

TrueMark™ Prostate Protein Assay

Catalog Number A46360

Pub. No. MAN0019000 Rev. A.0

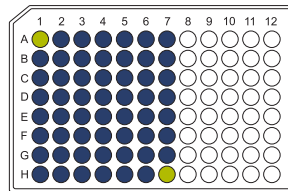
Note: For safety and biohazard guidelines, see the “Safety” appendix in the *TrueMark™ Prostate Protein Assay User Guide* (Pub. No. MAN0017766). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

This Quick Reference is intended as a benchtop reference for experienced users of the TrueMark™ Prostate Protein Assay. For detailed instructions, supplemental procedures, and troubleshooting, see the *TrueMark™ Prostate Protein Assay User Guide* (Pub. No. MAN0017766).

Assay procedure: Day 1

Prepare the Sample Stocks Plate

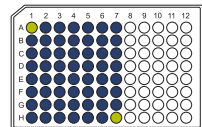
EDTA plasma samples
 Manually transfer aliquots of plasma test samples (●) plus the TrueMark™ Prostate Positive Control (●) to a Fast 96-well plate



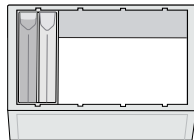
Sample Stocks Plate

Set up, then run the Plate 1 binding reactions on the Biomek™ FX^P instrument

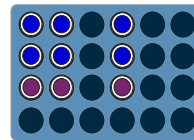
Plate 1 binding reactions (KLK2, PSP94)



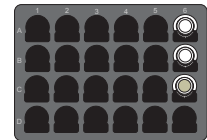
Sample Stocks Plate



Sample Buffer



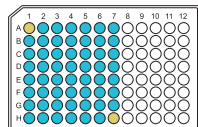
Probes (left) and Calibrators (right)
 (● KLK2, ● PSP94)



○ 2 x Probe Buffer, ● PSP94 Buffer

Ensure that all reagents, plates, samples, and tips are in the correct location on the instrument deck (see Figure 1)

Run Binding Reaction Setup Method—Plate 1 (~50 minutes)



Sample Dilutions Plate

Proceed to set up Plate 2 binding reactions

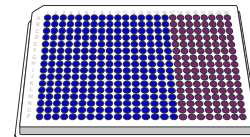


Plate 1 (samples + Probes + Calibrators)

>16 hours, 4°C (binding reaction)

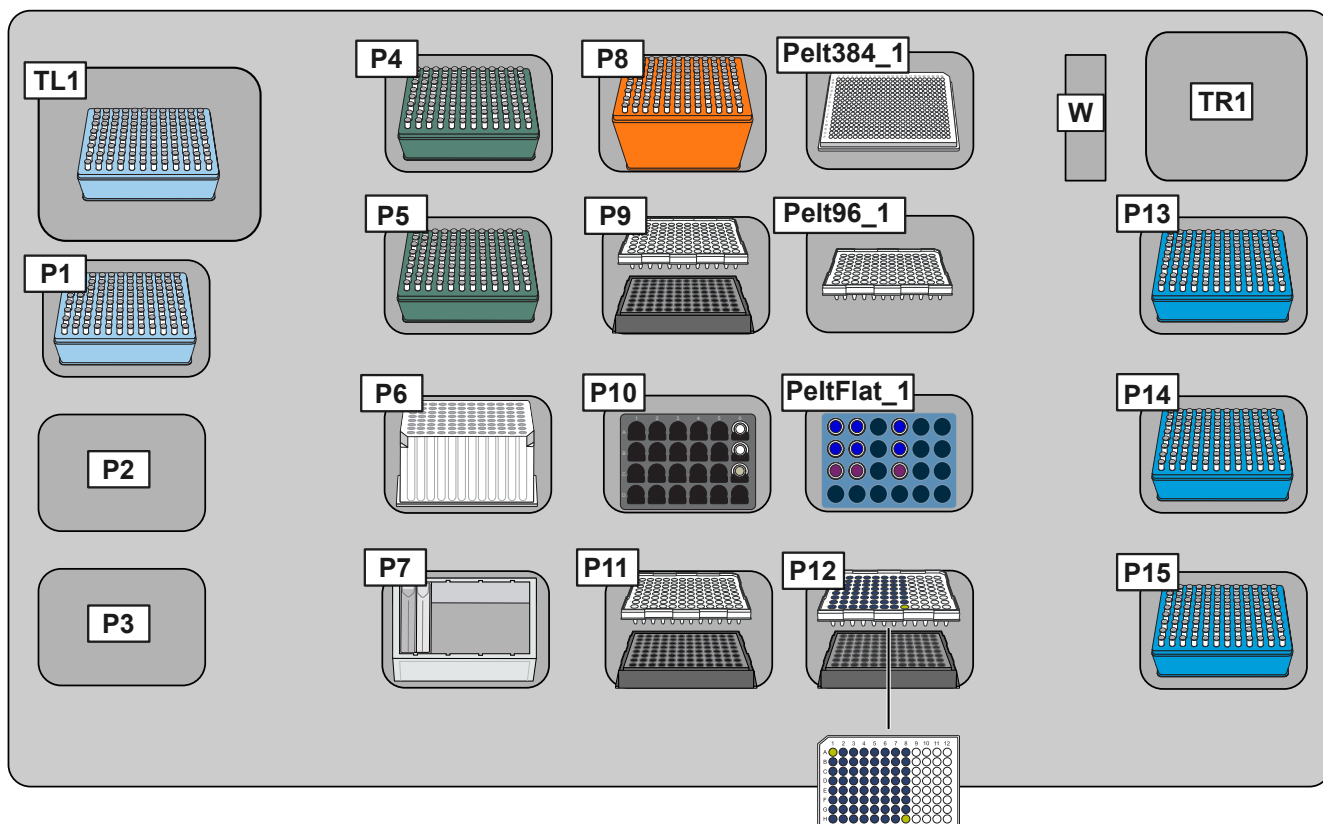
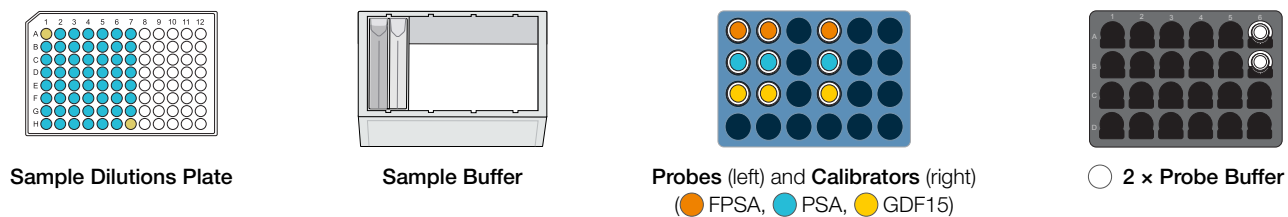


Fig. 1 Biomek™ FXP deck setup for Binding Reaction Setup Method—Plate 1

TL1	AP96 P20 Tips	Pelt384_1	Empty Plate 1 (384-well EnduraPlate™)
P1	AP96 P20 Tips	Pelt96_1	Fast 96-well plate
P4	Span-8 P250 Tips	PeltFlat_1	0.5-mL rack (Probes A and B, Calibrators for KLK2, PSP94)
P5	Span-8 P250 Tips	P12	Sample Stocks Plate on 96-well base
P6	1.0-mL deep-well plate	W	Liquid waste
P7	Sample Buffer (quarter reservoir in frame, liquid on the left)	TR1	Tip waste
P8	Span-8 P1000 Tips	P13	Span-8 P20 Tips
P9	Fast 96-well plate on 96-well base	P14	Span-8 P20 Tips
P10	2-mL rack (2 × Probe Buffer, PSP94 Buffer)	P15	Span-8 P20 Tips
P11	Empty Sample Dilutions Plate (Fast 96-well plate) on 96-well base		

Set up, then run the Plate 2 binding reactions on the Biomek™ FX^P instrument

Plate 2 binding reactions (FPSA, PSA, GDF15)
(start after Binding Reaction Setup Method—Plate 1 is complete)



Ensure that all reagents, plates, samples, and tips are in the correct location on the instrument deck (see Figure 2)

Run Binding Reaction Setup Method—Plate 2 (~45 minutes)

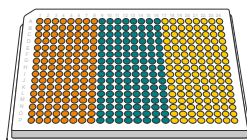


Plate 2 (samples + Probes + Calibrators)

>16 hours, 4°C (binding reaction)

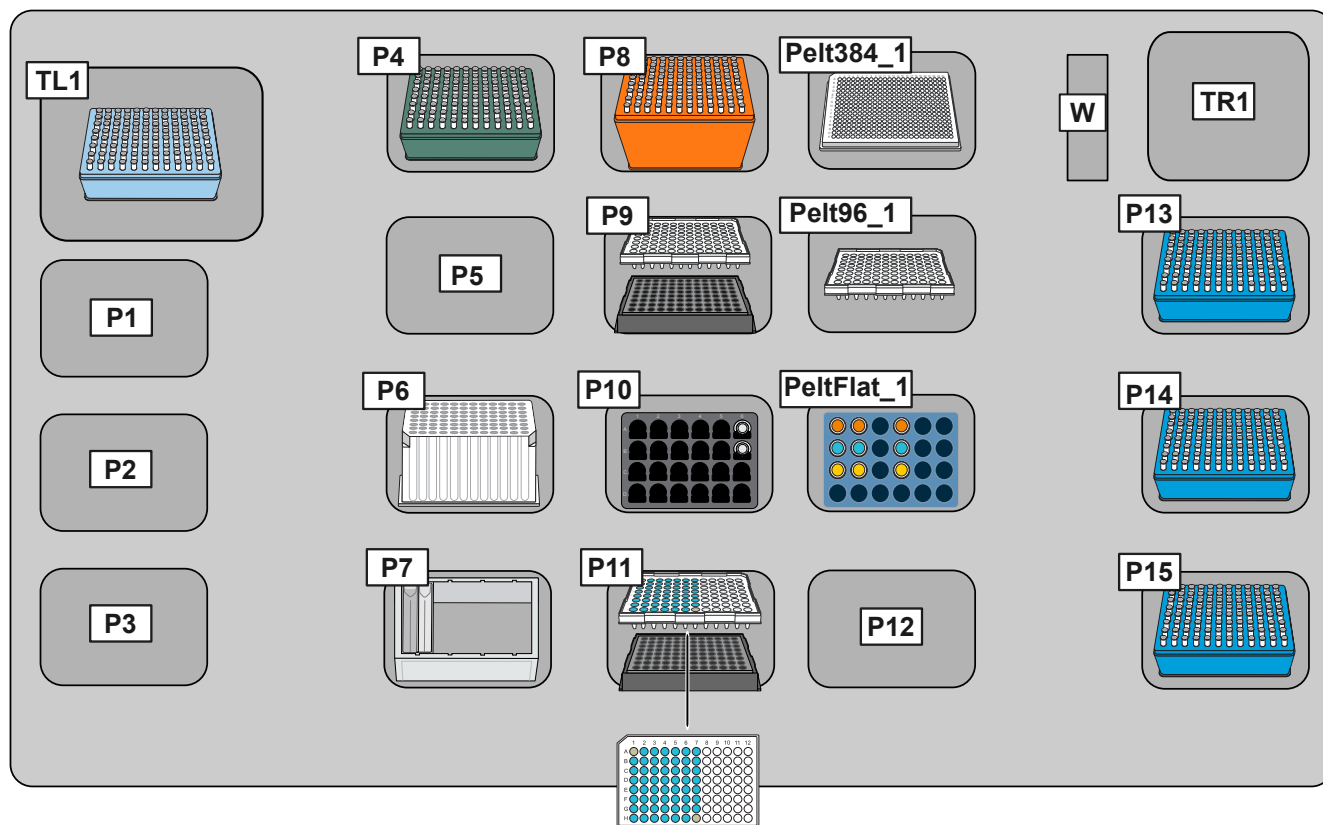


Fig. 2 Biomek™ FX^P deck setup for Binding Reaction Setup Method—Plate 2

TL1	AP96 P20 Tips	Pelt384_1	Plate 2 (<i>empty</i> 384-well EnduraPlate™)
P4	Span-8 P250 Tips	Pelt96_1	Fast 96-well plate
P6	1.0-mL deep-well plate	PeltFlat_1	0.5-mL rack (Probes A and B, Calibrators for FPSA, PSA, GDF15)
P7	Sample Buffer (quarter reservoir in frame, liquid on the left)	W	Liquid waste
P8	Span-8 P1000 Tips	TR1	Tip waste
P9	Fast 96-well plate on 96-well base	P13	Span-8 P20 Tips
P10	2-mL rack (2 x Probe Buffer)	P14	Span-8 P20 Tips
P11	Sample Dilutions Plate on 96-well base	P15	Span-8 P20 Tips

Assay procedure: Day 2

Set up, then run Ligation Reaction Setup Method on the Biomek™ FX^P instrument
Perform this procedure for Plate 1, then repeat for Plate 2.

Plate 1 (KLK2, PSP94)

IMPORTANT! Keep Plate 2 at 4°C during Plate 1 ligation and real-time PCR.

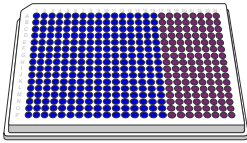
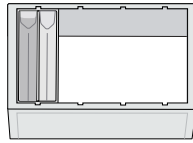
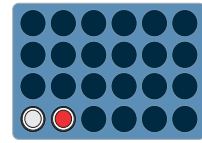


Plate 1
(probe binding to Plate 1 samples complete)



Ligation-qPCR Master Mix



DNA Ligase, Master Mix Additive

Ensure that all reagents, plates, samples, and tips are in the correct location on the instrument deck (see Figure 3)

Run Ligation Reaction Setup Method (~16 minutes)

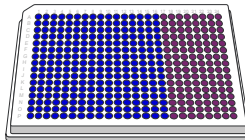


Plate 1 (ligation reaction complete)

Proceed to "Run the ligation reaction on the ProFlex™ thermal cycler" on page 6

Plate 2 (FPSA, PSA, GDF15)

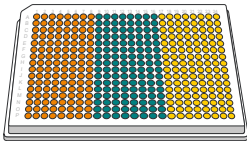
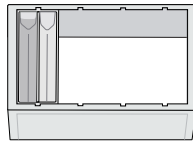
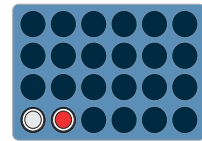


Plate 2
(probe binding to Plate 2 samples complete)



Ligation-qPCR Master Mix



DNA Ligase, Master Mix Additive

Ensure that all reagents, plates, samples, and tips are in the correct location on the instrument deck (see Figure 3)

Run Ligation Reaction Setup Method (~16 minutes)

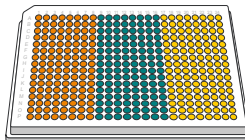


Plate 2 (ligation reaction complete)

Proceed to "Run the ligation reaction on the ProFlex™ thermal cycler" on page 6

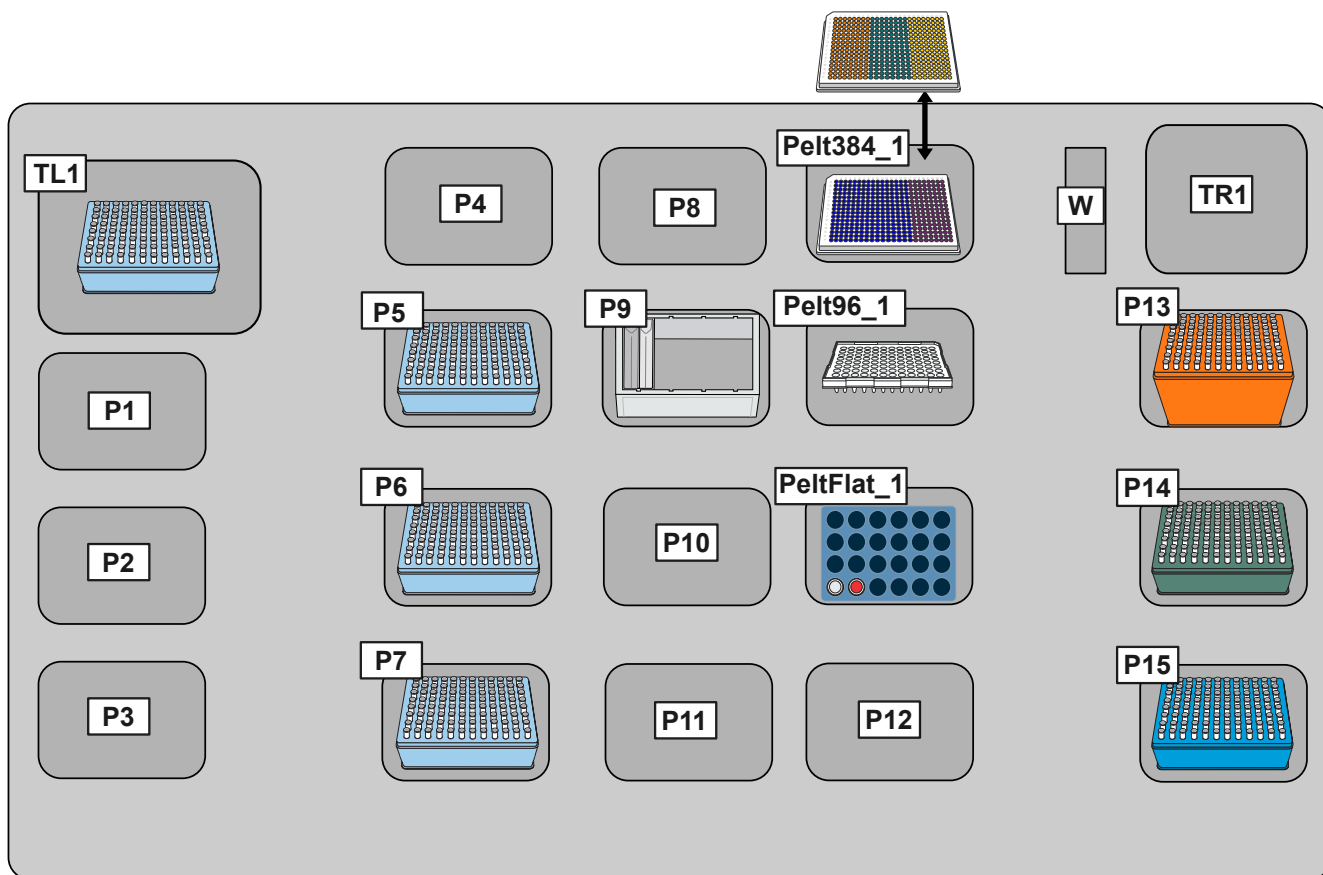


Fig. 3 Biomek™ FXP deck setup for Ligation Reaction Setup Method

TL1	AP96 P20 Tips	Pelt384_1	Plate 1/Plate 2
P5	AP96 P20 Tips	Pelt96_1	Fast 96-well plate
P6	AP96 P20 Tips	PeltFlat_1	0.5-mL rack (D1—DNA Ligase; D2—Master Mix Additive)
P7	AP96 P20 Tips	P13	Span-8 P1000 Tips
P9	Ligation-qPCR Master Mix (quarter reservoir in frame, liquid on the left)	P14	Span-8 P250 Tips
		P15	Span-8 P20 Tips

Run the ligation reaction on the ProFlex™ thermal cycler

Perform this procedure for Plate 1, then repeat for Plate 2.

1. Ensure that the heated cover setting is off.
2. Place the plate in the ProFlex™ thermal cycler, next to an empty 384-well EnduraPlate™.
3. Apply a compression pad over each plate, close the cover, then run the following program.

Temperature	Time
25°C	25 minutes
95°C	5 minutes
4°C	∞

4. Remove the plate from the thermal cycler, then place on ice or at 4°C for 5 minutes to ensure that the entire plate equilibrates to 4°C.
5. Tap the plate firmly on a benchtop 5–10 times to mix, centrifuge at 1,400 × g for 1 minute, then proceed to the next section.

Run the real-time PCR on the QuantStudio™ 12K Flex Instrument

1. In the QuantStudio™ 12K Flex Software, in the **Experiment** pane, click **Create**.
2. In the **Experiment Properties** screen, enter the experiment name, then select the experiment properties.

Property	Setting
Block	384-well
Experiment type	Standard Curve
Reagents	TaqMan™ Reagents
Instrument run	Fast

3. Click **Import** ▶ **Import Plate Setup**, navigate to, then select the plate setup file.
 4. Click **Start Import**.
 5. In the **Assign** screen, ensure that the assays and samples have been assigned to the appropriate wells.
 6. In the **Run Method** screen, enter the thermal cycling parameters.
 - Reaction volume per well—20 µL
- | Stage | Temperature | Time |
|-----------------|-------------|------------|
| Hold | 95°C | 20 seconds |
| PCR (40 cycles) | 95°C | 1 second |
| | 60°C | 20 seconds |
7. Save the experiment (EDS file), then load the plate into the instrument.
 8. Start the run.
 9. After the run is complete, click **Analysis**.
 10. In the **Analysis** screen, click **Analysis Settings**, ensure the following default C_t settings are selected, then click **Analyze** ▶ **Save**.
 - Baseline—Auto
 - Threshold—0.2
 11. In **Experiment Menu**, click **Export**, then select the format, data, and location for export.
 - Data format—QuantStudio 12K Flex
 - File type—*.txt
 - Data—select the checkboxes in the **Sample Setup** and **Results** tabs.
 12. Click **Start Export**.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Life Technologies Corporation | 6055 Sunol Blvd | Pleasanton, CA 94566

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. MAN0019000

Revision	Date	Description
A.0	27 February 2020	New document for new product launch.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.