



DiscoverRx Next Steps with Compound Preparation and Shipping Instructions

*****Please note: Projects will not be scheduled until receipt of both the compound(s) and P.O.*****

Step 1. Send in Payment Information

- **Purchase Order:** Include Payment information, Bill To and/or email address; reference DRX quote # (ex. DRX-11100)
- **Fax, or Email,** to DiscoverRx:
 - **Fax:** +44 121 260 6143
 - **Email:** EUservices@discoverx.com

Step 2. Complete Compound Shipping Document (CSD):

- **Quote #:** Enter Quote Number (ex. DRX-11100)
- **Compound I.D.:** Enter reference # of each compound
- **Final Concentration to be Tested:** Enter final highest screening concentration to be tested in μM for each compound
- **Company Originating From:** Enter your company information
- **Customer Name:** Enter your name and email address
- **Report Issued To:** Enter the email addresses to those receiving the data report
- **! Complete the Custom Declaration Form (Page 2 of the Compound Shipping Document)**

NOTE: Please complete only the columns that apply

For Solids (**up to 10*) Complete:

- **Weight Shipped**
- **Molecular Weight**
- **Recommended Solvent** (DMSO, PBS or water)
- **Recommended Stock Concentration**
- **Recommended Storage Temperature** for Dry and Reconstituted Compound
 - Provide solids in labeled tube(s)
 - If more than 10 compounds will be delivered in powder form, an additional cost of 50 USD per additional compound will apply

For Liquids (**up to 10*) Complete:

- **Vehicle:** For Liquids, enter the solvent used to reconstitute each compound
- **Stock Volume:** For Liquids enter Stock Volume (μL) for each compound
- **Stock Concentration:** For Liquids enter Stock Concentration for each compound
- **Liquid Storage Temp:** Enter Storage Temperature for each compound
 - Provide liquids in labeled vial(s)

For Plated Liquids (**more than 10*) Complete:

- ****Please inquire regarding optimal plate setup****
- **Well Location:** Enter Well Location of each compound
- **Plate ID #:** Enter Plate ID Number for each compound

Step 3. Email Excel copy of CSD to: EUservices@discoverx.com

Step 4. Attach the Customs Declaration Form to the outside of the shipping box and include a hardcopy of the Compound Shipping Document inside the package.

Compound Preparation and Shipping instructions:

- **Concentration:** Provide compounds in 100% DMSO, PBS or water at 1,000X the desired final top screening concentration for 10-pt dose response curves. For profiling in single point, provide compound at 1,000X the desired final screening concentration.
- **Minimum Amount:**
 - 20 μl for the first target and 2 μl for each additional target
 - or-
 - 1-2mg in labeled vial(s).

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

Dr. Neil W. Charter
DiscoverRx Corp.
42501 Albrae St.
Fremont, CA 94538
Tel: (510) 979-1415 x128
USA

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