



## Customer Registration Form

<b>Organisation Legal Entity Name</b>

<b>Invoice Address</b>	<b>Delivery Address</b>
VAT/EORI Number:	
Invoicing Currency: <input type="checkbox"/> GBP <input type="checkbox"/> USD <input type="checkbox"/> EUR	
Please tick the agreed currency for payment. Payment Method: Bank Transfer	

<b>Finance Contact Person</b>	<b>Delivery Contact Person</b>
Name:	Name:
Tel:	Tel:
Email for Invoices and Statements:	Email for Order Confirmations:

I confirm that the information provided is correct, that I have read and agree to Revvity’s Terms and Conditions of Sale and that I am an authorised signatory of this organisation:

Print Name: \_\_\_\_\_

Date: \_\_\_\_\_

Position within organisation: \_\_\_\_\_

Signature: \_\_\_\_\_

Please email signed form to [Customerexperience.oxford@revvity.com](mailto:Customerexperience.oxford@revvity.com)