

ASSAY NAME: STI2_BR
(Sexually transmitted infections panel 2 for BioRad)

Quantity: 100 x 20µL PCR reactions

5-plex assay: *Neisseria gonorrhoeae*, *Mycoplasma genitalium*, *Treponema pallidum*, *Ureaplasma urealyticum*, and human RPP30 DNA

SKU: PNP-STI2-D-BR-100 (Bio-Rad)

(RUO). Research Use Only. Not for use in Diagnostic Procedures.

SCOPE OF THIS PRODUCT INFORMATION SHEET (PIS):

The oligonucleotide recipes are optimized for each instrument (BioRad, QuantStudio, MIC). The verification data presented in this PIS were performed using PNP-STI2-D-BR-100 on a BioRad CFX96. The performance of the other SKUs on their corresponding instrument should be similar. Contact PCRassays.com if you need to use a different qPCR instrument.

CONTENTS

The STI2 kit contains a mixture of primers/probes targeting an intergenic region in the *Neisseria gonorrhoeae* genome, the genes of the RluA family pseudouridine synthase and M23 family metalloproteinase (Protein IDs: WGV92268.1, WGV91925.1) in the *Treponema pallidum* genome, the gene of the MgpB adhesin protein (Protein ID: ALN98184.1) in the *Mycoplasma genitalium* genome, and the gene of the iron ABC transporter permease protein (Protein ID: QDI63787.1) in the *Ureaplasma urealyticum* genome. The primers and probes in the STI2 assay are provided in Tube 1 as a 5X concentrated working solution that detects 4 pathogens and a human extraction control.

Table of Dyes used in this assay:

Pathogen/Target	Dyes	Quencher	Refs.
<i>N. gonorrhoeae</i>	FAM	BHQ-1	1,2
RPP30-DNA control	HEX	BHQ-1	7
<i>T. pallidum</i>	TEX615	BHQ-2	3,4
<i>M. genitalium</i>	Cy5	BHQ-2	5
<i>U. urealyticum</i>	Cy5.5	BHQ-2	6

The probes are designed as TaqMan[®] cleavage mechanism and thus the reaction requires a DNA polymerase with 5'-exonuclease activity.

Assay contents:

Tube 1: Primer/Probe mix (5X) for *Neisseria gonorrhoeae*, *Mycoplasma genitalium*, *Treponema pallidum*, *Ureaplasma urealyticum*, and hRPP30DNA.

Tube 2: (Do NOT add to specimen unknowns) Positive control: 5000

copies/µl positive controls of synthetic 500 bp DNA fragments of *Neisseria gonorrhoeae*, *Mycoplasma genitalium*, *Treponema pallidum*, *Ureaplasma urealyticum*, and hRPP30DNA.

Tube 3: InhibiTaQ Standard qPCR enzyme Mastermix (enough for 100 rxns. with 20 µL total volume). This is a custom formulation from Fortis Life Sciences to the specifications of PCRassays.com.



ASSAY HANDLING

The STI2 assay is shipped at ambient temperature, but should be stored at -20 °C. The tubes should be kept on ice once thawed. Do not subject the enzyme to multiple freeze-thaw cycles.

Any contamination should be avoided by using appropriate personal protective equipment (PPE), powder free gloves, aerosol barrier pipette tips, and a clean hood.

EXPERIMENTAL

Perform nucleic acid extraction/purification (recommended).

Set up your reaction (20 µL) as follows on ice:

Component	Volume (µL)
InhibiTaQ qPCR enzyme mastermix (2X)	10
Primer/Probe mix (5X)	4
Sample	2
Water	4

Notes: To improve assay sensitivity, up to 6 µL of sample can be added (water volume adjusted accordingly) for a total reaction volume of 20 µL. For positive control rxns., add 2 µL of the solution from Tube 2. Molecular biology grade water (NOT included) should be used to prepare the PCR reactions.

A PCR protocol was used for verification on a BioRad CFX96 system, with the following program:

Step	Thermocycling Protocol:
1	Incubate @ 95 °C for 2 minutes
2	Incubate @ 95 °C for 3 seconds
3	Incubate @ 55 °C for 15 seconds
4	Plate Read
5	Go to Step 2, repeat 44× more

RESULT INTERPRETATION

After running the qPCR reaction, perform a regression analysis on the data to determine the quantification cycle, C_q. (C_q is preferred over Ct). Each fluorescence channel with a C_q < 38 cycles and final RFU > 200 (BioRad) is considered “positive” or “+” in the Table below.

<i>N. gonorrhoeae</i> FAM TM	<i>T. pallidum</i> TEX615 TM	<i>M. genitalium</i> Cy5 TM	<i>U. urealyticum</i> Cy5.5 TM	hRPP30 HEX TM	Recommended Interpretation
—	—	—	—	—	The PCR reaction failed. Please repeat the experiment
—	—	—	—	+	The sample does not contain DNA of interest. The sample contains human RPP30 DNA.
+	—	—	—	—	The sample contains <i>N. gonorrhoeae</i> DNA. The sample may not contain human RPP30 DNA.
+	—	—	—	+	The sample contains <i>N. gonorrhoeae</i> DNA and human RPP30 DNA.
—	+	—	—	—	The sample contains <i>T. pallidum</i> DNA. The sample may not contain human RPP30 DNA.
—	+	—	—	+	The sample contains <i>T. pallidum</i> DNA and human RPP30 DNA.
—	—	+	—	—	The sample contains <i>M. genitalium</i> DNA. The sample may not contain human RPP30 DNA.
—	—	+	—	+	The sample contains <i>M. genitalium</i> DNA and human RPP30 DNA.
—	—	—	+	—	The sample contains <i>U. urealyticum</i> DNA. The sample may not contain human RPP30 DNA.
—	—	—	+	+	The sample contains <i>U. urealyticum</i> DNA and human RPP30 DNA.
+	+	+	+	—	The sample contains <i>N. gonorrhoeae</i> , <i>T. pallidum</i> , <i>M. genitalium</i> , <i>U. urealyticum</i> DNA. The sample may not contain human RPP30 DNA.
+	+	+	+	+	The sample contains <i>N. gonorrhoeae</i> , <i>T. pallidum</i> , <i>M. genitalium</i> , <i>U. urealyticum</i> DNA and human RPP30 DNA.

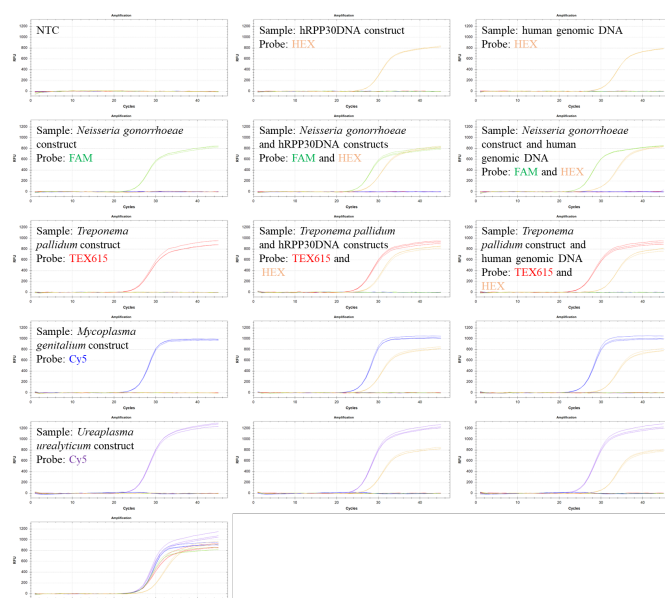


Figure 1: Verification experiments with single targets (given in text boxes for each panel). All sets of probes and primers are present in every reaction, but positive signal is only observed when the target(s) is present, indicating that the amplification is specific.

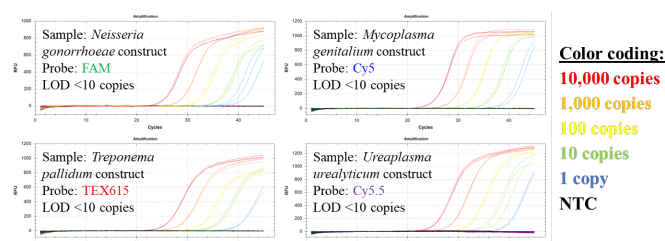


Figure 2: Serial dilution experiments show LOD <10 molecules for the synthetic DNA construct of each target.

VERIFICATION EXPERIMENTS

The STI2 assay verification was carried out as a 5-plex assay that simultaneously detects DNA from *Neisseria gonorrhoeae*, *Mycoplasma genitalium*, *Treponema pallidum*, *Ureaplasma urealyticum*, and human RPP30 DNA, which serves as a positive control assay.

Experiments were performed in triplicate using the experimental procedure given above, but with different samples added to each reaction. The samples used for the verification experiments contained 1×10^4 copies/reaction of synthetic 500 bp DNA constructs (from Twist Biosciences) harboring the regions of interest from the pathogen genomes, human RPP30 DNA gene, and human genomic DNA. **Figure 1** shows the results of these experiments, which indicate that the 5-plex specifically detects the different pathogens.

The limit of detection (LOD) was estimated by performing serial dilution experiments in triplicate (**Figure 2**). For dilution series only target construct was added. The results show a limit of detection (LOD) <10 copies/reaction.

NOTES

- ¹ FAMTM (Carboxyfluorescein), a trademark of Life Technologies Corporation.
- ² BHQ-1TM (Black Hole Quencher) is a trademark of Biosearch Technologies, Inc.
- ³ TEX615TM is a trademark of Thermo Fisher Scientific.
- ⁴ BHQ-2TM (Black Hole Quencher) is a trademark of Biosearch Technologies, Inc.
- ⁵ Cy5TM, a trademark of GE Healthcare.
- ⁶ Cy5.5TM is a trademark of GE Healthcare.
- ⁷ HEXTM (Hexachloro-fluorescein), a trademark of Thermo Fisher Scientific.
- ⁸ TaqManTM is a trademark of Roche Diagnostics, Inc.

Conclusion: The data in **Figure 1** indicate that the 5-plex primers and probes specifically detect and differentiate the pathogens and are also compatible with RPP30_DNA positive control primers. Human genomic DNA doesn't interfere with the detection of the pathogens.

CONTACT US

For assistance, please contact DNA Software using the link:
<https://www.pcrassays.com/contact/>

Address: Michigan Life Science and Innovation Center,
46701 Commerce Center Dr, Plymouth, MI 48170
Phone: (734) 222-9080