



Biotium Order Form

Please provide the information required in this form when placing an order. You may choose to send the completed form to us via fax (510) 265-1352 or E-mail (order@biotium.com).

Contact Name:		*Email:		
Organization:				
*Telephone:		Fax:		
Purchasing Method: <i>Purchase Order</i> <i>VISA</i> <i>MasterCard</i> <i>American Express</i>				
PO No. / CC No.				
CC Exp. Date		CC Holder Name:		
Shipping Address: <i>Commercial street address only.</i> <i>PO Box Number not acceptable.</i>		Billing Address:		
Catalog No.	Description	Unit Size	Unit Price(\$)	QTY
Special Instruction:				

Biotium, Inc.

46117 Landing Pkwy., Fremont, CA 94538
Tel: 1-510-265-1027; Fax: 1-510-265-1352

Quote or Order: order@biotium.com General Inquiry: btinfo@biotium.com
Technical Service: techsupport@biotium.com