

KINOMEScan™ Primary Screen: Next Steps with Compound Preparation and Shipping instructions

Step 1. Complete Project Quote:

- **Sign Page 2:** Include Payment information, Bill To mailing 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
- **Primary Screening Concentration:** Enter Screening Concentration in μM for each Compound (ex. "0.1, 1.0 or 10.0, etc.")

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
- **Comments:** Enter any Comments about the screen you may have
- **Send Data Report To:** Enter the email addresses to those receiving the data report
- **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:

- Less than 10 compounds:
 - 50 μL 10mM 100% DMSO in labeled tube(s) or labeled vial(s).
 - or -
 - 1-2 mg in labeled vial(s). Include volume required to make a 10mM stock solution in the Service Request Form
- More than 10 compounds:
 - 50 μL 10mM 100% DMSO in a labeled, **polypropylene**, 96 well plate (Costar 3363 or similar). Include well locations in the Service Request Form
 - *Do not send solids in plates*

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 compound(s)*****