

Shipment Instructions

Thank you for working with Reaction.

Our goal at Reaction's logistics department is to improve the service process, to facilitate customs clearance, to maximize the safety of your compound, and to reduce delivery timelines.

For this, we have prepared the following shipment instructions that may answer most of your questions. If you require any other information or advice, please contact us via email (RBE-logistics@reactionbiology.com) before initiating the shipment process. **We are happy to help!**

Where do I ship my parcel to?

Reaction Biology Europe GmbH

Logistics Department

RBE-logistics@reactionbiology.com

Phone: +49-761-769996-0

Project number PQxxxx _ yy department

Engesserstr.4

79108 Freiburg

GERMANY

We accept deliveries during business days

Monday – Friday,

8 am – 3 pm CET

Please fill in **yy department code:**

CP = Cellular Pharmacology

BP = Biochemical Pharmacology

IVP = In-Vivo Pharmacology

Which advice does Reaction provide for preparing my test articles?

- Use **vials** (with screw cap), **micronics** or **plates**, well closed and/or safely sealed.
- Deliver vials, micronics or plates in a **closed box**, please avoid bags.
- **Solid material** provided in vials needs to be **pre-weighed**.
- If using **pens** for labelling (better use **printed labels**), use **permanent** and **solvent proofed** versions.

What shall I include in the package?

- Your **items / compounds**
- A **print-out of the Assay Condition Form (ACF) filled in** (not necessary for in vivo studies)
- **Customs Invoice**

Generally, we prefer shipping with DHL Express

| For dry ice shipment | Amount of dry ice |
|----------------------|-------------------|
| NON-EU countries | 20 kg |
| EU countries | 5 kg |

What information shall I use to facilitate Customs' clearance?

- **TARIC Code** (Please state the correct code in the customs documents):

| Taric code | Items | |
|-------------|--|---|
| 294200 0000 | Reagents for diagnostic or laboratory purposes ("compounds") | If no Taric code seems to fit, please contact us via RBE-logistics@reactionbiology.com |
| 300259 0000 | Cell lines | |
| 05119985 00 | Tissue samples, tumors, sera | |
| 35079090 90 | Proteins, kinases, substrates | |
| 293499 0000 | DNA, plasmids | |

- Indicate:
"Only for laboratory testings, not for use in humans, only smallest amounts, no commercial value"
- Do not exceed a commercial value of 10 €//\$ per item and a total value of 150 €//\$.
- State number of items and estimated amounts (mg/µl), each.
- Use shipment Incoterm **"DDP"** (Delivered Duty Paid).

Who should I notify at Reaction?



Upon submission, please provide us with the **name of the selected carrier, and tracking number** by email to RBE-logistics@reactionbiology.com, indicating the project number PQxxxx in the subject line.