

# DOJINDO MOLECULAR TECHNOLOGIES, INC.

30 West Gude Drive, Suite 260, Rockville, MD 20850  
 Phone: 301-987-2667 Fax: 301-987-2687

Date	
Name	
e-mail	
Phone	

Billing Address	
How to receive an invoice	
e-mail / mail	
Attention	
Company/Institute	
Address	
Bldg. & Rm	
City	
State	
Zip Code	
e-mail	
Phone	

Shipping Address (If different)	
Attention	
Company/Institute	
Address	
Bldg. & Rm	
City	
State	
Zip Code	
Phone	
If you would like to use your account for the shipment, please provide your account number below.(UPS or Fedex only)	

Signature

Order Products	
Product Code	
Product Name	
Size / Quantity	/
Unit price	
Subtotal	
Product Code	
Product Name	
Size / Quantity	/
Unit price	
Subtotal	
Product Code	
Product Name	
Size / Quantity	/
Unit price	
Subtotal	

Payment Method	
<b>Purchase order</b>	
Purchase order #	
<b>Credit Card Order</b>	
Visa / Mastercard	
Name on Card	
Card Number	
Expiration Date	
Security Code	
Billing Zip Code	

Tax Exemption
Taxable or Tax Exempt
Taxexempt Certification#:

To place your order, please complete and info@dojindo.com or fax this form to 1-301-987-2687. Please indicate your PO# on the form. If you have any questions please feel free to call 1-877-987-2667 or email us at info@dojindo.com.