

**Molecular Devices 2010 Imaging Training**  
**Advanced MetaXpress® Software Training & In-Depth AcuityXpress™ Software Training**  
**REGISTRATION FORM**

Please select your events for and fax to 408.747.3562

<u>Downingtown, PA</u>			
<input type="checkbox"/>	Advanced MetaXpress® Software April 6 – 8, 2010 S9200-0126	\$1000	
<input type="checkbox"/>	AcuityXpress™ Software July 21-22, 2010 S9200-0127		\$1000
<input type="checkbox"/>	Advanced MetaXpress® Software October 5 – 7, 2010 S9200-0126	\$1000	

<u>Sunnyvale, CA</u>			
<input type="checkbox"/>	Advanced MetaXpress® Software April 27 – 29, 2010 S9200-0126	\$1000	
<input type="checkbox"/>	AcuityXpress™ Software May 12 – 13, 2010 S9200-0127		\$1000
<input type="checkbox"/>	Advanced MetaXpress® Software November 9 – 11, 2010 S9200-0126	\$1000	

Salutation	First name: *	Last name: *
Institution or Company: *		
Title:	Department:	
E-mail: *		Phone:
Address: *		
City: *		State or Province: *
Country: *		Zip/Postal code: *
*indicates a required field		

PAYMENT	
Card Number*	
Name as it appears on Card*	
Expiration Date*	
Signature*	

I would like to be put on a mailing list to receive updates about new products and events from Molecular Devices.

