



HAMBURG TRAINING CALENDAR 2021

LIFE SCIENCE AND DIAGNOSTICS

Applied Genomics	P/N	Site	Days	Jan	Feb	Mar	Apr	May	Jun	Jul	Aug	Sep	Oct	Nov	Dec
Janus® G3 User Training	EDU-AG1002HAM	Hamburg	3			9 - 11			15 - 17				5 - 7		
Maestro® Advanced Training	EDU-AG1004HAM	Hamburg	3			9 - 11			15 - 17				5 - 7		
LabChip® User Training	EDU-AG1001HAM	Hamburg	2						Organized upon request						
NGS Library Prep Training	EDU-AG1003HAM	Hamburg	1.5						Organized upon request						
Chemagen User Training	EDU-AG1005HAM	Hamburg	2						Organized upon request						
Application Consultation Day	EDU-AG1006HAM	Hamburg	1 - 5						Organized upon request						
plate::works™ User Training	N/A	Hamburg	1.5						Organized upon request						
High Content Screening	P/N	Site	Days	Jan	Feb	Mar	Apr	May	Jun	Jul	Aug	Sep	Oct	Nov	Dec
HCS University Stage 1	N0239824	Hamburg	3			23 - 25									
HCS University Stage 2	N0239824	Hamburg	3										19 - 21		
Radiometric Detection	P/N	Site	Days	Jan	Feb	Mar	Apr	May	Jun	Jul	Aug	Sep	Oct	Nov	Dec
RRD Basic Training - German	N0236520	Hamburg	3			1 - 3									
RRD Advanced Training - German	N0236522	Hamburg	2			4 - 5									
RRD Basic Training - English	N0236521	Hamburg	3											15 - 17	
RRD Advanced Training - English	N0236523	Hamburg	2											18 - 19	

Training Location

PerkinElmer Center of Excellence
 Schnackenburgallee 114
 22525 Hamburg
 Phone: +49 (0)40 307 090 476
 E-mail: Training.Hamburg@perkinelmer.com
<https://perkinelmer-appliedgenomics.com>

PerkinElmer, Inc.

Waltham, MA 02451 USA
 Phone: (800) 762-4000 or (+1) 203-925-4602
www.perkinelmer.com