

ORDERING INFORMATION

■ QUOTATION REQUEST

Email your request for a quotation to sales@cerep.com and copy our customer support teams:

- ▶ *In vitro* pharmacology requests . customer.support@cerep.fr
- ▶ ADME-Tox & PK requests . customersupportUS@cerep.com

Please detail your information including location, phone and fax #, the assays (and reference numbers) to be quoted, the number of compounds for each assay and the number of concentrations for each compound. **Standard quotations** for catalog assays will be returned to you by email within 48 hours maximum, and in most cases within 24 hours, of receipt at customer support. **Customized quotations** will be returned within 5 business days.

Your quotation number becomes your study number.

Cerep Order form.xls document can be used to facilitate the quotation process, but is not required. A copy can be emailed to you, request it from sales@cerep.com.

NOTE: Please review your quotation carefully to ensure that it includes the assays, the test concentrations and other specifications that concur with your request. We will try our best to interpret your inquiry, however, the quotation is what we will use to determine how and what we perform, and once the assays have been launched in production, it will be due as per the quotation.

■ MASTER SERVICE AGREEMENT

Simultaneous to quotation processing and if it is the first time you work with Cerep, we will prepare a Master Service Agreement for your company and send it in an e-mail. This also serves as your confidentiality agreement. Your legal counsel may have questions or propose changes. These must be reviewed by our in-house counsel at legal@cerep.fr. Once the content has been agreed upon by both parties, fax a signed copy, and send two signed originals to Cerep legal department who will have them signed, and will return one fully executed agreement to you.

- ▶ **Counsel - Legal department**
Cerep - Le Bois l'Evêque - 86600 Celle l'Evescault - FRANCE
Tel +33 (0)5 49 89 30 00 - Fax +33 (0)5 49 43 21 70 - e-mail: legal@cerep.fr

■ REQUESTED COMPOUND INFORMATION

To reduce the registration time and ensure that all the appropriate information is available to start the study in a shortest possible timeframe, please use **Cerep compound submission form**¹ or MS Excel file², and provide the following **compound information**:

- ▶ **Name (compound ID) / Batch # / Molecular weight**³ / **Formula weight**⁴ / **Stock concentration / Stock solvent / Quantity / Unit**⁵ / **Form / Storage conditions / Solubility**, as well as **Plate ID / plate position** for compounds delivered in plates, **Comments**⁶, and **Quotation number**.

NOTE: Impurity and colored compounds might affect the results (compound color information is mentioned in the study report).

- ▶ **General remarks:**
 - **If compound(s) are supplied as a stock solution in plate(s)** (preferred format for any submission of 10 or more compounds), please leave columns 1 and 12 empty in a 96VV plate. The 384WV plate format is also acceptable with columns 1, 2, 23 and 24 empty. For any other plate format, please inquire.
 - **If compound(s) are not soluble in 100% DMSO**, please provide any useful information concerning the solubility of the compound. The following solvents are compatible with most of our assays: DMSO (Cerep standard), H₂O, Methanol, Tris/HCl 10 mM pH 7.4.
 - **Organic solvents such as acetone, chloroform, ether, acetonitrile** (except for CYP assays), **tetrahydrofuran and trifluoroacetic acid are not recommended** as they will significantly affect the results from many *in vitro* assays, even at very low concentrations.

¹ Cerep compound submission form will be emailed to you with your quotation. A copy can be requested from sales@cerep.com, or downloaded from Cerep website: www.cerep.com/CatalogOnline

² Systematically required for studies of 10 compounds or more.

³ Molecular weight (MW) of free acid or base form.

⁴ Formula weight (FW) including salt form and/or hydrate form if applicable.

⁵ mg^g, mL^g

⁶ e.g. useful information such as sensitivity to light, stability or hygroscopicity issues.

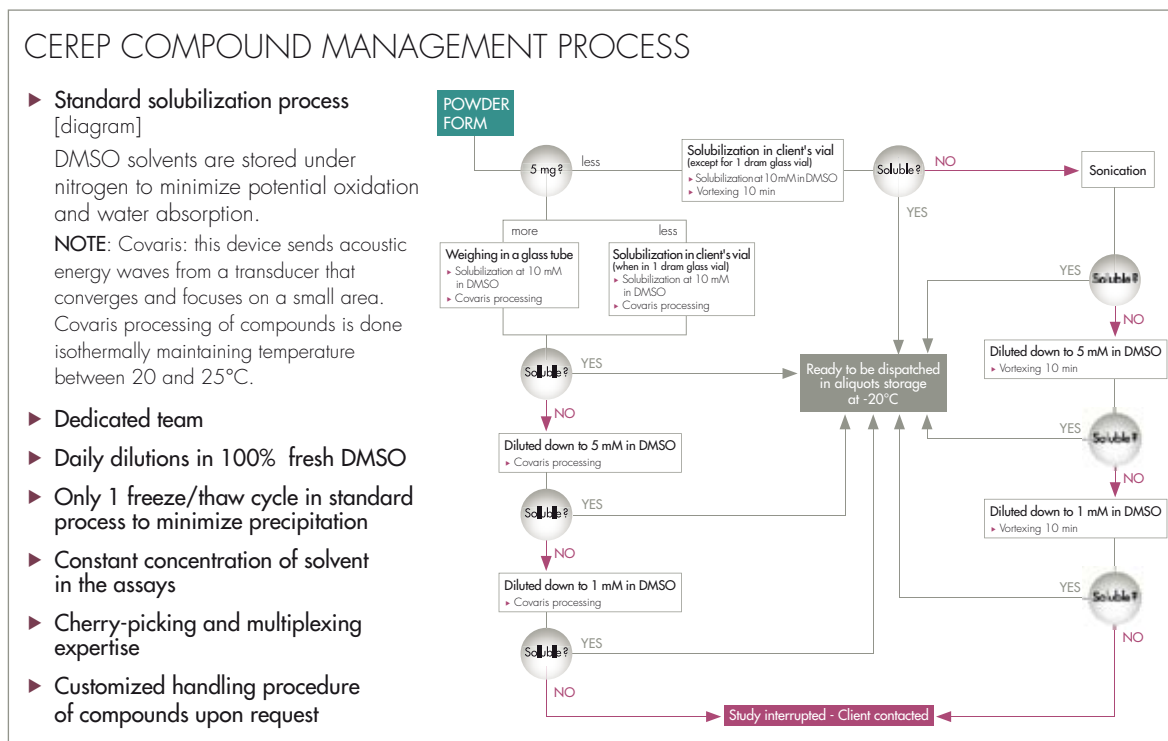
► ORDERING INFORMATION

► Remarks for ADME-Tox assays:

- Any study including mass spectrometry (MS) assays (such as stability, permeability, protein binding) requires both molecular weight (MW) and formula weight (FW) of each tested compound, even if they are the same. The FW is required to prepare the stock solution of each compound at the correct concentration. The MW gives indication on which expected molecular ion to look for in the LC-MS analysis of samples.
- Chemical structure or list of the ionizable groups is required for ionization constant (pKa) assay.
- Solubility information is useful and recommended for pKa and genotoxicity testing (due to high test concentrations).

WARNING: Cerep will apply the standard solubilization process described in the flow chart below when compounds are received at the testing site, unless special instructions are provided with the compounds.

Customized handling procedure of compounds can be accommodated, please inquire for pricing conditions.



■ SAMPLE SIZE [minimum compound amount - including IC₅₀/EC₅₀ follow-up studies]

■ BINDING, ENZYME AND CELLULAR ASSAYS

Assuming a molecular weight ≤ 500 g/mol and a testing concentration of 10 μ M in duplicate (including a possible retest).

	SCREENING		SCREENING + FOLLOW UP with highest concentration at 10 μ M ¹	
	WEIGHT (pre-weighed)	VOLUME (100% DMSO)	WEIGHT (pre-weighed)	VOLUME (100% DMSO)
INDIVIDUAL BINDING, ENZYME AND CELLULAR CATALOG ASSAYS				
1 to 3 assays	1 mg	25 μ L@ 10 mM	1 mg	60 μ L@ 10 mM
4 to 5 assays	1 mg	35 μ L@ 10 mM	1 mg	70 μ L@ 10 mM
6 to 10 assays	1 mg	50 μ L@ 10 mM	1 mg	100 μ L@ 10 mM
11 to 15 assays	1 mg	65 μ L@ 10 mM	1 mg	110 μ L@ 10 mM
16 to 20 assays	1 mg	75 μ L@ 10 mM	1 mg	125 μ L@ 10 mM
21 to 40 assays	1 mg	100 μ L@ 10 mM	1.5 mg	250 μ L@ 10 mM
41 to 50 assays	1 mg	150 μ L@ 10 mM	2 mg	275 μ L@ 10 mM
51 to 70 assays	1.5 mg	225 μ L@ 10 mM	2.5 mg	400 μ L@ 10 mM
71 to 100 assays	2 mg	300 μ L@ 10 mM	3 mg	550 μ L@ 10 mM
101 to 135 assays	2 mg	350 μ L@ 10 mM	4 mg	650 μ L@ 10 mM
136 to 150 assays	2.5 mg	400 μ L@ 10 mM	4 mg	750 μ L@ 10 mM
151 to 200 assays	3 mg	500 μ L@ 10 mM	5 mg	1000 μ L@ 10 mM
201 to 250 assays	3.5 mg	600 μ L@ 10 mM	inquire	inquire

¹ Assuming ~10% of test in IC₅₀. Usually, for 1 IC₅₀: 30 μ L@ 10 mM and + 25 μ L@ 10 mM by additional IC₅₀.

■ TISSUE ASSAYS

For screening at 3 concentrations in duplicate (2 tissues) or follow-up: sufficient amount to prepare 150 µL of a 1000-fold concentrated solution relative to the highest test concentration for each assay.

Maximum tolerable final DMSO concentration: 0.1%.

■ ADME-Tox ASSAYS

Assuming a molecular weight ≤500 g/mol and assays tested at default concentrations (including retest).

	WEIGHT (pre-weighed)	VOLUME (100% DMSO)
INDIVIDUAL ADME-Tox CATALOG ASSAYS		
1 to 20 assays	1 - 2 mg	200 µL@ 10 mM (50 µL min)
1 to 10 IC ₅₀ follow-up assays	1.5 - 2 mg	250 µL@ 10 mM (50 µL min)
Weights needed in addition to the general weight required for 1 to 20 ADME-Tox assays:		
. pKa	3 x 1.5 mg	-
. CYP1A, 2B6 & 3A induction (3 donors)	6 mg (2 mg/CYP)	120 µL@ 100 mM
. Ames (3 strains)	4 mg	800 µL@ 10 mM
. <i>in vitro</i> micronucleus	1 mg	40 µL@ 50 mM
. typical rat PK ¹	25 mg	-
. typical mouse PK (serial sampling) ¹	10 mg	-
. typical mouse PK (parallel sampling) ¹	20 mg	-
. typical rat BBB ²	15 mg	-
. typical mouse BBB ²	10 mg	-
. Plasma sample for quantitative bioanalysis	-	100 µL

¹ 2 routes, 1 dose: 1 mg/kg for IV and 5 mg/kg for PO, n=3

² 1 route, 1 dose: 1 mg/kg for IV, 3 time points, n=3

■ STANDARD PROFILES

Values are based on a molecular weight ≤500 g/mol and a typical testing concentration of 10 µM in duplicate (including a possible retest).

	SCREENING		SCREENING + FOLLOW UP with highest concentration at 10 µM ¹	
	WEIGHT (pre-weighed)	VOLUME (100% DMSO)	WEIGHT (pre-weighed)	VOLUME (100% DMSO)
STANDARD PROFILES				
ExpresSProfile	1 mg	125 µL@ 10 mM	1.1 mg	220 µL@ 10 mM
High-Throughput profile	1 mg	170 µL@ 10 mM	1.4 mg	270 µL@ 10 mM
ExpresSDiversity kinase profile	1 mg	140 µL@ 10 mM	1 mg	190 µL@ 10 mM
Cellular functional GPCR profile	3.25 mg	650 µL@ 10 mM	10 mg	2000 µL@ 10 mM
PDE high-throughput profile	1 mg	200 µL@ 10 mM	1.4 mg	275 µL@ 10 mM
Non-kinase enzyme profile	1 mg	185 µL@ 10 mM	1.5 mg	300 µL@ 10 mM
Diversity profile	1 mg	200 µL@ 10 mM	2 mg	400 µL@ 10 mM
BioPrint® profile (Full profile)	3.4 mg	475 µL@ 10 mM	6.8 mg	1350 µL@ 10 mM
BioPrint® <i>in vitro</i> pharmacology profile	1.4 mg	275 µL@ 10 mM	3.8 mg	750 µL@ 10 mM
Organ safety profile	9 mg	360 µL@ 50 mM	9 mg	360 µL@ 50 mM
Comprehensive kinase profile	2.5 mg	500 µL@ 10 mM	4.5 mg	850 µL@ 10 mM
ADME-Tox - Option I [Bioavailability]	1 - 2 mg	125 µL@ 10 mM	inquire	inquire
ADME-Tox - Option II [Lead selection/Prioritization]	1 - 2 mg	125 µL@ 10 mM	inquire	inquire
Genotoxicity profile	5 mg	200 µL@ 50 mM	inquire	inquire
CYP-based drug-drug interaction profile	10 mg	-	inquire	-
P-gp mediated drug-drug interaction profile	20 mg	-	inquire	-

¹ Assuming ~10% of test in IC₅₀. Usually, for 1 IC₅₀: 30 µL@ 10 mM and + 25 µL@ 10 mM by additional IC₅₀.

■ SHIPMENT INSTRUCTIONS

Please include **study number** (i.e. quotation number, see above Quotation request) on the shipping documents, and include the **Compound submission form** or MS Excel file (see Requested compound information, page 1) and attach it to the compounds.

► Form

- Dry powder, or
- 10 mM DMSO stock solution.

► Format

- Securely closed vials, or
- Individual 2D barcode tubes, or

► ORDERING INFORMATION

- Sealed 96-well plates (all SBS format standard plates are accepted): columns 1 and 12 empty and test compounds placed in A2-H11 column-wise, or
- Sealed 384-well plates (all SBS format standard plates are accepted): columns 1, 2, 23 and 24 empty and test compounds placed in A3-P22 column-wise.

For any other form and/or format: please inquire.

For frozen samples, please allow for enough dry ice to arrive frozen at Cerep site. In addition, it is very important to show on documents and mark on package that compound(s) must be kept in dry ice during all the shipping period (from its collection to its delivery at the Cerep site).

Compounds can be shipped⁷ to any of the two sites below (Cerep will manage the compound transfer to the other testing site when necessary):

- | | |
|--|--|
| <p>► Cerep
Attn: Compound receipt manager
Le Bois l'Evêque
86600 Celle l'Evescault - FRANCE
Tel. +33 (0)5 49 89 30 00</p> | <p>► Cerep
Attn: Compound receipt manager
15318 NE 95th St
Redmond, WA 98052 - USA
Tel. +1 425 895 8666</p> |
|--|--|

Every shipment of compounds going through customs must include a commercial invoice.

For shipments going through US customs: a **TSCA form** (Toxic Substance Control Act) is required (templates are available upon request).

For customers located in China, compounds can be shipped to:

- **Cerep (西海珀)**
Attn: Compound receipt manager
上海张江高科技区爱迪生路326号 302-1室 (326 Aidisheng Road, B 302-1)
Zhangjiang High-Tech Park
Shanghai 201203 - CHINA
tel. +86 21 5132 0568

To assist Cerep in efficiently processing compounds upon receipt, the following information is helpful to ensure the appropriate follow-up of shipment and streamlined initiation of the testing:

- Advance notice of shipment to Cerep – e-mail: compoundreceipt@cerep.com
- Shipment tracking number

■ STUDY INITIATION

To initiate a study (following compound shipment): Fax your signed quotation or attach it to the e-mail accepting it⁸ with purchase order number, to:

- For *in vitro* pharmacology assays: Customer support France: Fax +33 (0)5 49 43 21 70 – e-mail: customer.support@cerep.fr
- For ADME/PK-Tox assays: Customer support USA: Fax +1 (425) 895 8668 – e-mail: customersupportUS@cerep.com

An acknowledgement of shipment receipt will be emailed to you when the shipment has arrived at the laboratory location.

Your signed quotation is your commitment that you wish us to run exactly the assays and the concentrations that are included in the quotation. Please check it carefully.

■ TURNAROUND TIME

The turnaround time indicated in the quotation starts the day following receipt of compounds at the Cerep testing site for compounds received before 12:00 pm, unless otherwise specified with the client, and considering that the test compounds are received along with all the necessary information to initiate the study (quotation acceptance and requested compound information). If received after 12:00 pm, the study is initiated on the second day following receipt.

→ For majority of assays, **ExpresScreen option (5 business day turnaround time)** is available: please inquire.



■ RESULTS REPORTING

A **Data Online account** allowing visualization of the results as they are produced will be set-up when we receive your signed quotation to initiate the study. Your user name and password will be faxed to you. Please modify them to ensure confidentiality.

Log on to <https://www.cerep.com/secure> to view your results that are posted as soon as validated. Daily updates are highlighted with "NEW" information since your last log-in and are listed first. You will have the option to convert the data to MS Excel format and download it. Use the user name and password Demo to try out the site.

⁷ Preferred carriers: FedEx for room temperature and World Courier for dry ice shipment.

⁸ Your study will not begin until a signed quotation with purchase order # is received.

FRANCE
Le Bois l'Evêque
86600 CELLE L'EVESCAULT
tel. +33 (0)5 49 89 30 00

USA
15318 N.E. 95th Street
REDMOND, WA 98052
tel. +1 (425) 895 8666

CHINA
326 Aidisheng Road, B 302-1
Zhangjiang High-Tech Park
SHANGHAI 201203
tel. +86 21 5132 0568

■ **QUESTIONS OR CONCERNS?**
Please contact us: sales@cerep.com

sales@cerep.com
www.cerep.com