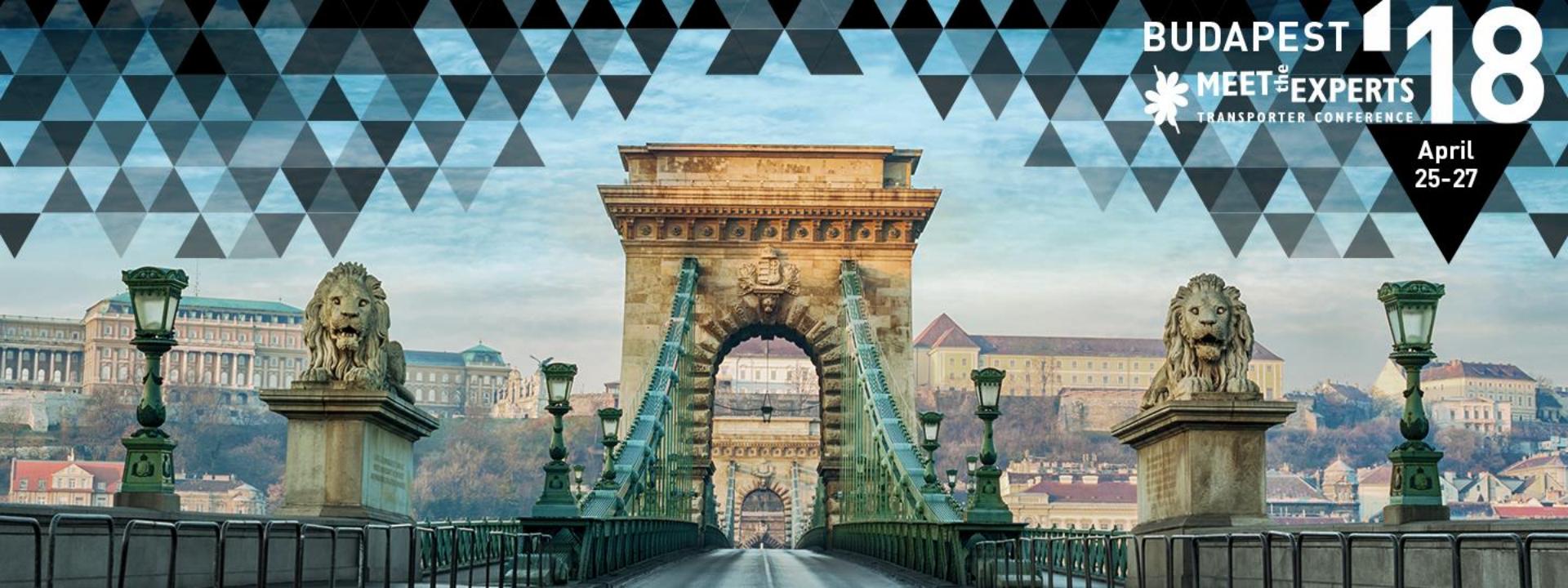


April  
25-27



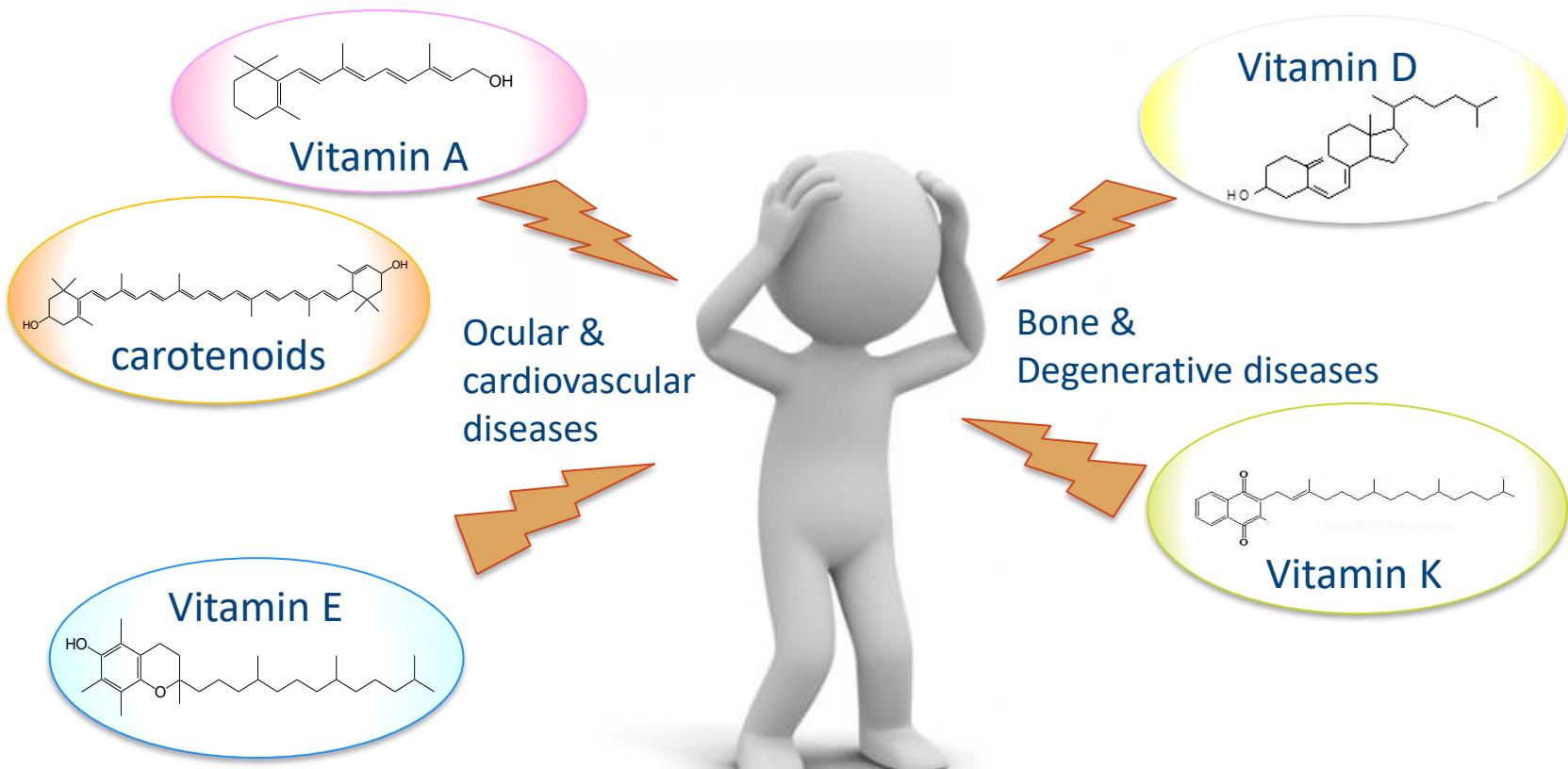
# Intestinal absorption of fat-soluble vitamins

Emmanuelle Reboul

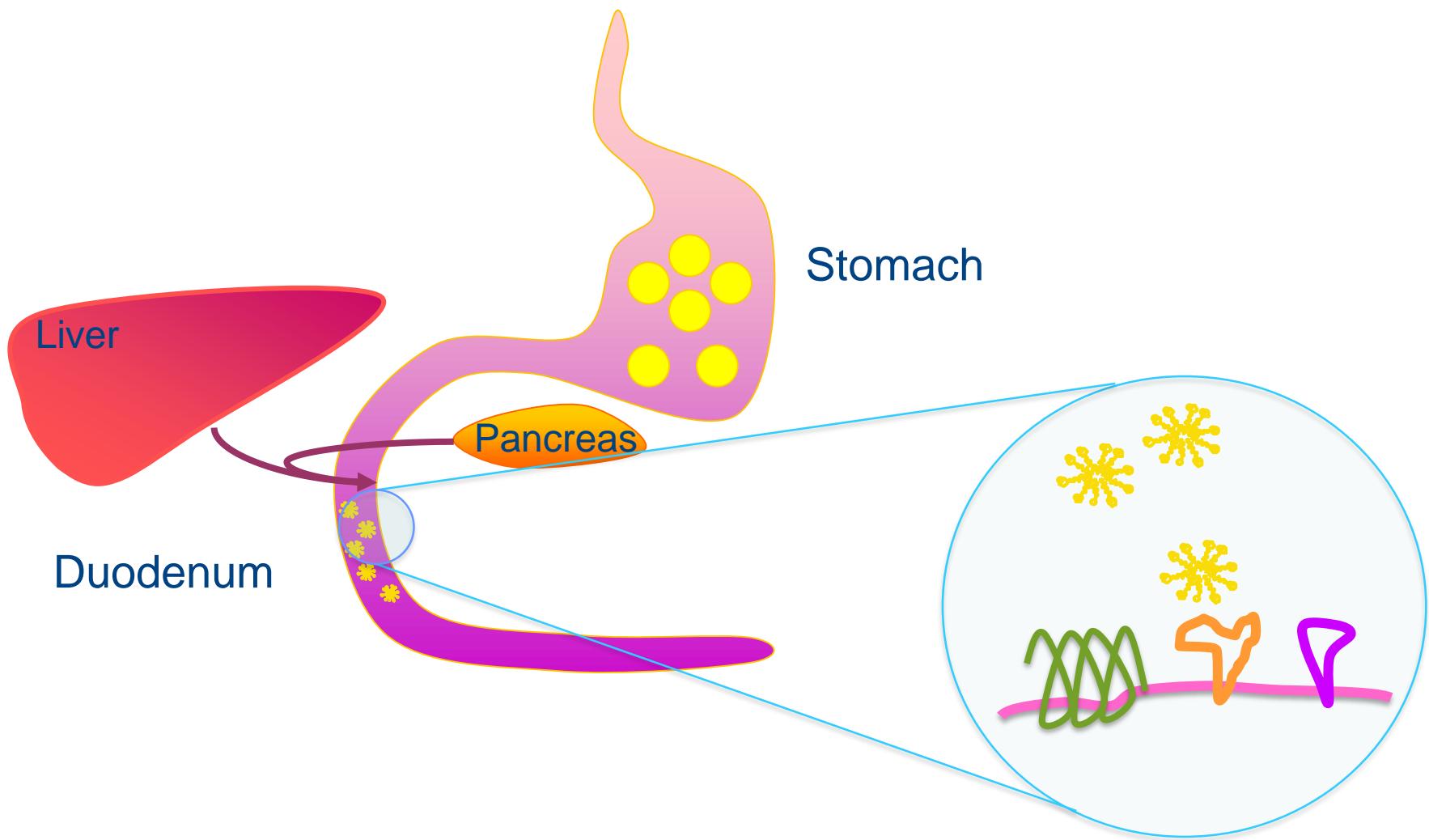
Human Micronutrition team  
Cardiovascular and Nutrition Center (C2VN)  
Marseille - FRANCE

# The lipid micronutrients

- Vitamins are non-energetic molecules present in small amounts in our diet (< 1g/d), necessary for our development and health. 4 fat-soluble vitamins: A, D, E, K.
- Carotenoids: not essential yet.
- ◆ Sub-deficiencies are very common

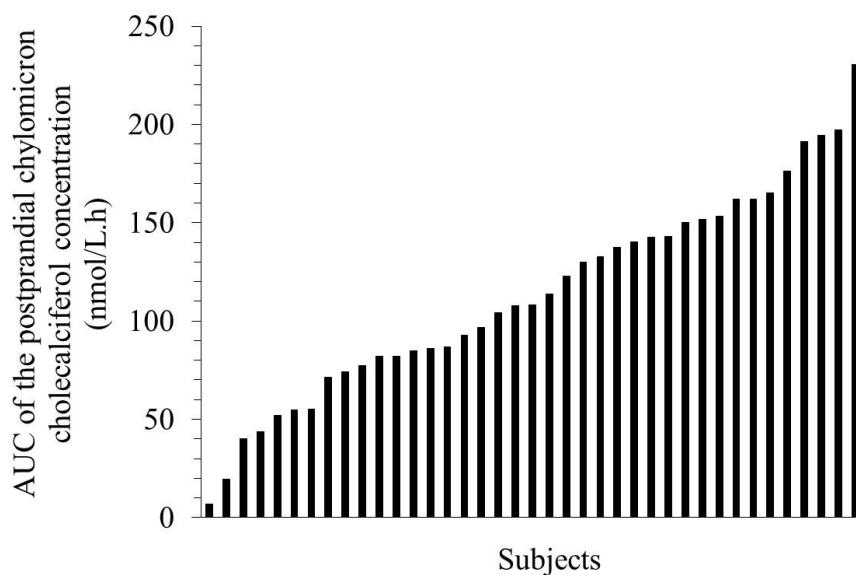


# Digestion of lipid micronutrients



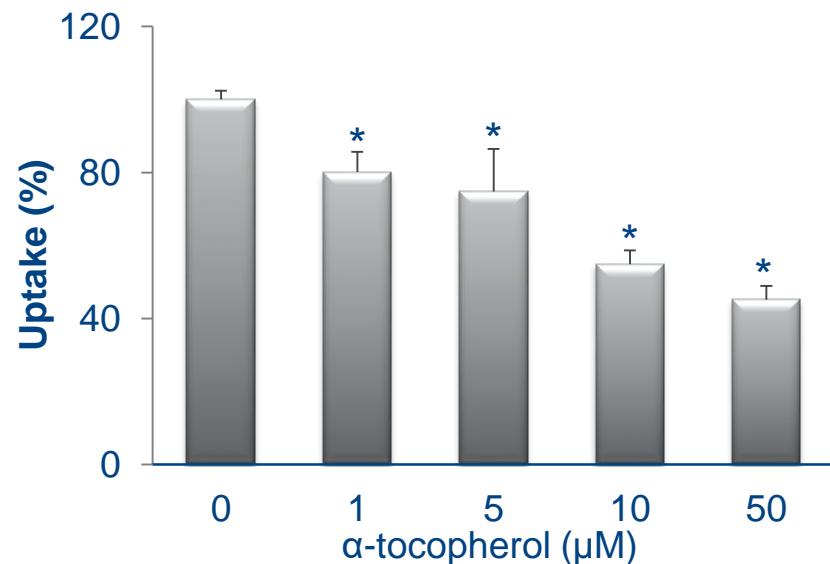
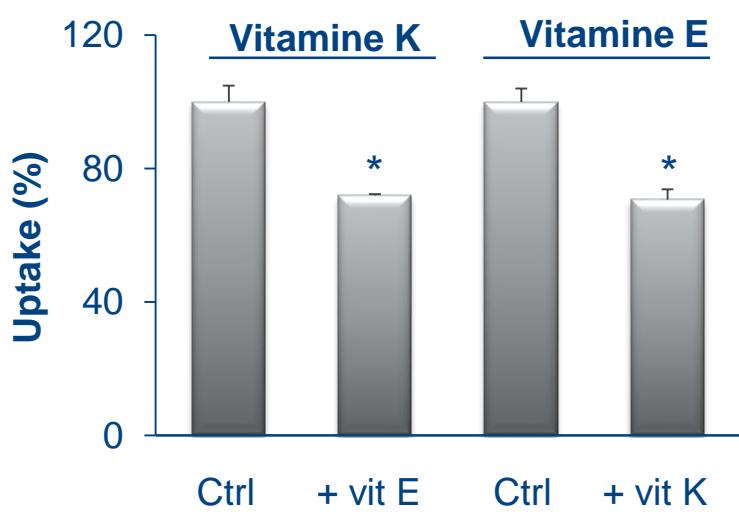
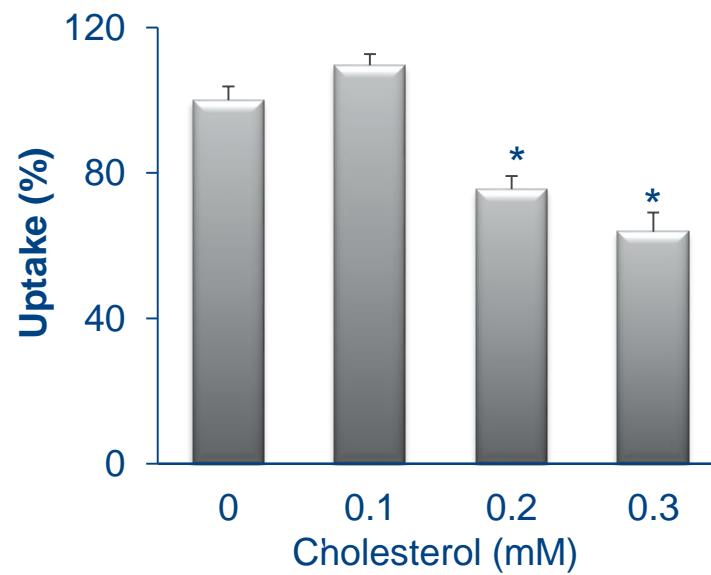
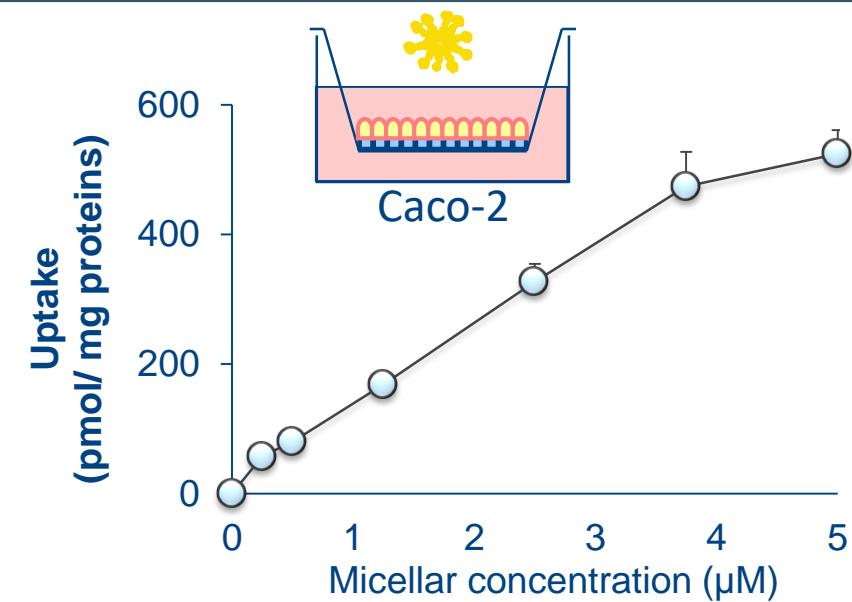
# Absorption by the enterocyte

- ◆ Absorption mechanisms initially defined by Hollander et al. in the 70s
  - Passive diffusion for vitamin D, E and carotenoids
  - Active process for vitamin A and K
- ◆ Inconsistent data
  - Competition for absorption
  - **Very high** interindividual variability

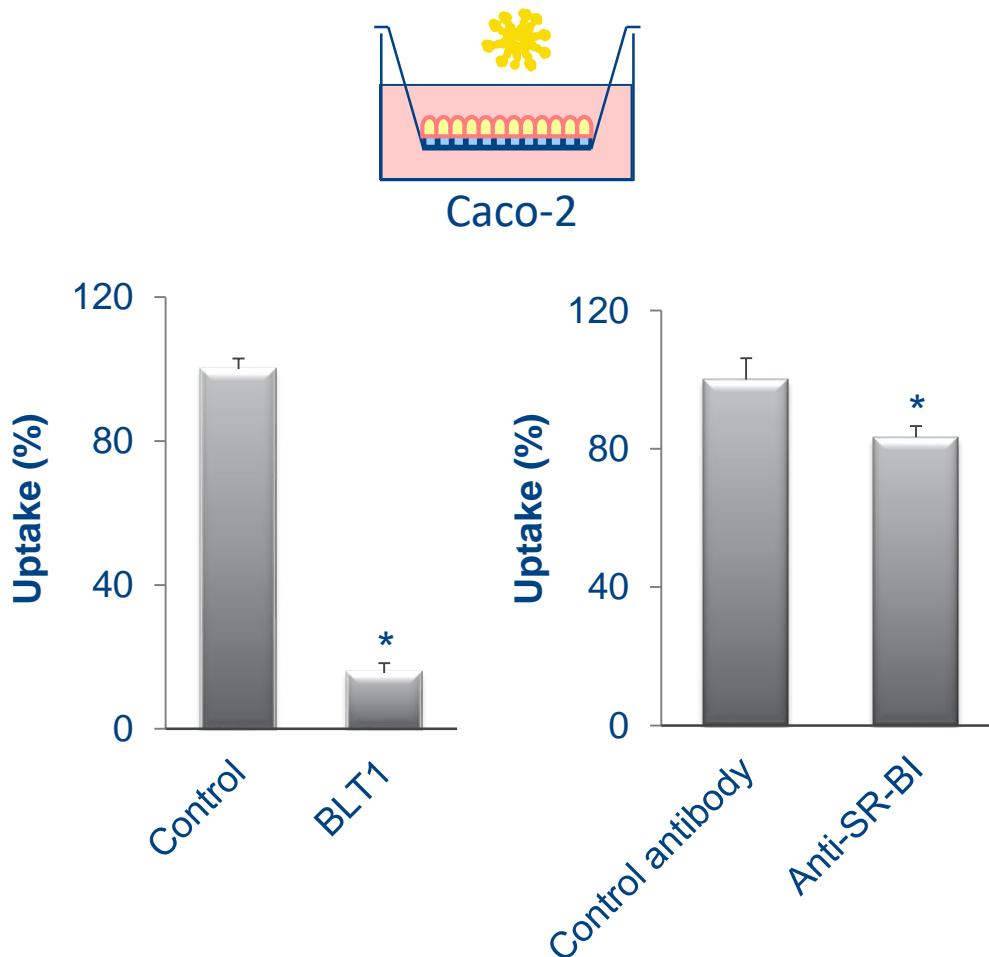


→Reinvestigation of lipid micronutrient absorption since 2002

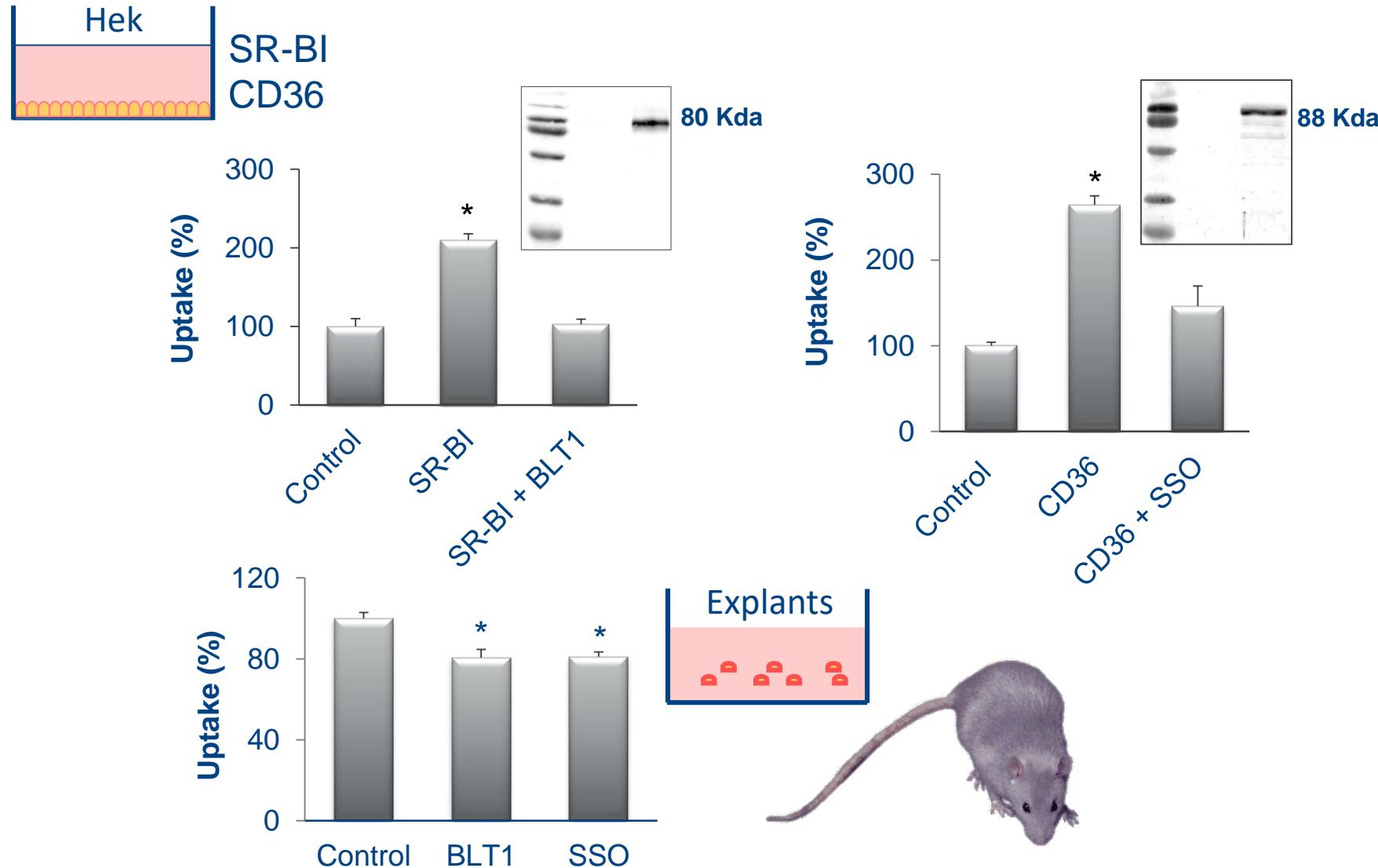
# Vitamin K absorption



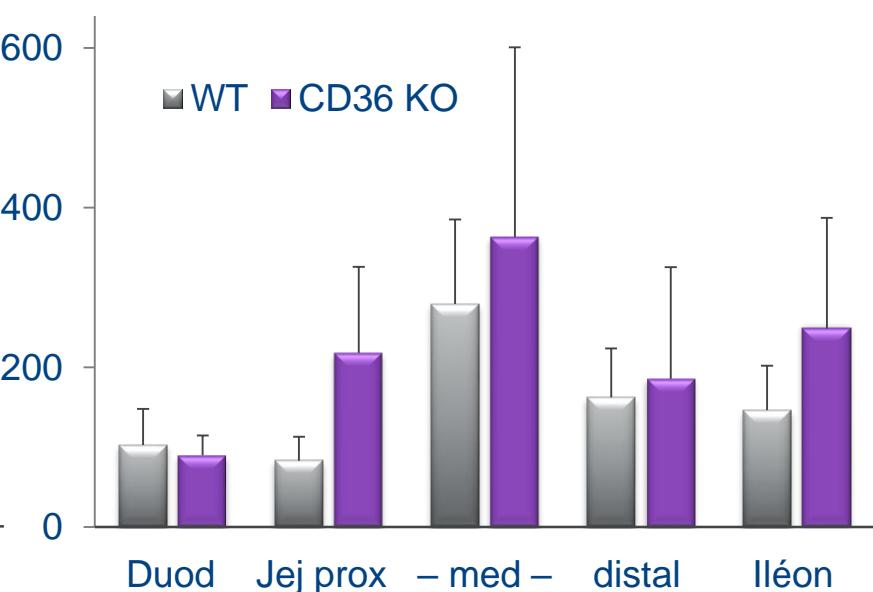
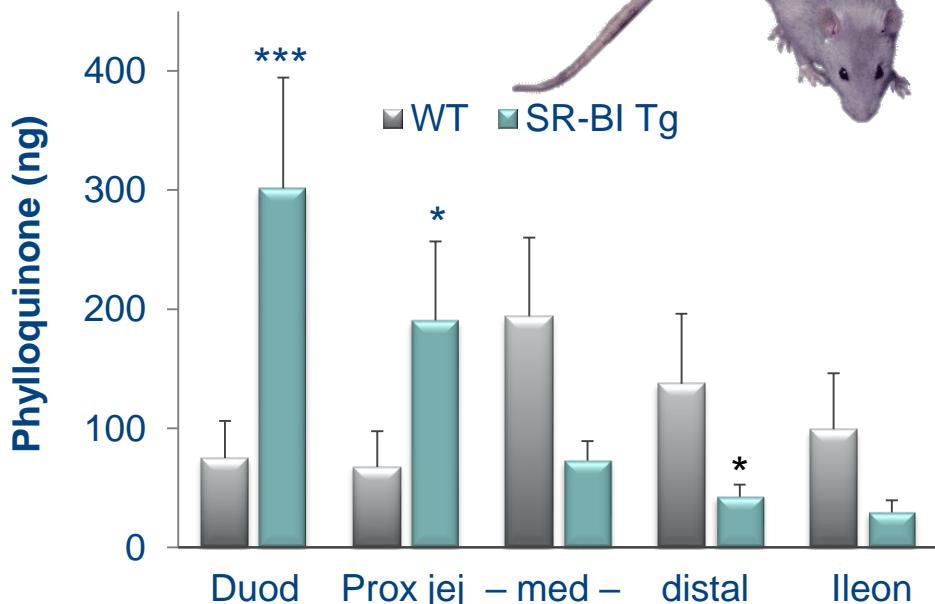
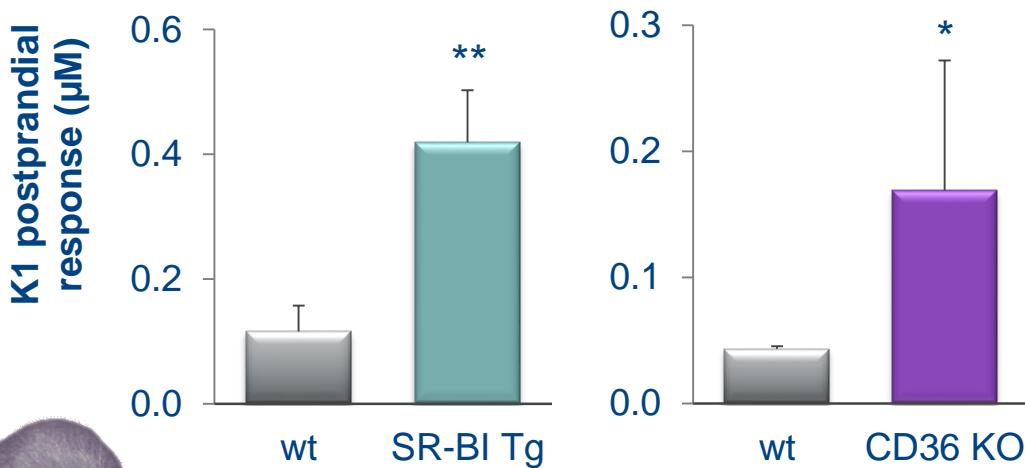
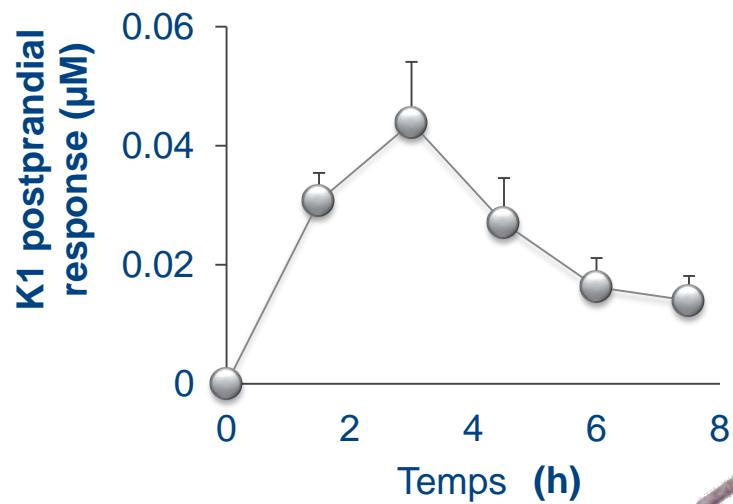
# Vit K and SR-BI in Caco-2 cells



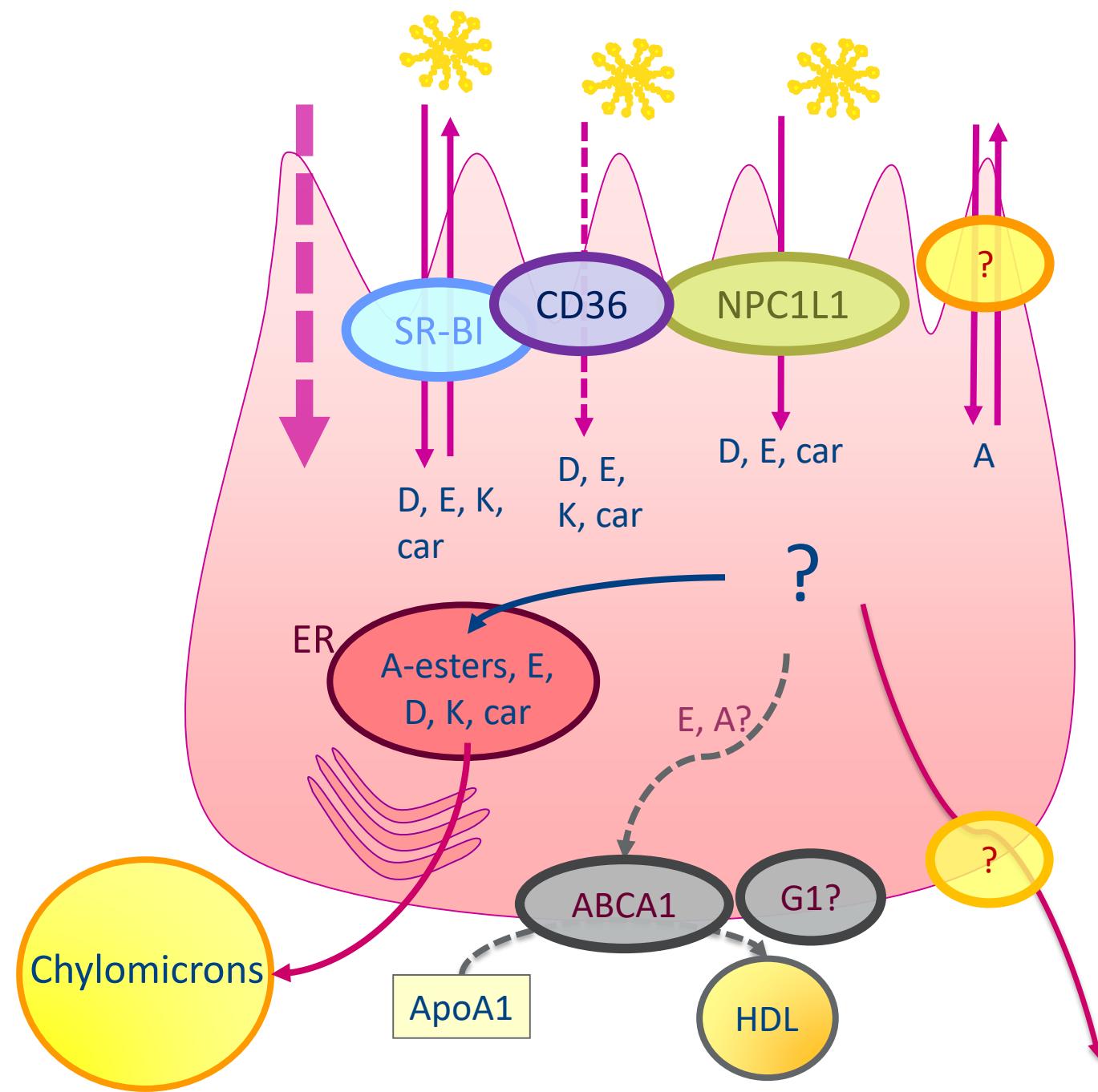
# Vit K & scavengers in transfected HEK cells and mouse intestinal explants



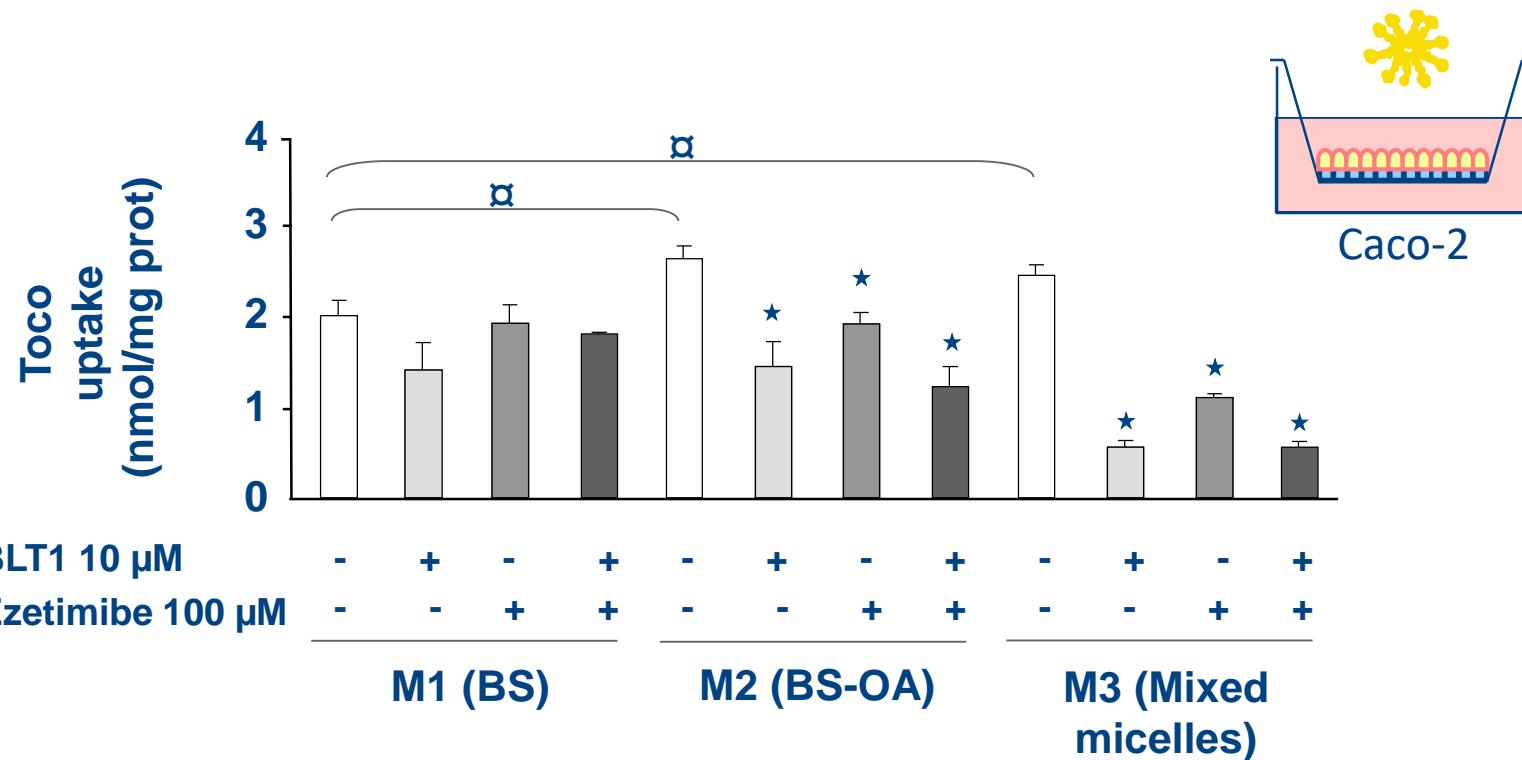
# Vit K postprandial response in mice



Reboul et al. 2005, 2006,  
2009, 2011; Moussa et al.  
2008; Borel et al. 2013;  
Goncalves et al. 2014 a, b.  
Van Benneckum et al. 2005;  
During et al. 2005, 2007;  
Anwar et al. 2006, 2007;  
Narushima et al. 2008; Nicod  
et al. 2013.

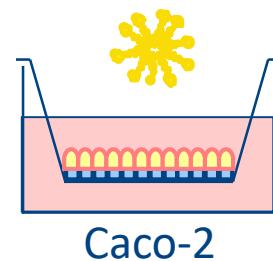
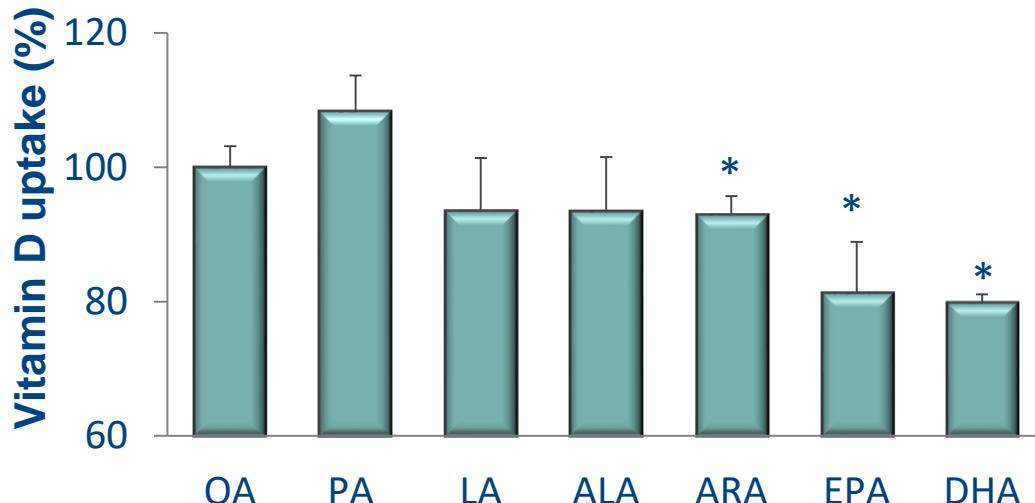


# Importance of the vehicle



- ◆ Presence of bile salts + fatty acids required for a transporter-dependent uptake

# Importance of the vehicle

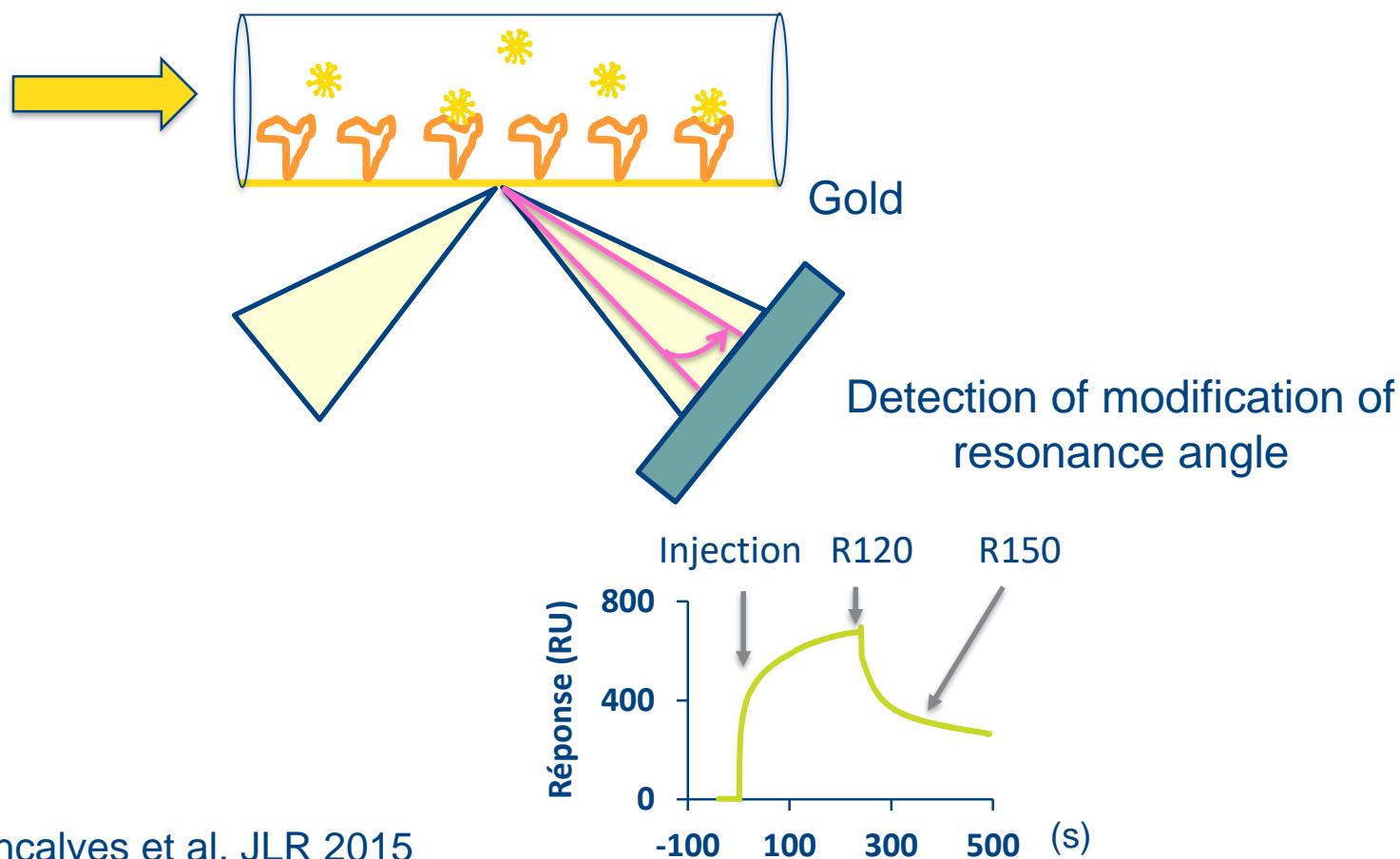


	OA	PA	LA	ALA	ARA	EPA	DHA
Zeta (mV)	-30,3 ± 2,4	-22,7* ±3,9	-34,5* ±2,7	-32,3* ±3,9	-31,0* ±1,9	-36,9* ±3,4	-38,8* ±3,9
Taille (r. nm)	5,4 ±0,2	5,1* ±0,4	3,6* ±0,6	3,5* ±0,5	3,4* ±0,1	3,5* ±0,2	3,6* ±0,1

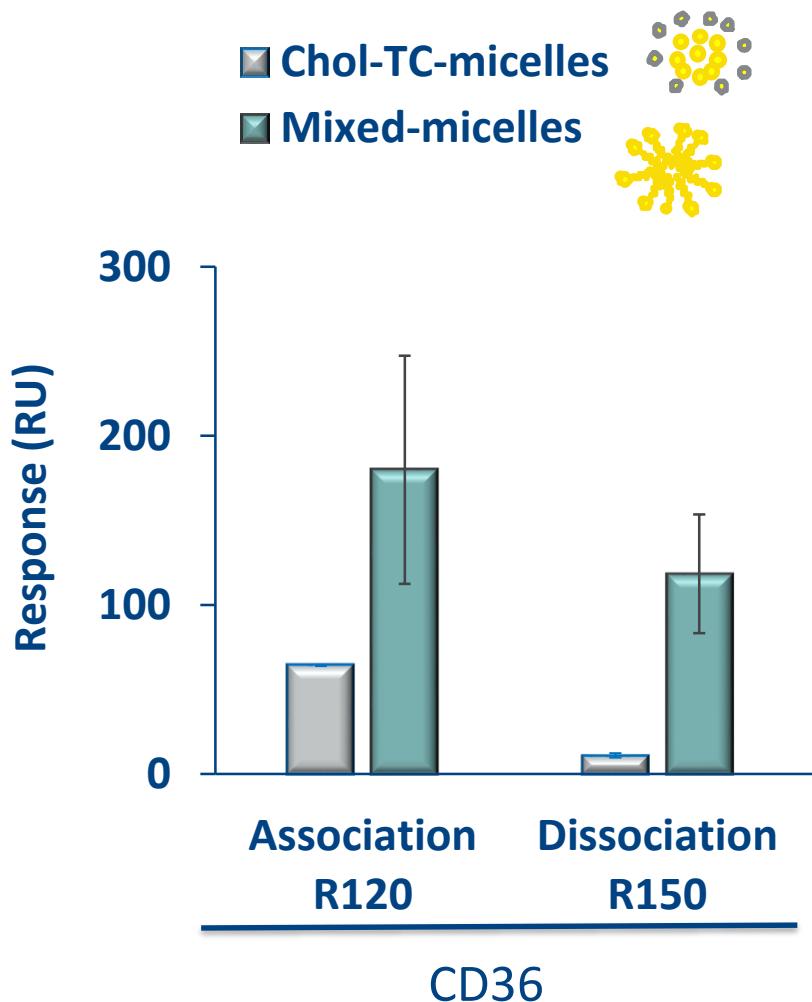
◆ Modification of micellar affinity with transporters?

# Intestinal scavenger receptor functioning

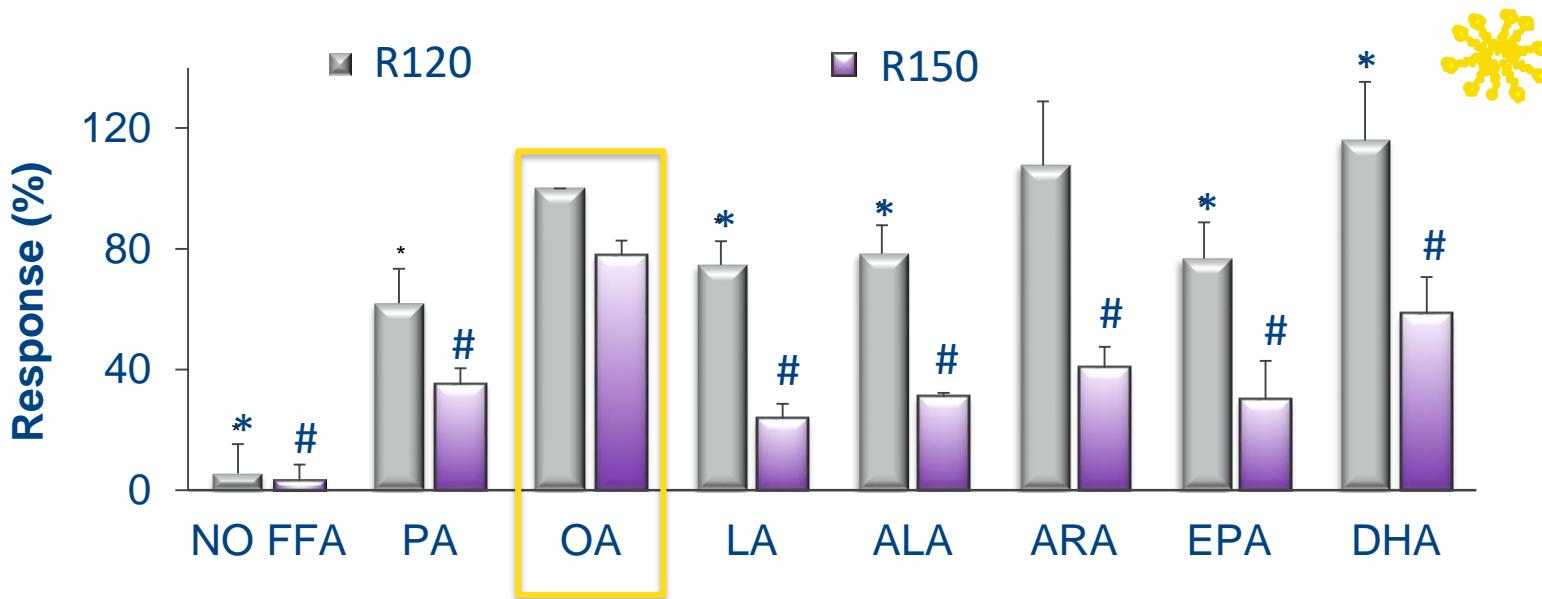
- ◆ Real-time analysis of protein-micelles interactions with Surface Plasmon Resonance
- ◆ Extracellular loop of SR-BI and CD36
- ◆ Synthetic mixed micelles



# Importance of micelle type

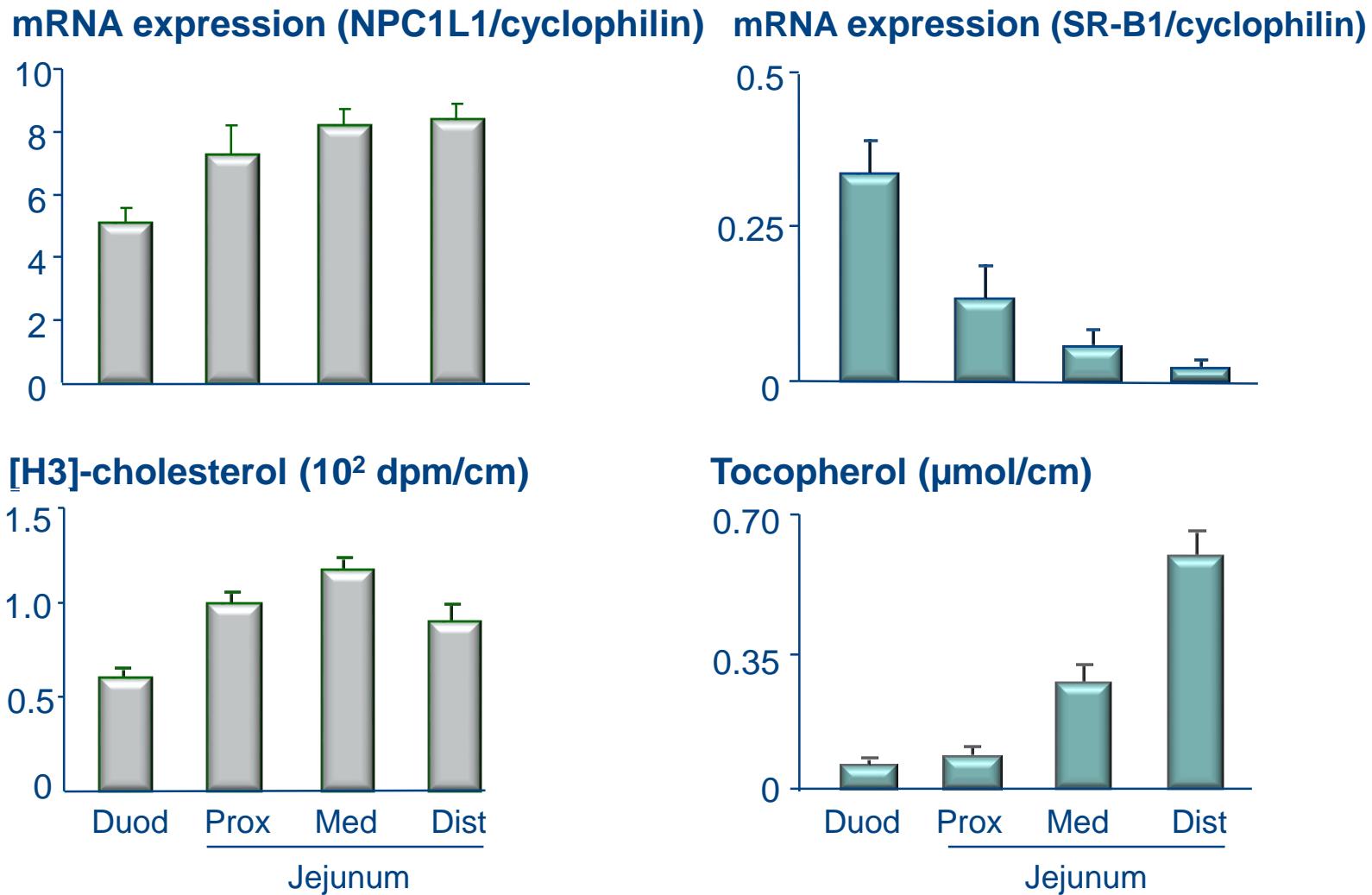


# Effect of micellar fatty acids



- ◆ Among the different micellar components (chol, PC), fatty acids are the most important regarding micelle-scavenger interactions

# Absorption sites of lipid micronutrients



# Take away messages

- ◆ Vitamin D, E, K and carotenoids share similar absorption pathways but are not absorbed at the same level of the intestine.
- ◆ Lipid micronutrient uptake through transporters require an adequate vehicle.
- ◆ The specificity of the interaction lipid micronutrient - transporter is not known yet.



Marseille  
Center for CardioVascular  
and Nutrition research



## ✓ Laboratory

Patrick Borel (PI)

Aurélie Goncalves (PhD)

Marielle Margier (PhD)

Marion Nowicki (RT)

Charlotte Halimi (RT)

Gabriel Masset (Postdoc)

MS-Nutrition (Start-up)



## ✓ Collaborations

Brigitte Gontero (BIP, CNRS, Marseille)

Catherine Caris (SQPOV, Avignon)

Isabelle Niot (Inserm 866, Dijon)

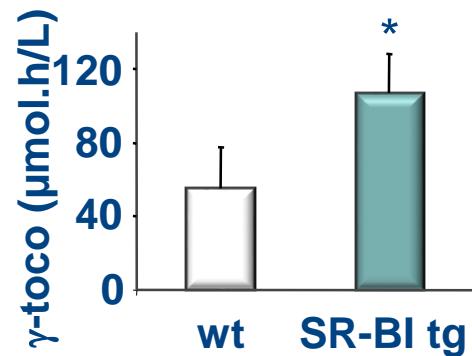
Xavier Collet (UMR 1048, Toulouse)

Wilfried le Goff (UPMC, Paris)

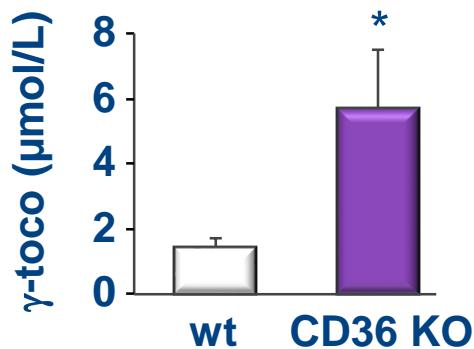
Anne Lespine (INRA, Toulouse)



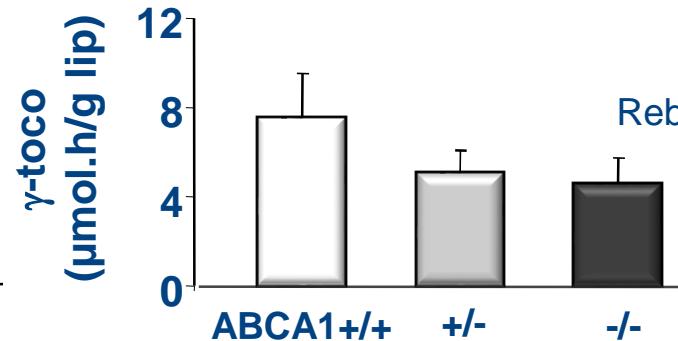
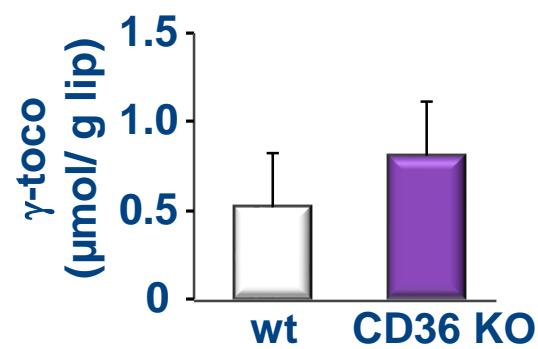
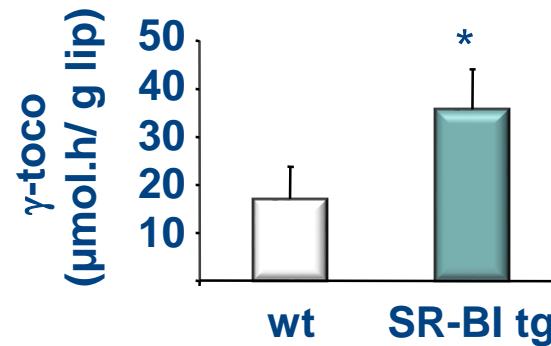
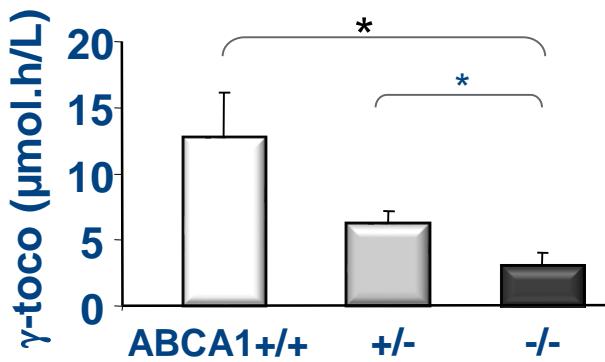
# Direct or indirect transport?



Reboul et al. JBC 2006



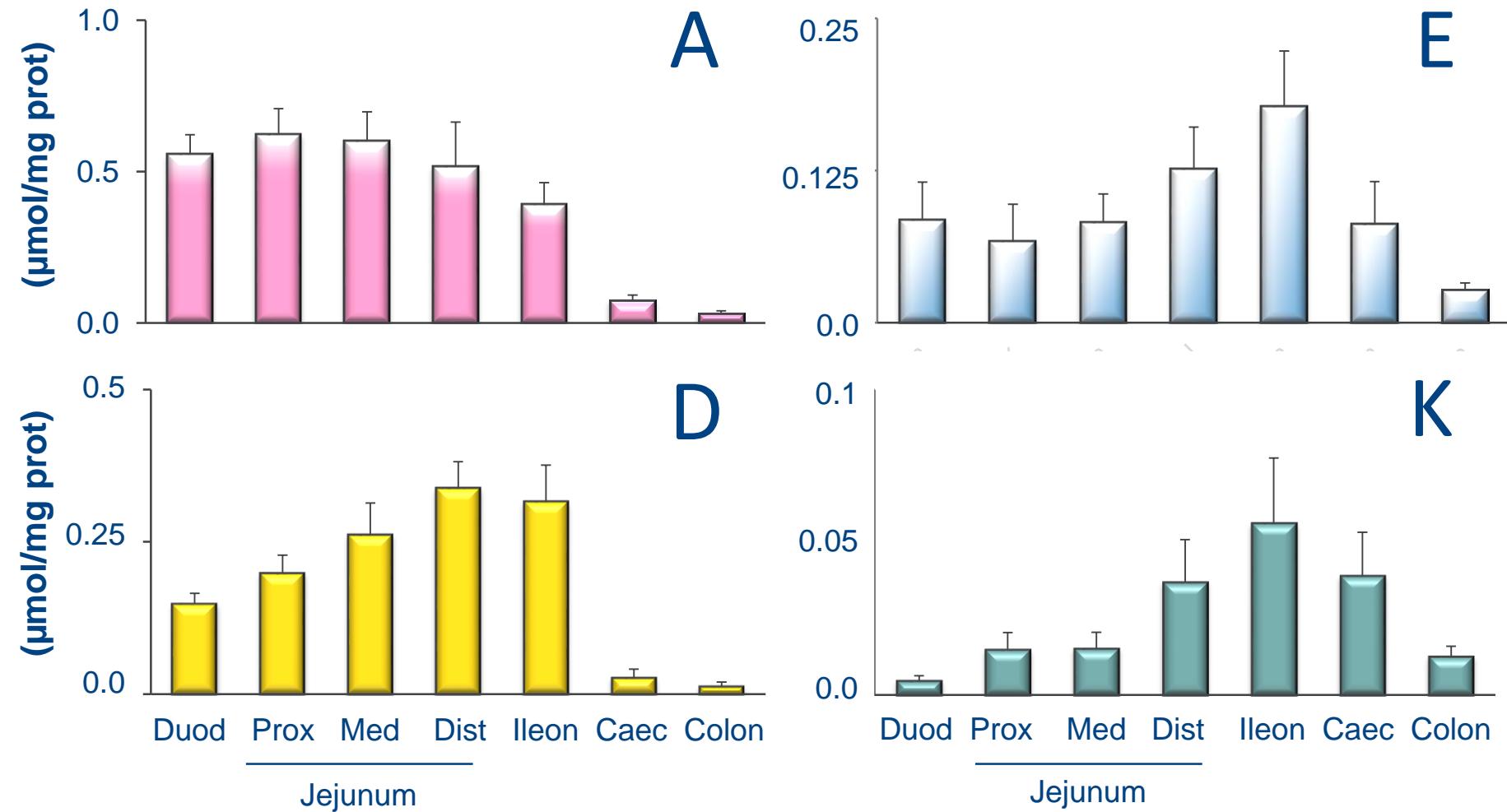
Goncalves et al. MNFR 2014



Reboul et al. Am J Clin Nutr 2009

- ◆ Direct: SR-BI?
- ◆ Indirect: CD36, ABCA1?

# Absorption sites of lipid micronutrients



◆ Different absorption sites → different absorption mechanisms?