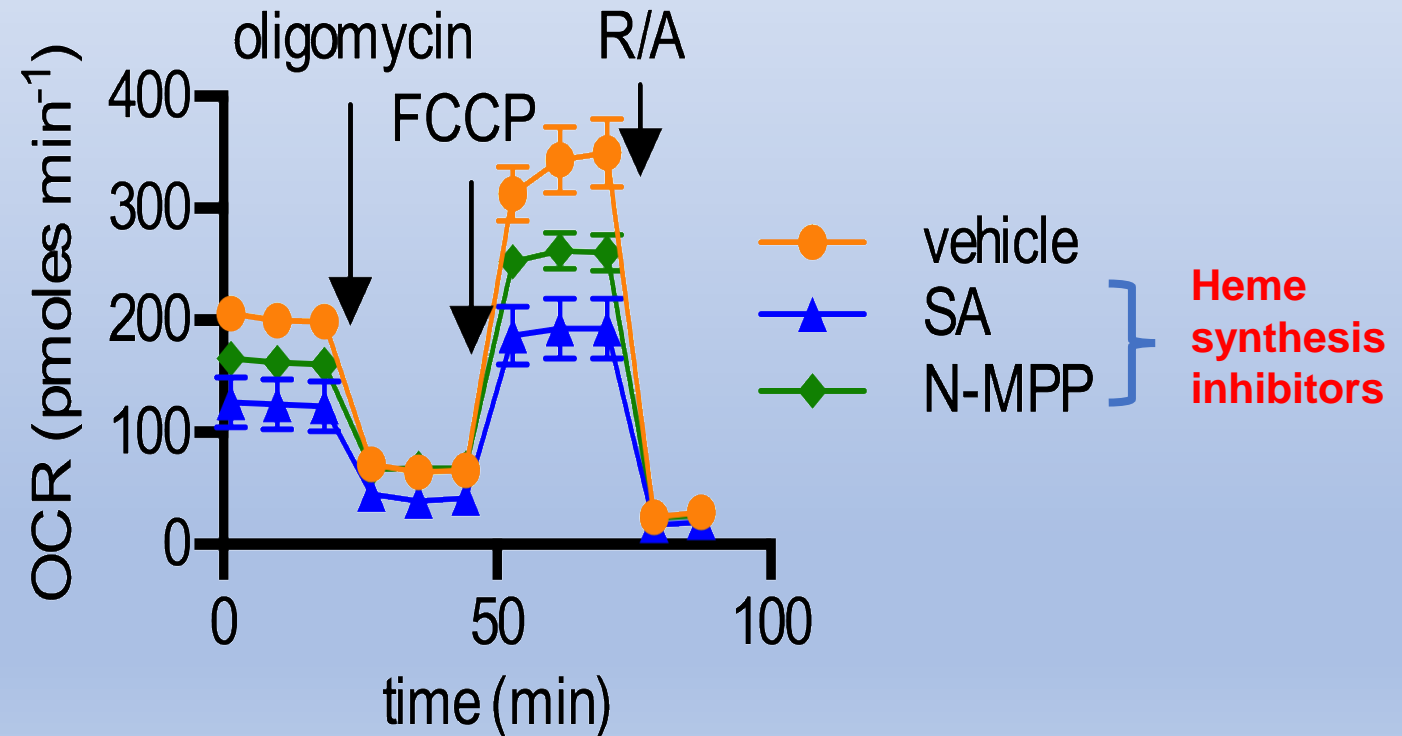
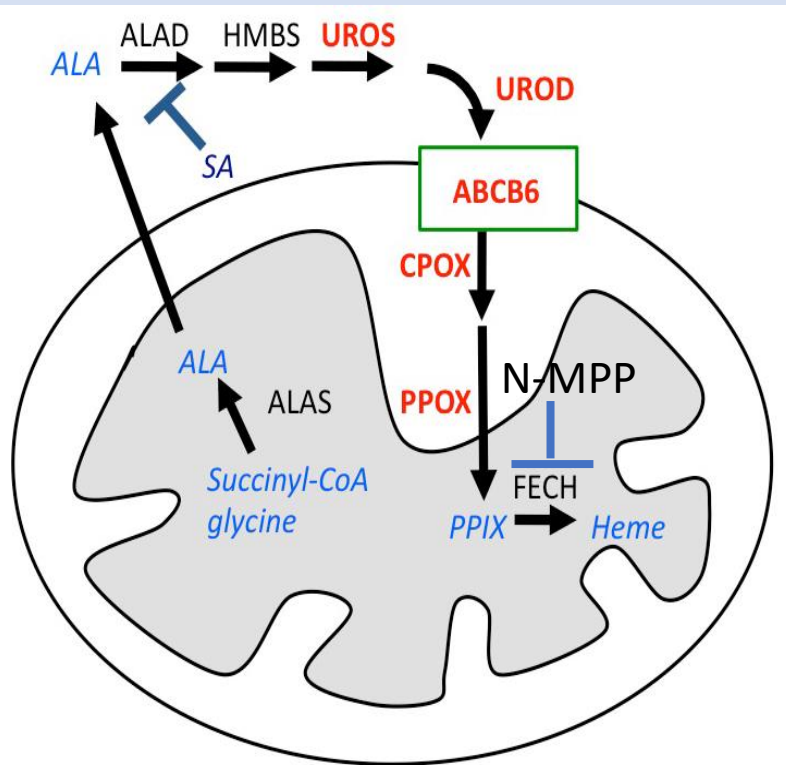


What is the purpose of increased heme biosynthesis ?

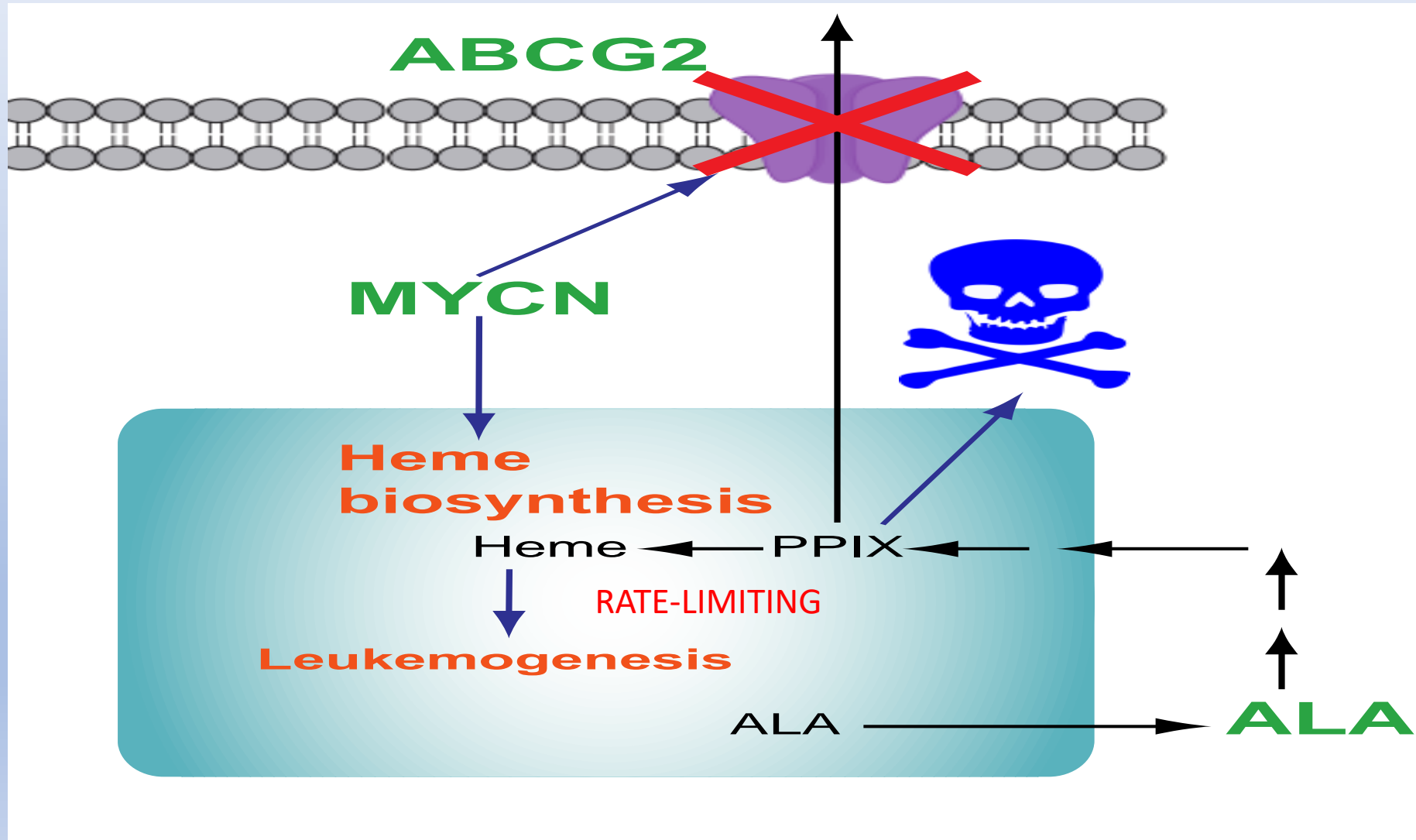
# In Leukemic Progenitors MYCN upregulates heme/porphyrin synthesis, UROD and ABCG2 expression



# Heme synthesis is required for maximal mitochondrial oxygen consumption (OCR) in MYCN-Leukemic Progenitors



Hypothesized that upregulated heme biosynthetic pathway would be a metabolic vulnerability in MYCN driven AML



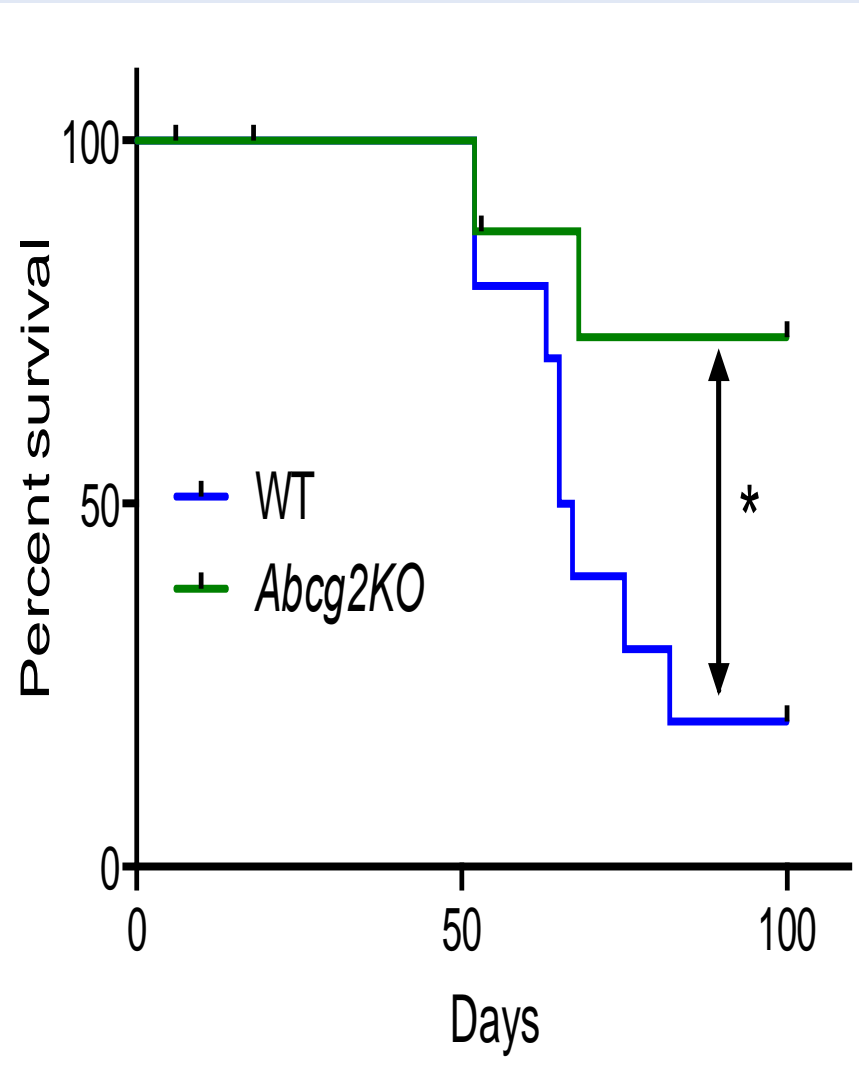
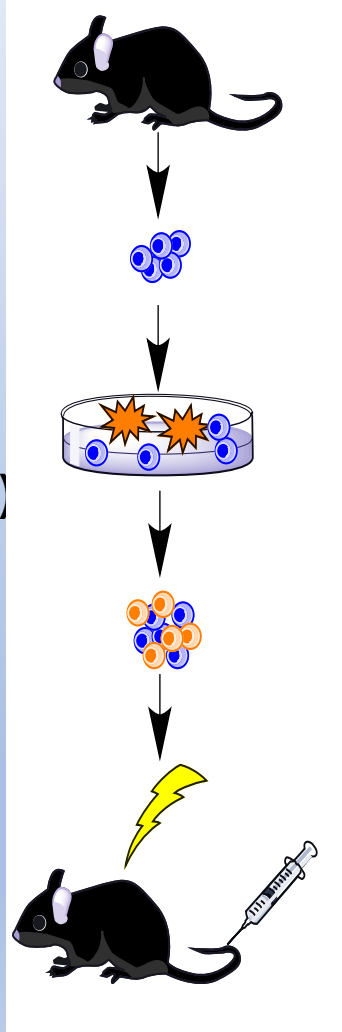
# A mouse model shows that loss of *Abcg2* increases MYCN driven myeloid leukemia survival

WT or *Abcg2*KO donor mouse

HPCs

MYCN transduction  
(Leukemic progenitors)

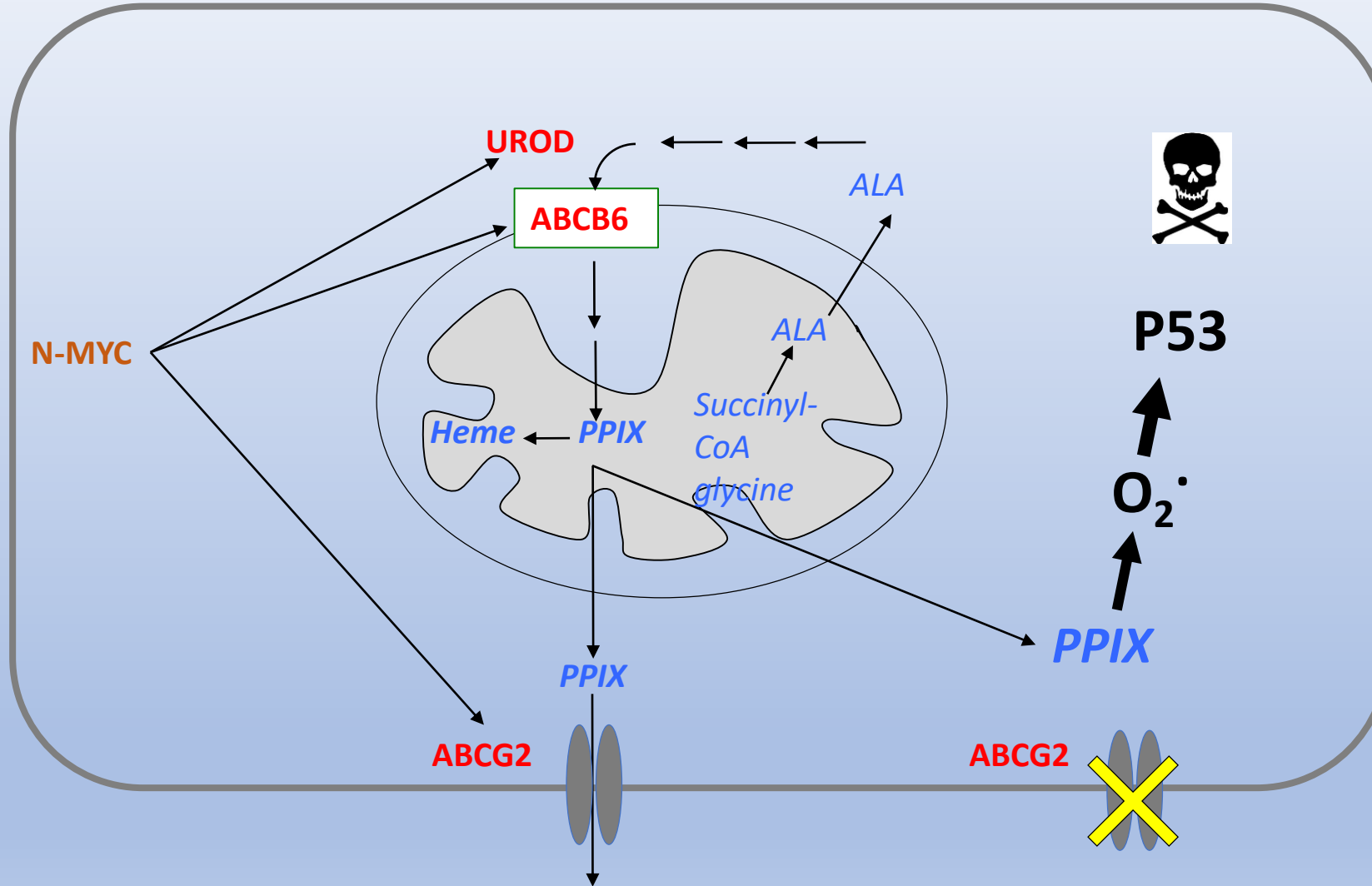
Transplant into lethally irradiated congenic recipient mice



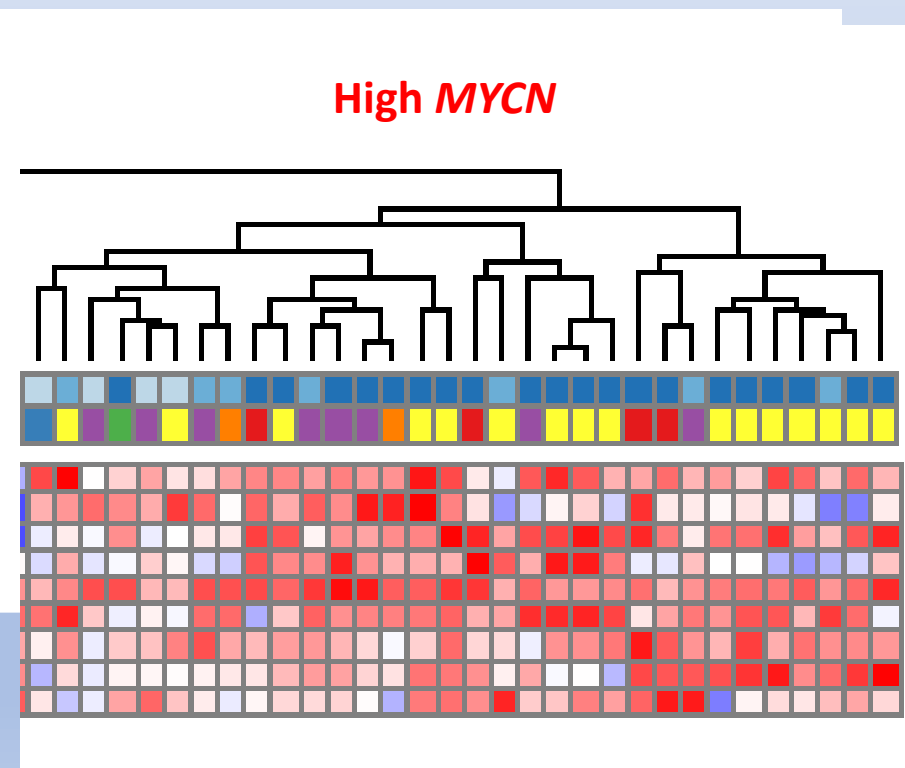
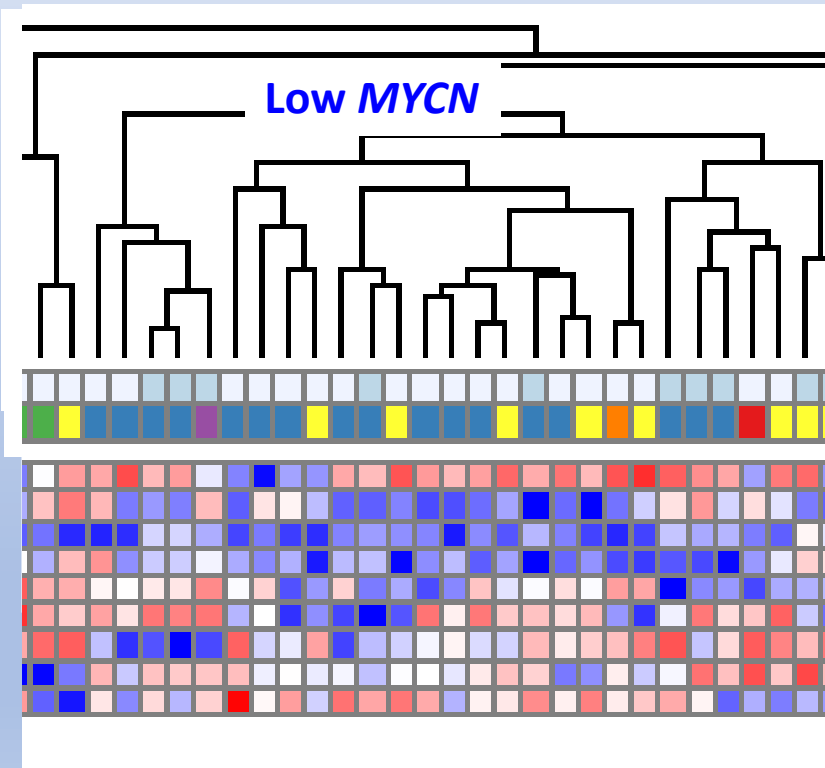
PPIX



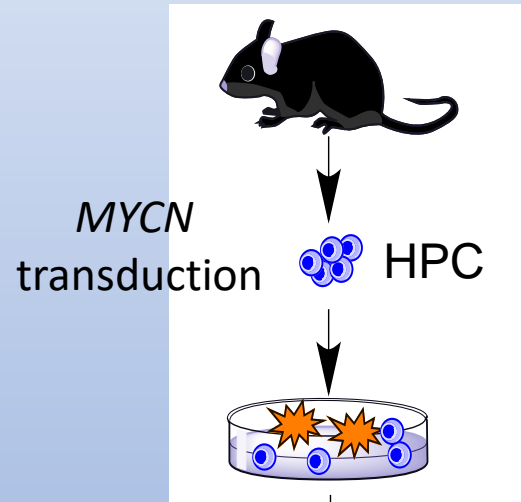
# Does PPIX activate P53?



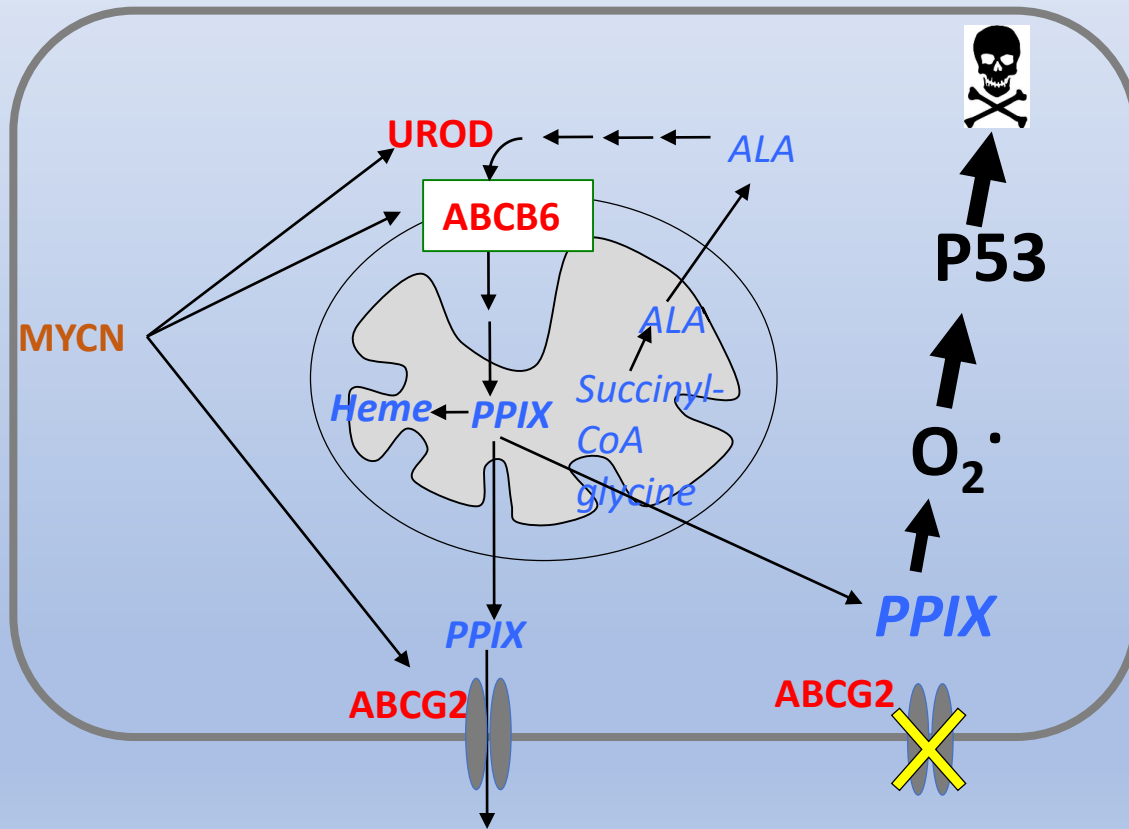
# Higher *MYCN* correlates with p53 activation in SJ AML patients



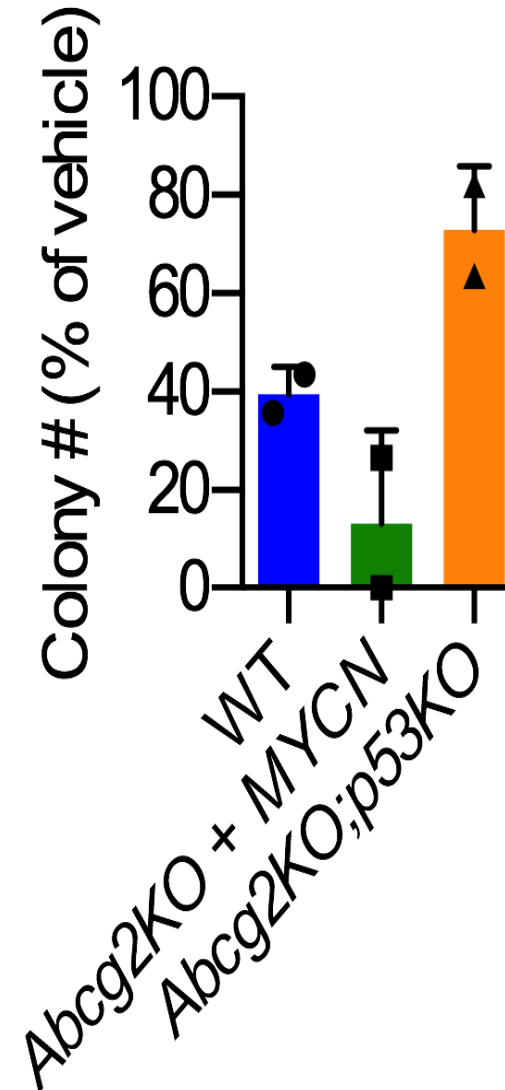
# P53 is activated in MYCN leukemic progenitors



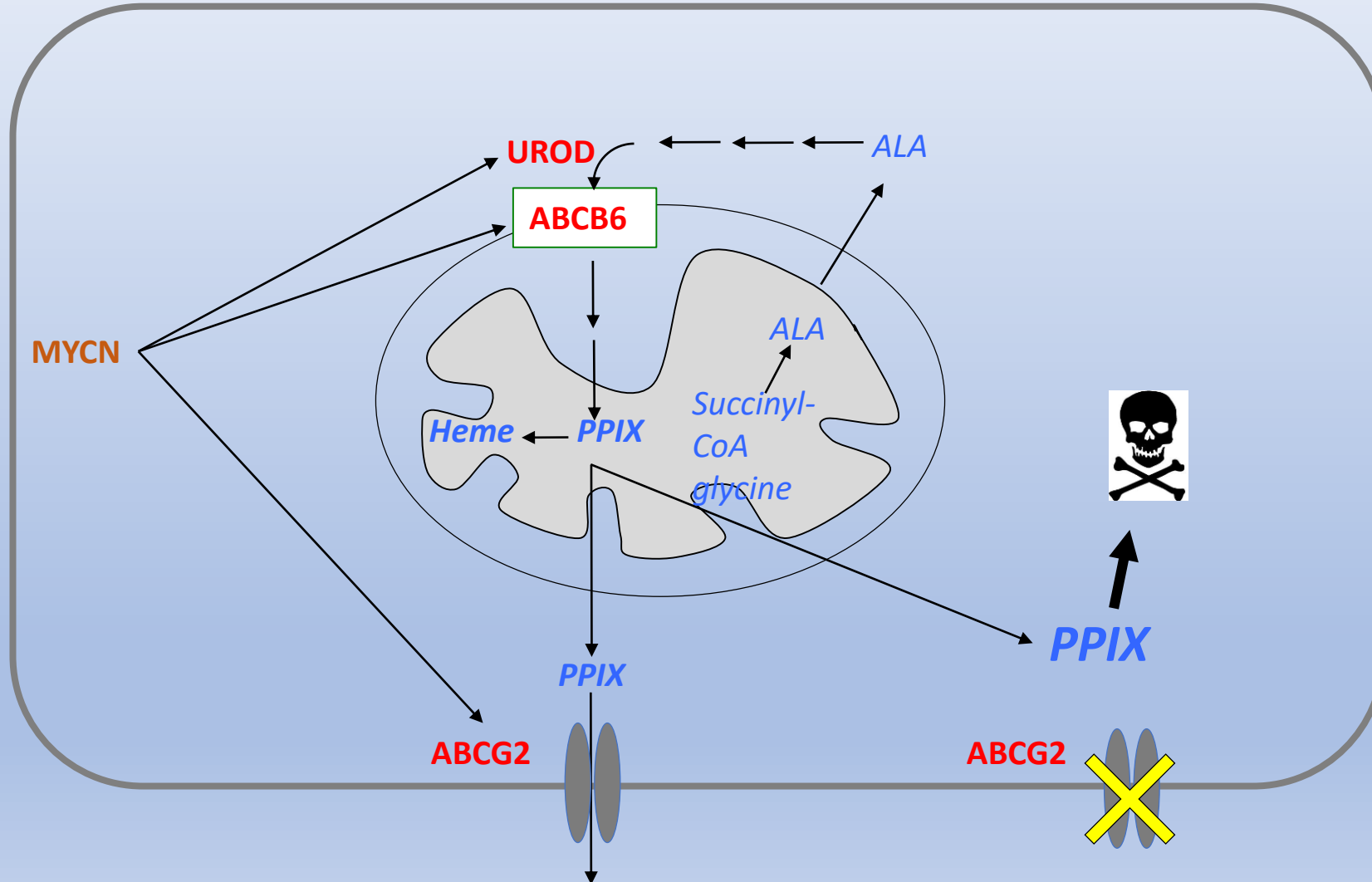
# P53 deletion mostly rescues self-renewal of porphyrin treated Abcg2KO MYCN cells



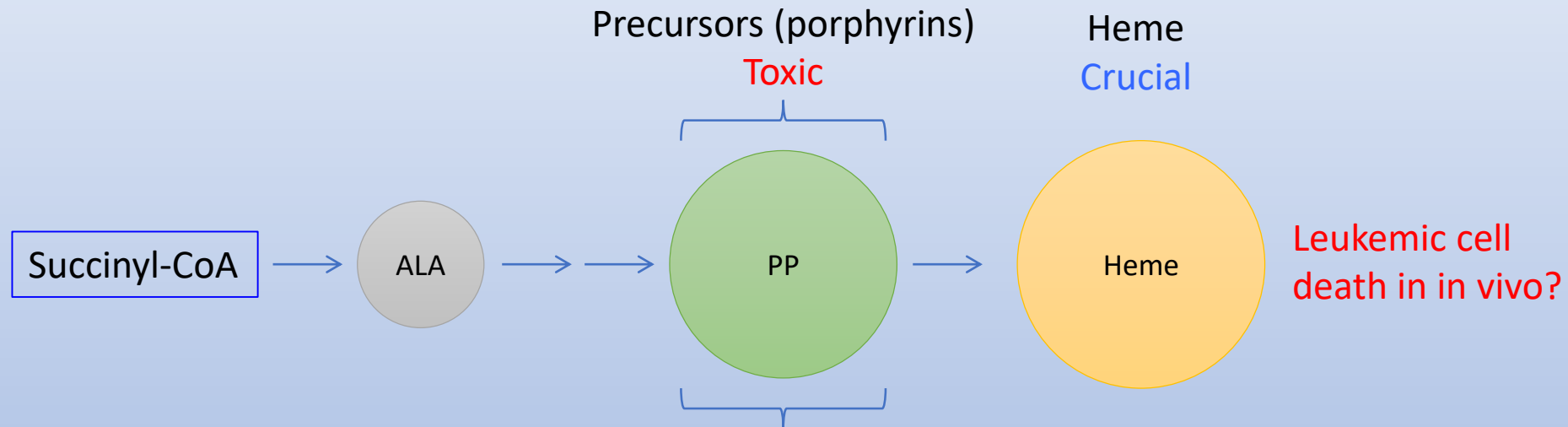
p53 null mice- Martine Roussel and Frederique Zindy



# Do additional PPIX-protein interactions produce cytotoxicity?

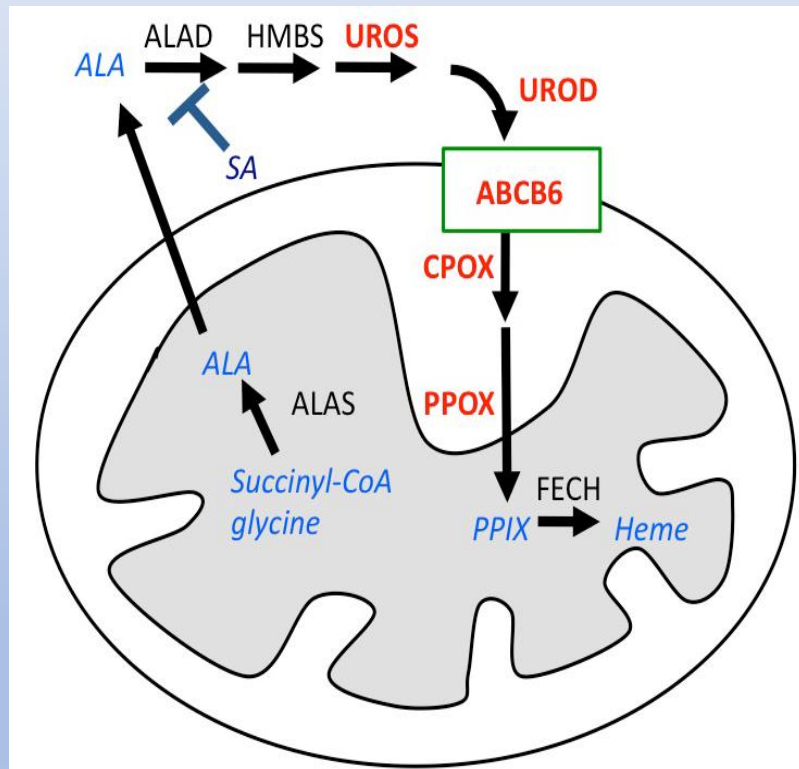


# Targeting the heme biosynthetic pathway in leukemic progenitors

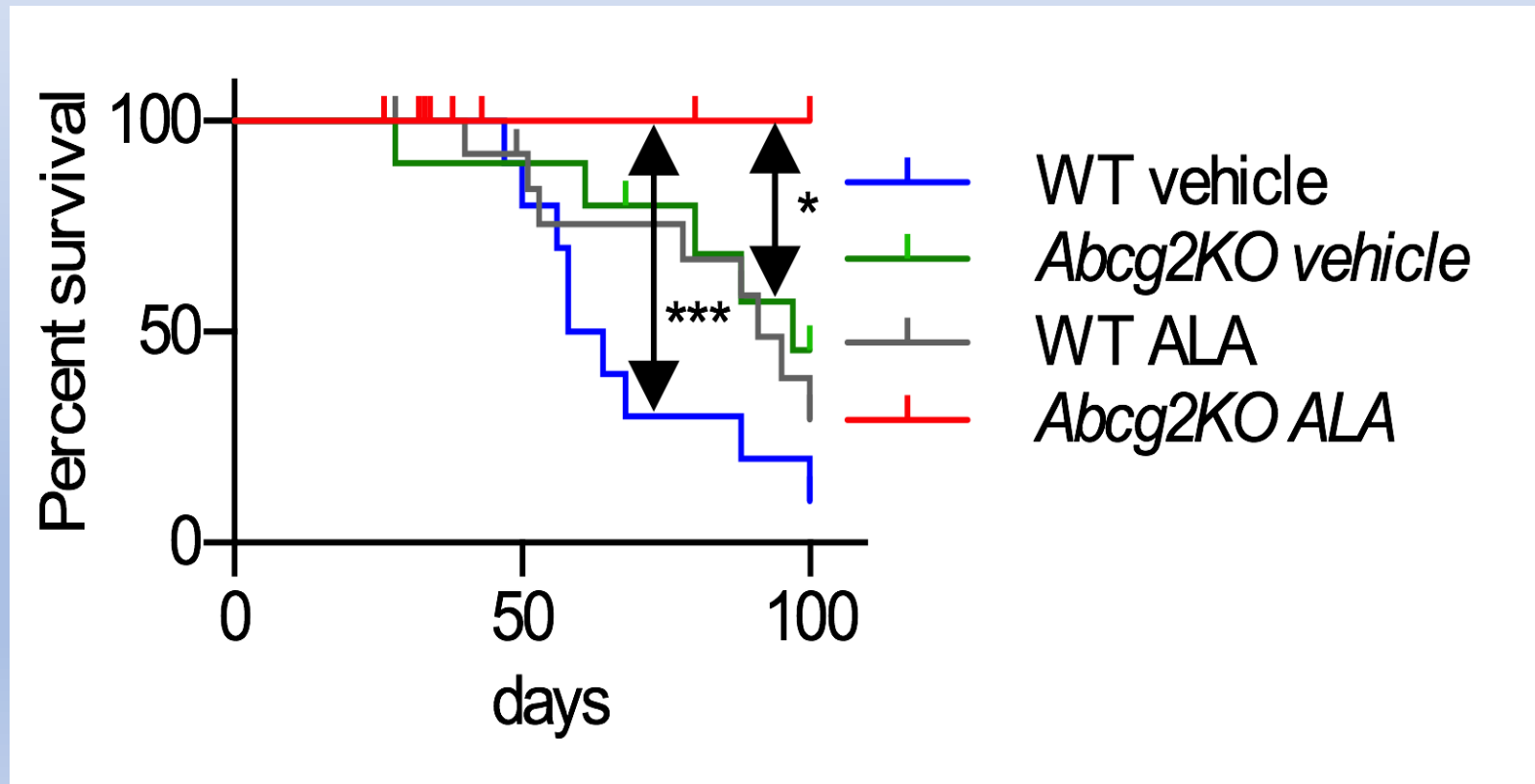


**Loss of Abcg2**  
**P53 activation**  
**Other targets?**

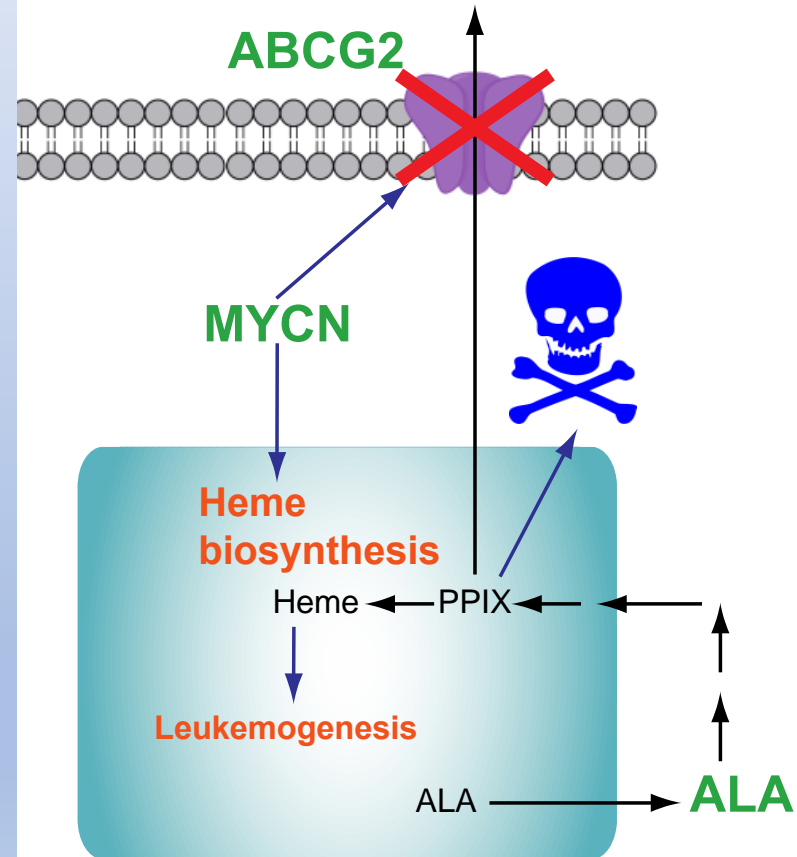
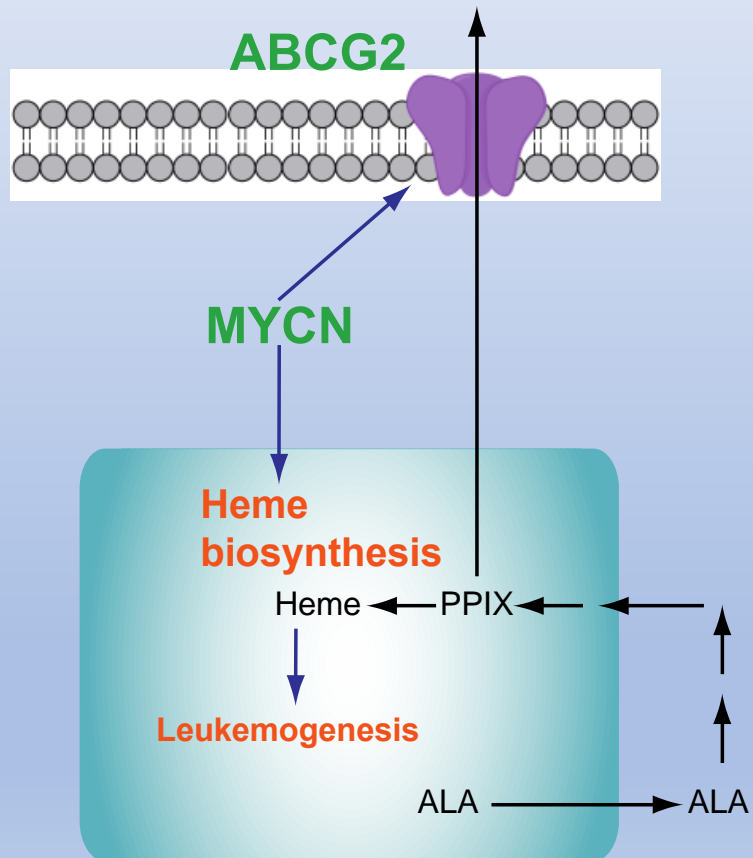
# Does increasing PPIX cure AML?



# ALA-treatment of MYCN Leukemic Progenitors cures ABCG2 KO AML



# Summary

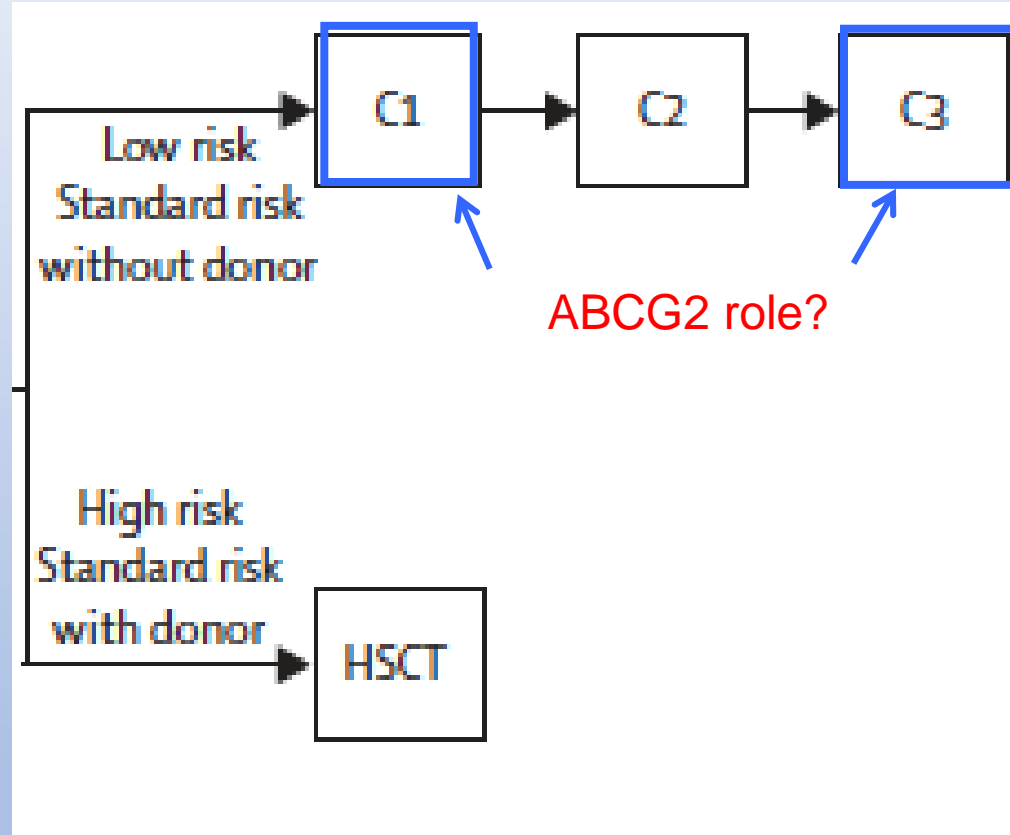


# ABCG2 substrates are used in conventional AML therapy

## AML02 treatment scheme

### Induction

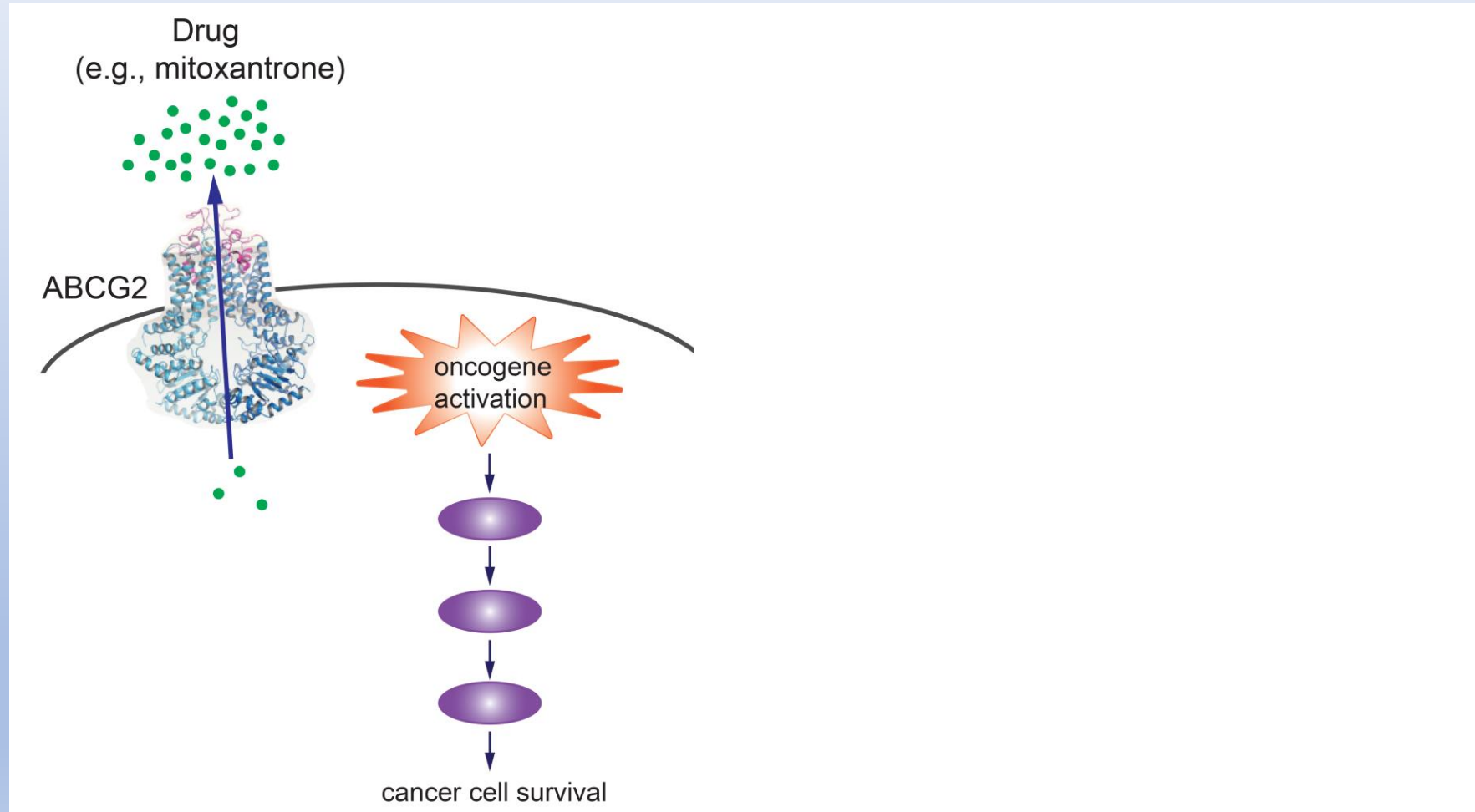
Daunorubicin  
etoposide  
cytarabine



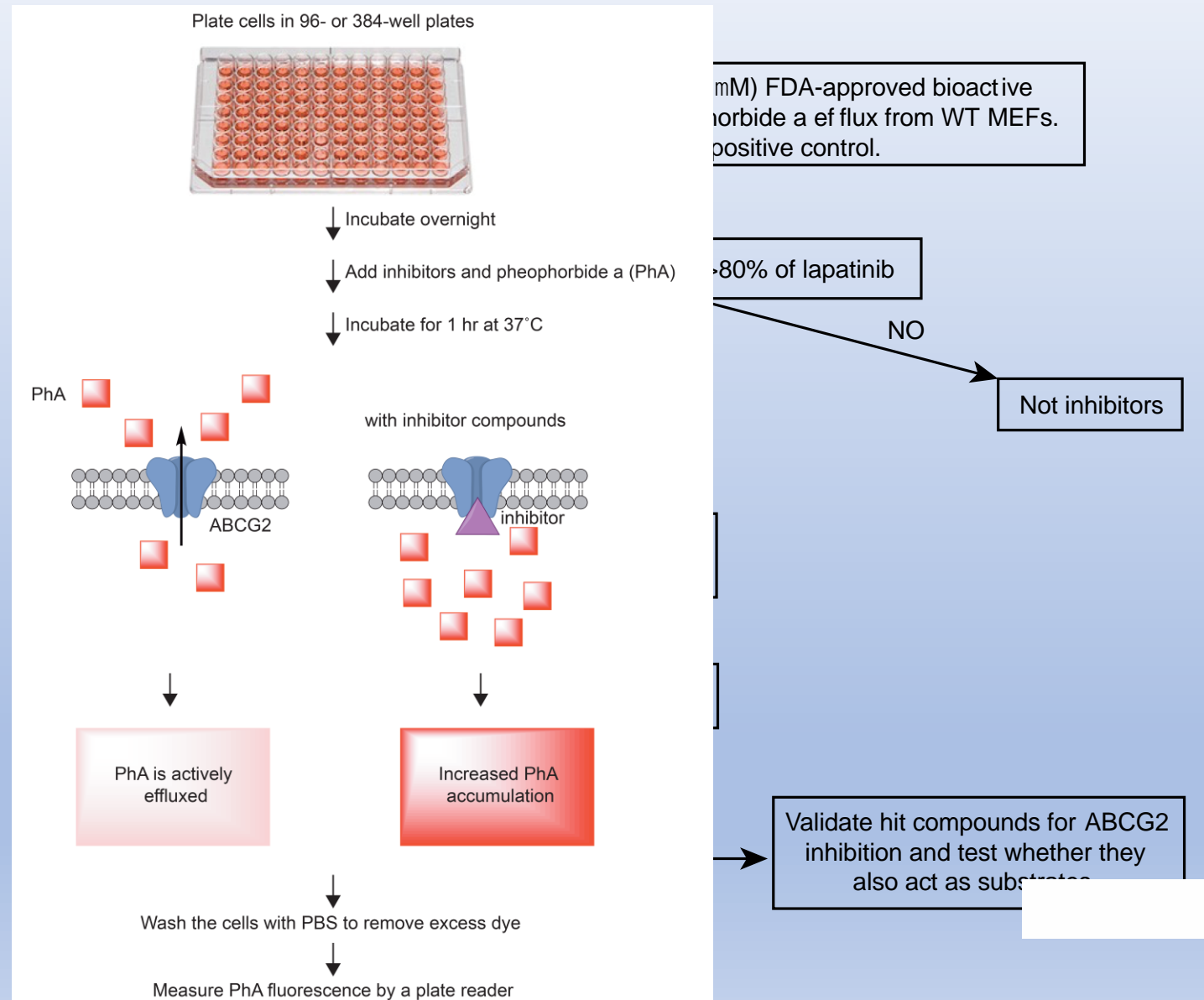
### Consolidation

C1: cytarabine & **cladribine**  
cytarabine & etoposide  
cytarabine & **mitoxantrone**  
C2: cytarabine & L-asparaginase  
C3: cytarabine & **mitoxantrone**

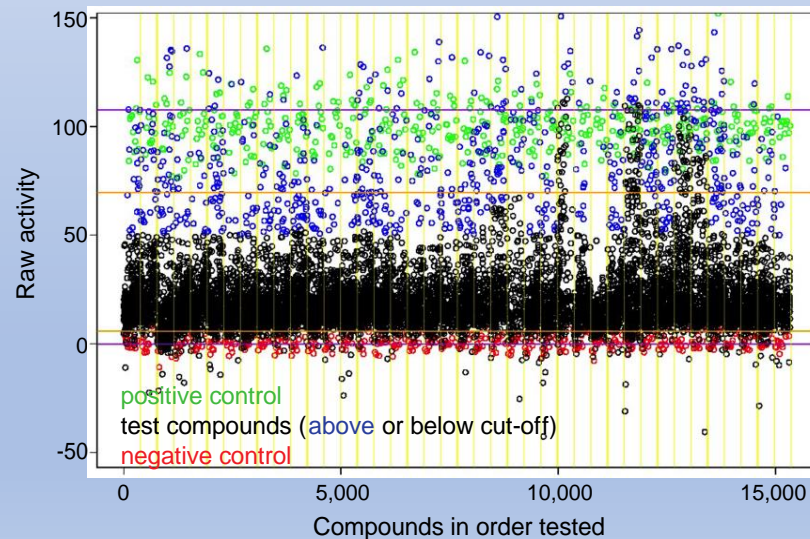
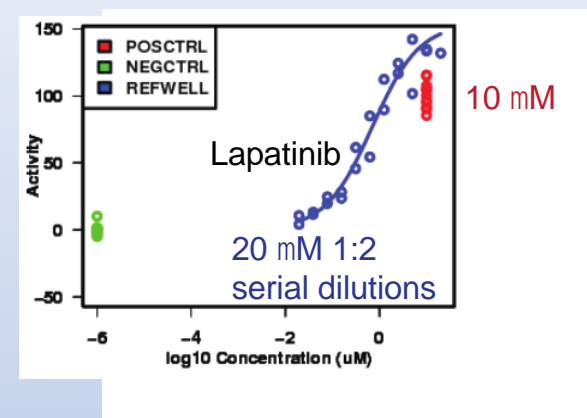
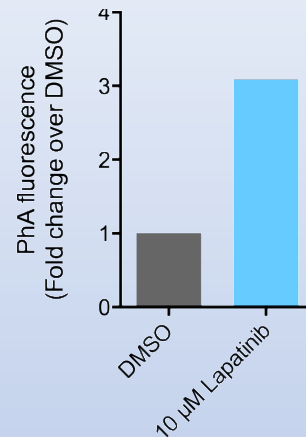
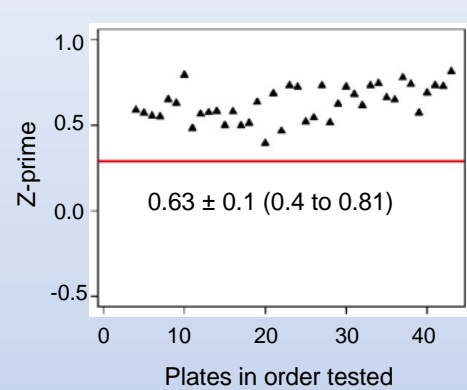
# Screening for ABCG2 and Cancer liability inhibitors



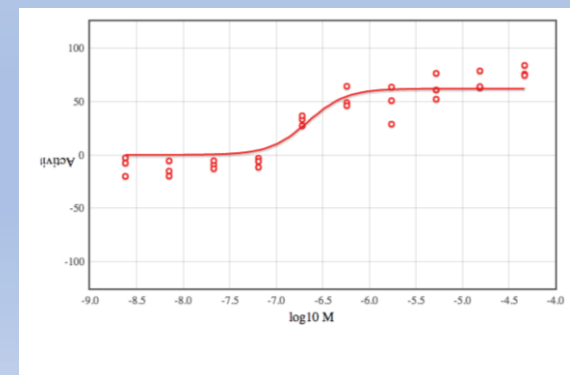
# High throughput screening to identify ABCG2 inhibitors



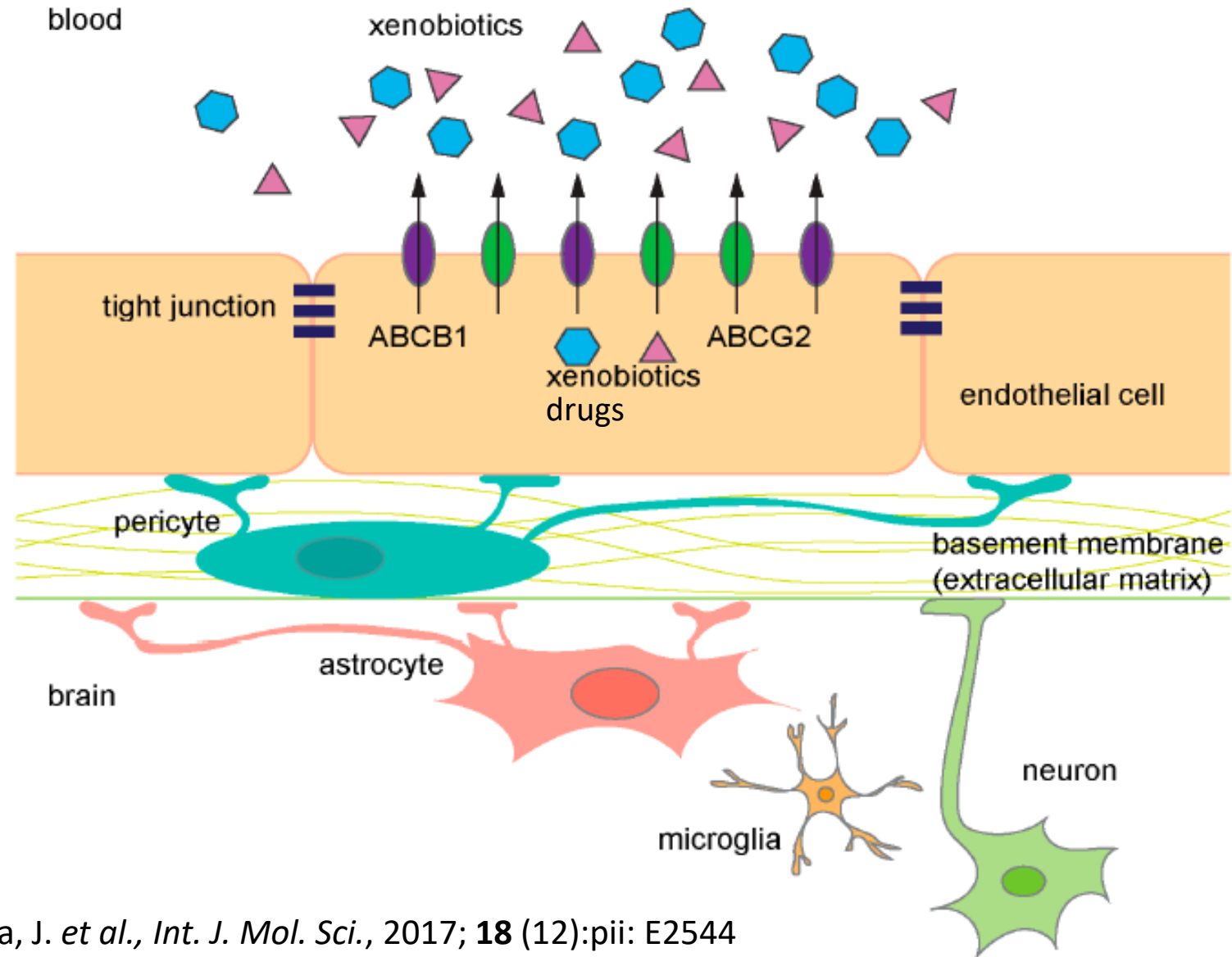
# High throughput screening for ABCG2 inhibitors identified novel inhibitors



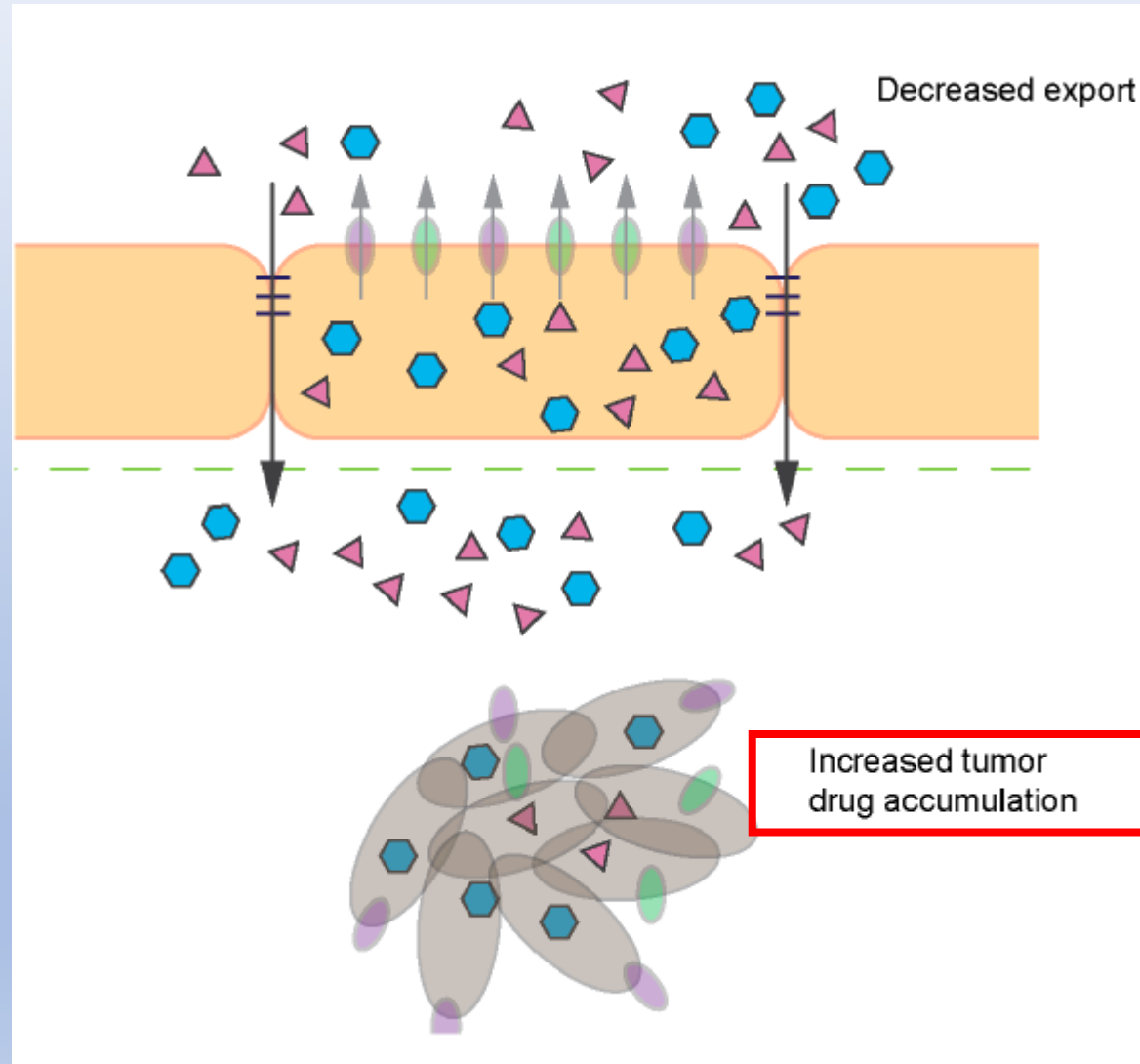
SJ000831433



# Inhibition of ABCG2 at the blood-brain barrier (BBB)



# ABCG2 inhibition may allow the drugs accumulation in the tumor



# “Teamwork to make the Dream work”



AP  
Naren  
CCHMC



Jin  
Zhang  
UCSD

## Funding

- P30 Cancer Center Support Grant
- ALSAC
- NIH R01s



**J Schuetz Lab-2018**

Yu  
Fukuda

## St Jude Collaborators



Martine  
Roussel  
Tumor Cell Biol



Jiyang  
Yu  
Comp. Bio



Taosheng  
Chen  
Chem. Bio.



Junmin  
Peng  
Struc.  
Biol