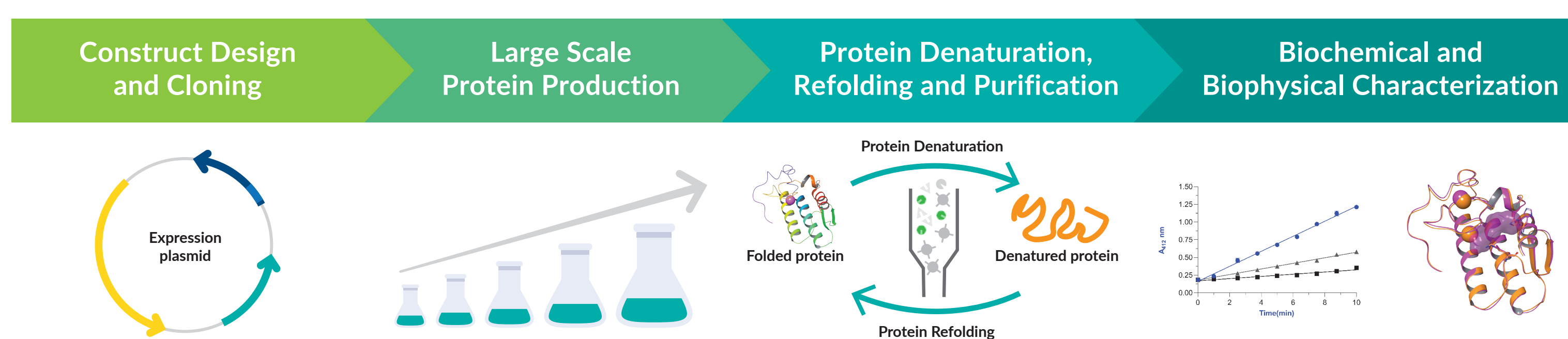


KEY FINDING sPLA₂ enzyme activity was compared in the presence and absence of inhibitors. sPLA₂-IIA apo and inhibitor-bound structures were solved with 1.72 Å and 1.68 Å resolutions, respectively.

INTRODUCTION

Secretory phospholipase A₂ (sPLA₂) catalyzes the hydrolysis of *sn*-2 acyl bond of membrane-bound phospholipids yielding a free fatty acid and a lysophospholipid. The sPLA₂ family contains 10 catalytically active (IB, IIA, IID, IIE, IIF, III, V, X, and XIIA) and one inactive isoform (XIIB) in mammals.^{1,2,3} This family of enzymes are disulfide-rich, low molecular weight, Ca²⁺-requiring lipolytic enzymes with a His-Asp catalytic dyad. Individual mammalian sPLA₂s have distinct enzymatic properties and display distinct tissue expression patterns, suggesting that each enzyme acts on a distinct phospholipid membrane.³ Here we describe the comparative activity of sPLA₂-IIA, sPLA₂-IID, sPLA₂-IIE, sPLA₂-V, and sPLA₂-X in the presence and absence of inhibitors. The closely related isoforms (sPLA₂-IIA, sPLA₂-IID, and sPLA₂-IIE) showed preference for diheptanoyl thio-PE (Thio-PE), whereas the other isoforms (sPLA₂-V and sPLA₂-X) preferred diheptanoyl thio-PC (Thio-PC) as a substrate. Inhibition assays showed that varespladib and AZD2716 inhibited enzyme activity. In the presence of 100 nM varespladib, Thio-PC conversion was abolished for all the sPLA₂s except for sPLA₂-V. However, 100 nM AZD2716 completely inhibited Thio-PC activity for sPLA₂-V, sPLA₂-IID and sPLA₂-IIE followed by residual (~<20%) activity for sPLA₂-IIA and sPLA₂-X. In contrast, 100 nM varespladib showed 100% inhibition for sPLA₂-X and sPLA₂-IIA in the presence of Thio-PE. In addition, we have solved the crystal structures of sPLA₂-IIA, apo and inhibitors bound with 1.72 Å and 1.68 Å resolutions, respectively.

EXPERIMENTAL WORKFLOW



RESULTS

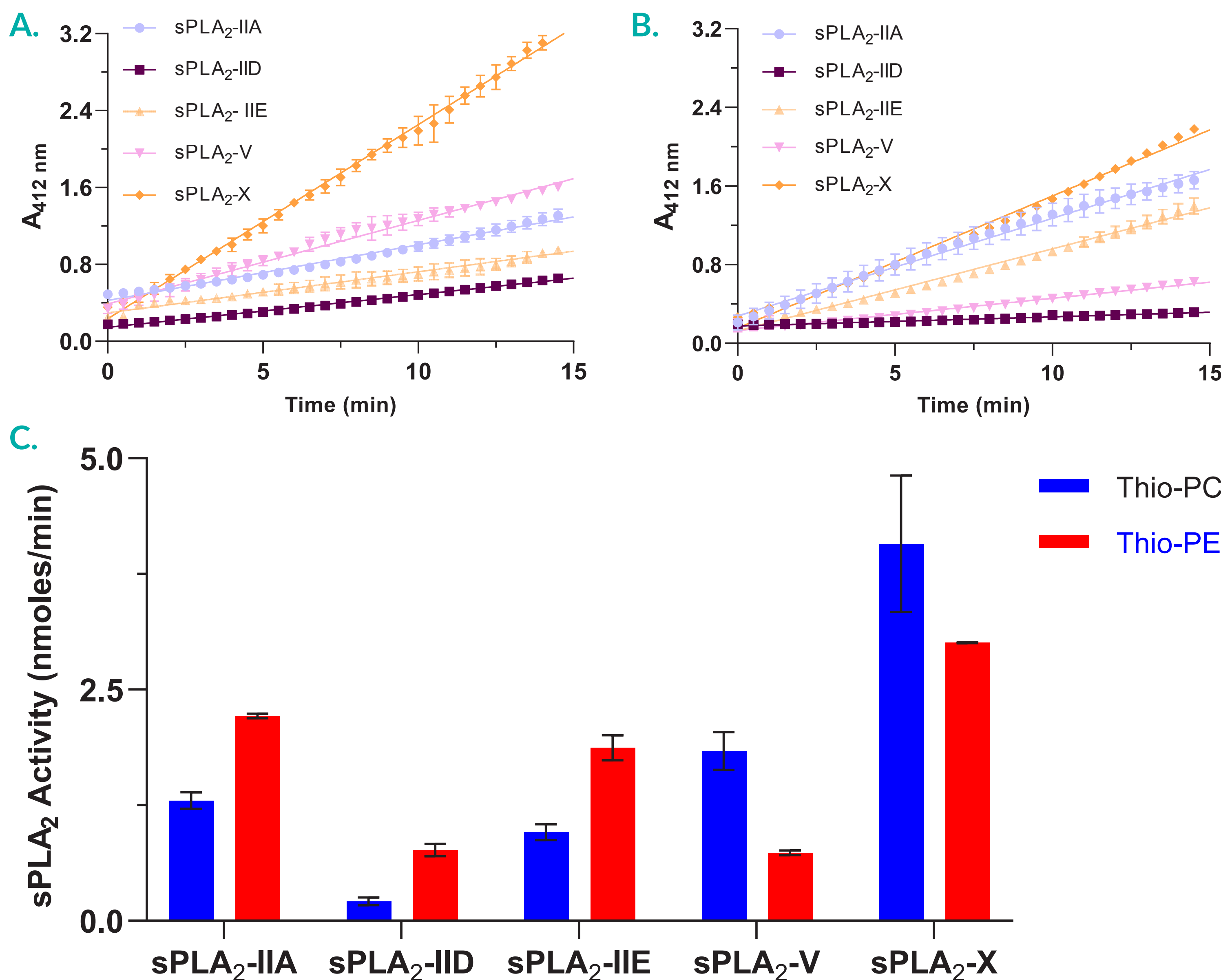


FIGURE 1 – sPLA₂ enzyme substrate specificity. sPLA₂ enzyme activity was determined using Cayman's sPLA₂ Assay Kit (Item No. 765001) with **A)** 1.48 mM Thio-PC (Cayman Item No. 62235) and **B)** 1.48 mM Thio-PE (Cayman Item No. 38245) as substrate. 200 ng of sPLA₂ enzyme was used in all reactions and assay was conducted for 15 minutes. **C)** Substrate preference for the sPLA₂ enzyme activity. sPLA₂-V and sPLA₂-X prefer Thio-PC as substrate, and sPLA₂-IIA, sPLA₂-IID, and sPLA₂-IIE prefer Thio-PE as substrate.

References

- Dennis, E.A. Diversity of group types, regulation and function of phospholipase A2. *J. Biol. Chem.* **269**, 13057-13060 (1994).
- Murakami *et al.* Secreted phospholipase A2 revisited. *J. Biol. Chem.* **150**, 233-255 (2011).
- Murakami *et al.* A new era of secreted phospholipase A2. **56**, 1248-1261 (2015).

RESULTS CONT.

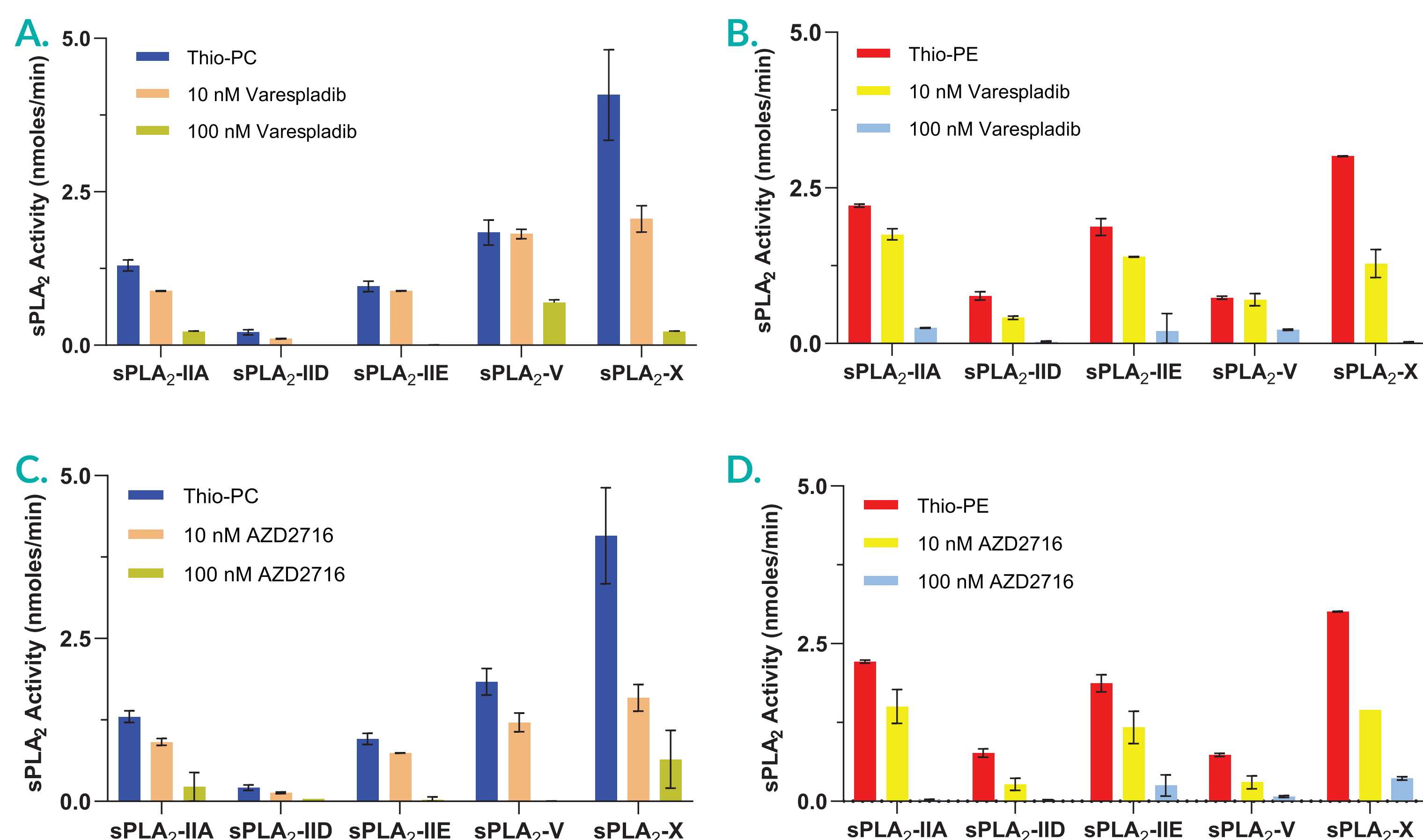


FIGURE 2 – sPLA₂ enzyme activity in the presence of the inhibitor. 200 ng of sPLA₂ was used in the assay with 10 nM and 100 nM of varespladib (Cayman Item No. 18267) and AZD2716 sPLA₂ inhibitors (Cayman Item No. 35067). **A)** and **B)** sPLA₂ enzyme activity inhibited by varespladib. sPLA₂ activity was measured in the presence of 10 nM and 100 nM varespladib in the presence of Thio-PC and Thio-PE, respectively. **C)** and **D)** sPLA₂ enzyme activity inhibited by AZD2716. sPLA₂ activity was measured in the presence of 10 nM and 100 nM AZD2716 in the presence of Thio-PC and Thio-PE, respectively.

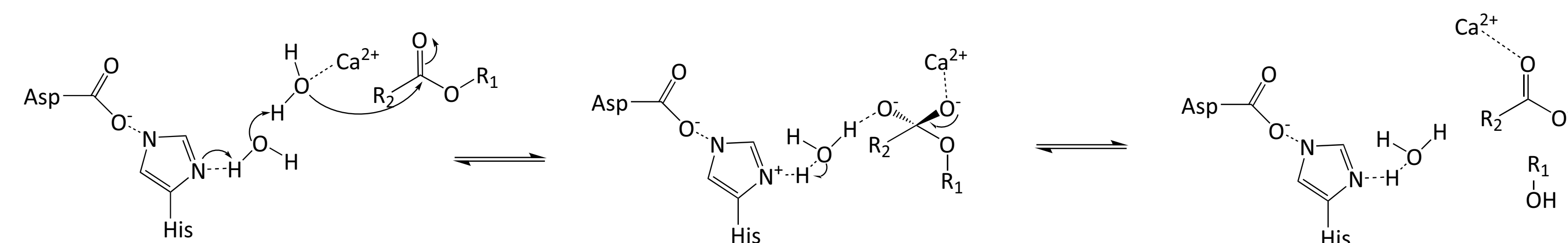


FIGURE 3 – Ca²⁺-dependent sPLA₂ enzyme reaction mechanism. The water molecule is activated by the presence of a histidine/aspartic acid dyad in a Ca²⁺-dependent manner.

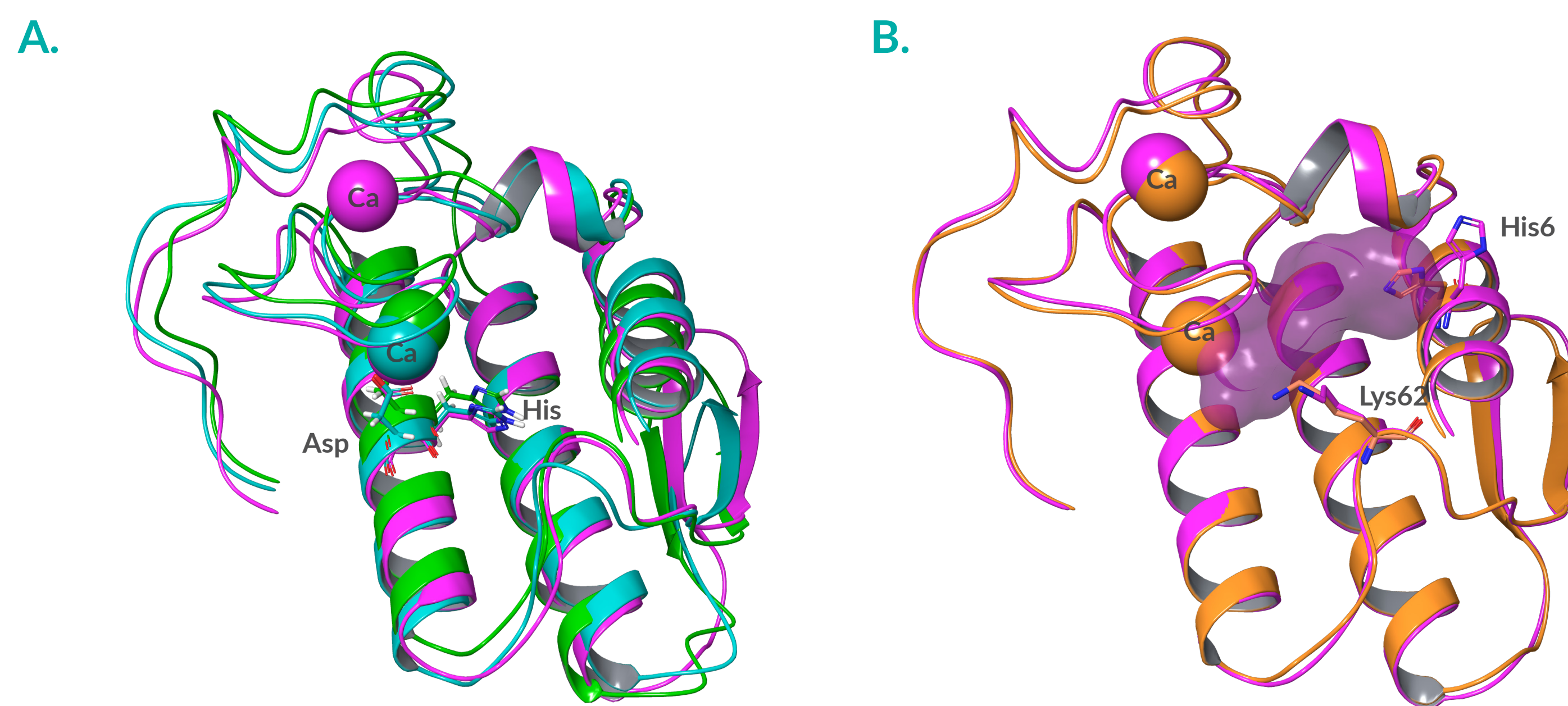


FIGURE 4 – Structure comparison of sPLA₂ enzymes. **A)** Overlay with ribbon model of sPLA₂-IIA (1.68 Å crystal structure solved by Cayman SBBC (unpublished) (magenta), sPLA₂-IIE (5WZW.pdb)(cyan), and sPLA₂-X (6G5J.pdb)(green). Calcium is shown as sphere in corresponding color. Catalytic His and Asp residues in active site shown in the corresponding colors as stick models. **B)** Overlay with ribbon models of sPLA₂-IIA complexed with inhibitor shown as a space filling model (1.68 Å crystal structure solved by Cayman SBBC (unpublished) (magenta).

TABLE 1 – Variables obtained from SPR analysis using Biacore™ Insight software using varespladib inhibitor.

	sPLA ₂ -IIA	sPLA ₂ -X
1:1 binding k_b (M ⁻¹ s ⁻¹)	2.05e+06	8.21e+05
k_d (s ⁻¹)	2.01e-02	1.65e-03
K_D (M)	9.77e-09	2.02e-09

CONCLUSIONS

- sPLA₂-IIA, sPLA₂-IID, sPLA₂-IIE, sPLA₂-V, and sPLA₂-X enzymes were purified from inclusion bodies and refolded.
- sPLA₂-IIA, sPLA₂-IID, sPLA₂-IIE preferred Thio-PE, and sPLA₂-V, and sPLA₂-X preferred Thio-PC as substrate.
- Varespladib and AZD2716 inhibited sPLA₂ enzyme activity.
- Crystal structures of sPLA₂-IIA, apo and inhibitor bound were solved with 1.72 Å and 1.68 Å resolution, respectively.



Download this poster & learn more about our protein services