

RAPID' *Salmonella* Capsule / Supplement

356-4710

356-4709

356-4712

SCOPE OF APPLICATION

The RAPID' *Salmonella* capsule and RAPID' *Salmonella* Supplement are a selective supplement for addition to Buffered Peptone Water as part of the RAPID' *Salmonella* method – a short, NF VALIDATION and Nordval-certified protocol (see RAPID' *Salmonella* technical data sheet), for detection of *Salmonella* spp. in human and animal food products, and in environmental samples.

FORMATS

- **100 RAPID' *Salmonella* capsules**
100 x QSP 250 ml **code 356-4710**
- **100 RAPID' *Salmonella* capsules (10 times concentrated)**
100 x QSP 2.5 Litres **code 356-4709**
- **RAPID' *Salmonella* Supplement box**
1 x QSP 100 analyses **code 356-4712**

STORAGE / SHELF LIFE / BATCH

- At +2 – 8°C away from light.
- The expiry date and batch number are indicated on the packaging.

PRINCIPLE

The RAPID' *Salmonella* capsule and RAPID' *Salmonella* Supplement are a selective supplement for addition to Buffered Peptone Water for carrying out selective salmonella enrichment. The formulation enables highly effective *Salmonellae* growth, even when stressed, while limiting the growth of interfering flora.

TYPICAL FORMULA

RAPID' *Salmonella* Capsule **code 356-4710**

QSP 250 ml

Selective mixture 15 mg
Excipient qsp 575 mg

RAPID' *Salmonella* Capsule (10 times concentrated) **code 356-4709**

QSP 2.5L

Selective mixture 150 mg
Dye 3.75 mg
Excipient qsp 575 mg

RAPID' *Salmonella* Supplement box

code 356-4712

QSP 100 analyses

Selective mixture 1500 mg
Dye 37.5 mg
Excipient qsp 5,75 g

PRODUCT(S) REQUIRED (NOT PROVIDED)

• Buffered peptone water

6 x 225 ml bottles (ex. code **355-4170**)
500 g (ex. code **356-4684**)
5 x 2.3 l bags (ex. code **355-5789**)
2 x 5 l bags (ex. code **355-5790**)

• RAPID' *Salmonella* medium

90 mm x 20 boxes (code **356-3961**)
90 mm x 100 boxes (code **356-3962**)
500 g (code **356-4705**)

Preparation of samples by the alternative RAPID' *Salmonella* method – short protocol

Sample preparation with direct addition of the capsule supplement in the enrichment broth

Note: In the context of NF VALIDATION mark, no samples of over 25 grams were tested.

Dilute 7 g or 7 mL of the sample in 9 x 7 mL Buffered Peptone Water.

Example: dilute 25 g or 25 mL of the sample in 225 mL Buffered Peptone Water and dilute to 1/10th.

ISO standard 6579 specifies how to prepare the stock suspension (cocoa, acid foods etc.).

Homogenise in a Stomacher type blender.

Open a RAPID' *Salmonella*, QSP 250 ml capsule (code **356-4710**) and pour the contents directly into the broth.

Homogenise by agitating vigorously.

Note: Either the whole capsule or its contents only can be added before the stomacher phase. In order to make handling easier, we recommend opening the capsule and pouring out the contents (see PRECAUTIONS FOR USE).

RAPID' *Salmonella* Capsule / Supplement

Sample preparation with addition of the capsule supplement as a concentrated solution, to the enrichment broth

The capsule contents can be diluted in Buffered Peptone Water or Sterile Distilled Water first, for incorporation in liquid form.

- Dilute η g or η mL of the sample in 9 x η ml of Buffered Peptone Water.

- Homogenise in a Stomacher blender.

- Where RAPID' *Salmonella* QSP 250 ml capsules (code **356-4710**) are used: Open n capsules and pour the contents directly into n x 10 ml Buffered Peptone Water to obtain a concentrated supplement solution.

Add η x 0.4ml of the concentrated supplement solution to the sample to be analysed.

Homogenise by agitating vigorously.

- Where RAPID' *Salmonella* QSP 2.5 Litre capsules are used (code **356-4709**): Open n capsules and pour their contents directly into an empty recipient. Fill with n x 10 ml Buffered Peptone Water or n x 10 ml of Sterile Distilled Water.

Homogenise by agitating vigorously to obtain a red concentrated solution.

Add η x 0.04ml of the concentrated supplement solution to the sample diluted in the Buffered Peptone Water.

Homogenise by agitating vigorously.

- Where RAPID' *Salmonella* Supplement (code 356-4712), 1 x QSP 100 analyses are used: Open the box and fill with 100 ml of Buffered Peptone Water or Sterile Distilled Water. Homogenise by agitating vigorously to obtain a red concentrated solution.

Add η x 0.04ml of the concentrated supplement solution to the sample diluted in the Buffered Peptone Water.

Homogenise by agitating vigorously.

Example for a 10g sample:

- Dilute the 10g sample in 90 ml Buffered Peptone Water.

- Homogenise using a Stomacher type blender.

- Dilute 1 RAPID' *Salmonella* QSP 250 ml capsule (code 356-4710) in 10 ml Buffered Peptone Water to obtain a concentrated supplement solution

- Add 4 ml of concentrated supplement solution to the 90 ml Buffered Peptone Water diluent + sample in order to achieve the correct capsule dilution ratio.

NB.: The concentrated solution, once reconstituted with Buffered Peptone Water or Sterile Distilled Water can be stored for 1 week at ambient temperature, or at +2-8°C.

PRECAUTIONS FOR USE

- Good Laboratory Practice must be observed. (EN ISO 7218)

- If the whole capsule is added to the Buffered Peptone Water, sterile tweezers must be used to add the capsule to the bag. We recommend that you check that the capsule actually opened during the stomaching stage.

- If the capsule contents are handled with the fingers, it cannot be added to the enrichment broth due to the risk of contamination.

- The RAPID' *Salmonella* capsule and RAPID' *Salmonella* Supplement contain selective agents and an excipient. Selective agents dissolve very well. The excipient however, remains in suspension and may create a deposit when the contents of the capsule are diluted in a small quantity of Buffered Peptone Water or Sterile Distilled Water. Always shake well therefore before using the concentrated solution.

QUALITY CONTROL

All products manufactured and sold by Bio-Rad are subject to a quality assurance system, from receipt of raw materials to commercialisation of the finished products.

Each finished product batch is subject to quality control and is only marketed if it is compliant with the applicable acceptance criteria.

Documentation relating to batch production and control is kept on file.

MICRO-ORGANISMES	Culture des micro-organismes en 24 heures à 37°C
<i>Salmonella</i> Enteritidis ATCC 13076	Colonies magenta
<i>Salmonella</i> Typhimurium ATCC 14028	Colonies magenta
<i>Escherichia coli</i> ATCC 25922	Inhibition totale ou partielle Colonies non colorées
<i>Enterococcus faecalis</i> ATCC 19433	Inhibition totale

KEYWORDS

RAPID' *Salmonella*/Capsule/*Salmonella*/Food products/Detection/Chromogenic/Medium.

RAPID' *Salmonella* Capsule / Supplement

BIBLIOGRAPHY

- **PERRY J.D., FORD M., TAYLOR J., JONES A.L., FREEMAN R., GOULD F.K.** 1999: ABC medium, a new chromogenic agar for selective isolation of *Salmonella* spp. J Clin Microbiol, 37 (3): 766-8.
- **LAUER W., MARTINEZ F.** 2009: RAPID' *Salmonella* chromogenic medium. Performance Tested Method 050701. J AOAC Int. 2009 Nov-Dec;92(6): 1871-5