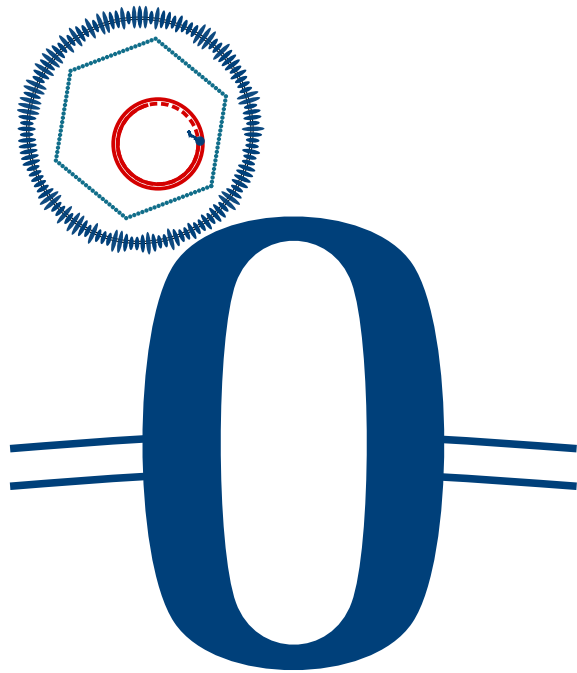


Optimization of an *in vitro* assay for identification of NTCP inhibitors



Pieter Van Brantegem
KU Leuven

Meet the Experts Transporter
Conference Budapest
26 April 2018

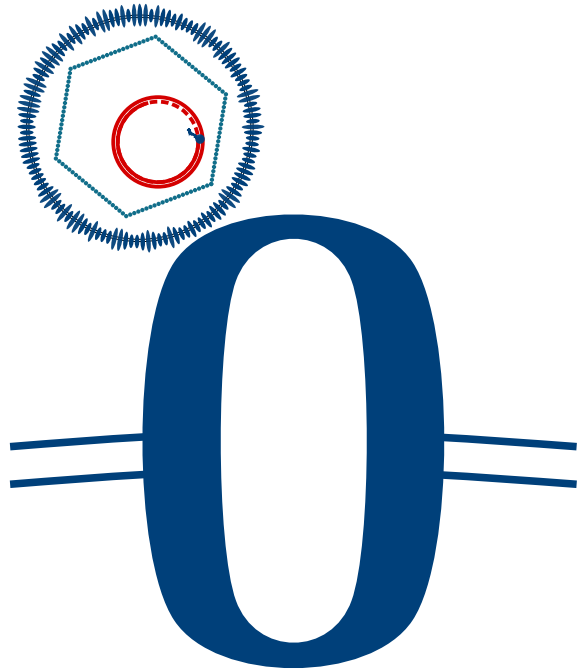
Chronic hepatitis B virus infection

Affects 250 million people world wide

Is a common cause of cirrhosis and hepatocellular carcinoma

Cannot be eradicated yet

Optimization of an *in vitro* assay for identification of NTCP inhibitors



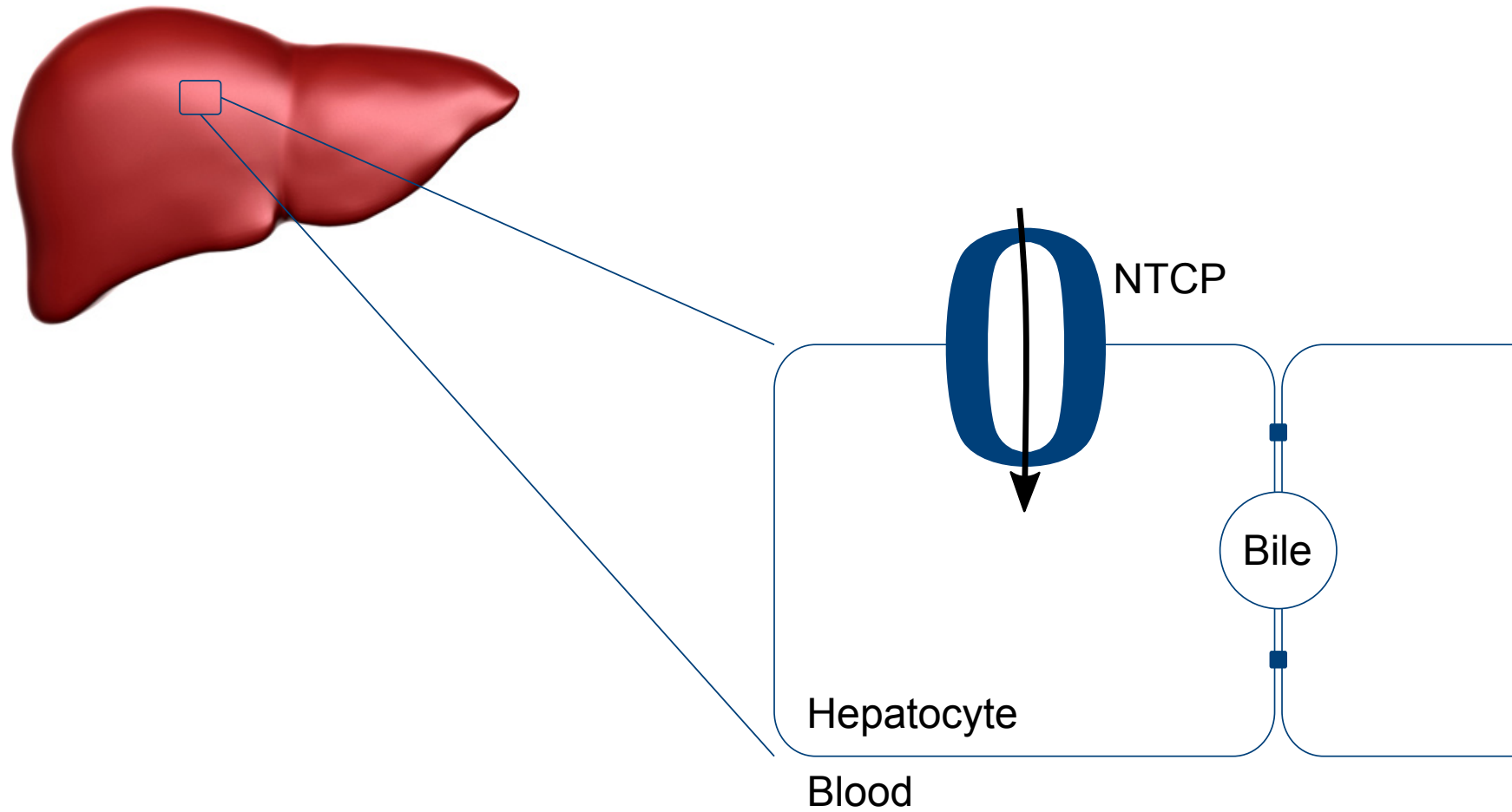
Rationale

Methods

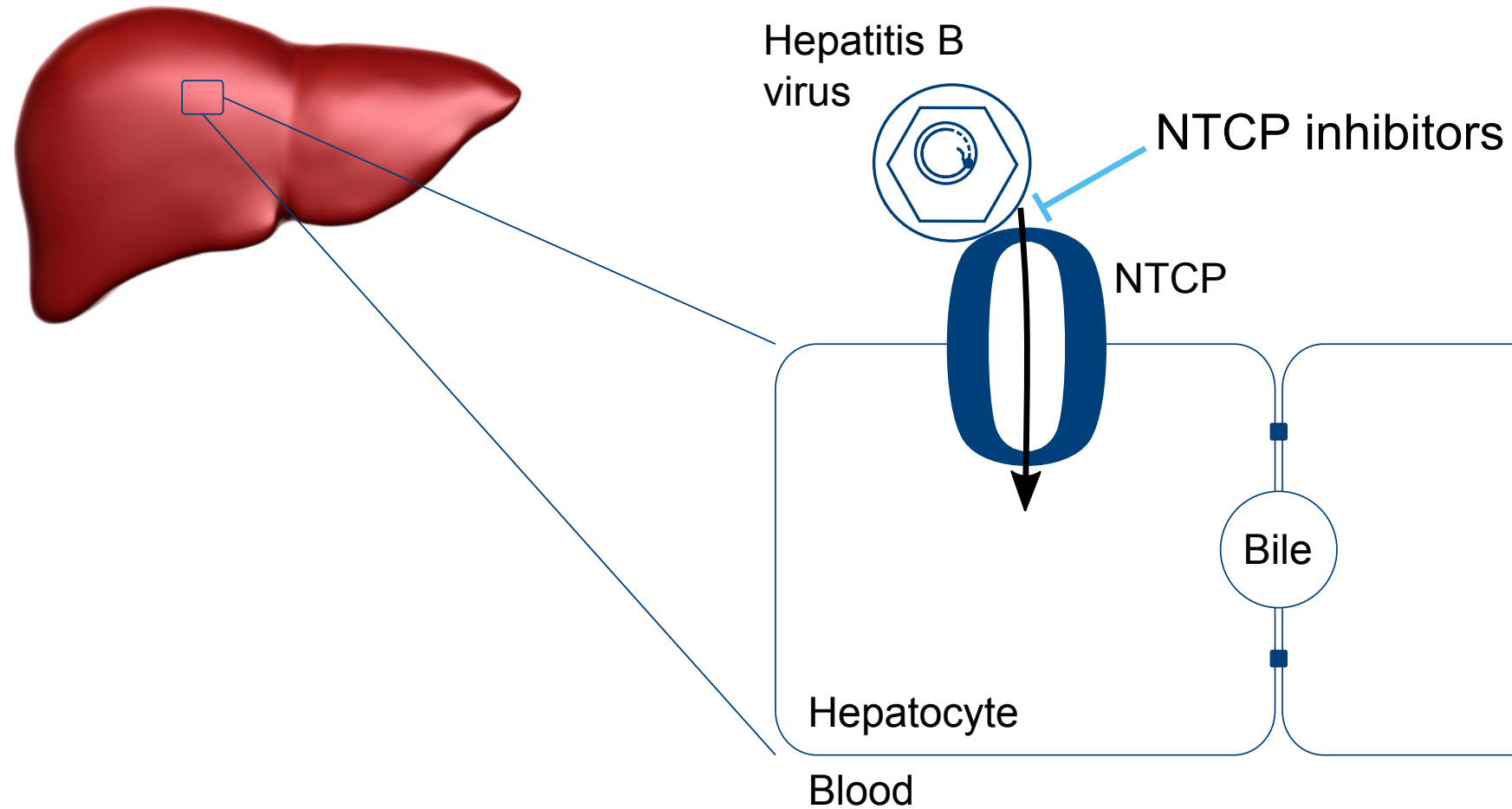
Results

Conclusion

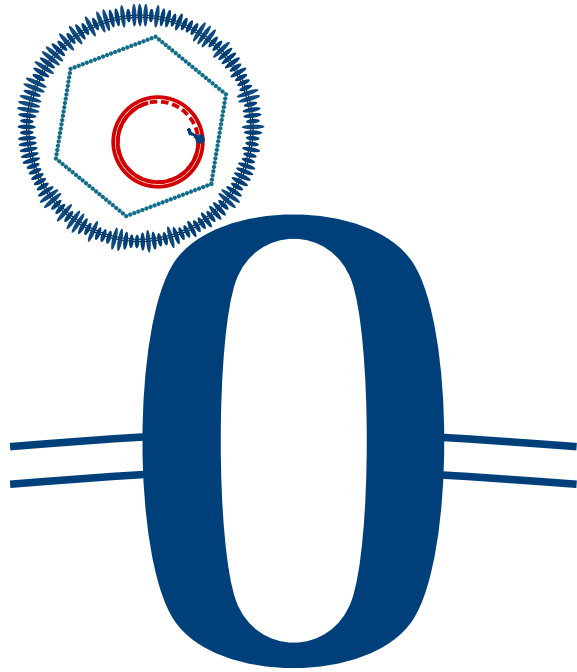
NTCP inhibition prevents HBV entry into hepatocytes



NTCP inhibition prevents HBV entry into hepatocytes



Optimization of an *in vitro* assay for identification of NTCP inhibitors



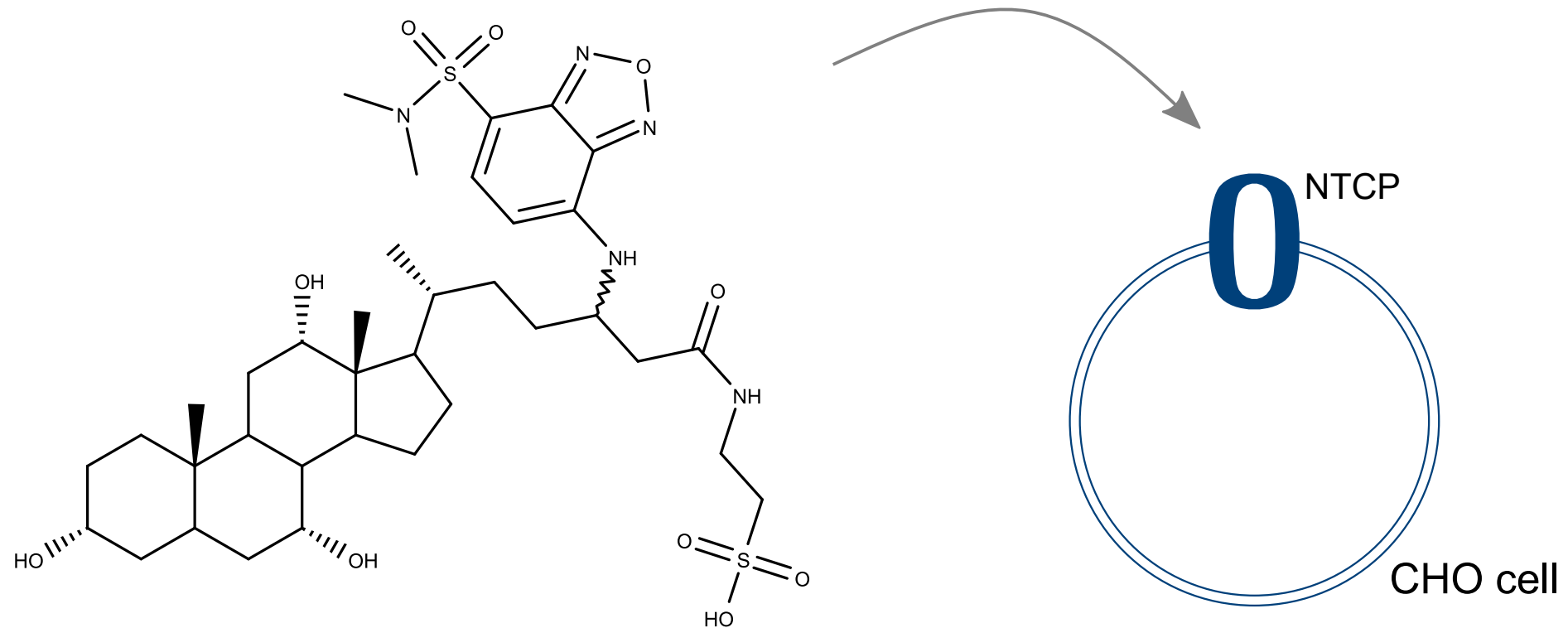
Rationale

Methods

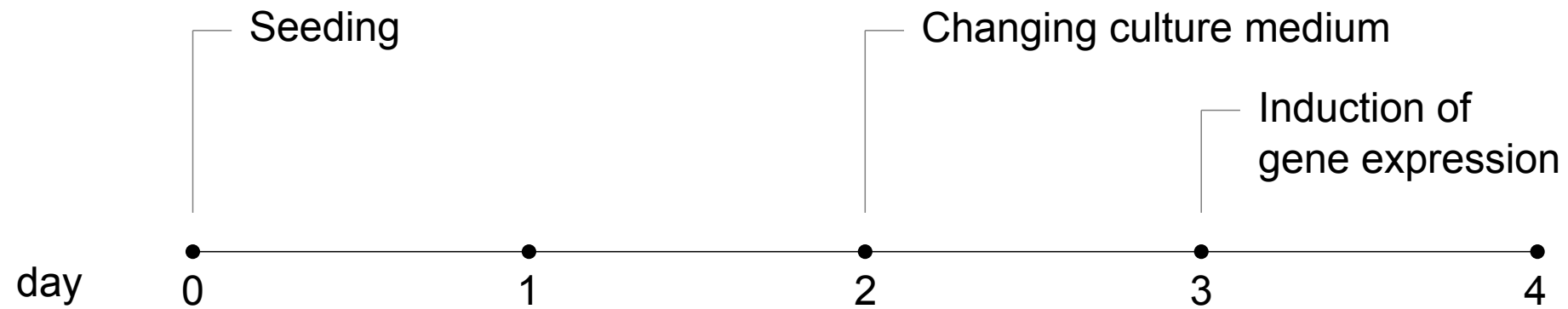
Results

Conclusion

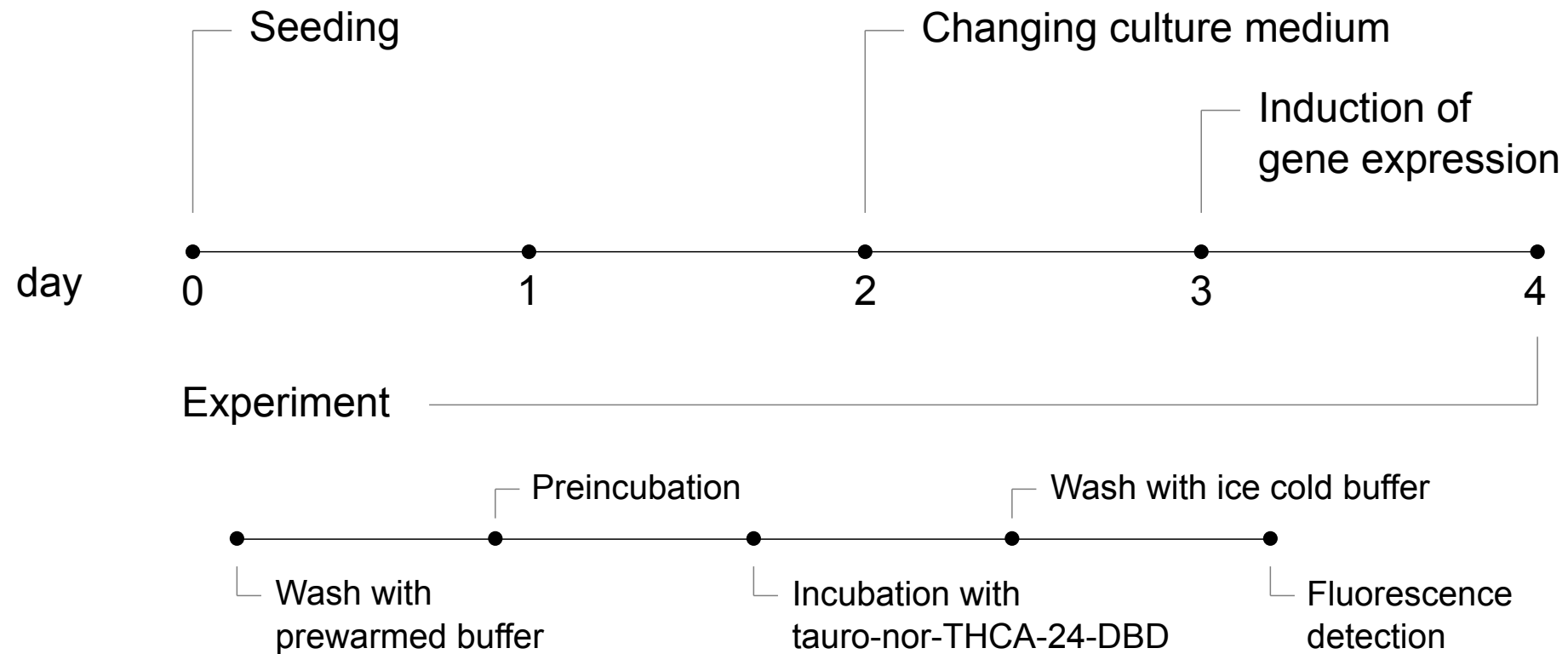
Uptake of the fluorescent NTCP substrate tauro-nor-THCA-24-DBD was determined in NTCP transfected CHO cells



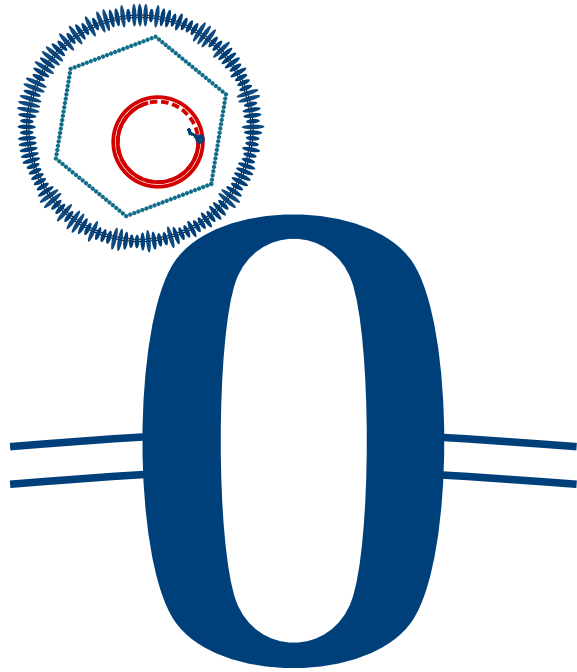
Uptake of the fluorescent NTCP substrate tauro-nor-THCA-24-DBD was determined in NTCP transfected CHO cells



Uptake of the fluorescent NTCP substrate tauro-nor-THCA-24-DBD was determined in NTCP transfected CHO cells



Optimization of an *in vitro* assay for identification of NTCP inhibitors



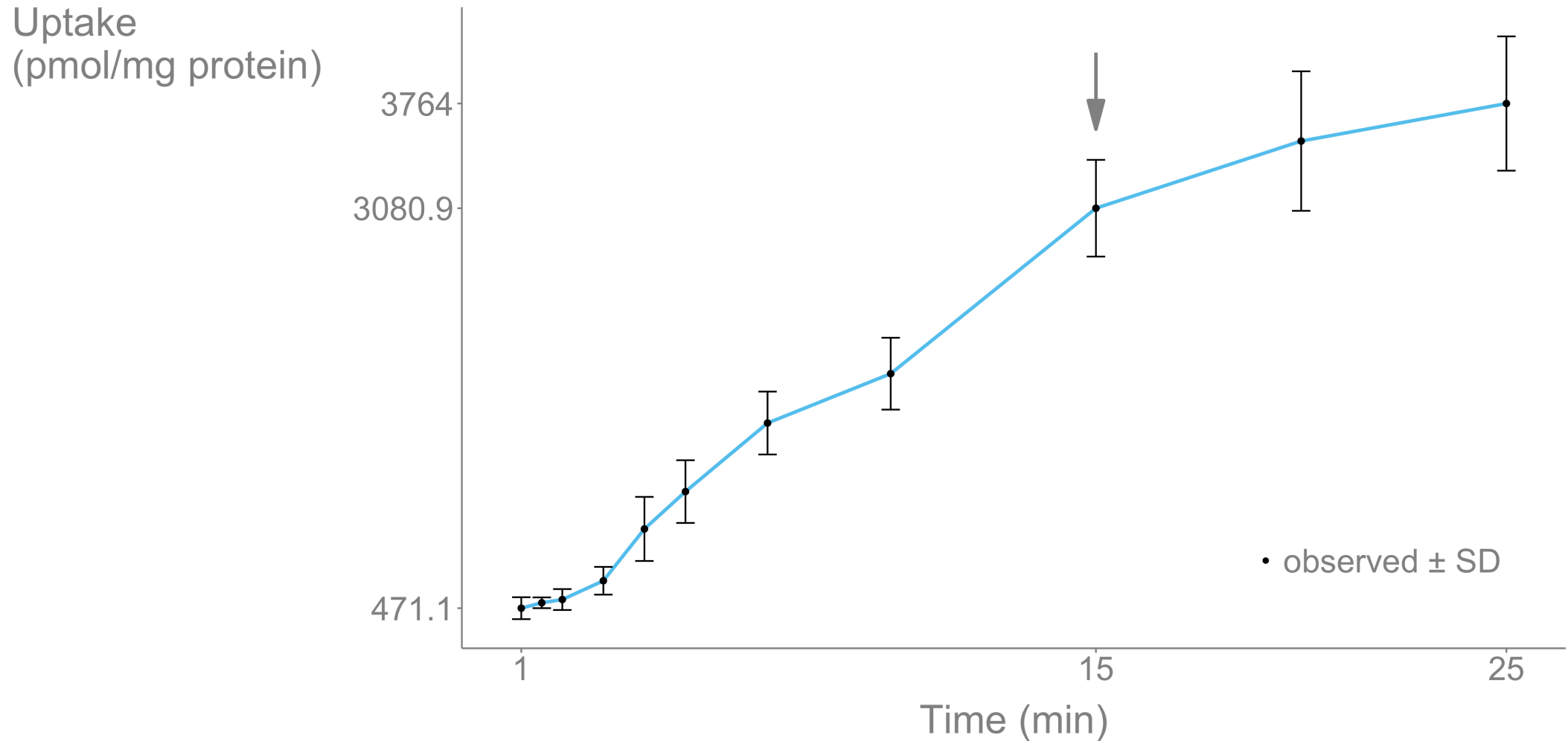
Rationale

Methods

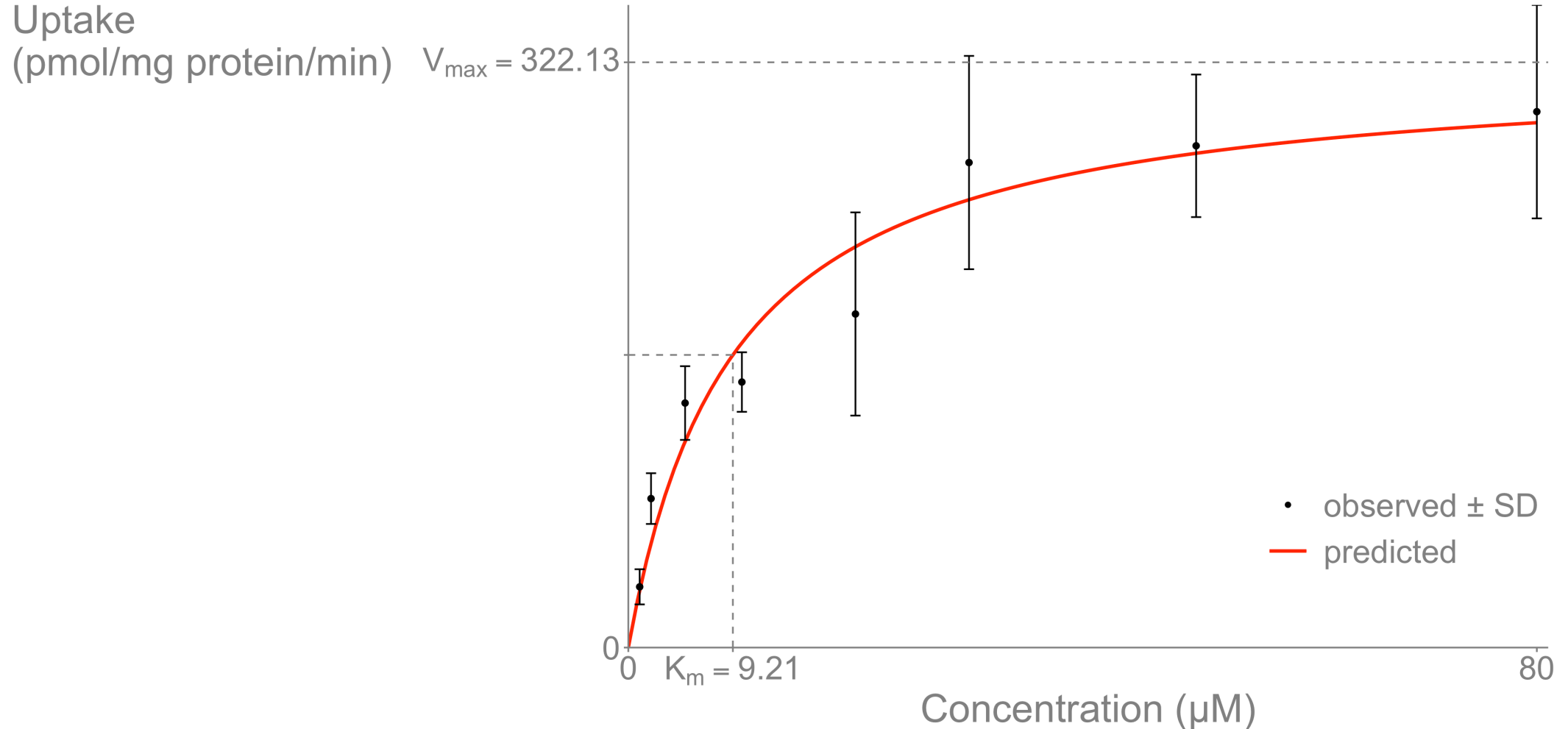
Results

Conclusion

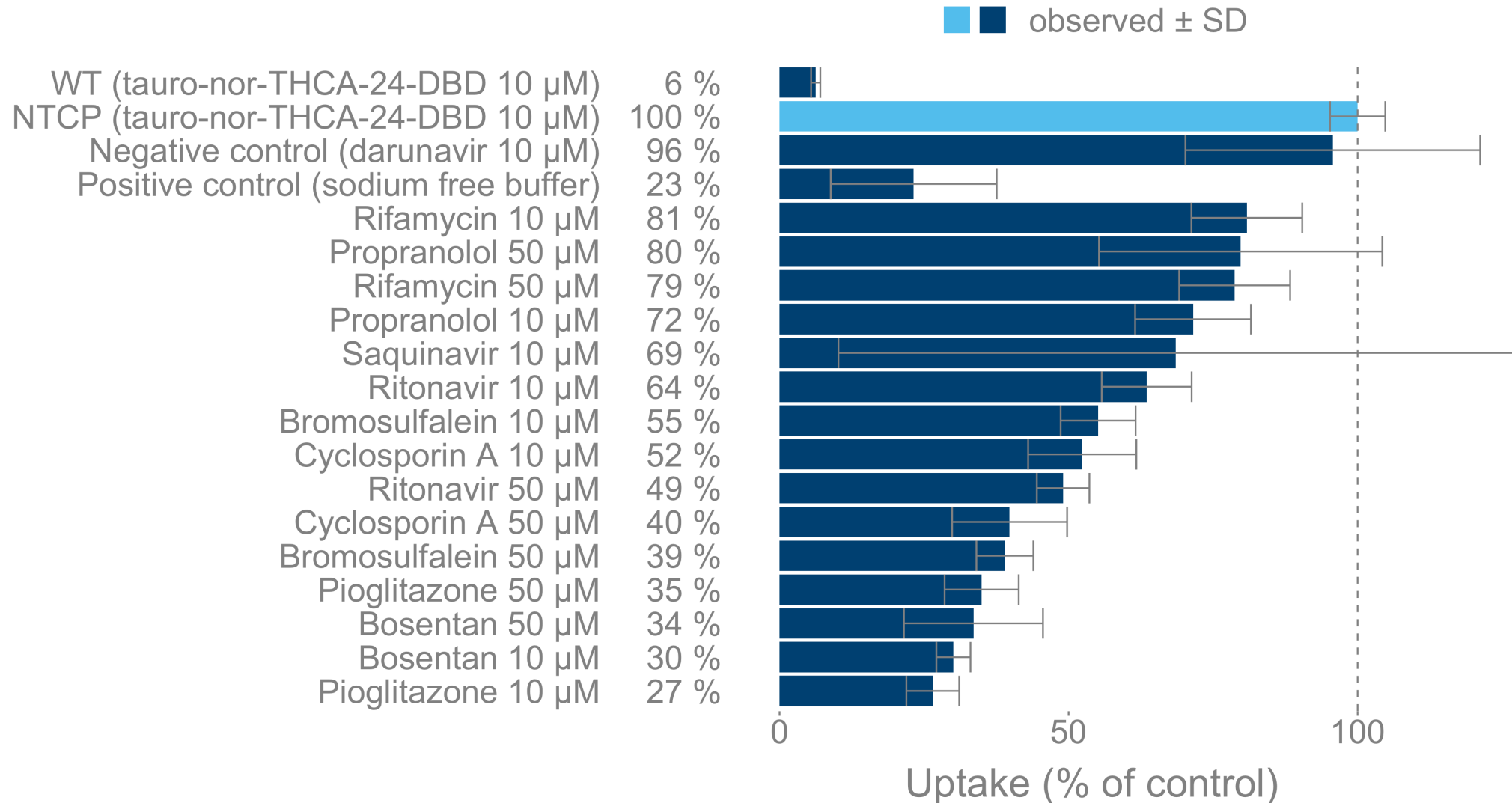
Uptake of 10 μM tauro-nor-THCA-24-DBD increased up to 20 minutes incubation



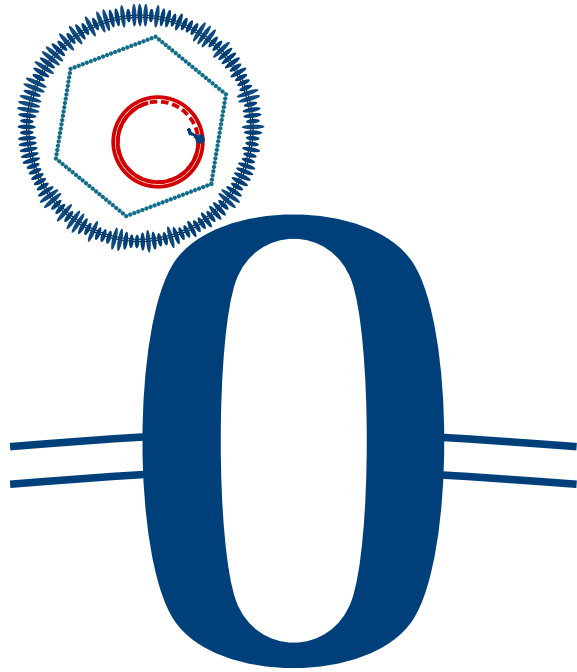
Concentration dependent uptake of tauro-nor-THCA-24-DBD (15 minutes) shows Michaelis-Menten kinetics



The assay was validated using known NTCP inhibitors



Optimization of an *in vitro* assay for identification of NTCP inhibitors



Rationale

Methods

Results

Conclusion

The assay was optimized and will be used for identification of novel NTCP inhibitors

Uptake in wild-type cells remained below the LOQ at all time points

Inhibitory potential of 2000 compounds will be explored

Uptake will expressed to the control of each batch

A (proteo)chemometric model will be developed

Acknowledgement



Laboratory of Liver Infectious Diseases



Prof. Bruno Stieger
University Hospital Zürich

Optimization of an *in vitro* assay for identification of NTCP inhibitors

Pieter Van Brantegem, Philip Meuleman and Pieter Annaert

fwo

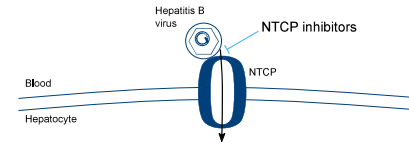
KU LEUVEN

Chronic hepatitis B virus (HBV) infection

affects up to 250 million people world wide
is a common cause of cirrhosis and hepatocellular carcinoma
cannot be eradicated yet

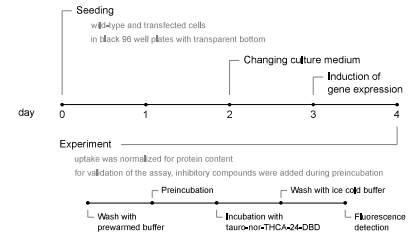
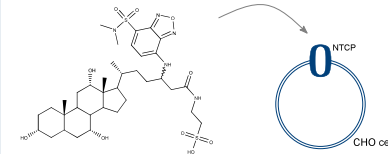
Sodium taurocholate co-transporting polypeptide (NTCP)

is a hepatic transport protein
is located at the sinusoidal membrane
transports bile acids from the blood into hepatocytes
has been identified as receptor for entry of the HBV into hepatocytes

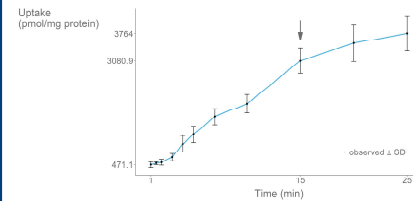


NTCP inhibition to prevent HBV entry has been shown to be successful → We need to find small molecules that inhibit NTCP
Therefore, we developed an *in vitro* assay that allows high throughput screening

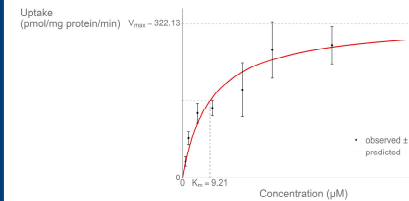
Uptake of the fluorescent NTCP substrate tauro-nor-THCA-24-DBD was determined in NTCP transfected CHO cells



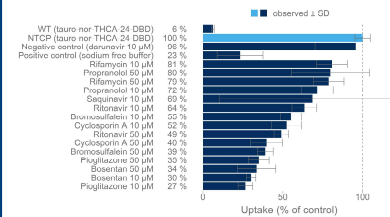
Uptake of 10 μ M tauro-nor-THCA-24-DBD increased up to 20 minutes incubation



Concentration dependent uptake of tauro-nor-THCA-24-DBD (15 minutes) shows Michaelis-Menten kinetics



The assay was validated using known NTCP inhibitors



The assay was optimized and will be used for identification of novel NTCP inhibitors

Uptake of tauro-nor-THCA-24-DBD in *wild-type* cells remained below the limit of quantification at all time points as a consequence, no correction for uptake in *wild-type* cells will be made

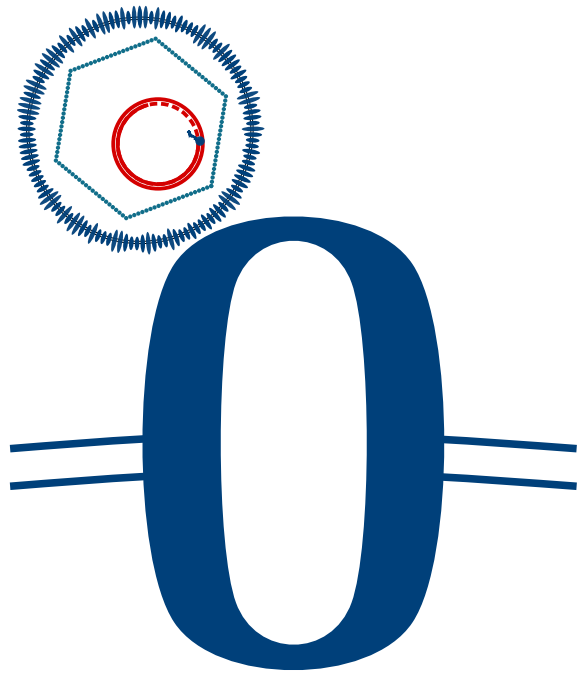
The inhibitory potential of 2000 compounds will be explored at equimolar concentrations of 10 μ M and 15 minutes incubation time

Uptake will be expressed relative to the control for each batch

Compounds will be considered strong inhibitors if at least 50% uptake inhibition can be obtained

Cyclosporin A 50 μ M will be used as positive control

Optimization of an *in vitro* assay for identification of NTCP inhibitors



Pieter Van Brantegem
KU Leuven

Meet the Experts Transporter
Conference Budapest
26 April 2018