

# 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](mailto: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](mailto: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 <a href="mailto:RBE-logistics@reactionbiology.com">RBE-logistics@reactionbiology.com</a>
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](mailto:RBE-logistics@reactionbiology.com), indicating the project number PQxxxx in the subject line.