

# The manganese efflux transporter SLC30A10 is essential for gastrointestinal manganese excretion

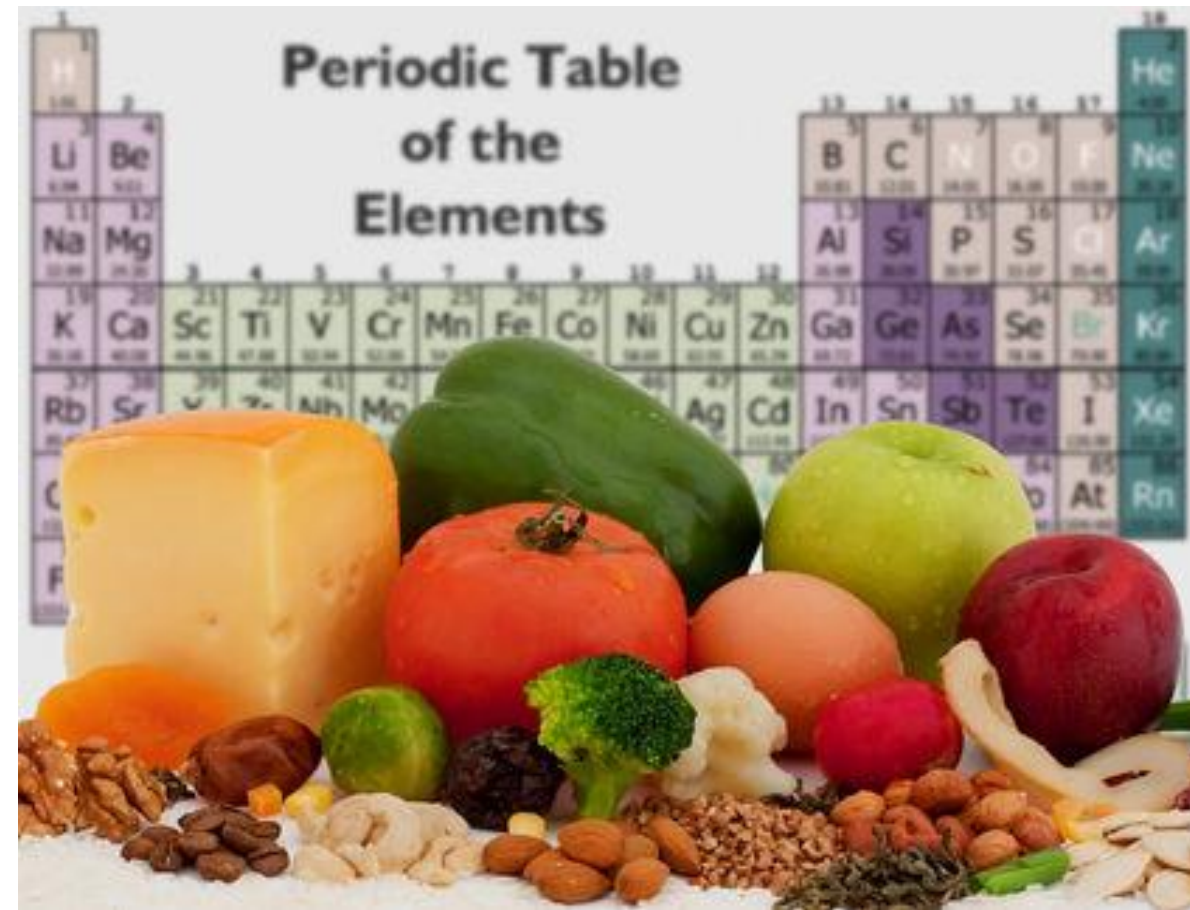
Courtney Mercadante, Ph.D.

Bartnikas Laboratory

Brown University, Providence, RI

September 5, 2019

# Acquisition of Manganese (Mn)

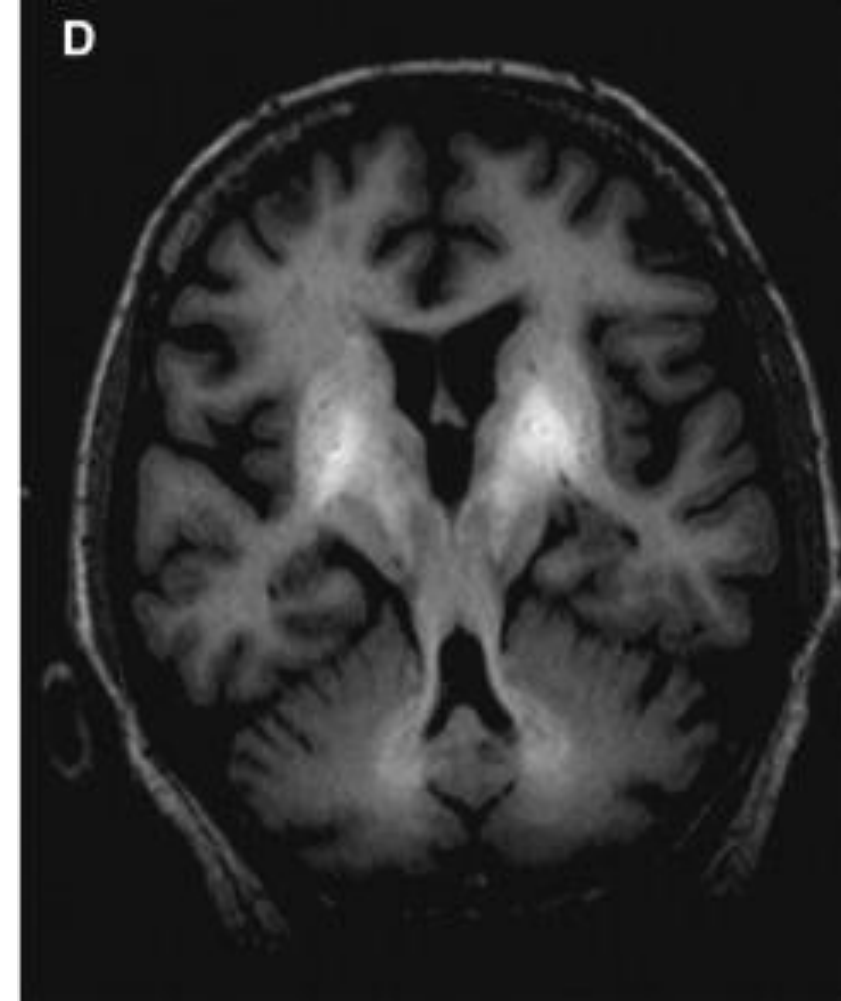


- Cofactor of essential proteins
- Oxidant defense, bone growth, immunity
- Men and women require 2.3 and 1.8 mg/day, respectively
- Plentiful dietary sources
- Only 3-5% of dietary Mn is absorbed
- Mn deficiency is very rare
- Mn related diseases are often diseases of excess

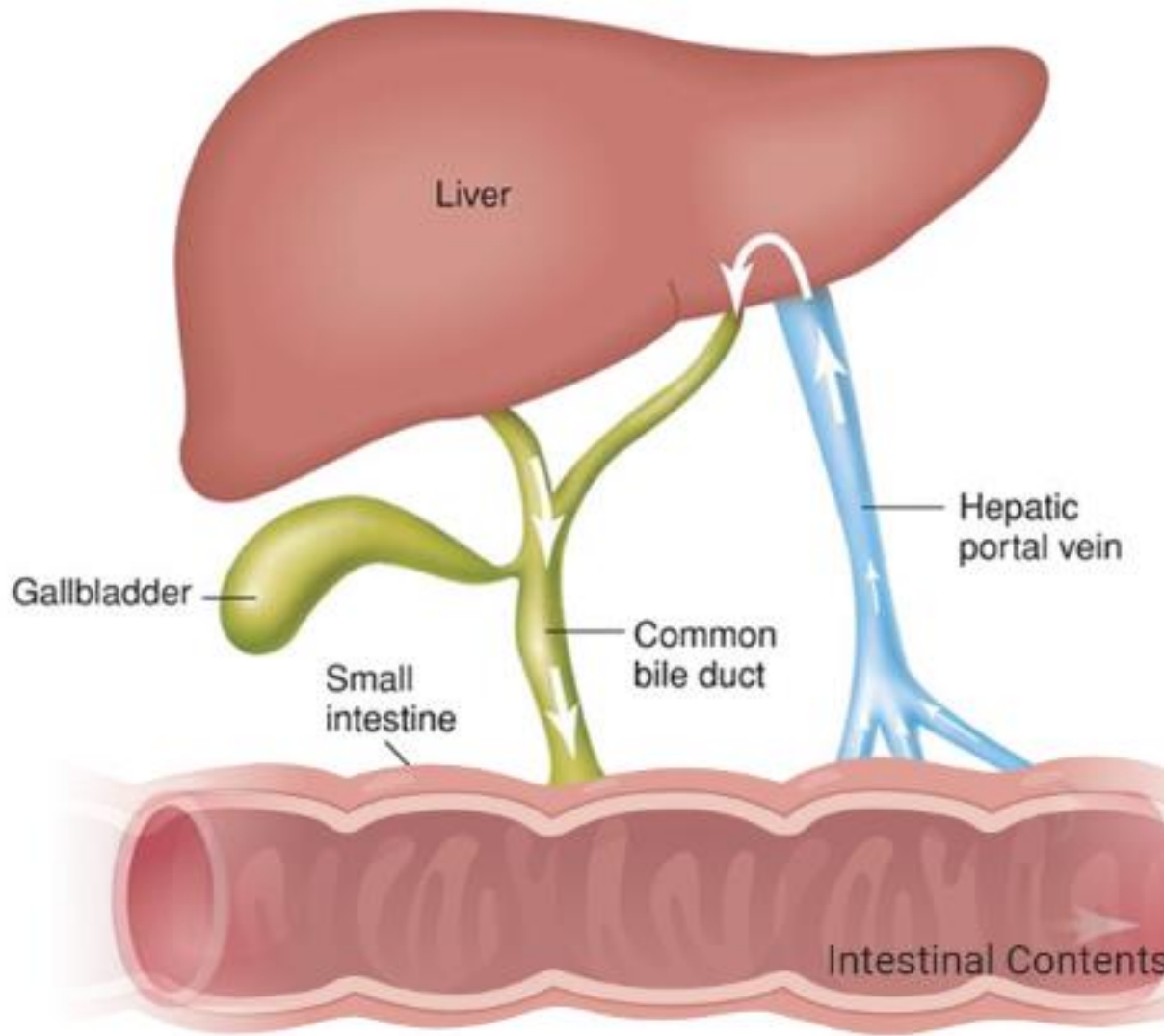
<http://www.chemistryexplained.com/elements/L-P/Manganese.html>

# Manganese in excess is toxic

- Mitochondrial dysfunction, compromised antioxidant defense mechanisms
- Manganism: Parkinson's- like disorder
  - Behavioral changes, slow and clumsy movements, tremors, cock-walk gait

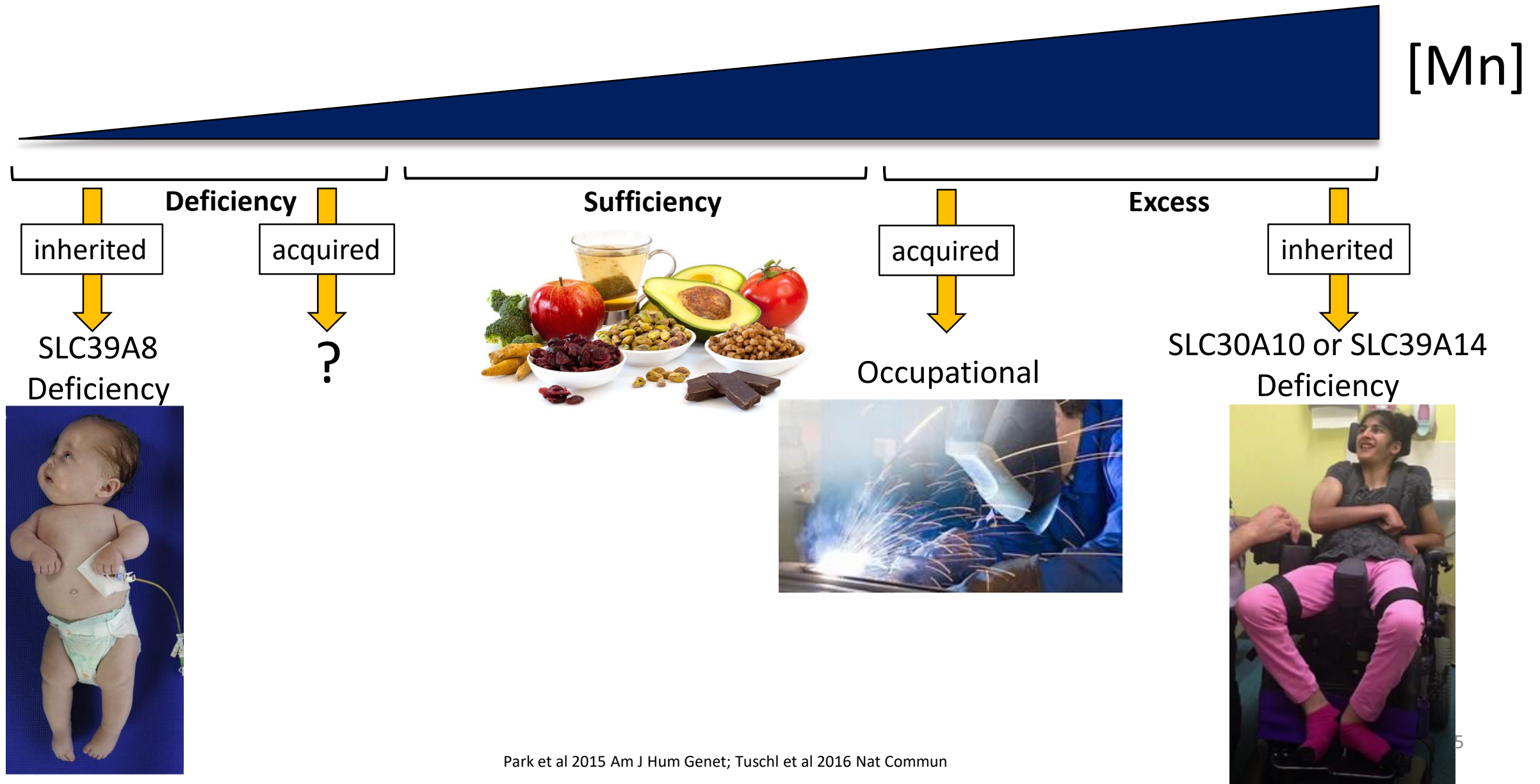


# Physiological Mechanisms of Mn Homeostasis

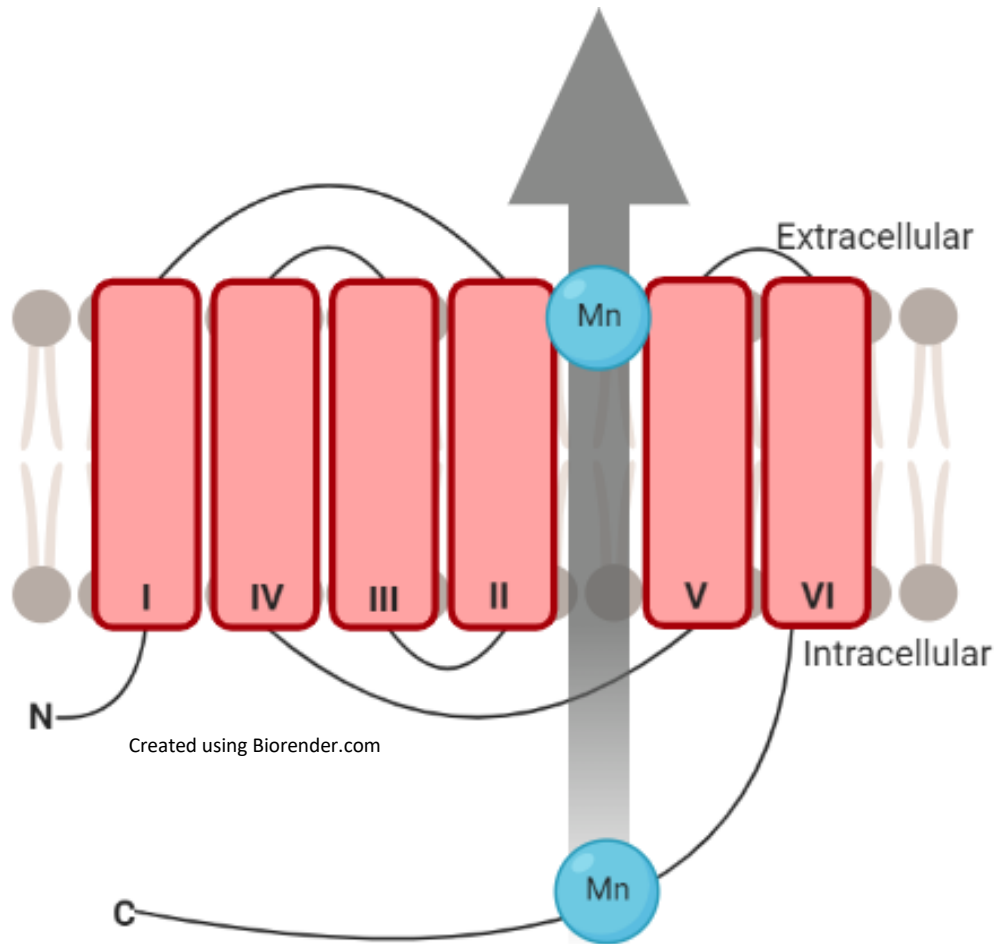


Modified from: [https://www.78stepshealth.us/human-physiology/images/3204\\_738\\_974-liver-gallbladder-portal-vein-intestines.jpg](https://www.78stepshealth.us/human-physiology/images/3204_738_974-liver-gallbladder-portal-vein-intestines.jpg), Created using Biorender.com

# Manganese excess: Acquired vs. Inherited



# Manganese excess: Acquired vs. Inherited

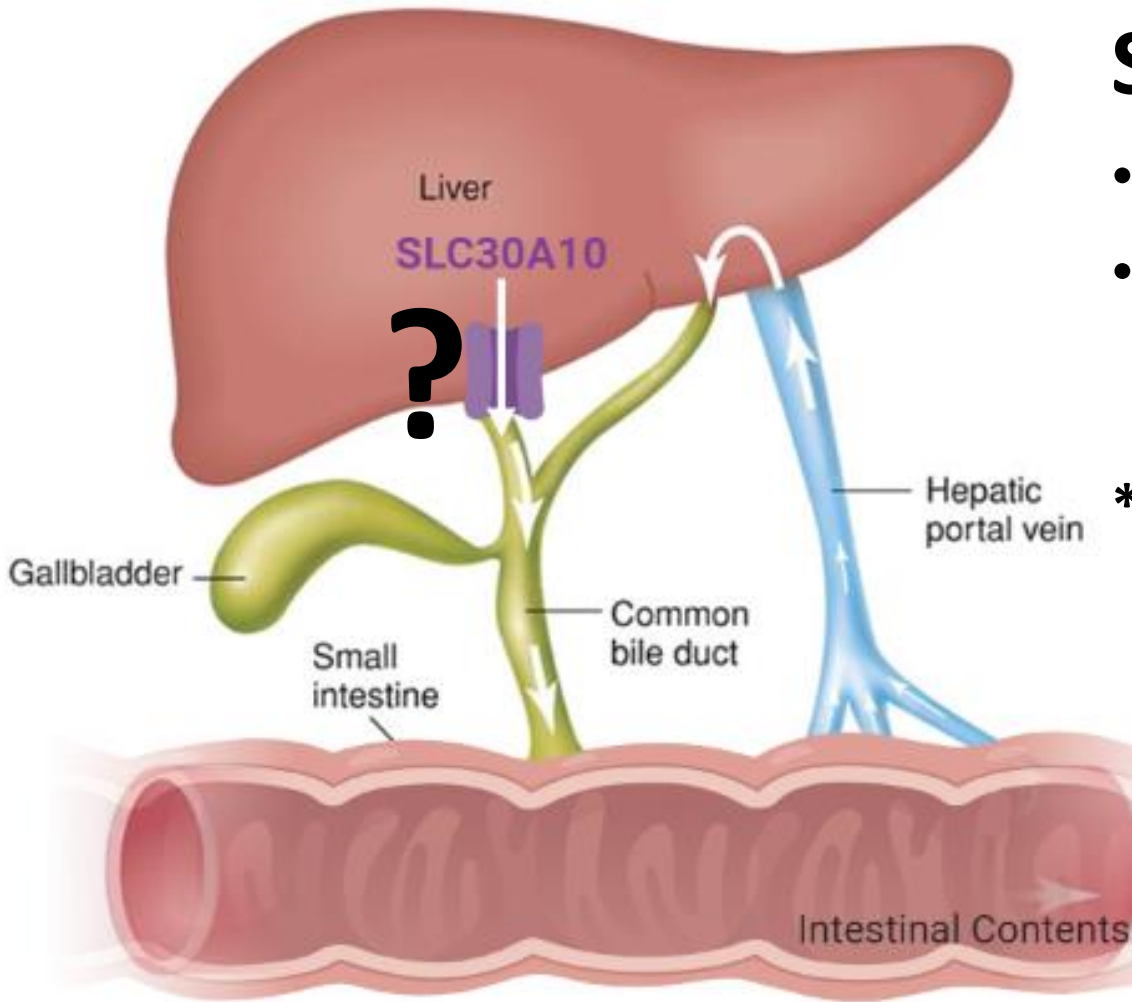


## SLC30 family of proteins

- Originally thought to be a zinc transporter (Znt10)
- SLC30 proteins- cation diffusion facilitator superfamily of metal transporters
  - Transport divalent metal cations (Fe, Zn, Cu, Ni, Co, Cd, Mn)
  - Efflux metals from cytosol to extracellular space
- 6 transmembrane domains with a cytoplasmic N and C termini



# Manganese excess: Acquired vs. Inherited



## SLC30A10 deficiency

- Autosomal recessive
- Mutations in *SLC30A10* identified in patients with hepatic cirrhosis, polycythemia, manganism, and high Mn levels

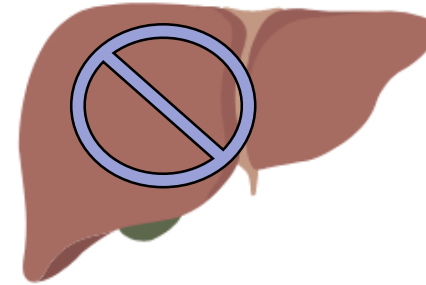
\* **No history of environmental Mn exposure**

# Generating Slc30a10-deficient mice

Whole body Slc30a10 deficiency  
*Slc30a10<sup>KO/KO</sup>*

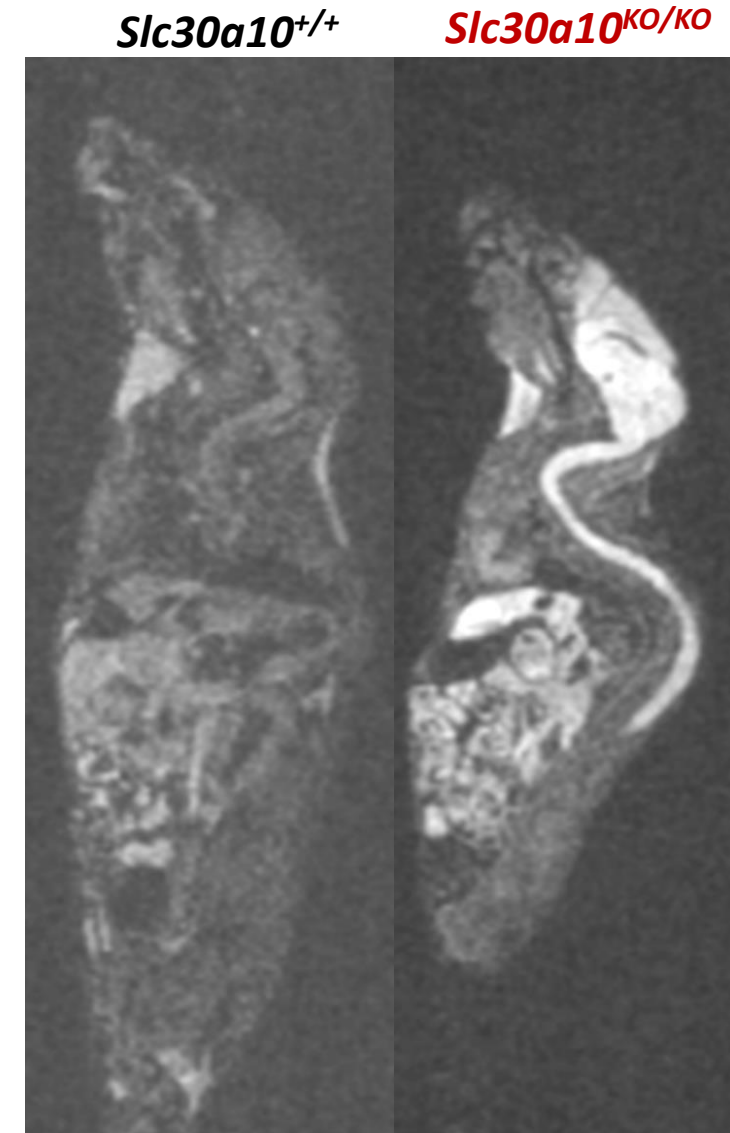
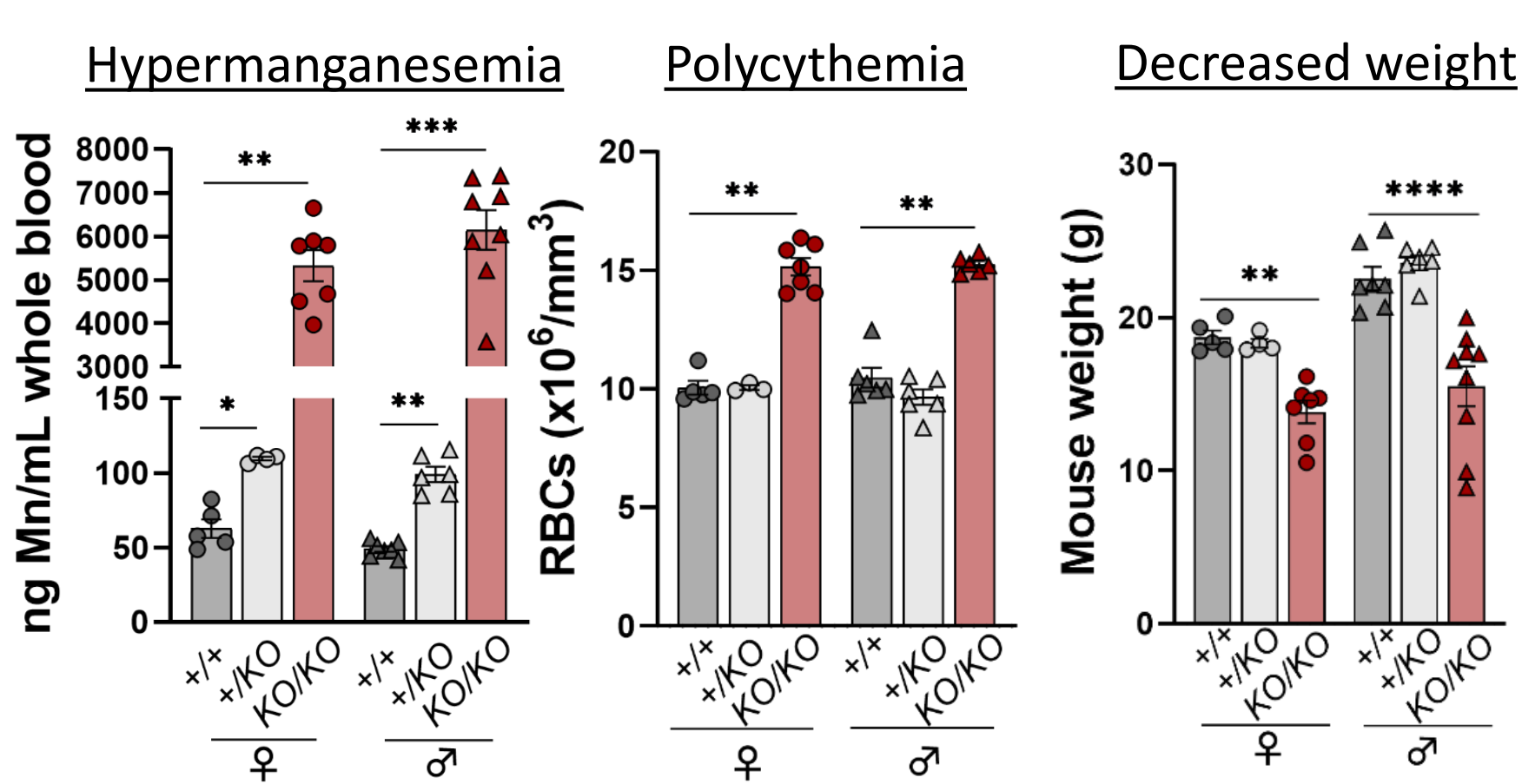


Hepatic  
Slc30a10 deficiency  
*Slc30a10<sup>lox/lox</sup> Alb*

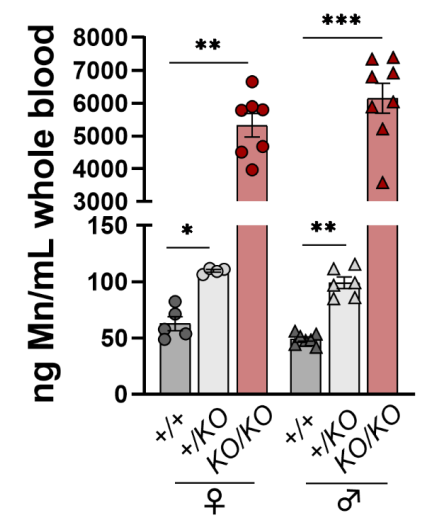
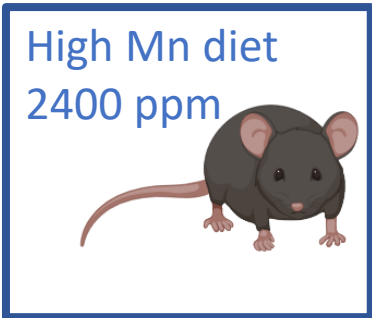
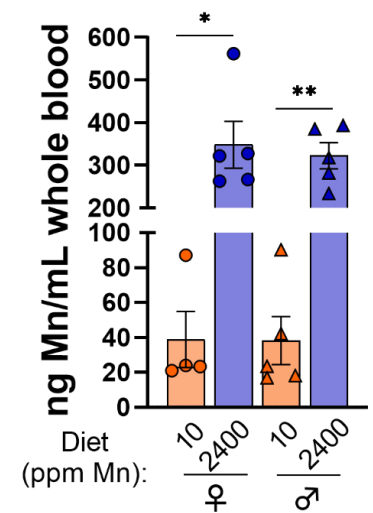




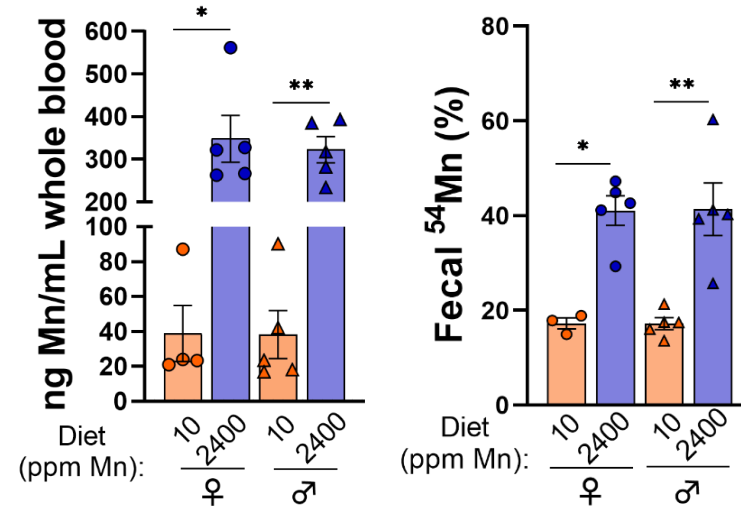
# Slc30a10-deficient mice recapitulate human disease phenotypes



# Wild-type mice on high Mn diet accumulated high Mn levels



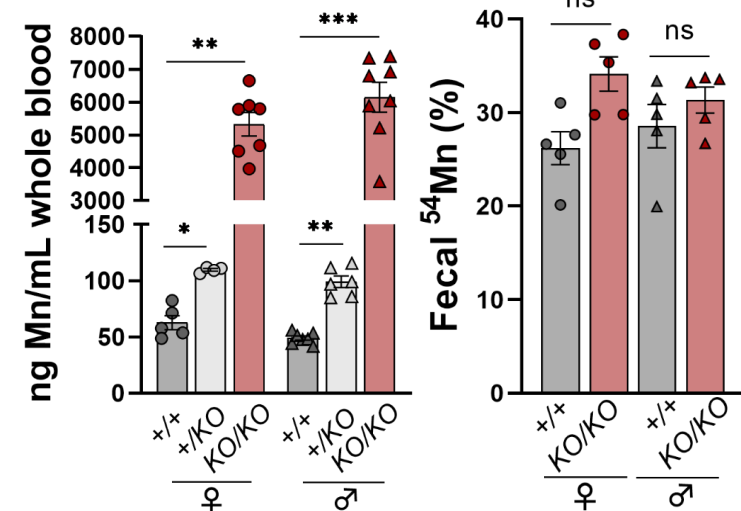
# Slc30a10-deficient mice have impaired Mn excretion



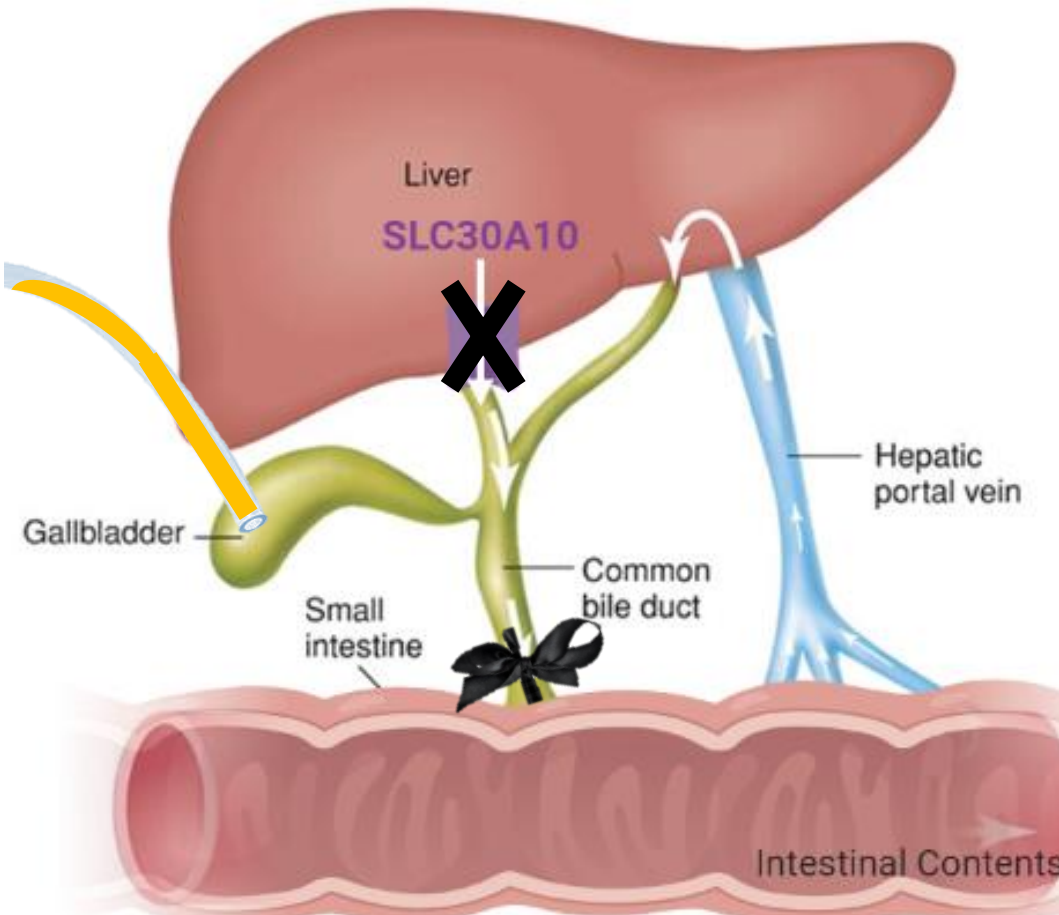
Control Mn diet  
10 ppm



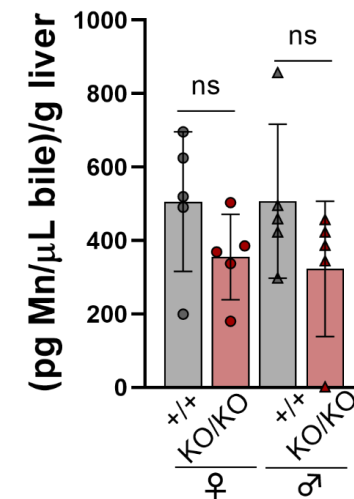
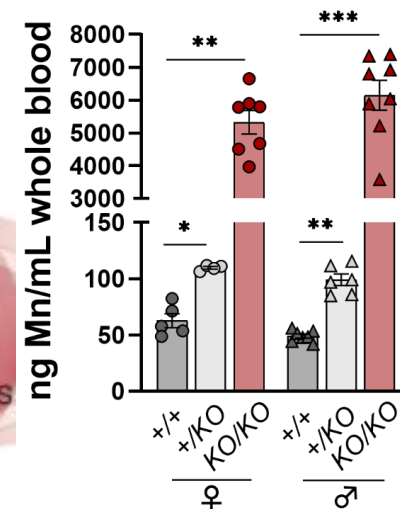
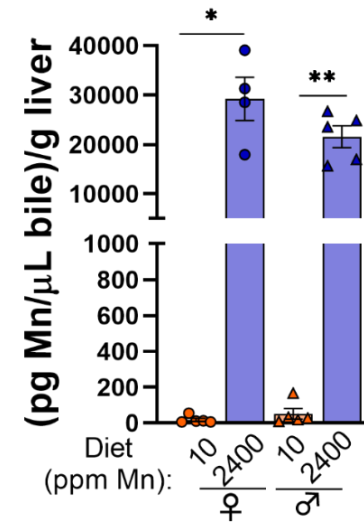
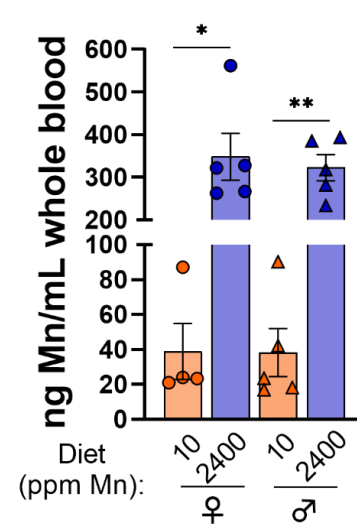
High Mn diet  
2400 ppm



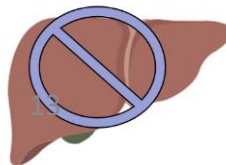
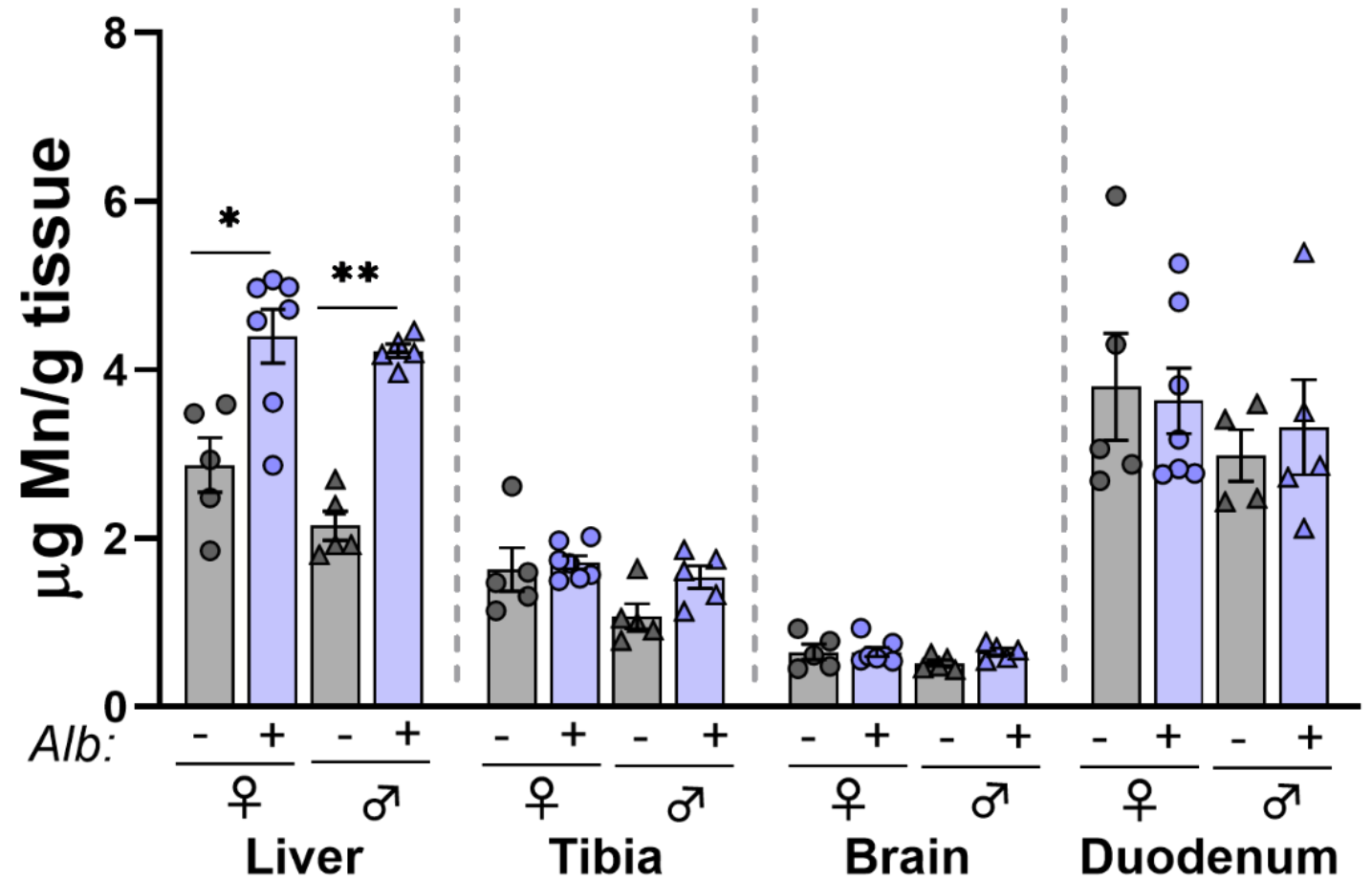
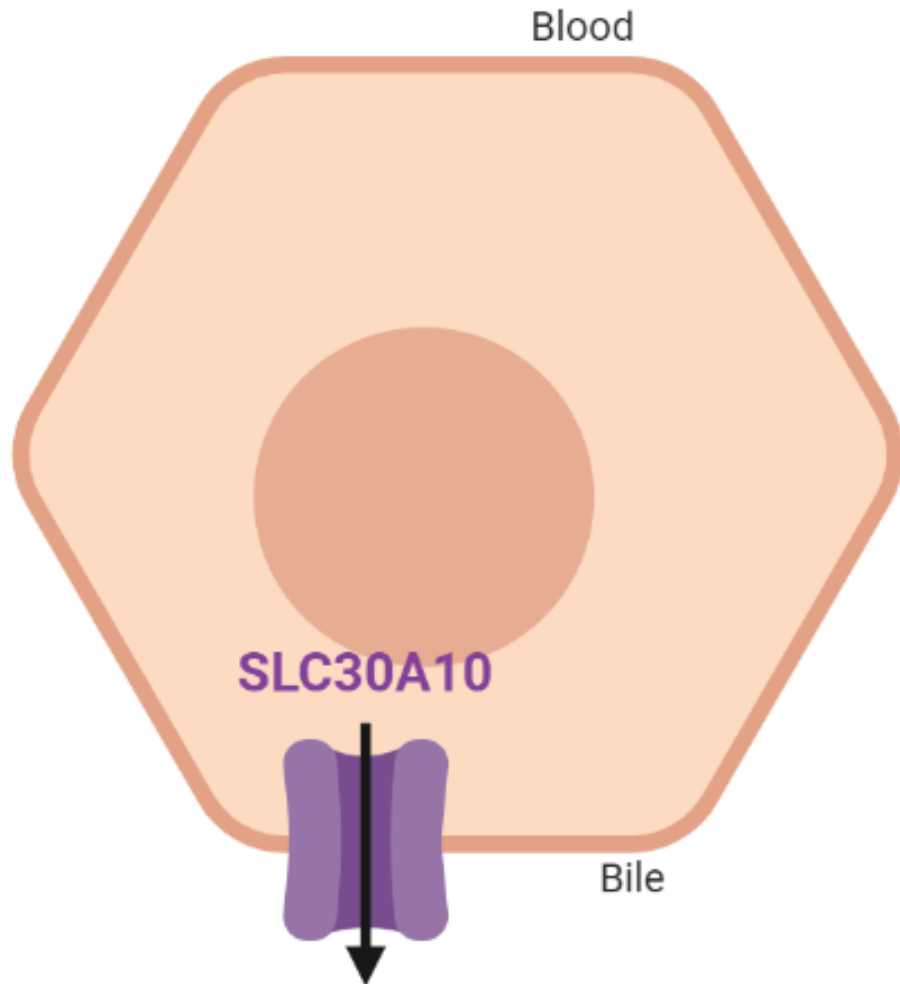
# Slc30a10-deficient mice have impaired Mn excretion



Modified from: [https://www.78stepshealth.us/human-physiology/images/3204\\_738\\_974-liver-gallbladder-portal-vein-intestines.jpg](https://www.78stepshealth.us/human-physiology/images/3204_738_974-liver-gallbladder-portal-vein-intestines.jpg). Created using Biorender.com

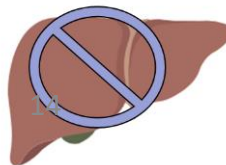
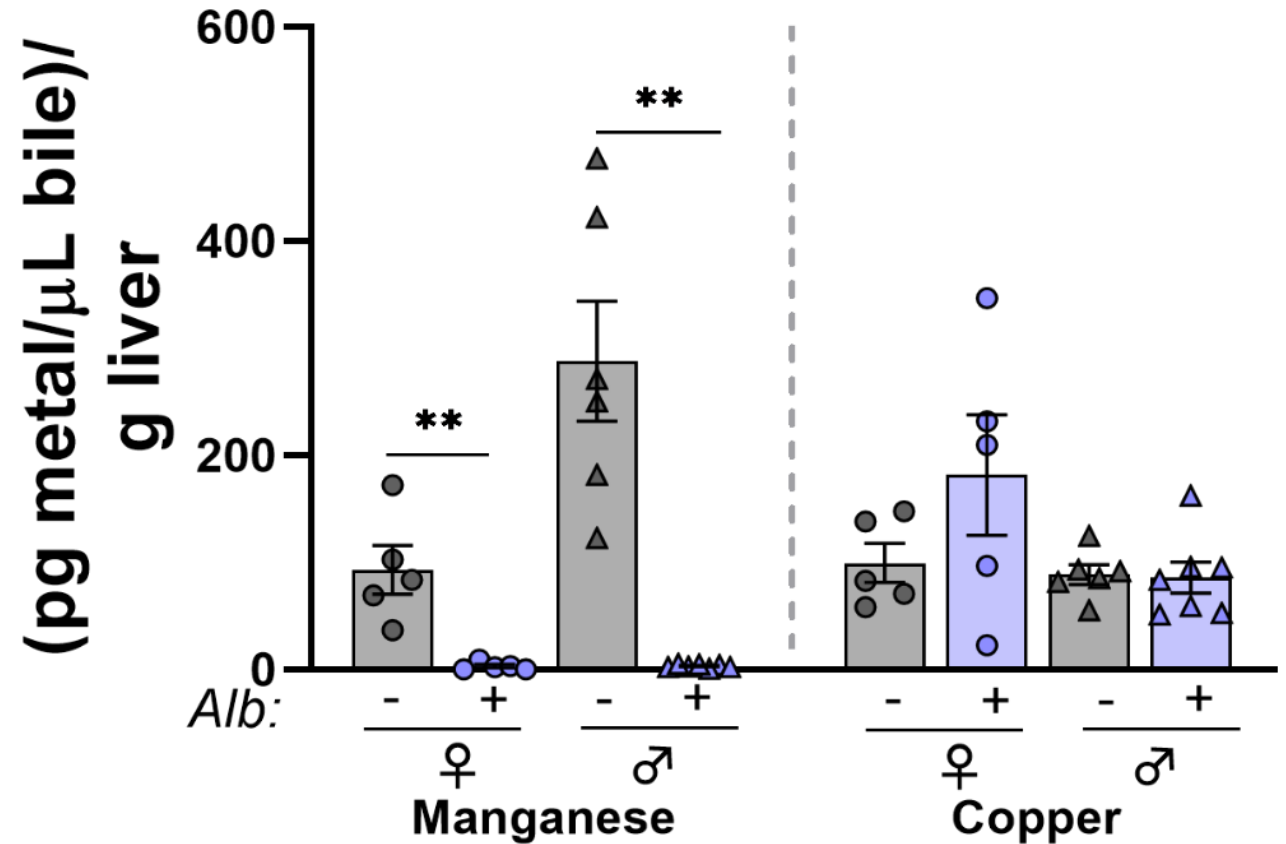
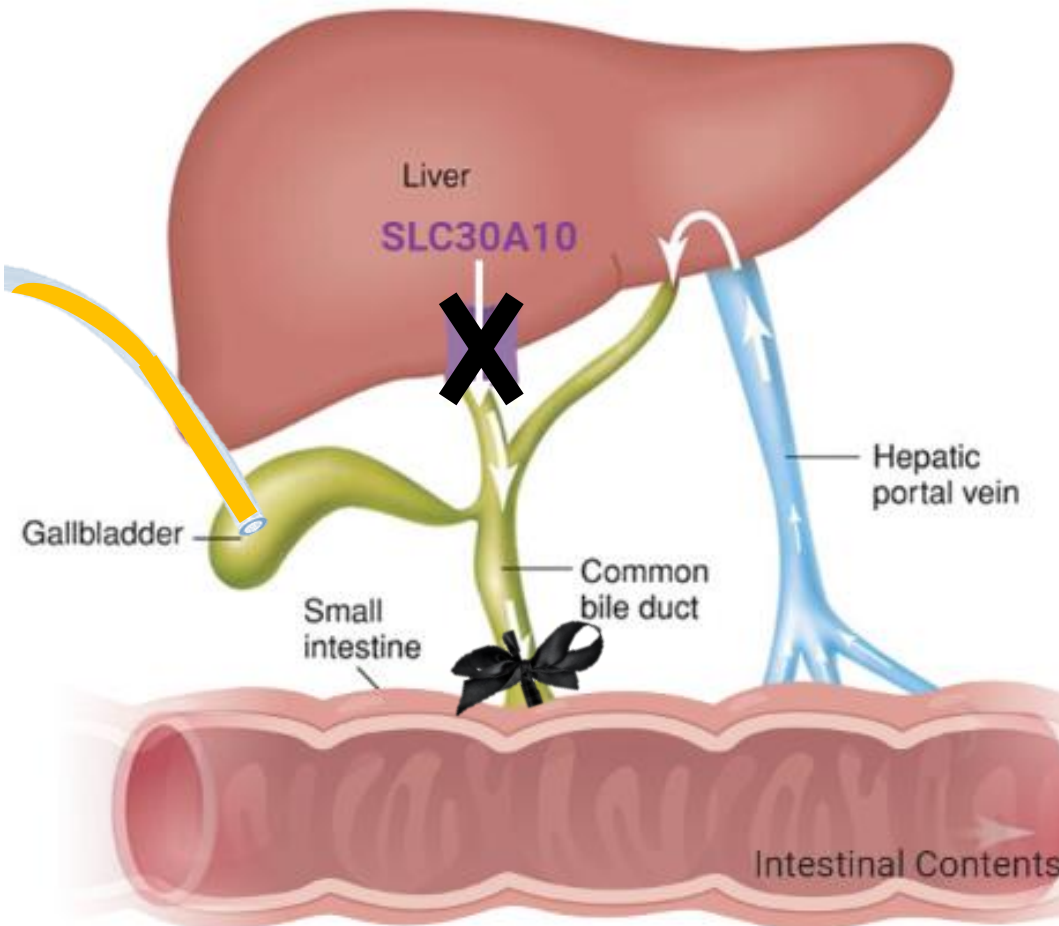


# Hepatic-Slc30a10 deficiency leads to minimal Mn excess





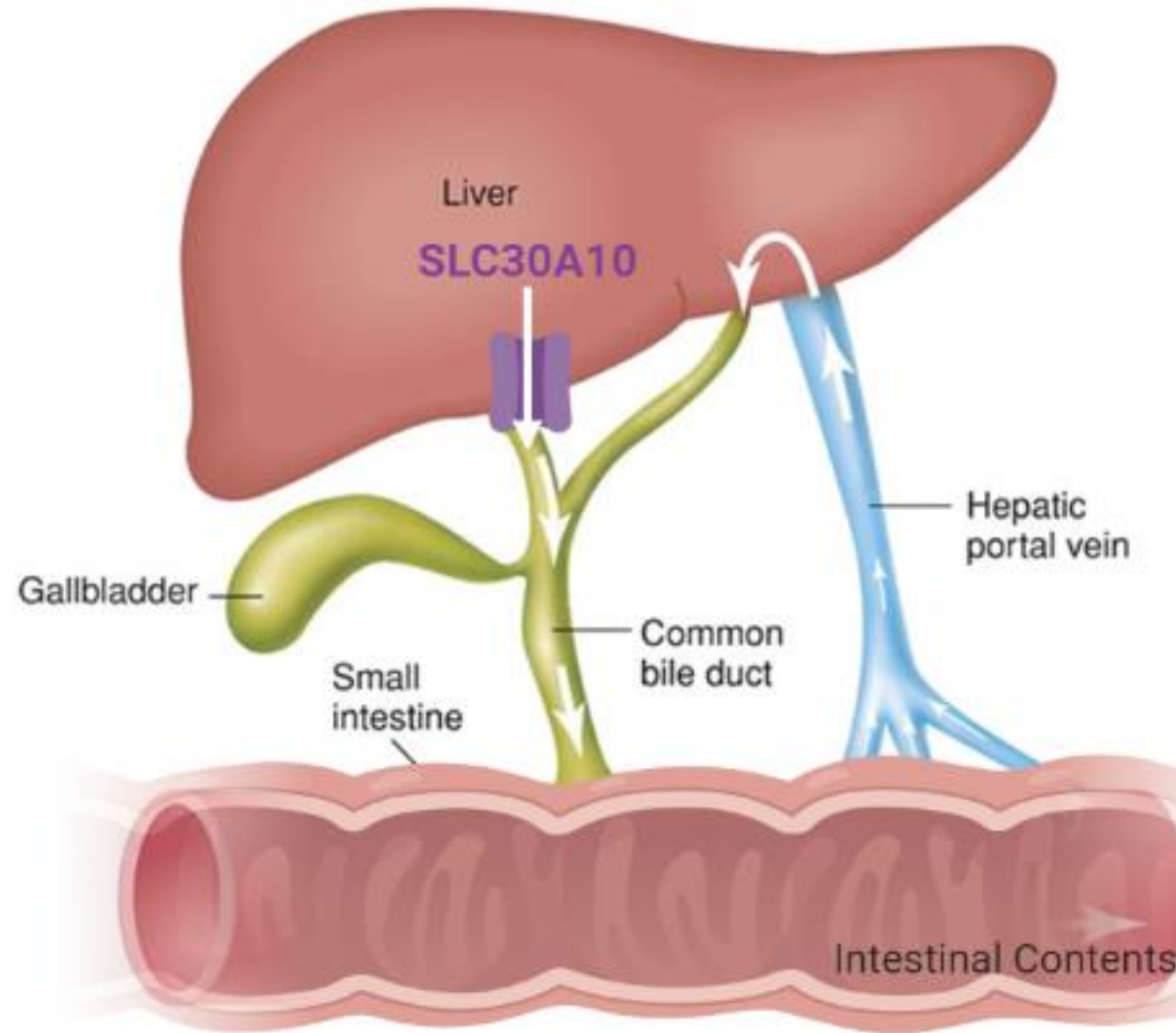
# Hepatic Slc30a10-deficient mice have impaired biliary Mn excretion, but minimal Mn excess



# Why do hepatic Slc30a10-deficient mice have minimal Mn excess?

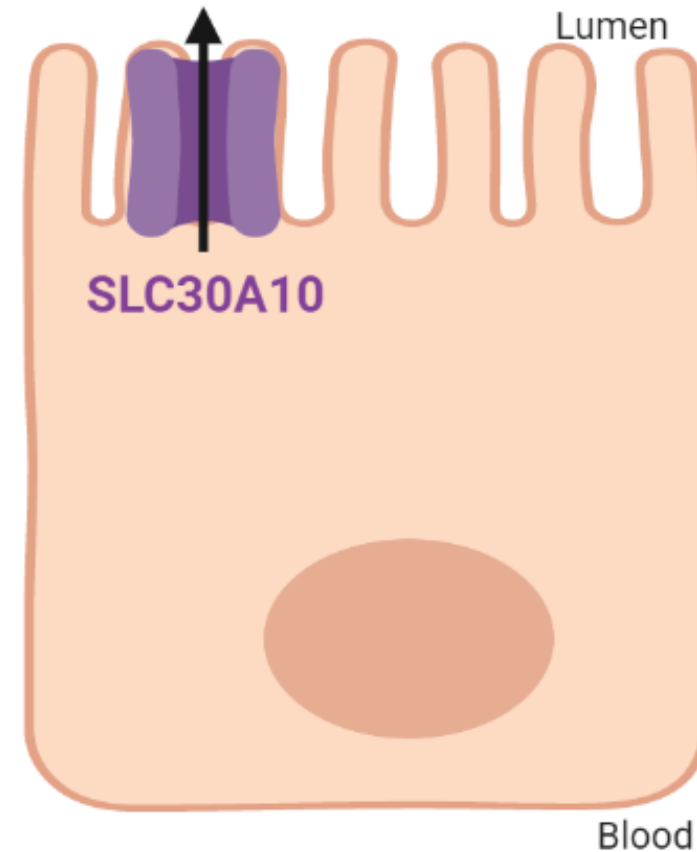
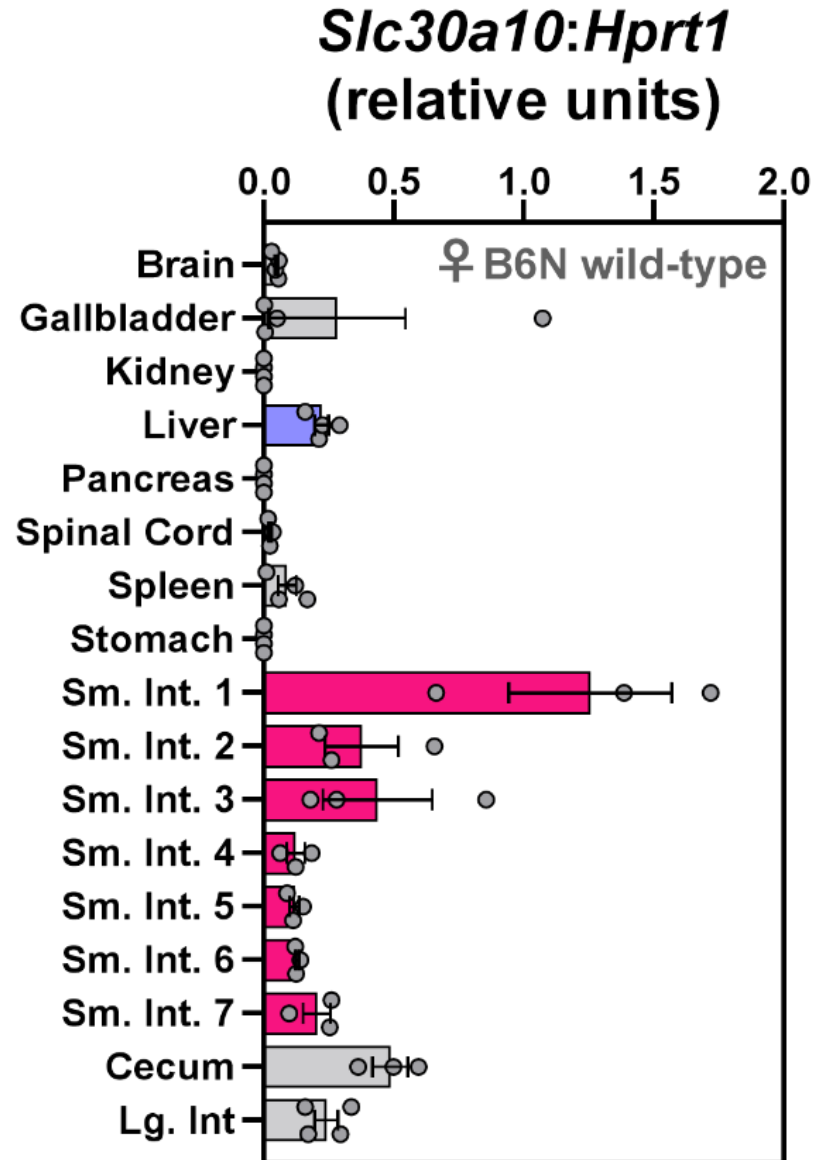
Hepatic Slc30a10 is essential for hepatobiliary Mn excretion, but hepatic Slc30a10 deficiency only results in a minimal Mn excess?

- Late onset of Albumin-Cre transgene (P21)
- Another tissue contributes to Mn excretion





# Slc30a10 is highly expressed in intestines



# Generating Slc30a10-deficient mice

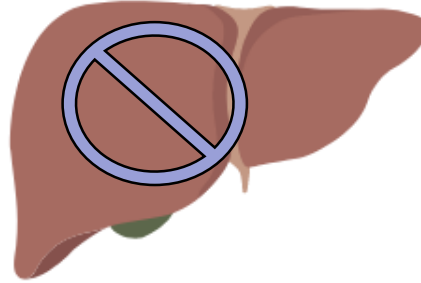
Whole body Slc30a10 deficiency

***Slc30a10<sup>KO/KO</sup>***



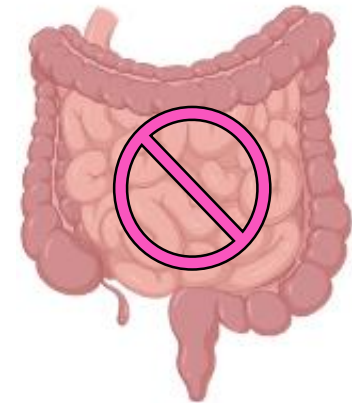
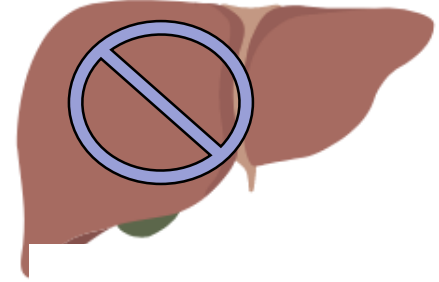
Hepatic  
Slc30a10 deficiency

***Slc30a10<sup>lox/lox</sup> Alb***



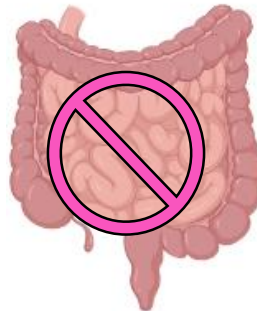
Hepatic and Intestinal  
Slc30a10 deficiency

***Slc30a10<sup>lox/lox</sup> Alb Vil***

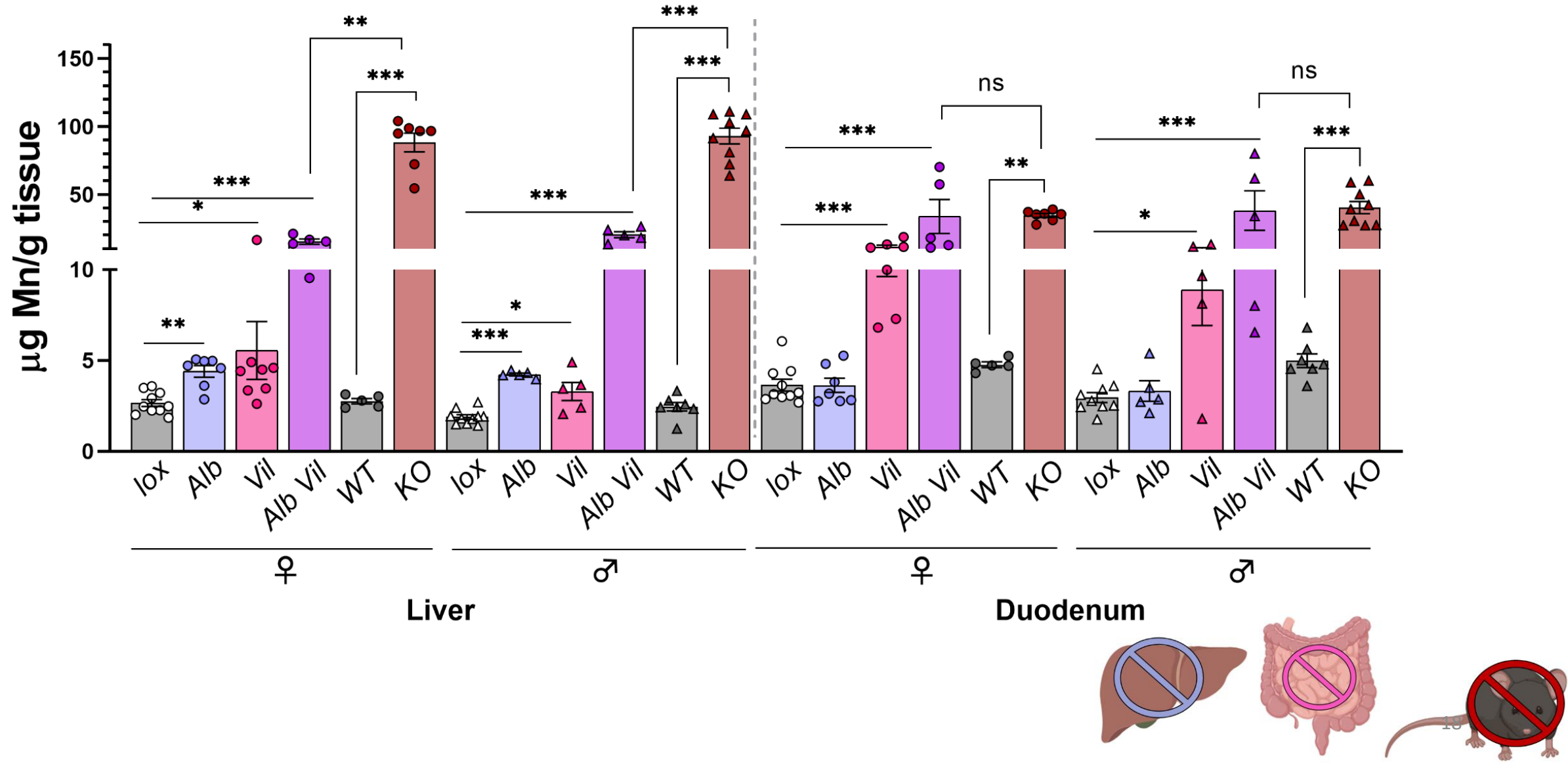


Intestinal  
Slc30a10 deficiency

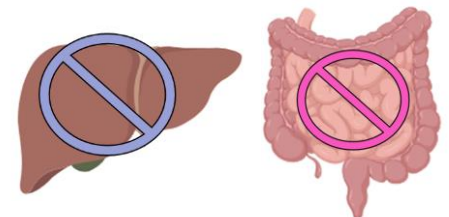
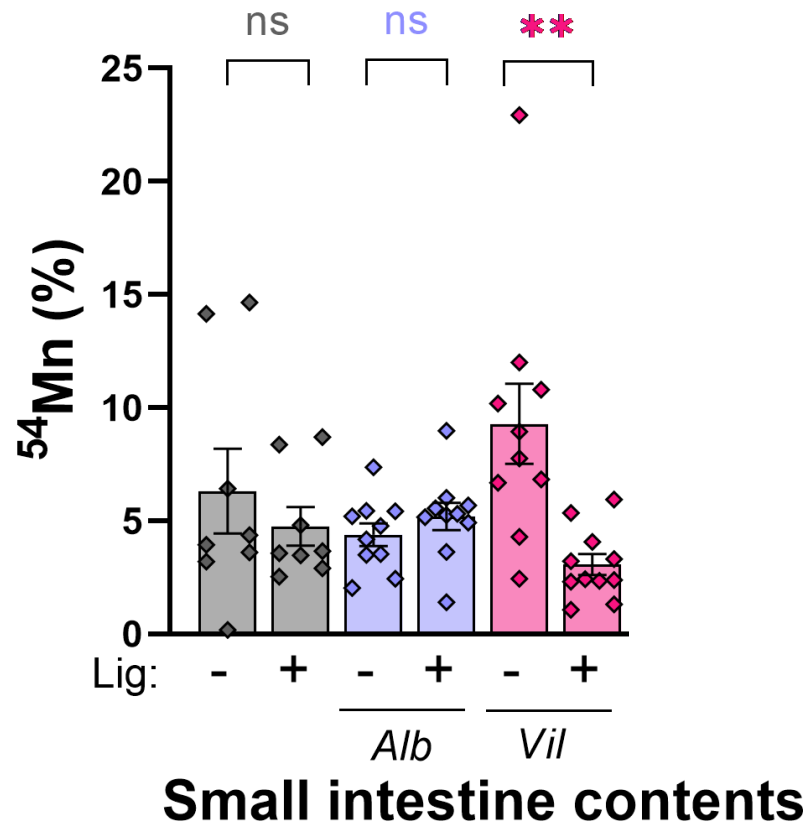
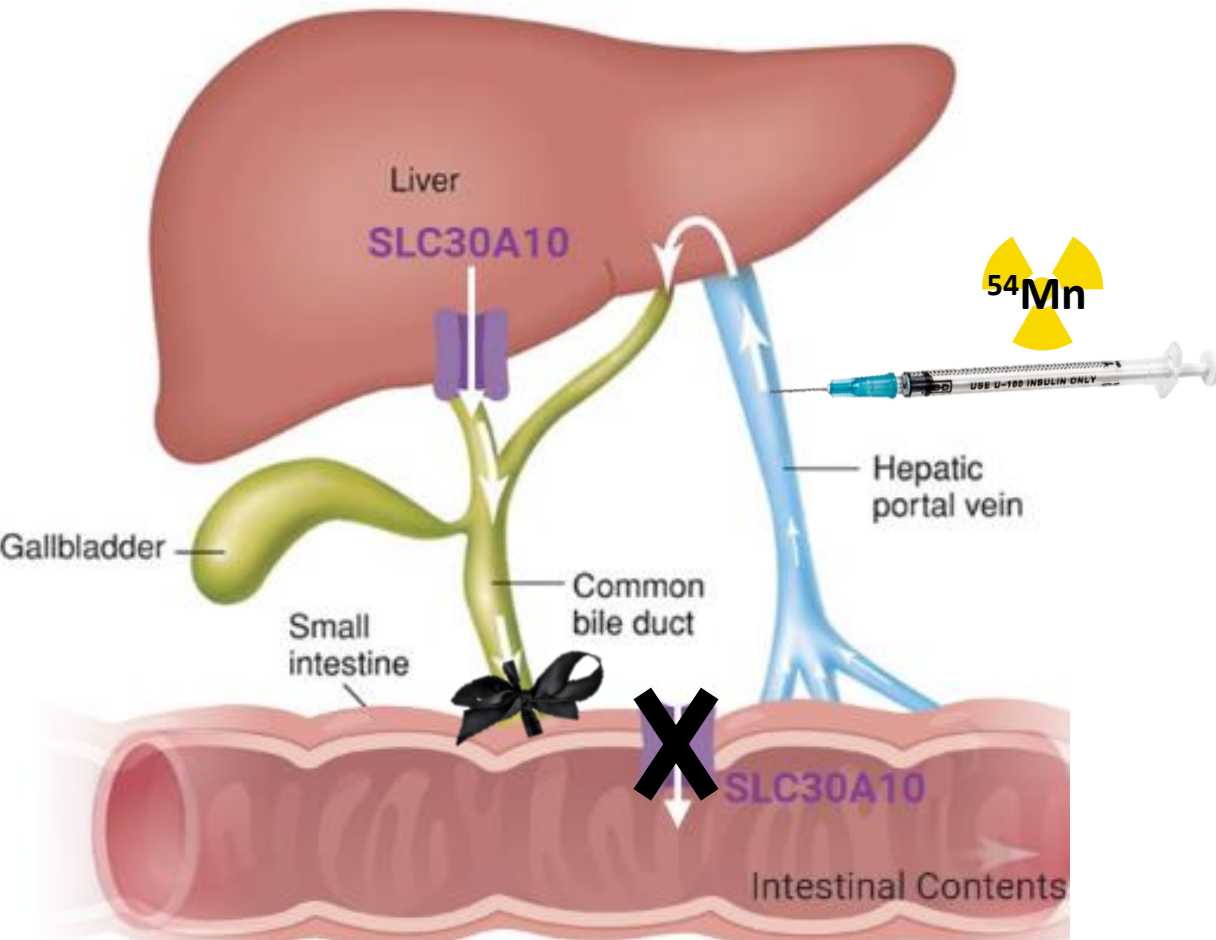
***Slc30a10<sup>lox/lox</sup> Vil***



# Intestinal and hepatic Slc30a10 contribute to Mn homeostasis

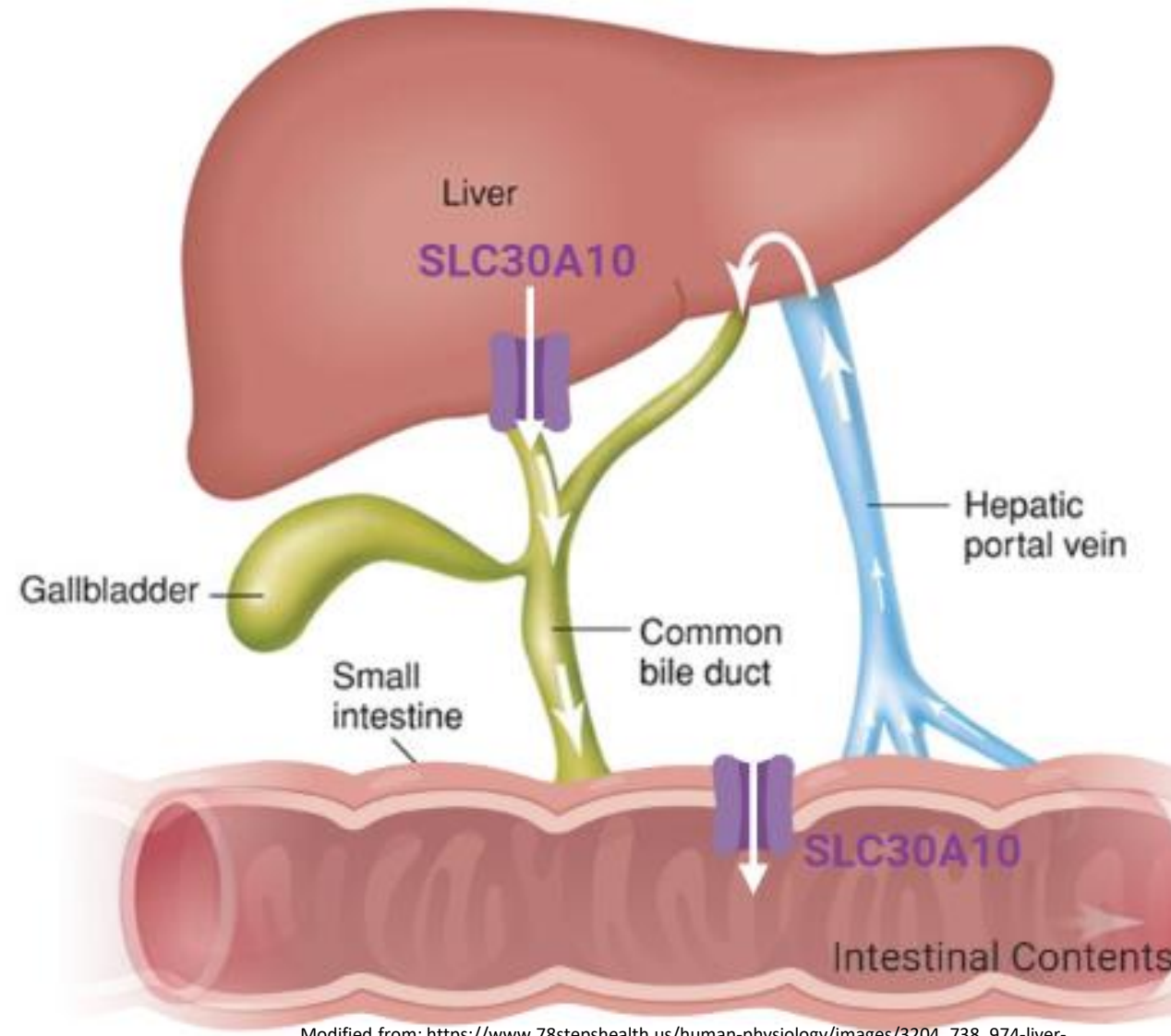


# Intestinal Slc30a10 contributes to intestinal Mn excretion



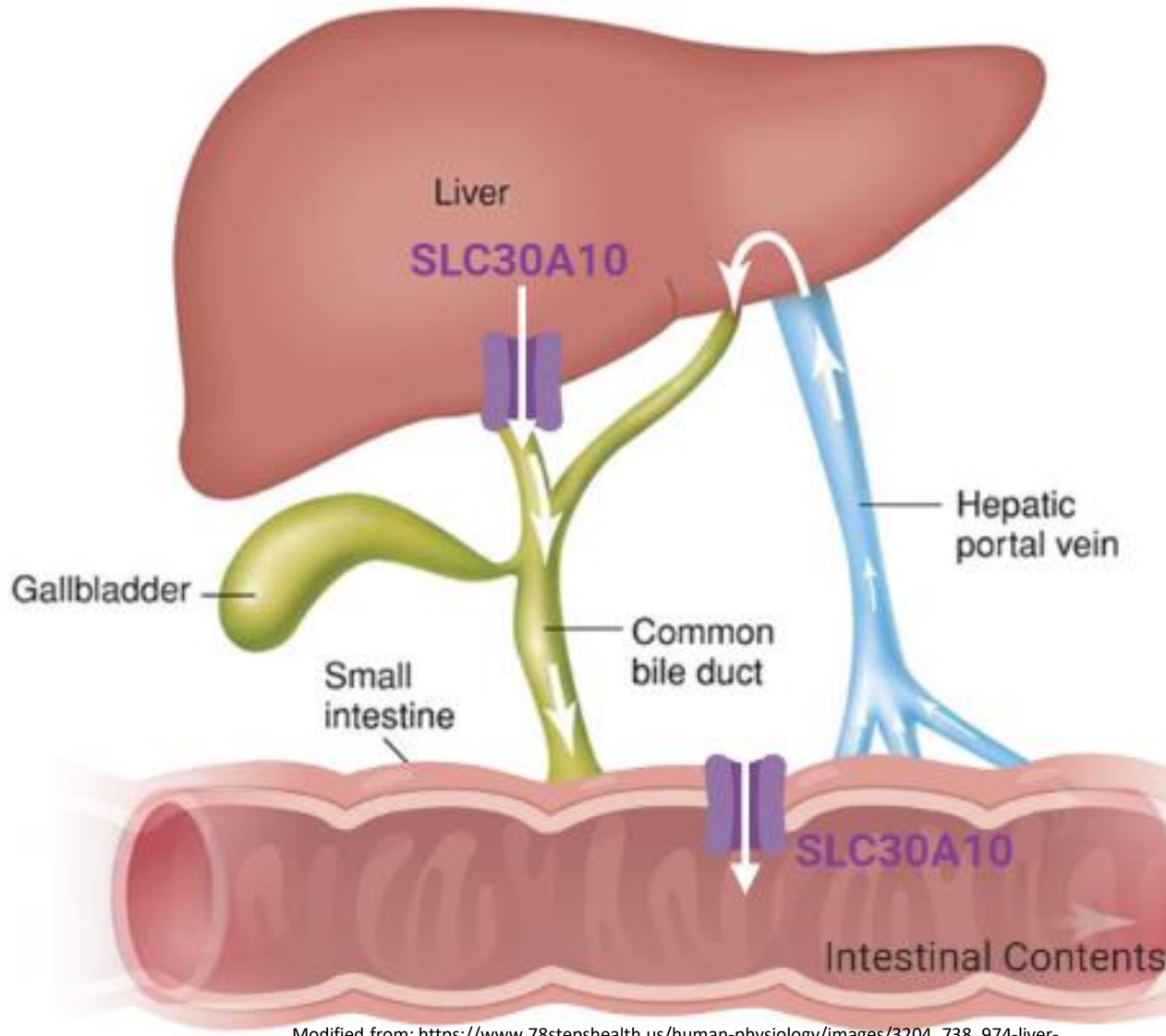
# Conclusions

- Slc30a10-deficient mice have impaired Mn excretion
- Hepatic Slc30a10 is essential for hepatobiliary Mn excretion
- Hepatic and intestinal Slc30a10 contribute to manganese homeostasis
- Intestinal Slc30a10 contributes to excretion of Mn into intestinal lumen



# Future Directions

- Determine the contribution of Slc30a10 to Mn homeostasis in developing mice
- Determine the role of Slc30a10 in cecum and large intestine to Mn homeostasis
- What about Mn absorption?



Modified from: [https://www.78stepshealth.us/human-physiology/images/3204\\_738\\_974-liver-gallbladder-portal-vein-intestines.jpg](https://www.78stepshealth.us/human-physiology/images/3204_738_974-liver-gallbladder-portal-vein-intestines.jpg), Created using Biorender.com



# Acknowledgments



## Bartnikas laboratory

Dr. Tom Bartnikas

Dr. Milan Prajapati

Heather Conboy

Miriam Dash

Carolina Herrera

Mike Pettiglio

## Genomics Core

Dr. Christoph Schorl

## Metal analysis

Dr. Joseph Orchard

## Funding

NIDDK- R01 DK110049, F31 DK117524

NIEHS- T32 ES007272

