



Ordering Instructions

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 at orders@cellbiolabs.com

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



Order Form

Contact Information	Name:		Phone:	
	Company or Institution Name:			
	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: