

Lysosomal Phospholipase A2 (LPLA2) Inhibitor Screen

K-7000I (384 tests)

Support: echelon@echelon-inc.com

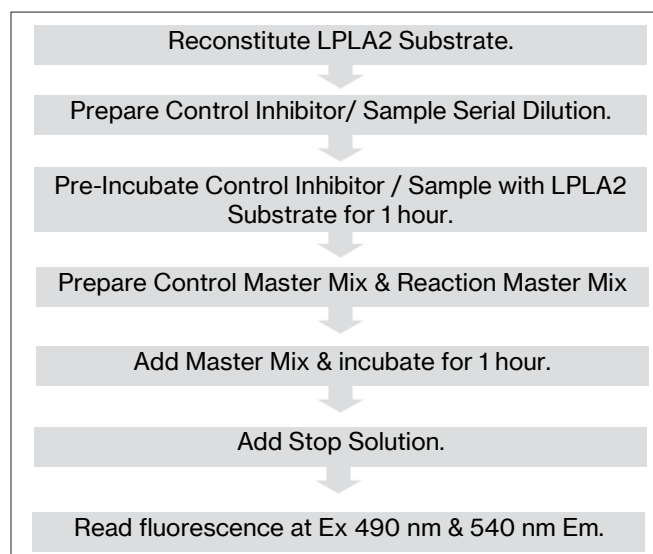
Description: Description: LPLA2 Inhibitor Screen (K-7000I) is a homogeneous assay designed to screen molecules that potentially trigger phospholipidosis (PL) through inhibition of LPLA2.

Storage: Upon receipt, store assay kit at -20 °C. Opened and reconstituted reagents are less stable. Please see assay notes for additional information.

Materials Provided

| Catalog # | Description | Quantity |
|-----------|---|------------|
| K-7001I | Human LPLA2 (Lyophilized) | 1 vial |
| K-7002I | LPLA2 Substrate (Dried) | 1 vial |
| K-7003I | 4X Reaction Buffer | 1 bottle |
| K-7004I | Control Inhibitor – Amiodarone (1 µmol) | 1 vial |
| K-7005I | 5X Stop Solution | 1 bottle |
| K-GS01 | Protein Stabilizer | 2 x 600 mL |
| K-DIL7 | Diluent | 1 bottle |
| --- | 384-well Low Volume Black Plate | 1 plate |
| --- | Plate Sealers | 2 seals |

Quick Protocol



Additional Materials Provided by User

- Pipettes (capable of delivering between 5 and 1,000 µL solution with appropriate tips)
- Reagent grade water
- Dimethyl sulfoxide (DMSO)
- (optional) plate shaker
- Microplate reader capable of reading with 490 nm excitation & 540 nm emission (see assay note #1).

Echelon Biosciences products are sold for research and development purposes only and are not for diagnostic use or to be incorporated into products for resale without written permission from Echelon Biosciences. Materials in this publication, as well as applications and methods and use, may be covered by one or more U.S. or foreign patents or patents pending. We welcome inquiries about licensing the use of our trademarks and technologies at echelon@echelon-inc.com

Background

The human lysosomal phospholipase A2 (LPLA2) enzyme is responsible for normal lipid metabolism. It is unique from other known PLA2s in that LPLA2 is only active in an acidic environment such as the lysosome (~pH 4.5). Furthermore, it demonstrates negligible activity at pH 7.4 which is normally found for the cytosol. In addition, LPLA2 acts as a transacylase with C2-ceramide as a preferred lipophilic acceptor.¹ Research has shown LPLA2 to be involved in drug-induced phospholipidosis (DIPL).² Phospholipidosis (PL) is a condition resulting from the excessive accumulation of intracellular phospholipids, causing tissue inflammation and organ damage. PL commonly manifests in patients taking cationic amphiphilic drugs (CADs) such as fluoxetine (Prozac[™], Sarafem) and Amiodarone. The FDA has determined DIPL a serious drug safety issue.³ Although the complete cause of DIPL is unknown, evidence is accumulating that DIPL is the result of 3. certain CADs directly inhibiting the actions of the lysosomal phospholipase A2 (LPLA2). Thus, LPLA2 inhibition is a potential predictor of drug-induced phospholipidosis.

Assay Design

Echelon's LPLA2 Inhibitor Screen (K-7000I) is designed to screen a drug's ability to inhibit LPLA2 activity in vitro, a potential predictor of drug-induced phospholipidosis, in a high throughput format (HTS) using "smart probe" technology. A quenched fluorogenic substrate is liberated by LPLA2 activity, resulting in a bright fluorescent product (Figure 1). Therefore, the fluorescent signal is significantly reduced in the presence of an LPLA2 inhibitor. This direct biochemical approach provides a quantitative measurement in a robust and simple to use in vitro plate-based assay, providing a greatly improved throughput when compared to traditional microscopy methods of tissue cultured cells. The assay has been validated with a group of known PL-inducing and non-PL inducing drugs. A known PL-inducing cationic amphiphilic drug (CAD), amiodarone, is included in the assay as a control.

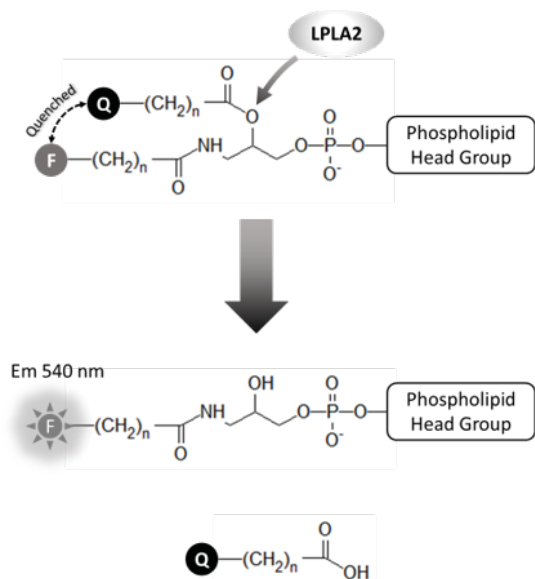


Figure 1. Assay Design
Q = Quencher, F = Fluorophore

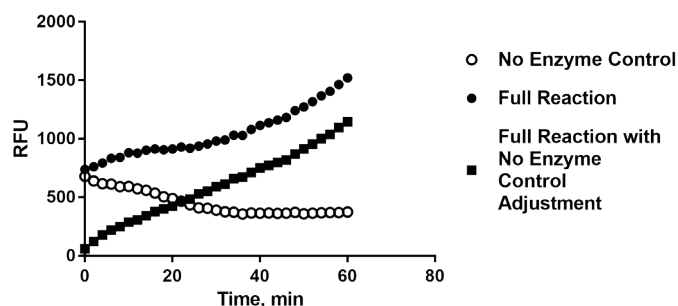
Assay Performance

For best results, please follow the protocols provided. Not following instructions may result in suboptimal performance of the kit and failure to produce accurate data.

Assay Notes

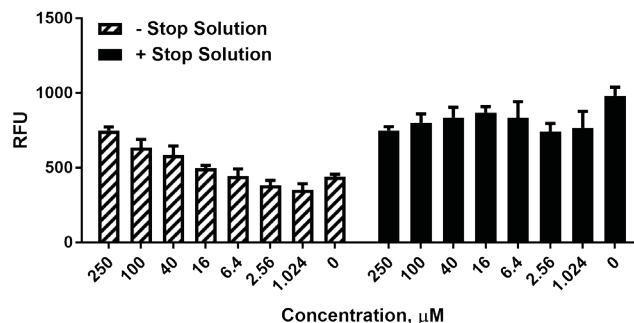
1. The LPLA2 Inhibitor Screen is designed to be run as a one-time use only assay. If the inhibitor screen is to be run on separate occasions, the remaining kit reagents can be stored at -20 °C. However, a reduction in the strength of the fluorescent signal has been observed after freeze-thaw. Always include the Control Inhibitor as a standard to ensure accurate results. DO NOT COMBINE different vials of reconstituted LPLA2 Substrate (K-7002I).
2. If an alternative solvent to DMSO is used in preparing the test compounds, include a "solvent only" control without the test compound to detect potential solvent effects. To ensure assay reproducibility, it is advised to use consistent solvent concentrations between experiments.
3. The LPLA2 Inhibitor Screen can also be read in kinetic mode, reading every 5 minutes for 1 hour, using the same excitation/emission settings as step 17. However, photobleaching might occur as demonstrated by the "No Enzyme Control" (Figure 2). Including this "No Enzyme Control" allows one to account for photobleaching and is recommended for data analysis.

Figure 2. Photobleaching



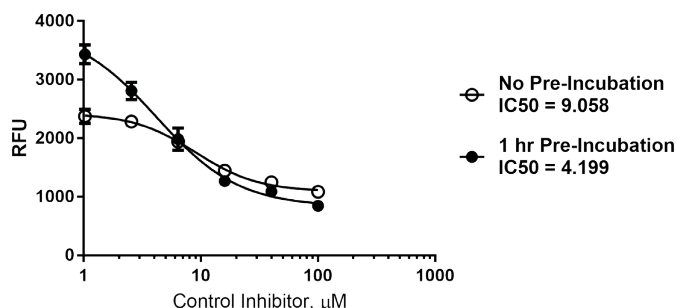
4. Fluorescent interference, including self-quenching and inhibitor related auto-fluorescence, may result in false positives or false negatives when screening. If fluorescent interference is a concern, screening the inhibitor without the LPLA2 enzyme will serve as control. We have observed that the addition of Stop Solution (K-7005I) significantly improves this self-quenching or auto-fluorescence effect when screening drugs (Figure 3).

Figure 3. Clomipramine Background (No Enzyme Controls)



- For optimal results, we suggest using a consistent pre-incubation time (step 8). Varying the pre-incubation time significantly affects the IC₅₀ of the Control Inhibitor (K-7004I, Figure 4).

Figure 4. Drug-Substrate Pre-Incubation



Assay Protocol

Reagent volumes can be adjusted when not using the entire assay, please see assay note #1 for more details.

- Place LPLA2 Enzyme (K-7001I) on ice. Bring the remaining reagents to room temperature (RT).
- Reconstitute the Control Inhibitor (K-7004I) with 20 μL of DMSO for a 50 mM Stock. Vortex to mix. Keep at RT.
- Add 2 mL of reagent grade water to the LPLA2 Substrate (K-7002I). VORTEX VIGOROUSLY FOR 5 MINUTES to form liposomes. Keep at RT.
- Prepare the 7-point Control Inhibitor serial dilutions in microcentrifuge tubes. See Table 1. Keep at RT and protect from light.

Table 1, Control Inhibitor Dilutions

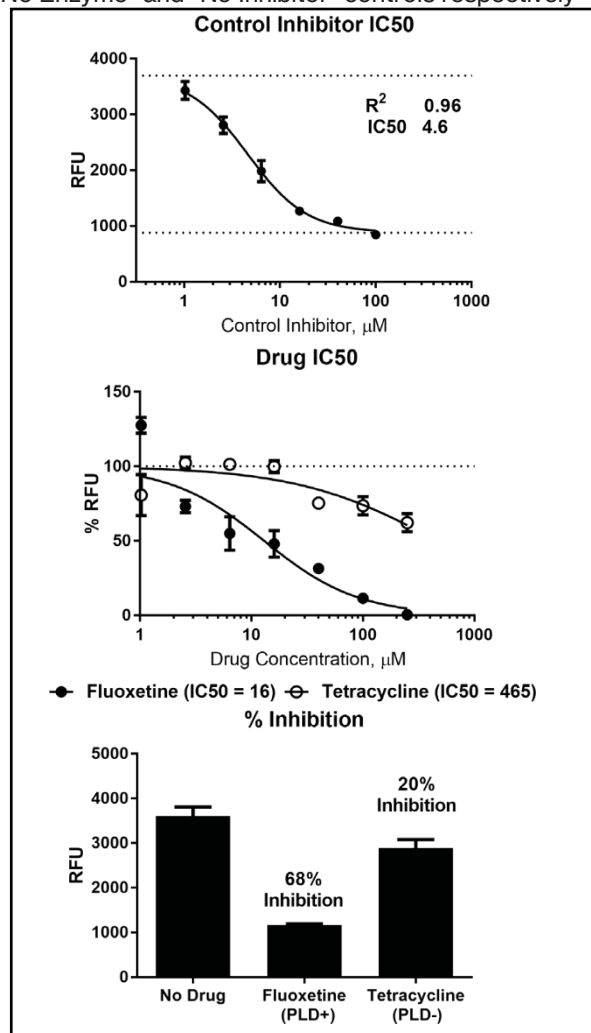
| Standard | 1X Control Inhibitor | 16X Control Inhibitor | Previous Dilution Needs | DMSO to add |
|----------|----------------------|-----------------------|-------------------------|-------------|
| A | 100 μM | 1600 μM | 5 μL of K-7004I | 151 μL |
| B | 40 μM | 640 μM | 30 μL of 1600 μM | 45 μL |
| C | 16 μM | 256 μM | 30 μL of 640 μM | 45 μL |
| D | 6.4 μM | 102.4 μM | 30 μL of 256 μM | 45 μL |
| E | 2.56 μM | 40.96 μM | 30 μL of 102.4 μM | 45 μL |
| F | 1.024 μM | 16.384 μM | 30 μL of 40.96 μM | 45 μL |
| G | 0 μM | 0 μM | None | 45 μL |

- Prepare the compounds to be tested in DMSO at 16X the desired final concentration. See assay note #2 if you use a solvent other than DMSO.

- Add 1 μL/well of the 16X Control Inhibitor serial dilutions (step 4.) and 1 μL/well of the 16X compounds to be tested. For the “No LPLA2” control, add 1 μL/well of DMSO or the solvent of your choice. It is suggested that samples be run in triplicate. A suggested plate layout is shown below (Table 2). See assay note #4 for information on drug auto-fluorescence.
- Add 4 μL of the LPLA2 Substrate (K-7002I), prepared in step 3, to all wells.
- Cover the plate with a plate seal and pre-incubate the control inhibitors or test compounds with the LPLA2 substrate for 1 hour at RT with shaking and protected from light. See assay note #5 for more information on pre-incubation.
- Reconstitute the Human LPLA2 (K-7001I) with reagent grade water with the volume labeled on the vial. GENTLY PIPETTE UP & DOWN AND INCUBATE ON ICE FOR 10 MINUTES to ensure full reconstitution. Keep on ice.
- Control Master Mix: Prepare 137.5 μL of “Control Master Mix” solution by combining the following in a Microcentrifuge tube:
 - 50 μL 4X Reaction Buffer (K-7003I)
 - 50 μL Diluent (K-DIL7)
 - 37.5 μL Reagent Grade Water
 - Vortex to mix. Keep at RT.
- Prepare 1,700 μL of the human LPLA2 by further diluting the reconstituted human LPLA2 (step 9) with Diluent (K-DIL7) according to the dilution factor labeled on the vial. Vortex for 3 seconds to mix. Keep at RT.
- Reaction Master Mix: Prepare 4,675 μL of “Reaction Master Mix” by combining the following in an appropriate container.
 - 1,700 μL 4X Reaction Buffer (K-7003I)
 - 1,700 μL LPLA2 Enzyme diluted in Diluent (step 11.)
 - 1,275 μL Reagent Grade Water
 - Vortex to mix. Keep at RT.
- Following the 1-hour LPLA2 Substrate / Inhibitor pre-incubation step, add 11 μL of the Control Master Mix into the “No LPLA2” control wells. Add 11 μL of the Reaction Master Mix into the remaining wells to start the reactions.
- Cover the plate with a plate seal and incubate at RT for 1 hour with shaking. Protect from light. If the plate shaker is not available, gently tap plate to mix before incubation. See assay note #3 for additional information regarding kinetic analysis.
- After 1 hour, add 4 μL of the 5X Stop Buffer (K-7005I) to each well to stop the reactions. Incubate at RT in the dark for 15 minutes. Stop Solution must be added prior to fluorescence recording for end-point assay. See assay note #4 for additional information.
- Record the fluorescence using 490 nm excitation and 540 nm emission settings. See assay note #3 for additional information. additional information regarding kinetic analysis.

Data Analysis

Examples of the Control Inhibitor's IC₅₀ curve (Top), an end-point inhibitor screen (Middle), and the analysis of the drugs, fluoxetine and tetracycline, and their respective IC₅₀ curves (Bottom). IC₅₀ is analyzed using the log(agonist) vs. response – variable slope curve fit from GraphPad Software with top & bottom constrained by “No Enzyme” and “No Inhibitor” controls respectively



References

1. Abe, A.; Gregory, S.; Lee, L.; Shayman, J. A., Use of sulfobutyl ether beta-cyclodextrin as a vehicle for D-threo-1-phe- nyl-2-decanoylamino-3-morpholinopropanol- related glucosylceramide synthase inhibitors. Anal Biochem 2000, 287 (2), 344-7.
2. Halliwell, W. H., Cationic amphiphilic drug-induced phospholipidosis. Toxicol Pathol 1997, 25 (1), 53-60.
3. Shayman, J. A.; Kelly, R.; Kollmeyer, J.; He, Y.; Abe, A., Group XV phospholipase A(2), a lysosomal phospholipase A(2). Prog Lipid Res 2011, 50 (1), 1-13.

Related Products

| Products | Catalog Number |
|----------------------|----------------|
| Enzyme | |
| Human LPLA2 | E-7000 |
| Assay | |
| LPLA2 Activity Assay | K-7000A |

Please visit our website at www.echelon-inc.com for more enzyme and lipid products.

Table 2, Suggested Plate Layout

| | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 | 13 | 14 | 15 | 16 | 17 | 18 | 19 | 20 | 21 | 22 | 23 | 24 |
|---|------------|---|---|-------------|---|---|-------------|---|---|-------------|----|----|-------------|----|----|-------------|----|----|--------------|----|----|--------------|----|----|
| A | No LPLA2 | | | Compound 9 | | | Compound 25 | | | Compound 41 | | | Compound 57 | | | Compound 73 | | | Compound 89 | | | Compound 105 | | |
| B | Standard A | | | Compound 10 | | | Compound 26 | | | Compound 42 | | | Compound 58 | | | Compound 74 | | | Compound 90 | | | Compound 106 | | |
| C | Standard B | | | Compound 11 | | | Compound 27 | | | Compound 43 | | | Compound 59 | | | Compound 75 | | | Compound 91 | | | Compound 107 | | |
| D | Standard C | | | Compound 12 | | | Compound 28 | | | Compound 44 | | | Compound 60 | | | Compound 76 | | | Compound 92 | | | Compound 108 | | |
| E | Standard D | | | Compound 13 | | | Compound 29 | | | Compound 45 | | | Compound 61 | | | Compound 77 | | | Compound 93 | | | Compound 109 | | |
| F | Standard E | | | Compound 14 | | | Compound 30 | | | Compound 46 | | | Compound 62 | | | Compound 78 | | | Compound 94 | | | Compound 110 | | |
| G | Standard F | | | Compound 15 | | | Compound 31 | | | Compound 47 | | | Compound 63 | | | Compound 79 | | | Compound 95 | | | Compound 111 | | |
| H | Standard G | | | Compound 16 | | | Compound 32 | | | Compound 48 | | | Compound 64 | | | Compound 80 | | | Compound 96 | | | Compound 112 | | |
| I | Compound 1 | | | Compound 17 | | | Compound 33 | | | Compound 49 | | | Compound 65 | | | Compound 81 | | | Compound 97 | | | Compound 113 | | |
| J | Compound 2 | | | Compound 18 | | | Compound 34 | | | Compound 50 | | | Compound 66 | | | Compound 82 | | | Compound 98 | | | Compound 114 | | |
| K | Compound 3 | | | Compound 19 | | | Compound 35 | | | Compound 51 | | | Compound 67 | | | Compound 83 | | | Compound 99 | | | Compound 115 | | |
| L | Compound 4 | | | Compound 20 | | | Compound 36 | | | Compound 52 | | | Compound 68 | | | Compound 84 | | | Compound 100 | | | Compound 116 | | |
| M | Compound 5 | | | Compound 21 | | | Compound 37 | | | Compound 53 | | | Compound 69 | | | Compound 85 | | | Compound 101 | | | Compound 117 | | |
| N | Compound 6 | | | Compound 22 | | | Compound 38 | | | Compound 54 | | | Compound 70 | | | Compound 86 | | | Compound 102 | | | Compound 118 | | |
| O | Compound 7 | | | Compound 23 | | | Compound 39 | | | Compound 55 | | | Compound 71 | | | Compound 87 | | | Compound 103 | | | Compound 119 | | |
| P | Compound 8 | | | Compound 24 | | | Compound 40 | | | Compound 56 | | | Compound 72 | | | Compound 88 | | | Compound 104 | | | Compound 120 | | |