



# TURKU TRAINING CALENDAR 2021

REPRODUCTIVE HEALTH	SITE	JAN	FEB	MAR	APR	MAY	JUN	JUL	AUG	SEP	OCT	NOV	DEC
<b>MATERNAL &amp; FETAL HEALTH COURSES</b>													
Vanadis® NIPT System	Turku (Finland)			15-19		17-21	31.5.-4.6.			13-17		8-12	
LifeCycle™ 7.0 Software	Turku (Finland)			15-19		17-21	31.5.-4.6.			13-17		8-12	
Prenatal Screening with AutoDELFIA®/DELFIAXpress / VICTOR™ 2D System	Turku (Finland)			15-19		17-21	31.5.-4.6.			13-17		8-12	
<b>NEONATAL SCREENING COURSES</b>													
GSP® System	Turku (Finland)			15-19		17-21	31.5.-4.6.			13-17		8-12	
Neonatal Screening with AutoDELFIA®/VICTOR™ 2D System	Turku (Finland)			15-19		17-21	31.5.-4.6.			13-17		8-12	
EnLite™ and Neonatal TREC Kit	Turku (Finland)			15-19		17-21	31.5.-4.6.			13-17		8-12	
Migele™ System & RESOLVE™ hemoglobin kit	Turku (Finland)			15-19		17-21	31.5.-4.6.			13-17		8-12	
QSight® MD System & NeoLSD MSMS Kit	Turku (Finland)			15-19		17-21	31.5.-4.6.			13-17		8-12	
QSight® MD System & NeoBase 2 MSMS Kit	Turku (Finland)			15-19		17-21	31.5.-4.6.			13-17		8-12	
Specimen Gate® Essentials	Turku (Finland)			15-19		17-21	31.5.-4.6.			13-17		8-12	
Eonis™ System & SMA, SCID & XLA Assays	Turku (Finland)			15-19		17-21	31.5.-4.6.			13-17		8-12	

#### Training Location

PerkinElmer Center of Excellence  
 Mustionkatu 6  
 20750 Turku Finland  
 +358 2 2678 111  
 training.turku@perkinelmer.com

#### PerkinElmer, Inc.

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