

Phospholipid Substrates for Lysosomal Phospholipase A2 (LPLA2)

Fok Vun Chan[§], James A. Shayman[‡], Paul O. Neilsen[§], L. Jeff Johnson[§]

[§]Echelon Biosciences, Inc. Salt Lake City, Utah, United States

[‡]Internal Medicine – Nephrology, University of Michigan, Ann Arbor, MI, United States



1. OVERVIEW

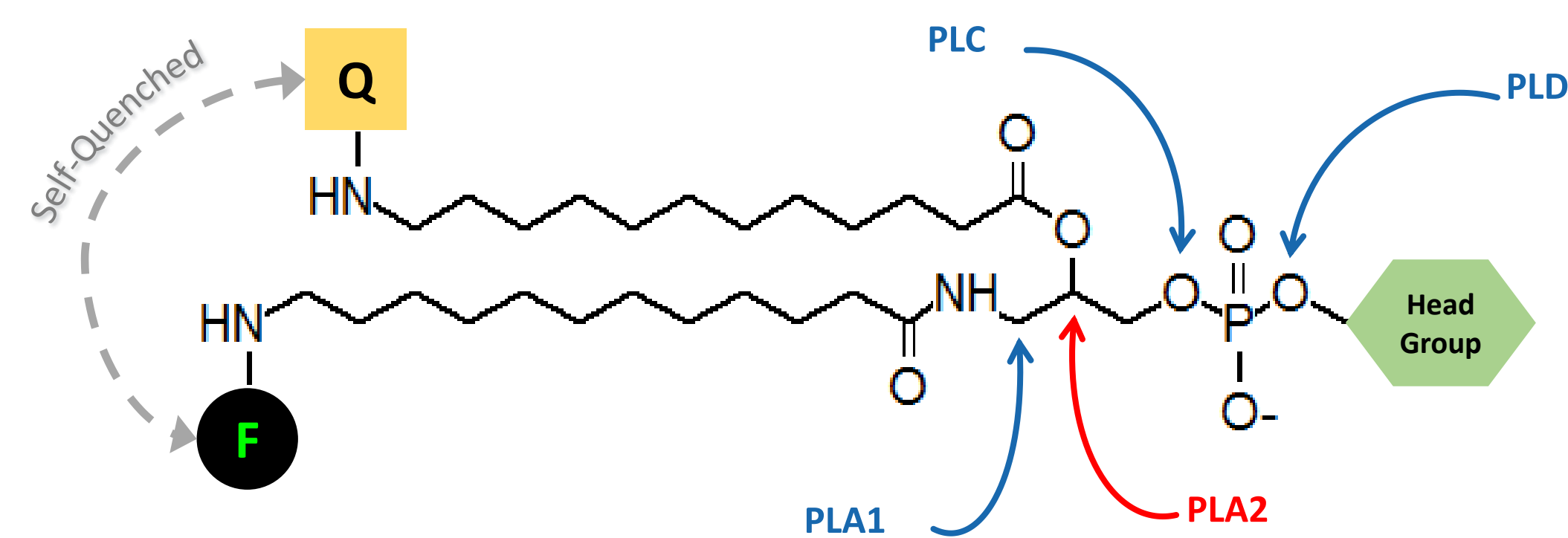
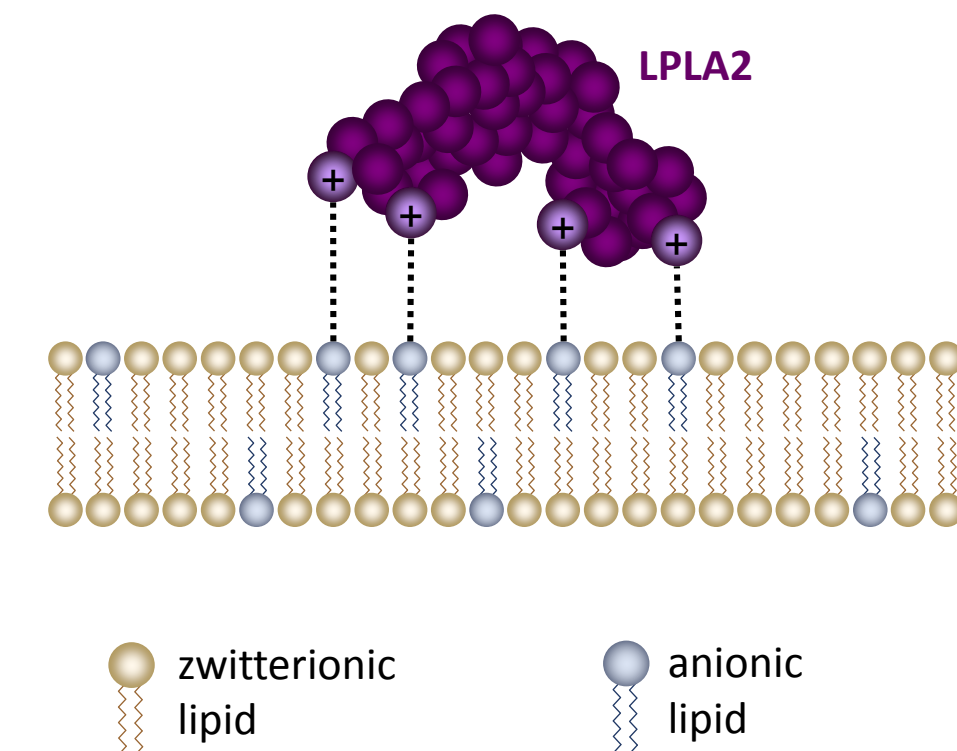
➤ Lysosomal phospholipase A2 (LPLA2) is involved in both drug-induced phospholipidosis (DIPL) & drug-induced lupus (DIL)

➤ Evaluated phospholipid charge & structure in relation to LPLA2 activity

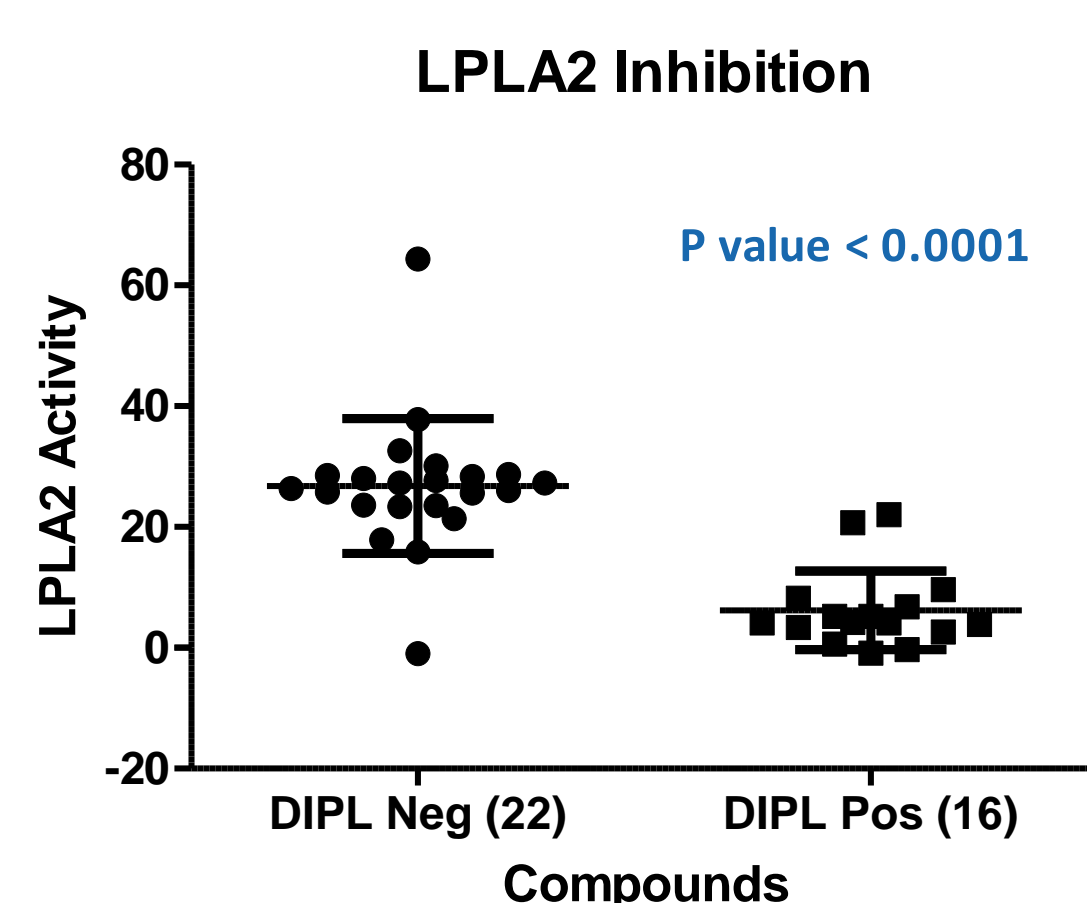
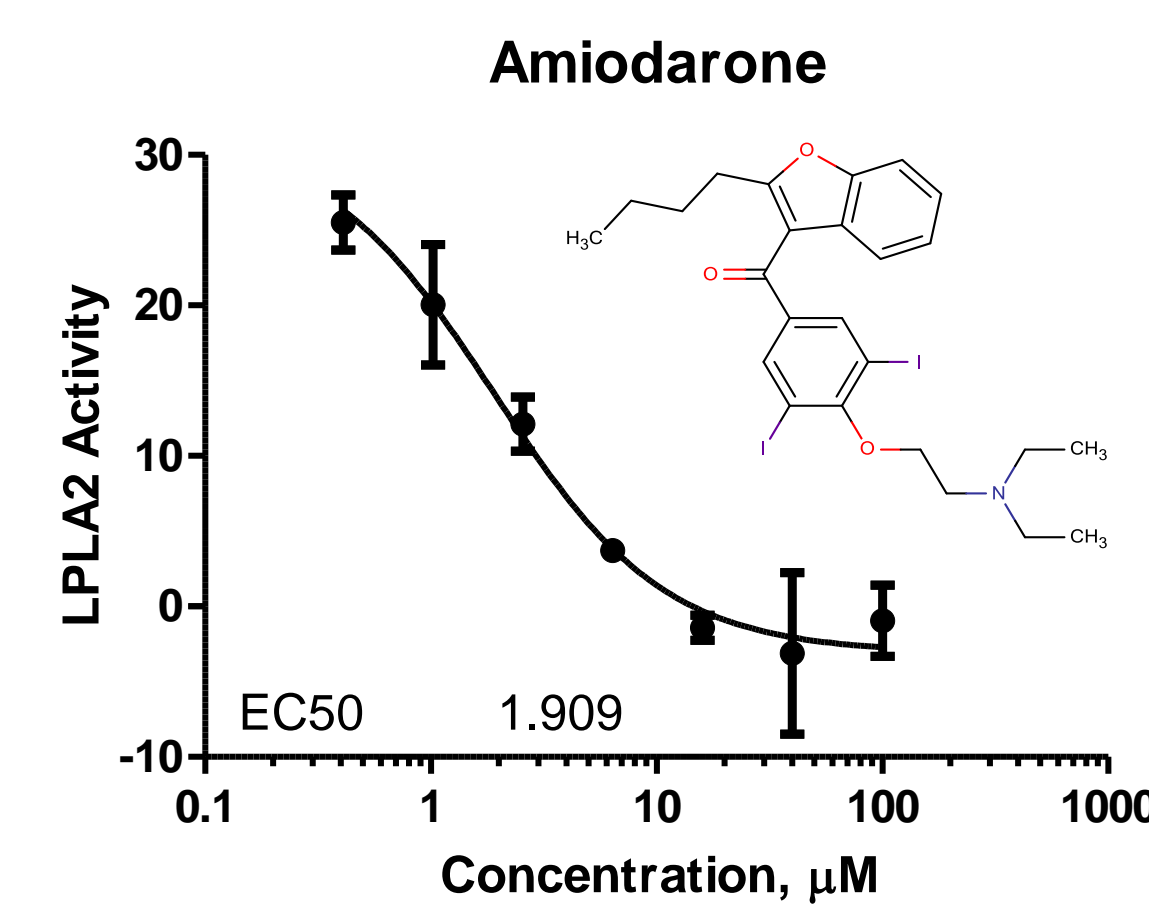
➤ Studied plasma LPLA2 substrate profile using a self-quenched fluorogenic probe specifically designed for LPLA2 in acidic environment

2. LYSOSOMAL PHOSPHOLIPASE A2 (LPLA2)

- LPLA2:
 - Localized in lysosome with optimal pH at 4.5
 - Has both PLA1 and PLA2 activity
 - Also has transacylase activity when an acceptor, such as N-acetyl-sphingosine (NAS), is present
- Detection method:
 - A self-quenched fluorogenic probe is synthesized specifically for PLA2 activity
 - The self-quenched fluorogenic probe is incorporated into liposomes and used as a substrate for LPLA2
 - Reaction is performed under acidic condition

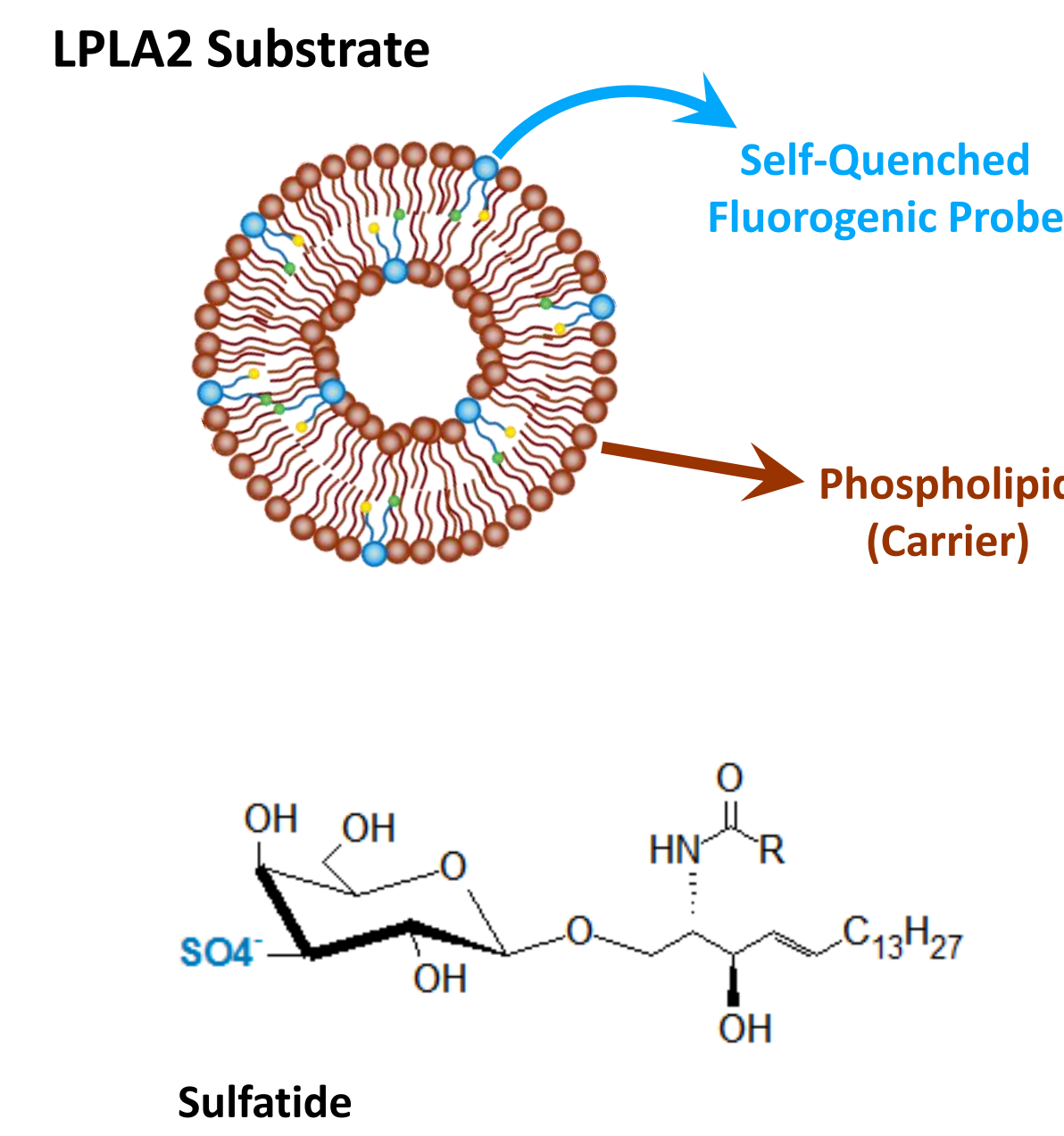


3. LPLA2 & DISEASE

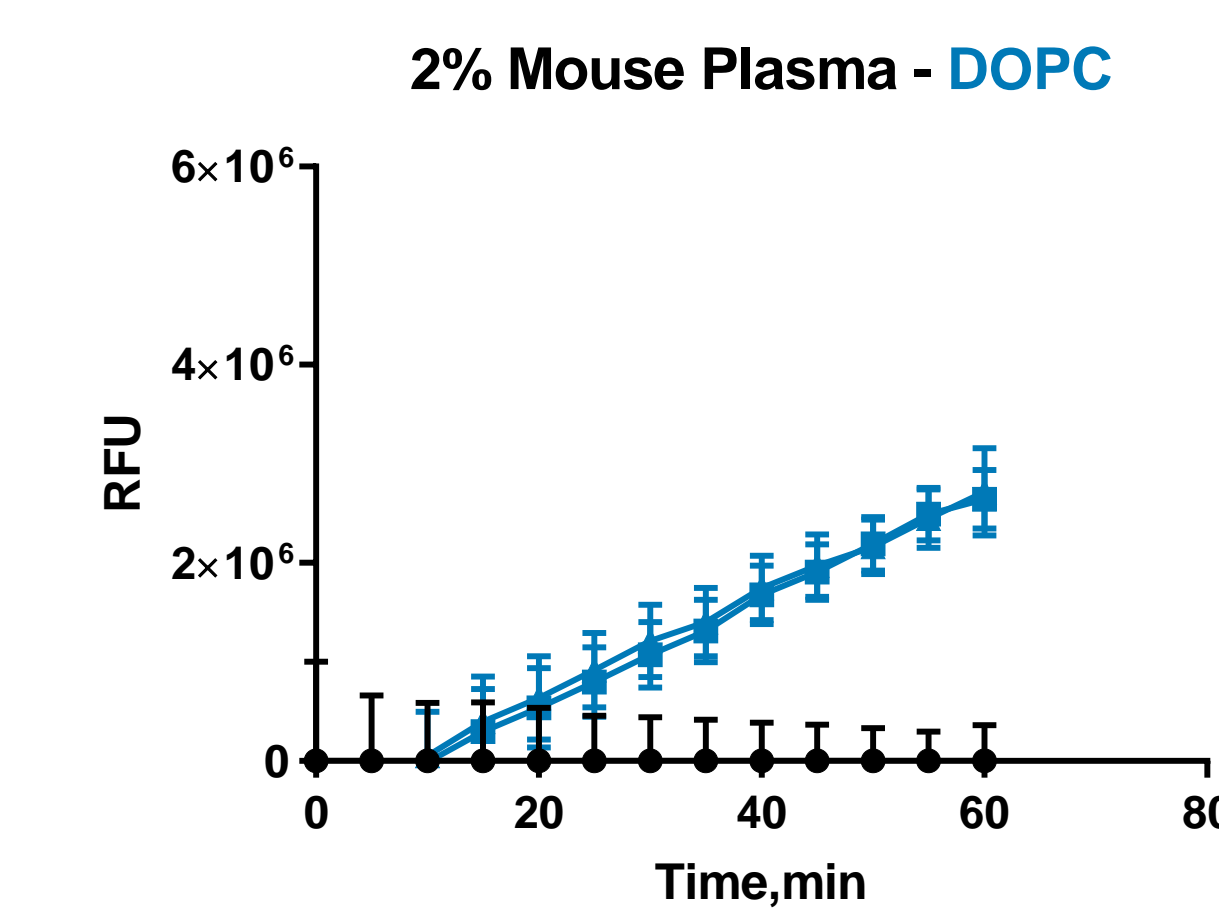
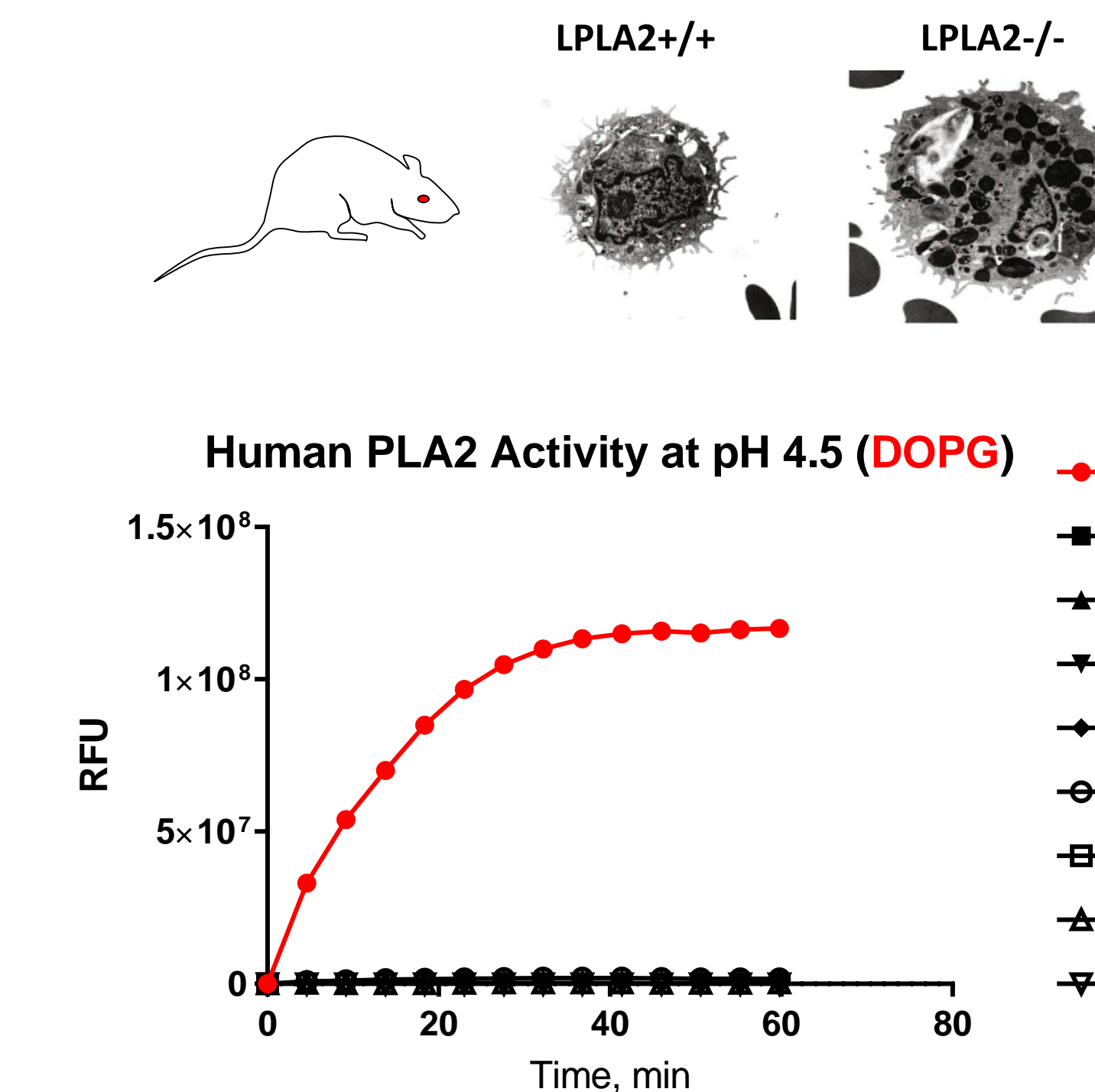
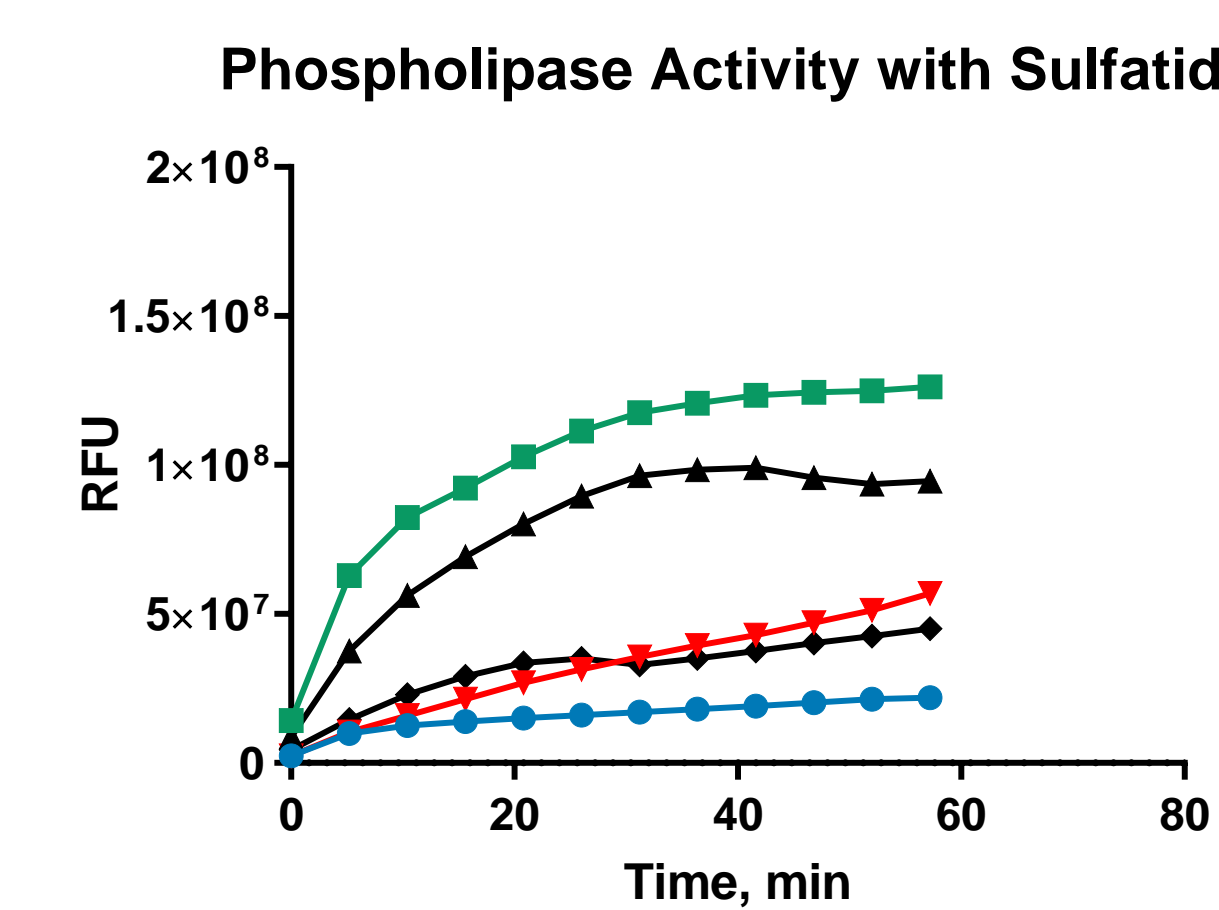
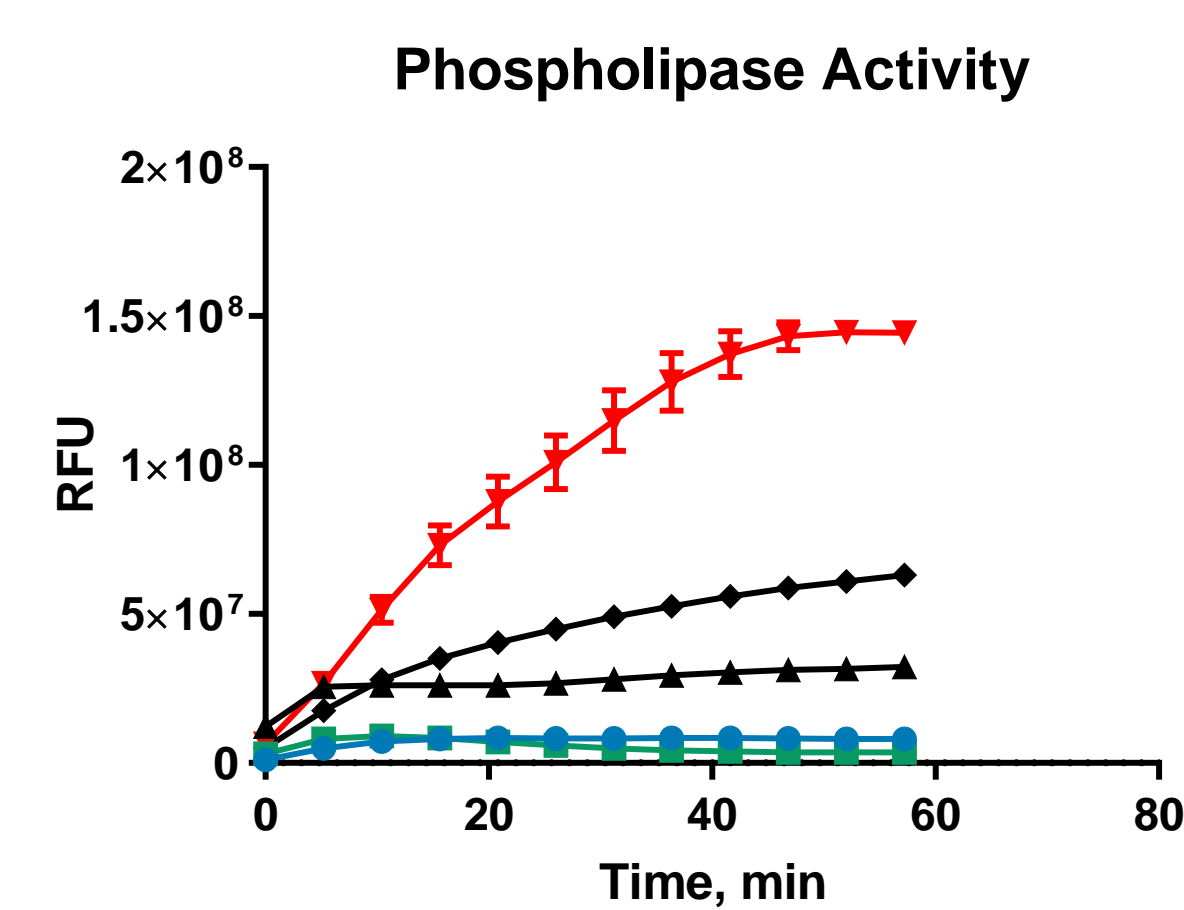


- Drug-induced phospholipidosis (DIPL)
 - Condition of excessive accumulation of intracellular phospholipids caused by common cationic amphiphilic drugs (CADs) on the market
 - CADs significantly inhibit LPLA2 activity in vitro
- Drug-induced lupus (DIPL)
 - Systemic autoimmune disease when immune system attacks own tissues and organs
 - 5-10% lupus is triggered by long-term drug use
 - LPLA2 KO mice express phenotypes similar to lupus

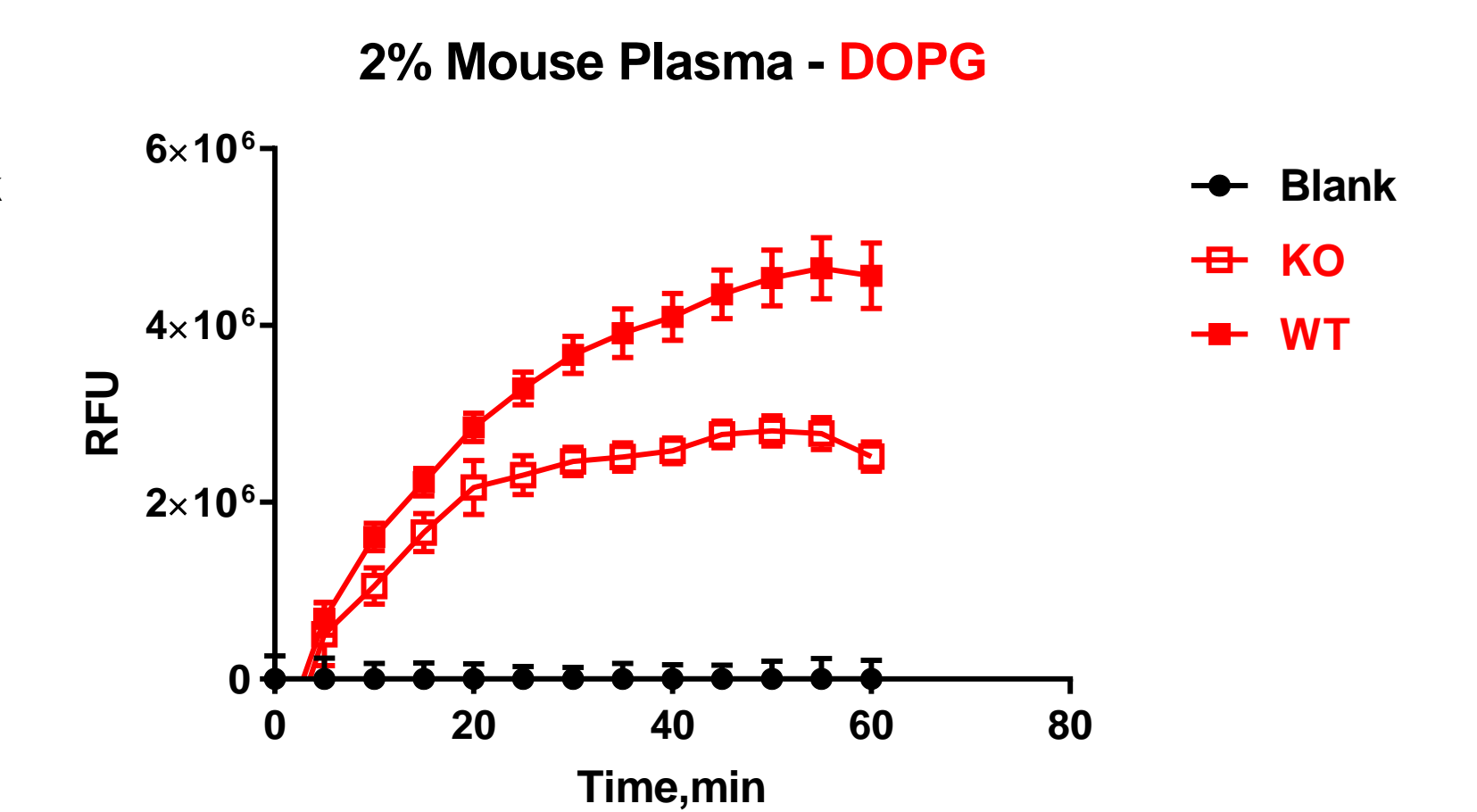
4. LPLA2 ACTIVITY & MEMBRANE PROPERTIES



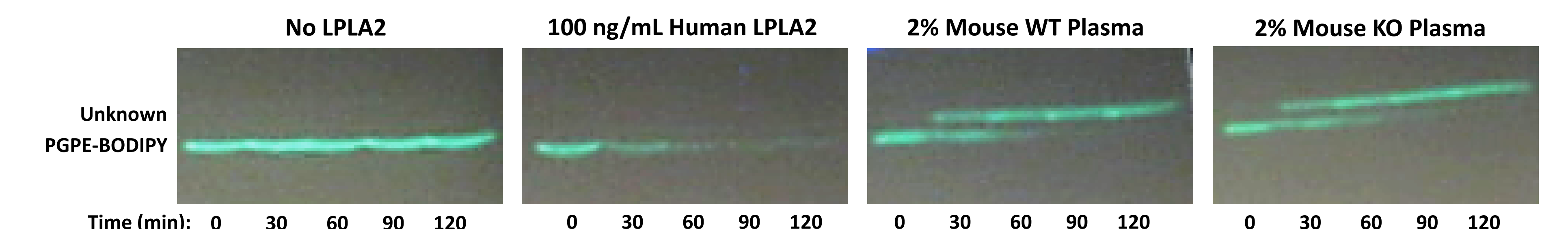
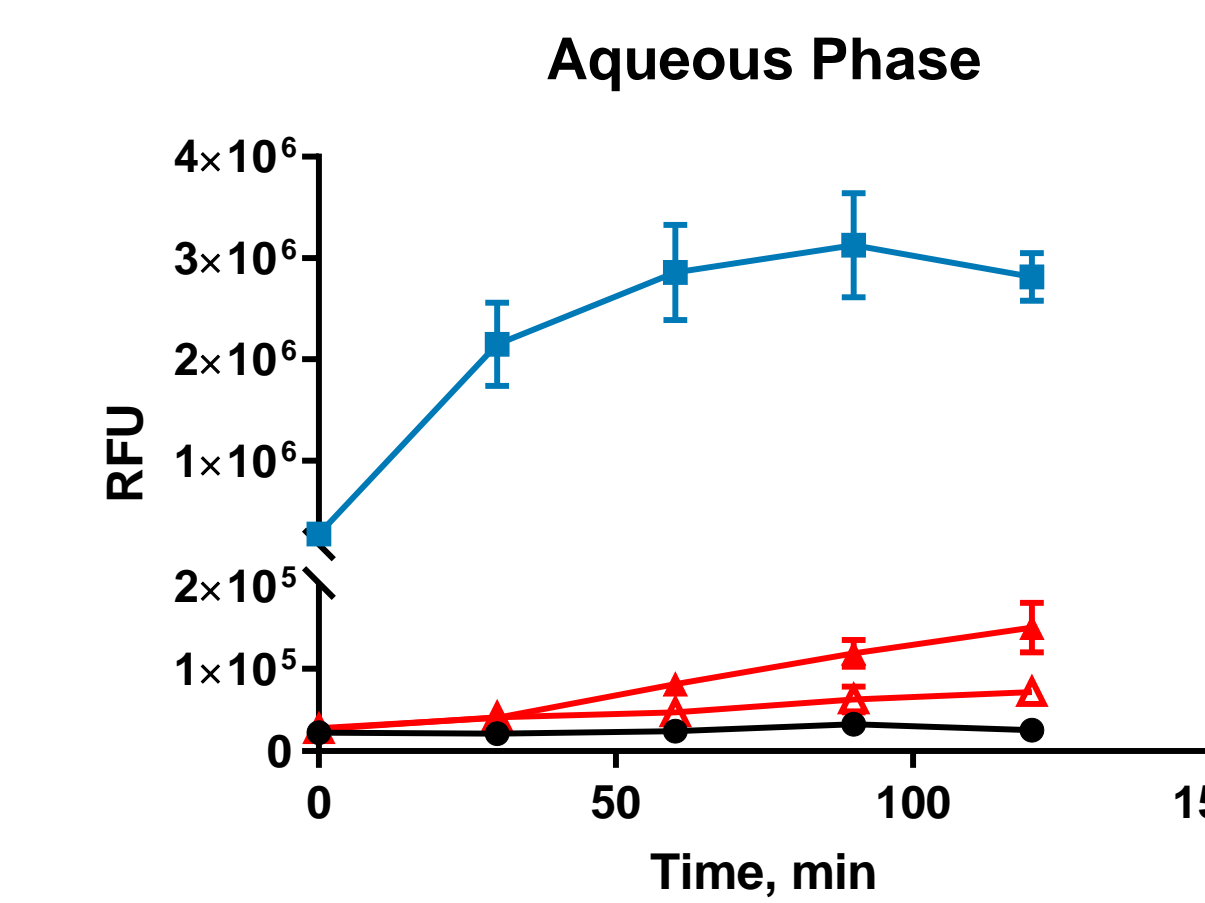
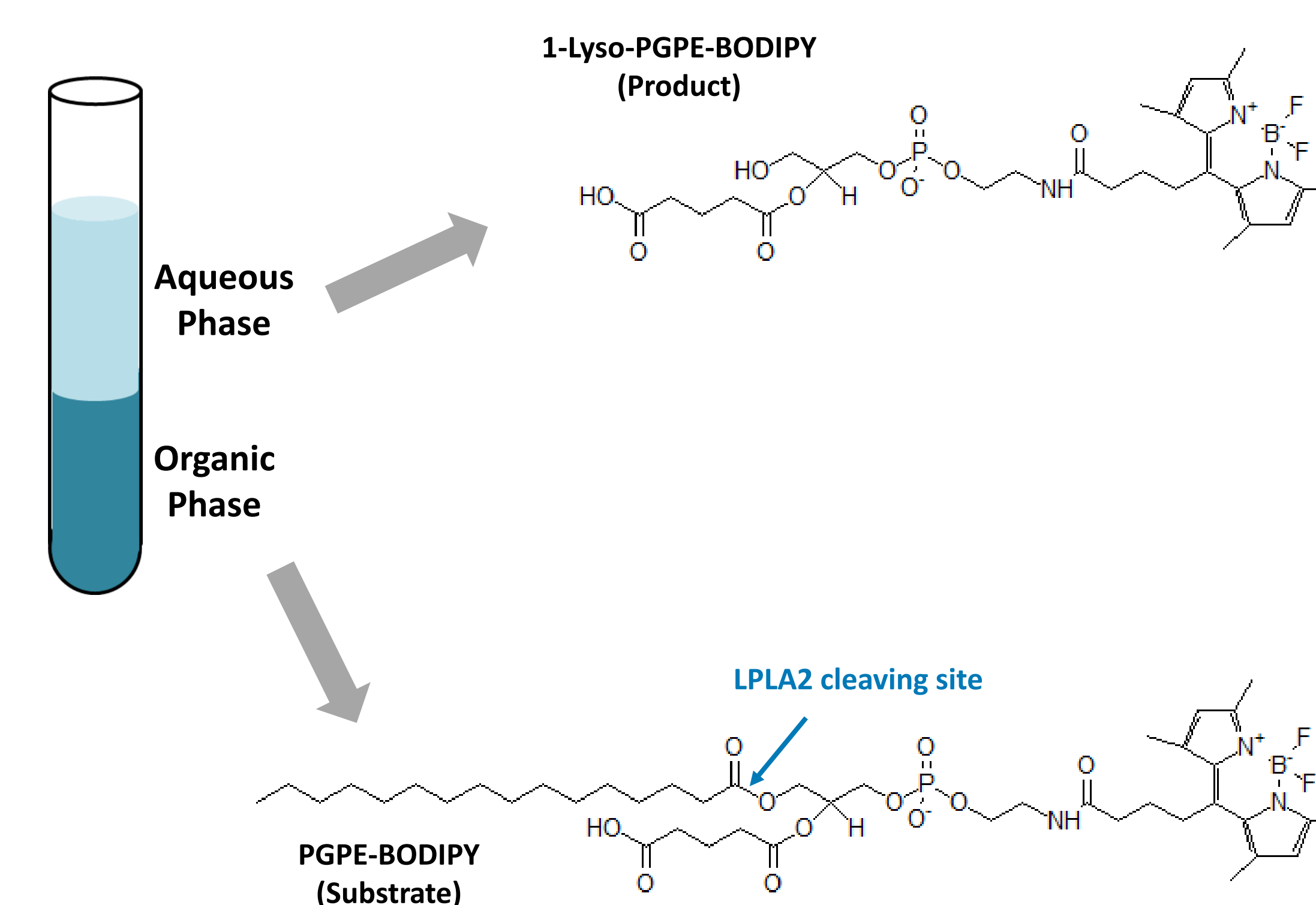
- The self-quenched fluorogenic probe incorporated into negative charged phospholipids such as DOPG & DOPEth results in significantly higher LPLA2 activity
- No significant LPLA2 activity when the self-quenched fluorogenic probe incorporated into neutral phospholipids such as DOPC & DOPE
- The self-quenched fluorogenic probe also shows strong LPLA2 activity when incorporated in the bis(monoacylglycerol)phosphate (BMP), a special late endosome lipid
- The highly negatively charged lipid, sulfatide, significantly enhances the LPLA2 activity towards the self-quenched fluorogenic probe incorporated into neutral phospholipids such as DOPE



- LPLA2 KO mice were generated by systemic deletion of exon 5 (catalytic site) of the LPLA2 gene
- Some non-specific PLA2 activity was detected in both mouse WT & KO plasma
- A negatively charged phospholipid carrier in the self-quenched fluorogenic probe liposome substrate is critical for specific LPLA2 activity detection
- No other PLA2s except LPLA2 metabolize the self-quenched fluorogenic probe incorporated in negatively charged phospholipid (DOPG)



5. LPLA2 sn-1 PHOSPHOLIPASE ACTIVITY



- Substrate (PGPE-BODIPY incorporated in DODPC) depletion observed in both recombinant human LPLA2 and mouse plasma
- Human recombinant LPLA2 cleaves only at the sn-1 position on truncated & oxidized phospholipid fluorescent probe
- Unknown product (possibly by PAF-AH cleaving at the sn-2 position) is formed in mouse plasma
- Mouse WT plasma shows higher aqueous product formation than mouse LPLA2 KO plasma

6. CONCLUSIONS & ACKNOWLEDGEMENTS

- CAD interfering with LPLA2 activity is a promising mechanism for DIPL
- Membrane charge & structure are critical for LPLA2 substrate engagement
- Mouse WT and KO plasma LPLA2 activity presents different substrate profile

- We appreciate Dr. James M Willard from FDA in sharing the Phospholipidosis Working Group database
- We thank our collaborators Dr. Piotr W. Rzepecki for advice
- This study is supported by FDA, SBIR Contract, 5R44FD004052

Presenter Contact Information:
Fok Vun Chan
schan@echelon-inc.com
(801) 588-0455 ext. 378