

xMAP University 2008

Teaching the xPERT

Dates for currently scheduled training courses are located on the calendar. All courses are trained in the Austin, TX training facility unless otherwise noted.

Additional course dates and locations may be scheduled upon request. Methods for course delivery include web-based, in-house at a Luminex facility, on-site at a customer facility and at regional locations.

Find out more about Luminex Training Courses by visiting the Luminex Training site at www.luminexcorp.com/training.

For more information or to register for a training course, please contact Luminex at 1-512-219-8020, ext. 4386 or via email at training@luminexcorp.com.

IS 2.3 1-Day Basic Training (1-day)

xPONENT 1-Day Basic Training (1-day)

Fundamental Assay Techniques Training – Protein (2-days)

Fundamental Assay Techniques Training – Nucleic Acid (2-days)

Targeted Applications Workshop – Protein (1-day)

Targeted Applications Workshop – Nucleic Acid (1-day)

Technical Representative Training (2 days)

Field Service Training (5-days)

Luminex Corporation Internal Courses

Other Scheduled Course

* Training in Oosterhout, The Netherlands
 **xMAP University Regional Events: March 6th – Boston, MA, August 20th – San Francisco, Sept. 17th – Philadelphia

Q1							Q2							Q3							Q4							
JANUARY							APRIL							JULY							OCTOBER							
S	M	T	W	T	F	S	S	M	T	W	T	F	S	S	M	T	W	T	F	S	S	M	T	W	T	F	S	
		1	2	3	4	5			1	2	3	4	5			1	2	3	4	5			1	2	3	4		
6	7	8	9	10	11	12	6	7	8	9	10	11	12	6	7	8	9	10	11	12	5	6	7	8	9	10	11	
13	14	15	16	17	18	19	13	14	15	16	17	18	19	13	14	15	16	17	18	19	12	13	14	15	16	17	18	
20	21	22	23	24	25	26	20	21	22	23	24	25	26	20	21	22	23	24	25	26	19	20	21	22	23	24	25	
27	28	29	30	31			27	28	29	30				27	28	29	30	31			26	27	28	29	30	31		
FEBRUARY							MAY							AUGUST							NOVEMBER							
S	M	T	W	T	F	S	S	M	T	W	T	F	S	S	M	T	W	T	F	S	S	M	T	W	T	F	S	
					1	2				1	2	3						1	2							1		
3	4	5	6	7	8	9	4	5	6	7	8	9	10	3	4	5	6	7	8	9	2	3	4	5	6	7	8	
10	11	12	13	14	15	16	11	12	13	14	15	16	17	10	11	12	13	14	15	16	9	10	11	12	13	14	15	
17	18	19	20	21	22	23	18	*19	*20	*21	*22	23	24	17	18	19	**20	21	22	23	16	17	18	19	20	21	22	
24	25	26	27	28	29		25	26	27	28	29	30	31	24	25	26	27	28	29	30	23	24	25	26	27	28	29	
MARCH							JUNE							SEPTEMBER							DECEMBER							
S	M	T	W	T	F	S	S	M	T	W	T	F	S	S	M	T	W	T	F	S	S	M	T	W	T	F	S	
						1																						
2	3	4	5	**6	7	8	8	9	10	11	12	13	14	7	*8	*9	*10	*11	12	13	7	8	9	10	11	12	13	
9	10	11	12	**	13	14	15	15	16	17	18	19	20	21	14	*15	*16	**17	18	19	20	14	15	16	17	18	19	20
16	17	18	19	20	21	22	22	23	24	25	26	27	28	21	22	23	24	25	26	27	21	22	23	24	25	26	27	
23	24	25	26	27	28	29	29	30						28	29	30					28	29	30	31				
30	31																											