

Ordering Instructions

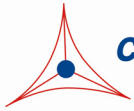
Shipments within the USA:

1. Please complete all information on the following Order Form, including payment information (Credit Card or Purchase Order number).
2. Return the completed form to Cell Biolabs:
 - By Fax: (858) 271 6514
 - By E-mail: orders@cellbiolabs.com
 - By Postal mail: Cell Biolabs Inc., Attn: Customer Service, 7758 Arjons Drive, San Diego, CA 92126
3. Our Customer Service department will apply any discounts to your order and add shipping and handling (S&H) charges. Sales tax will be added to shipments within California. Upon receipt of your order, we will contact you as soon as possible either by phone or by email to confirm your order and advise you on product availability and expected delivery time.

Shipments outside the USA:

1. Orders may be placed with our international distributors (see www.cellbiolabs.com for country specific contact information). International orders placed directly with Cell Biolabs must be pre-paid by credit card or wire transfer. Contact us at orders@cellbiolabs.com for wire transfer information.
2. Complete the following form, along with your FedEx account number for shipping, and return the form to Cell Biolabs:
 - By Fax: +1 (858) 271 6514
 - By E-mail: orders@cellbiolabs.com
 - By Postal mail: Cell Biolabs Inc., Attn: Customer Service, 7758 Arjons Drive, San Diego, CA 92126 USA
 - Our Customer Service department will apply any discounts to your order and add handling charges. Upon receipt of your order, we will contact you as soon as possible either by phone or by email to confirm your order and advise you on product availability and expected delivery time.

Thank you for ordering from Cell Biolabs. We appreciate your business!



Order Form

Contact Information	Name:		Phone:	
	Company or Institution Name:		Fax:	
	Account Number (if known):		E-mail:	

Address Information		Billing Address	Shipping Address (if different)
	Company or Institution Name:		
	Attention Line:		
	Bldg/Room #:		
	Street Address or Box Number:		
	City/State:		
	Zip/Postal Code:		
	Country:		

Payment Information	Payment Method:	<input type="checkbox"/> Credit Card	<input type="checkbox"/> Purchase Order	
	Name as Shown on Card:		Purchase Order Number:	
	Card Number:			
	Expiration Date:			
	FedEx Account Number (required outside USA only):			

Products Ordered	Cat. No.	Description	Qty	Unit Price	Subtotal Price
	Total				

Authorized Signature: _____ **Order Date:** _____

Print Name: _____ **Promotion or Quote #:** _____

Notes or Special Instructions: