

# MultiPROBE II Normalization Workstation

PerkinElmer Life Sciences has developed an integrated MultiPROBE II Normalization Workstation for quantification and normalization of nucleic acid samples. This workstation is built around the PerkinElmer's MultiPROBE® II System and Molecular Devices' SPECTRAmax® Microplate Readers.



## System Components:

### PerkinElmer

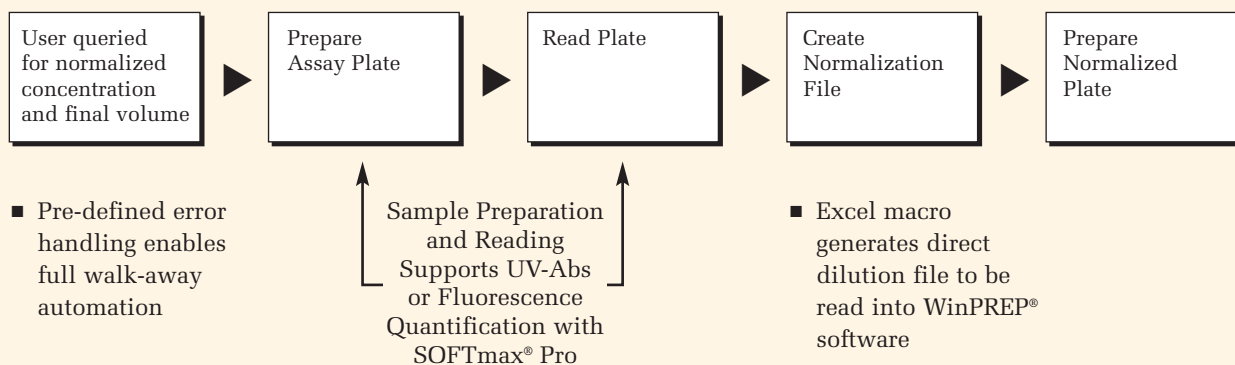
- MultiPROBE II (4-Tip or 8-Tip)
- Gripper™ Integration Platform
- Nucleic Acid Normalization Software Kit

### Molecular Devices

#### Select from

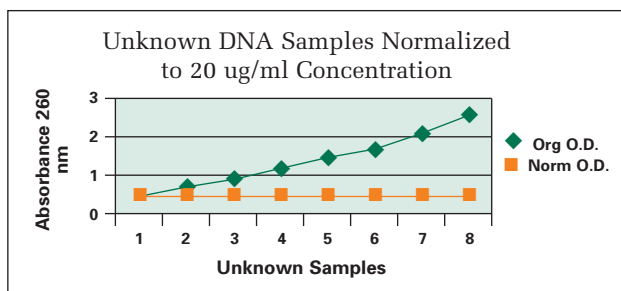
- SPECTRAmax 190 (Absorbance)
- SPECTRAmax PLUS 384 (Absorbance 96/384)
- SPECTRAmax GEMINI XS (Fluorescence)

## The Process:

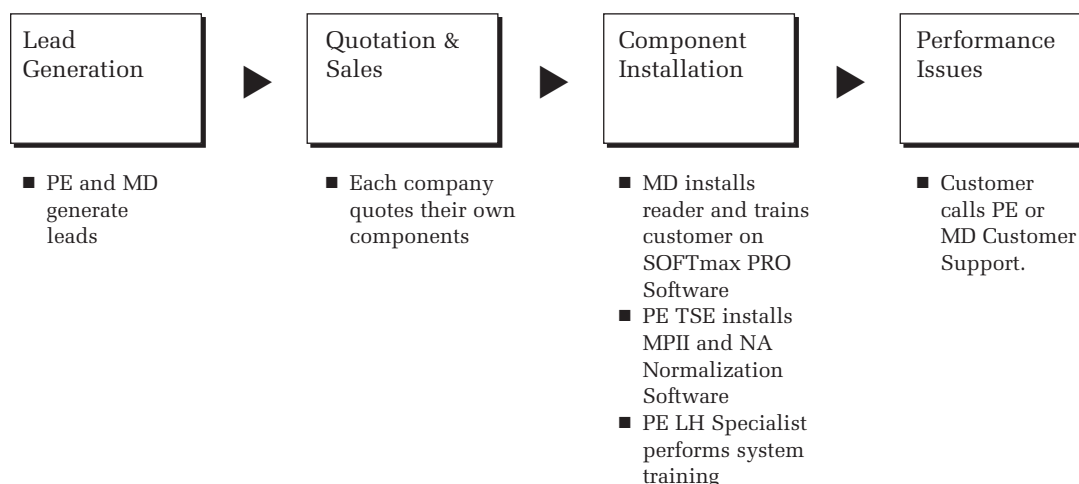


## System Performance:

- Run-time generation of normalization worklist enables full walk-away automation of quantification and normalization for one or many plates.
- Excellent accuracy and precision even at 1 uL volumes.
- Flexible platform supports the choice of:
  - Fixed or disposable tips to control carryover
  - 96- or 384- well plate formats
  - Quantification via UV absorbance or fluorescence
- A 96-well plate can be setup for quantitation in less than 7 minutes while normalization can be achieved in as little as 8 minutes.



## Workstation Support:



## Ordering Information:

Liquid Handling System		Reader
Domestic	International	
<b>MPII4W0 MP II Workstation (4-Tip) includes:</b> <ul style="list-style-type: none"> <li>■ Computer</li> <li>■ MP II EX</li> <li>■ Left &amp; Right Expansion modules</li> <li>■ VersaTip Plus Kit</li> <li>■ 500 uL syringes</li> <li>■ Gripper Integ. Platform</li> <li>■ Double reagent trough holder</li> <li>■ Trough holder, 60 ml 4 pos.</li> <li>■ 60 ml troughs</li> <li>■ 20 uL disposable tips</li> <li>■ 200 uL disposable tips</li> </ul>	<b>MPII8W1 MP II Workstation (4-Tip) includes:</b> <ul style="list-style-type: none"> <li>■ Computer</li> <li>■ MP II EX</li> <li>■ Left &amp; Right Expansion modules</li> <li>■ VersaTip Plus Kit</li> <li>■ 500 uL syringes</li> <li>■ Gripper Integ. Platform</li> <li>■ Double reagent trough holder</li> <li>■ Trough holder, 60 ml 4 pos.</li> <li>■ 60 ml troughs</li> <li>■ 20 uL disposable tips</li> <li>■ 200 uL disposable tips</li> </ul>	<b>0200-3850 - SpectraMax Plus 384</b> with SOFTmax Pro  <b>0200-5025 - SpectraMax 190</b> with SOFTmax Pro  <b>0200-3940 - Gemini XS</b> with SOFTmax Pro  <b>Optional:</b> 0200-2405 - SpectraTest Validation plate for absorbance plate readers  0200-5008 - SOFTmax Pro Software v4.0 Validation pkg
or	**Must Specify Voltage** 7001516 WinPREP Software Kit 8950000 Disk Media Kit	
<b>MPII8W0 MP II Workstation (8-Tip)</b> <ul style="list-style-type: none"> <li>■ Same parts as 4-Tip</li> </ul>	or	
and	MPII8W1 MP II Workstation (8-Tip)	
<b>7001957 Kit, NA Norm Software includes:</b> <ul style="list-style-type: none"> <li>■ MD Reader Integ. Kit</li> <li>■ NA Norm Software</li> </ul>	<ul style="list-style-type: none"> <li>■ Same parts as 4-Tip</li> </ul>	
	**Must Specify Voltage**	
	7001717 WinPREP Software Kit (HT) 8950000 Disk Media Kit	
	and	
	<b>7001957 Kit, NA Norm Software includes:</b> <ul style="list-style-type: none"> <li>■ MD Reader Integ. Kit</li> <li>■ NA Norm Software</li> </ul>	