



MEET ^{the} EXPERTS

TRANSPORTER CONFERENCE

BOSTON '19
CAMBRIDGE

SEPTEMBER 3-5



Closing remarks

Peter Krajcsi



LADME

| | Key message | Note |
|----------|--|--|
| LA | Understanding all LA parameters (pH dependent solubility, in stomach, precipitation in intestine, effect of micronization, efflux) is needed for right formulation strategy. | Betain HCL transiently restores pH and helps stomach pH dependent solubility |
| A | Precision-cut intestinal slices are good models to study P-gp | |
| D | Spheroid model > applicable for screening for CPPs and stapling enhances brain delivery | Calibration is needed for prediction of brain penetration |
| D | Good concordance between in vitro and in vivo approaches to predict brain exposure | Not applicable for zwitterions and acids (yet) |
| D | MDCKII-MDRI, MDCKII-BCRP monolayers are widely applied tools to predict brain penetration of drugs | |
| D | Specific inhibitors (Valspodar, Ko143) can be used to set-up chemical knock-out models for P-gp and BCRP | Ko143 is unstable in plasma so samples should be collected into chilled tubes containing NaF |
| (M) E | Good understanding of role of sinusoidal uptake by OATPs | Role of OAT2 and OCT1 is emerging |

DDI and Tox

| | Key messages | Notes |
|-----|---|--|
| DDI | Extensive understanding of P-gp-based tDDI | The ideal test system for DE is still investigated |
| DDI | Biomarkers > early DDI assesement | Full complexity should be looked at |
| DDI | Improved protocols (preincubation, calibration) provide more relevant data | Preincubation has an effect on IC50 data of multiple trps |
| DDI | Automation, high-speed bioanalysis and LIMS makes early application of ADME profiling possible | |
| DDI | HEK-BCRP membranes are superior to other BCRP-overexpressing membranes | |
| DDI | Accurate prediction of low fu values is possible | Current guidelines lead to overprediction of clinical DDI |
| Tox | High pos pred value (PPV) of hPTEC and high NPV of animal testing is best combination for nephrotox | Primary PTC cultures preserves expression of trps, including OATs |
| Tox | rBsep KD > best to monitor T3-BA and TCA-d4 in plasma | rBsep KD model > complex phenotype due to comp mech |
| Tox | ABCG2 inhibition (by SJ000831433?) may improve prognosis of high MYCN pediatric AML | Porhyrin inducing drugs > toxicity in patients with impaired ABCB6 |
| Tox | DILISym predicts liver injury > increasingly used in | |

Highlighted topics

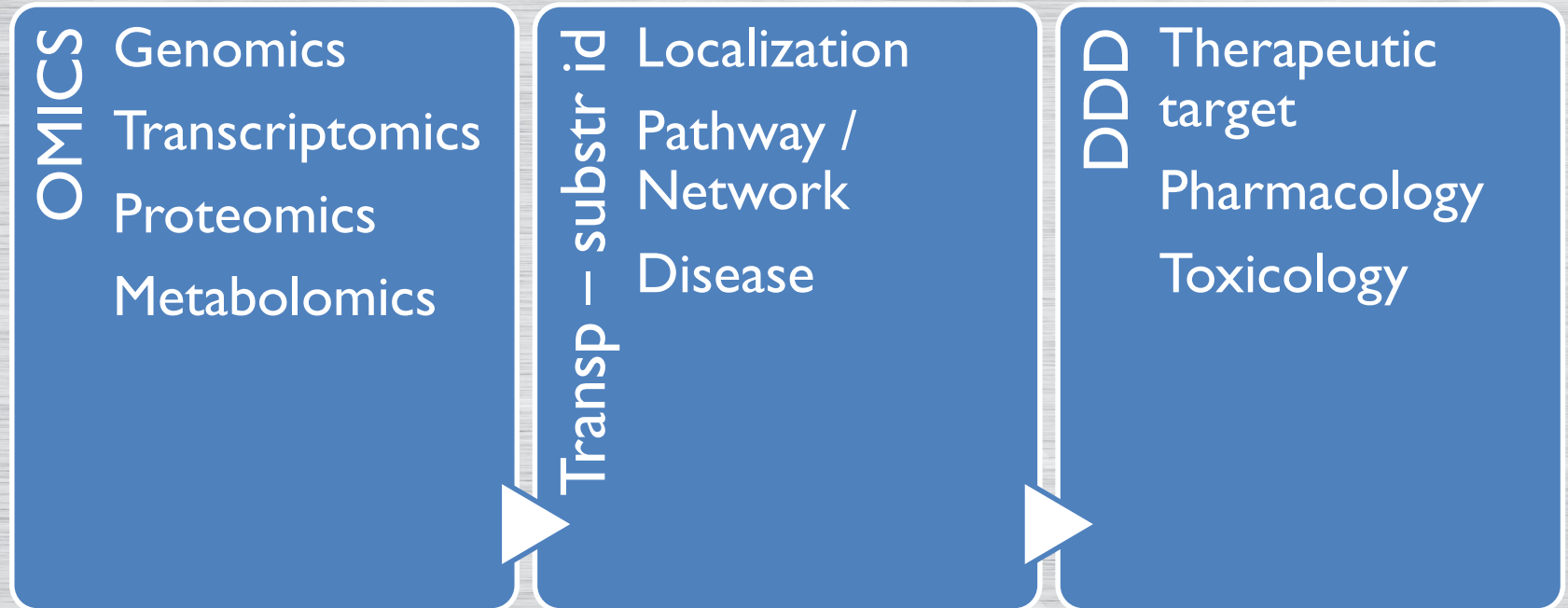
Microbiota

- Role in producing new metabolites (EPAC MII)
- Role in deconjugation of glucuronide conjugates of EPAC/metabolites
- Role in deconjugation (deamidation) and reduction of bile salts
- MoA of antibiotics induced DDI

Exosomes

- Communication between cells / organs
- Rich pool of biomarkers (represent protein, miRNA, etc profile of donor organs)
- Transfer of drug resistance is possible
- Way of tissue sampling

Technology and applications



Therapeutic target

Influx

- BBB OATPs transport statins through BBB > reduce ischemic damage
- LAT1 and ASCT2 > BBB penetration, cancer
- SLC30A10 > ameliorates excess of Mn
- Updated version of BDDCS illuminates druggable targets

Efflux

- ABCG2 inhibition potential cure for MYCN AML

Pharmacology – regulation of transporter activity

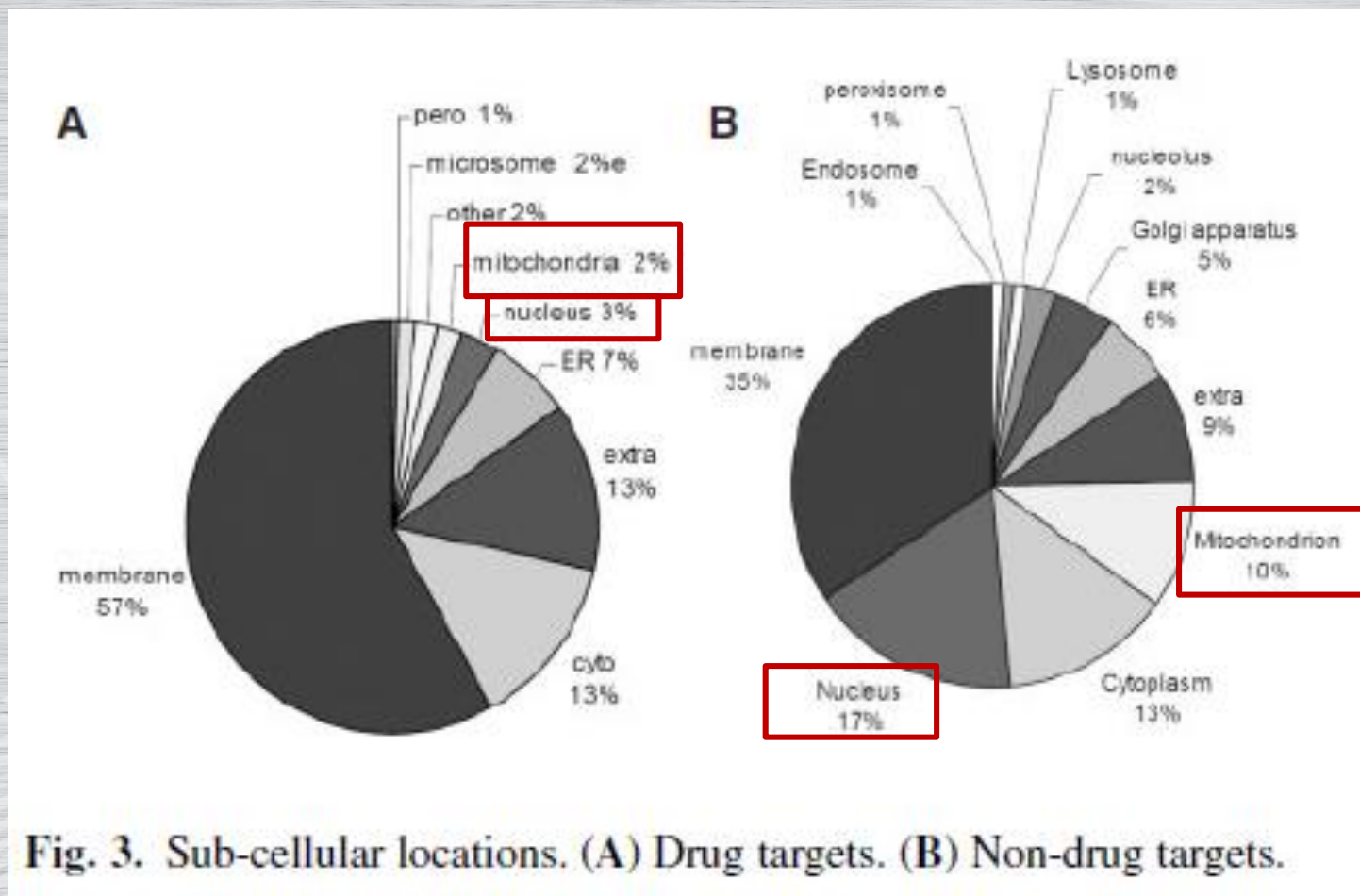
Inhibition

- **Non-covalent inhibitors (glifozins, urads, etc.)**
- Modulators of gene expression / activity (Crawford 2018 DMD)

Activation

- **Correctors (lumacraftor)**
- **Potentiators (ivacaftor)**
- Modulators of gene expression (ALK5 antagonists, ALK1 agonists > increased Oatp1a4 exp)
- Pathway modulators
 - GalNAc antisense to Tmprss6 > increased hepcidin > decreased Mn

Subcellular localization of drug targets



Toxicity /Adverse effects

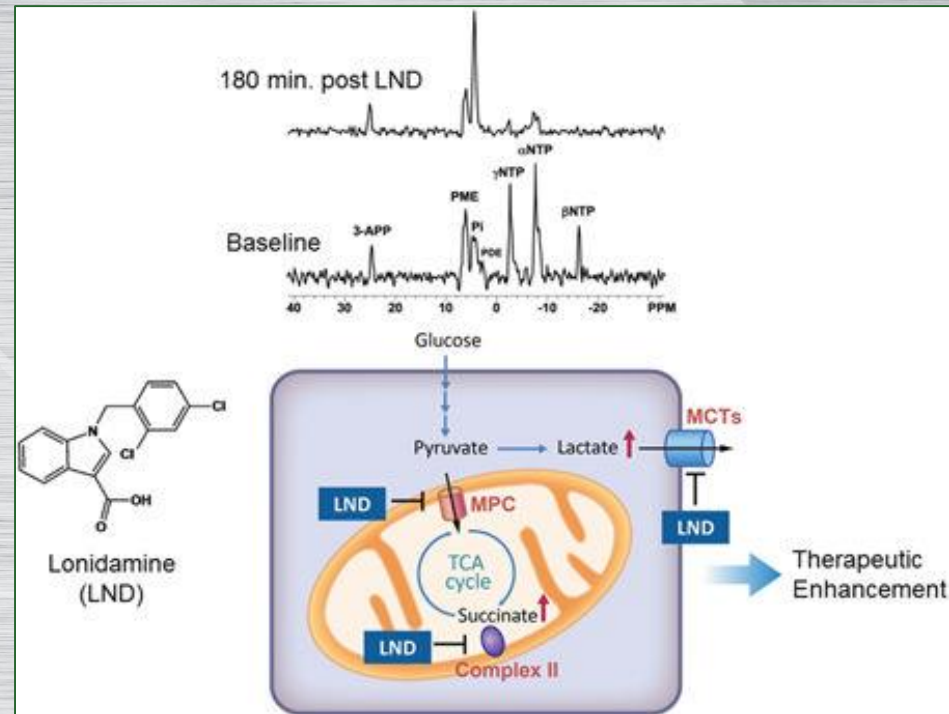
Physiological substrate

Modulation of transport
Modulation of metabolism

Impaired homeostasis

Toxicity
Adverse events

- Greater availability of data for plasma membrane transporters > plasma levels go up (bile acids/salts, urate, bilirubin, etc.)
- Fewer examples for intracellular transporters
 - Pb^{2+} , Mn^{2+} , Cd^{2+} enter mitochondria via Ca^{2+} channels > replace Ca^{2+} > toxicity
 - Lonidamine toxicity – not toxic to normal cells



Nath 2016 BBA

Transporter issues

Disease

- Transporters as therapeutic targets (BCRP, LAT1, ASCT2, Slc30a10)
- Transporters as determinants of PK of drugs (CNS drugs (P-gp, BCRP, Oatp1a4))
- Transporters as tissue/cell targeting of drugs (OCTN2, LAT1)

Preclinical testing

- in vitro / in vivo with reference to PK/PD properties (BBB models (transfectants, spheroids, in vivo), absorption (transfectants, precision cutslices)) excretion (kidney)
- tDDI tests (P-gp, OATPs, OCTs), biomarkers (cynos)
- Tox profile (BSEP, ABCB6, various kidney transporters, serotonin reuptake)
- Modeling / simulations (docking studies (LAT1, ASCT2), IVIVE (BBB), PBPK

Clinical testing

- tDDI (P-gp) / PGx / biomarkers
- Toxicity (hepatotoxicity, nephrotoxicity)
- PK / PD (plasma vs tissue levels)

Thanks for attending
Special thanks to presenters

Have a safe trip home

SW

Strength

- Understanding LADME parameters > right formulation strategy
- Extensive understanding of P-gp with regard DDI
- High positive predictive value (PPV) of hPTEC and high NPV of animal testing
- rBsep KD > best to monitor T3-BA and TCA-d4 in plasma
- Spheroid model > applicable for screening for CPPs
- Good concordance between in vitro and in vivo approaches to predict brain exposure
- Understanding role of sinusoidal uptake
- Improved protocols (preincubation, calibration) provide more relevant data

Weakness

- rBsep KD model > complex phenotype due to comp mech
- BBB spheroids > low dynamic range for efflux trp effects
- Not applicable for zwitterions and acids
- Preincubation has an effect on IC50 data of multiple trps

Opportunities

- Tools available to study molecular details of ADMETox (EPAC)
- Biomarkers > early DDI assessment
- Potentiation of Oatp-mediated uptake of statins > stroke treatment
- A selection of IBI/IB3 probes / inhibitors offer greater relevance of DDI studies
- Accurate prediction of fu values is possible
- ABCG2 inhibition > potential to cure MYCN driven AML
- In silico LAT1 and ASCT2 work tools to study reprogramming metabolic networks and developing leads
- Targeting NPs to ATB^{0,+} and OCTN2 increased uptake of nanoparticles
- DILI Sym predicts liver injury > increasingly used in reg submissions
- Automation, high-speed bioanalysis and LIMS makes early application of ADME profiling possible

Threats

- Potential interference with homeostasis of endogenous substrate (was not the case for EPAC)
- Current guidelines tend to overpredict clinical DDI
- Porphyrin inducing drugs > toxicity in patients with reduced/inhibited ABCB6 function