



Immunochemistry Technologies, LLC
 9401 James Avenue South, Suite #155
 Bloomington, MN 55431 USA
 952-888-8788; fax 952-888-8988
 1-800-829-3194
 www.immunochemistry.com

ORDER

1-800-829-3194; FAX 952-888-8988

Billina Address:	Shipping Address:
Organization:	Organization:
Department:	Department:
Name:	Name:
Title:	Title:
Address:	Address:
City/State/Zip/Country:	City/State/Zip/Country:
Phone:	Phone:
Fax:	Fax:
Payment method:	FedEx/shipping account #:
Tax ID#:	Other:

PURCHASER		P.O. NUMBER		SALES REP	DELIVERY	DATE
PART NUMBER	PRODUCT NAME	SIZE	QUANTITY	UNIT PRICE	TOTAL PRICE	
					SUBTOTAL	
					SALES TAX / DUTY	
					SHIPPING & HANDLING	
					TOTAL	

For credit card orders only, please fill out the information below:

circle one Visa/MasterCard # _____ - _____ - _____ - _____ Exp ____ / ____ mo/yr
Name on the card: _____ Signature: _____
Address to which the bill is sent: _____

If you have any questions concerning this order, call Sally Hed at 952-888-8788,
 or send an email to sally@immunochemistry.com.

THANK YOU FOR YOUR BUSINESS!