

KINOMEScan™ KdELECT: Next Steps with Compound Preparation and Shipping Instructions (if necessary)

Step 1. Complete Project Quote:

- **Sign Page 2:** Include payment information: Bill To address and/or email address
- **Fax, or Email,** pages 1 and 2 to:
 - **Fax:** (858)-630-4600 *ATTN: Aimee Allen*
 - **Email:** orders@KINOMEScan.com

Step 2. Complete Service Request Form (SRF):

- **Project Quote #:** Enter Project Quote Number from signed Quote (ex. "MAXXP99999A")
- **Compound Name:** Enter Name of each compound

NOTE: Please complete only the columns that apply

For Solids Complete:

- **µL of DMSO to add to make a 10mM stock solution:** Enter microliters (µL) of DMSO required to make a 10 millimolar (mM) stock solution for each Compound
- **Compound Weight:** Enter Weight in milligrams (mgs) for each Compound - (optional)

For Liquids Complete:

- **Stock Concentration:** For Liquids enter Stock Concentration (mM) for each Compound
- **Stock Volume:** For Liquids enter Stock Volume (µL) for each Compound

For Plated Liquids Complete:

- **Well Location:** Enter Well Location of each Compound
- **Plate ID #:** Enter Plate ID Number for each compound
- **KINOMEScan Gene Symbol:** Enter KINOMEScan Gene Symbol as shown on the Primary screen data report
- **Compound Onsite at KINOMEScan:** Enter "Yes" if KINOMEScan's Compound Management Team confirmed there is sufficient quantity of each compound on site to perform the requested number of Kd determinations. Enter "No" if there is not a sufficient amount of each compound to perform the requested number of Kd determinations.
- **Comments:** Enter any Comments about the screen you may have
- **Send Data Report To:** Enter the email addresses to those receiving the data report

NOTE: Only Complete Shipping and Tracking Information if Sending Additional Compound(s) or this is a first time screen:

- **Shipped by:** Enter the email address of the contact that shipped the compound(s)
- **Tracking Number:** Enter tracking number for KINOMEScan's Compound Management Team

Step 3. Email Excel copy of SRF to: orders@KINOMEScan.com

Step 4. Include a Hardcopy of SRF with Compound Shipment

Compound Preparation and Shipping instructions: (if necessary)

	For Liquids	For Solids
Number of Kinases per Compound	Minimum Volume (µL)	Minimum Weight (mgs)
≤ 6 kinases	50µL	1-2mgs
> 6 kinases	50µL + 6µL per each additional kinase*	2-3mgs

*EXAMPLE: Volume required for 10 kinases per compound	
first 6 kinases	50µL
Additional 4 kinases	6µL* 4 kinases = 24µL
50µL + 24µL = 74µL of 10mM in 100%DMSO	

Step 5. Ship compound(s) to the following address:*NEW ADDRESS:**

KINOMEScan, a division of DiscoverX
Compound Management
11180 Roselle Street, Suite D
San Diego, CA 92121

Tel: (800)-644-5687

*****An email confirmation will be sent upon receipt of the compound(s)*****