SmartNotes



Yes. The new Thermo Scientific™ Oxoid™ Columbia CNA Agar/Thermo Scientific™ Brilliance™ GBS Agar Biplate detects Group B Streptococci (GBS), including non-hemolytic GBS, as well as clinically-relevant Gram positives, like enterococci, staphylococci and streptococci. Results are available in 18-24 hours, without using a pre-enrichment step for GBS detection, and no need for reincubation.

Why is it a benefit to use a biplate combination with Oxoid Columbia CNA Agar and *Brilliance* GBS Agar?

The biplate combination of Oxoid Columbia CNA Agar/*Brilliance* GBS Agar enables the simultaneous detection of clinically-relevant Gram positives, like enterococci, staphylococci and streptococci on Oxoid Columbia CNA and also GBS on *Brilliance* GBS Agar, including non-hemolytic GBS which could be missed on CNA alone.

As a lab automation user, what is the benefit of using a biplate combination of Oxoid Columbia CNA Agar/ *Brilliance* GBS Agar?

By using this biplate you can use two samples instead of one at the same time with your lab automation for HVS (high vaginal swabs) and increase your output by 50%. You can also save space by using a biplate rather than two single plates in the incubator and plate stacking. Furthermore, by having both media available on one image you will have a more comprehensive picture of Gram positive bacteria present in the sample.

Thermo Fisher

thermoscientific

Do I need to use an enrichment step to detect GBS?

No. An enrichment step is not necessary to detect GBS on *Brilliance* GBS Agar since the medium incorporates patented Inhibigen™ technology, making the medium highly selective to background flora and very sensitive to GBS.

Why should I use an extra chromogenic media for GBS when I can detect these on a simple blood agar?

Whilst the significant proportion of GBS infections are hemolytic, approximately 4% of all GBS infections are non-hemolytic and therefore not identified on a standard blood plate. *Brilliance* GBS is able to also detect non-hemolytical GBS as bright pink colonies which are distinct and easy to read.

What kind of sample can I use with *Brilliance* GBS Agar?

Low vaginal swabs (LVS), high vaginal swabs, (HVS), vaginal or recto-vaginal samples can be used to detect GBS. The studies are available at thermofisher.com/brilliance-gbs-cna for further review.

How do I incubate the Oxoid Columbia CNA Agar/*Brilliance* GBS Agar Biplate?

Aerobically. No additional kits or jars are needed to create any specific atmosphere and only 24 hours of incubation are required, instead of 48 hours with traditional media.

Is it possible to perform automated antimicrobial susceptibility testing (AST) analysis directly with colonies from *Brilliance* GBS Agar?

Yes. It is possible to use the colonies from *Brilliance* GBS Agar directly to confirm GBS using automated AST systems.

International:

+44 (0) 1256 841144 microbiology@thermofisher.com

