

Sheep and Horse Blood/Additive

355-6641
355-6652

- Meat-yeast agar (fresh blood)

DEFINITION

Sheep and horse blood is included in the composition of certain culture media:

- either as an enrichment factor for the preparation of agars used for the culture of fastidious bacteria.
- or for investigating the hemolytic response of some of them, streptococci in particular.

PRESENTATION

• Sterile horse blood

5 ml x 1 vial

code 355-6641

• Sterile sheep blood

50 ml x 1 bottle

code 355-6652

STORAGE

- Ready-to-use: + 2°C to 8°C in a dry place.
- Expiration date and batch number are shown on the package.

N.B.: These products are delivered on subscription.

UTILISATION

1. Fresh blood agar

(study of hemolytic response)

Add 5-10% of blood to the sterile base agar maintained at a temperature of 44°C - 47°C, then pour into Petri dishes.

2. "Chocolate" agar (cooked blood agar)

After the addition of 10% of blood to the base agar, it is distributed in bottles and placed in a water-bath at 80°C for 10-15 minutes until a chocolate shade is obtained. Agitate and pour into Petri dishes.

Chocolate agar releases factors (in particular hemin) conducive to the growth of *Haemophilus* and *Neisseria*.

3. Agar with the addition of blood

(See preparation)

- Columbia agar (fresh or cooked blood)
- Blood agar (fresh or cooked blood)
- Blood-enriched agar (fresh or cooked blood)
- Blood agar + Nalidixic acid (fresh blood)
- Schaedler agar (fresh blood)
- Trypto-Casein-Soy (fresh blood)

QUALITY CONTROL OF MANUFACTURER

Every product manufactured and marketed by Bio-Rad is subject to a quality-assurance procedure at all stages, from the reception of raw materials to the marketing of the end-product. Each batch of finished product undergoes quality control and is marketed only if it satisfies the acceptability criteria.

Documentation relative to the production and control of each batch is kept on file.

KEY WORDS

Blood / Sheep / Horse / Chocolate Agar.