

## Bacterial Serotyping Guide for *Salmonella*



## Minimizing Risk

The monitoring of veterinary diseases and quality control of industrial products are public health issues. Microbial populations which cause infection vary over time depending on manufacturing and transport conditions. Consequently, to prevent such risks, it is important to take an active role in the overall food chain monitoring process. As part of such an approach, studying bacteria from an epidemiological standpoint is essential for monitoring and anticipating dynamic changes in microbial structures. An antiserum is a high performance reagent enabling accurate identification of bacteria. It thus makes classification and evaluation of changes in such microbial populations possible.

## Serotyping as an Identification Tool

Serotyping (serological typing) is based on the long-standing observation that microorganisms from the same species can differ in the antigenic determinants expressed on the cell surface. Serotyping is one of the classic tools for epidemiological study and is applied to numerous species that express different serotypes, such as: *Escherichia coli*, *Pseudomonas aeruginosa*, *Salmonella* species, *Shigella* species, *Yersinia* and *Vibrio cholerae*.

## Expertise in Serotyping

With an extensive range of immune sera for serotyping, Bio-Rad offers a wide range of reagents serving the needs of veterinary, food and service laboratories. The method's ease of use enables every laboratory to perform identification of bacterial strains. Data collected promotes enhanced prevention of veterinary, human, and industrial risks.

# Identification of *Salmonella* by Serotyping

Salmonellae are Gram-negative, flagellated, facultative anaerobic bacilli possessing three major antigens: H or flagellar antigen, O or somatic antigen, and Vi (capsular) antigen (possessed by only a few serovars). The different species are serotyped according to these three different antigens.

- **H (flagellar) antigen** may occur in either or both of two forms, phase 1 and phase 2. There are over 1800 known serovars which current classification considers being separate species. The organisms tend to change from one phase to the other.
- **O (somatic) antigens** occur on the surface of the outer membrane and are determined by specific sugar sequences on the cell surface.
- **Vi (capsular) antigen** is a superficial antigen overlying the O antigen (additional surface antigen). It is present in a few serovars, the most important being *Salmonella* Typhi, but also present in *Salmonella* Paratyphi C and *Salmonella* Dublin.

Once the O, H-phase 1 and H-phase 2 are identified, the antigenic formula can be used to identify the serotype by referring to a **Kauffman-White reference catalog**.

- The formula gives the O antigen(s) first followed by the H antigen(s). O antigens, Vi (when present), H antigens phase 1, H antigens phase 2 (when present).
- Colons separate the major antigens and commas separate the components of the antigens.
- Further conventions:
  - Underlined O factor is encoded by a bacteriophage (lysogenic strain)
  - [ ]: square bracketed factor may or may not be present (not phage-encoded)
  - { }: curly bracketed factor never coexists with others (exclusive)
  - ( ): parenthesis around a factor indicate weakly agglutinable factor

## Examples

### ***Salmonella enterica* serotype Typhimurium: 1,4,[5],12:i:1,2**

This strain has the O antigen factors 1, 4, [5], and 12, the flagellar H antigen i (1st phase) and the flagellar H antigens 1 and 2 (2nd phase).

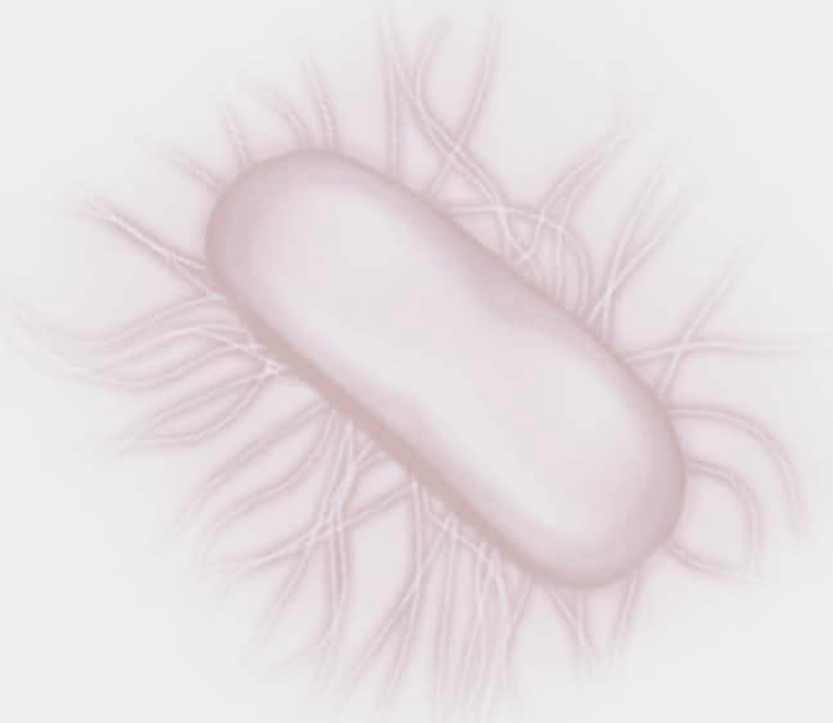
### ***Salmonella enterica* serotype Lagos: 1,4,[5],12:i:1,5**

This strain has the O antigen factors 1, 4, [5], and 12, the flagellar H antigen i (1st phase) and the flagellar H antigens 1 and 5 (2nd phase).

### ***Salmonella enterica* serotype Virchow: 6,7,14:r:1,2**


This strain has the O antigen factors 6, 7 and 14; the flagellar H antigen r (1st phase) and the flagellar H antigens 1, 2 (2nd phase).


For more examples, please refer to our Quick Guide (#xxx) including a table with some typical *Salmonella* serotypes isolated in food products.





# How to Perform *Salmonella* Serotyping

## A Simple Protocol

- 

1 After confirmation, isolate colonies on TSI slant (or recommended agar) for identification by antisera. It is important to use pure cultures.
- 

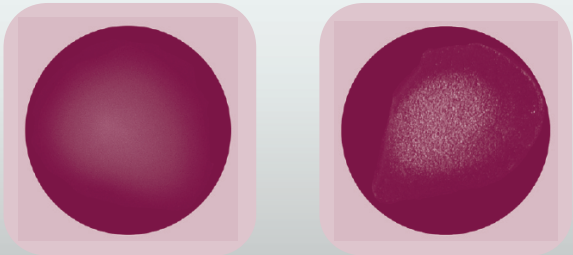
2 Place 1 free falling drop of antisera onto a slide for agglutination and add a pure colony.
- 

3 Mix reagents thoroughly.
- 

4 Rock the slide in the circular motion for 30 seconds and observe for agglutination.

- Serotyping is performed after identification of the species on a fresh, pure culture of *Salmonella* isolated on a non-selective agar medium. There are several media recommended for use, including Müller-Hinton or Nutrient Agar, TSI (Triple Sugar Iron) and/or LIA (Lysine Iron Agar) slants or TSA (Tryptic Soy Agar).
- Where polyvalent and monovalent antisera are available, start by testing agglutination with polyvalent sera, then with the specific monovalent sera corresponding to the mixture giving marked agglutination.
- Agglutination should appear between 1–10 seconds. If agglutination occurs > 60 sec, the antigens can not be identified correctly.

## Easy Interpretation

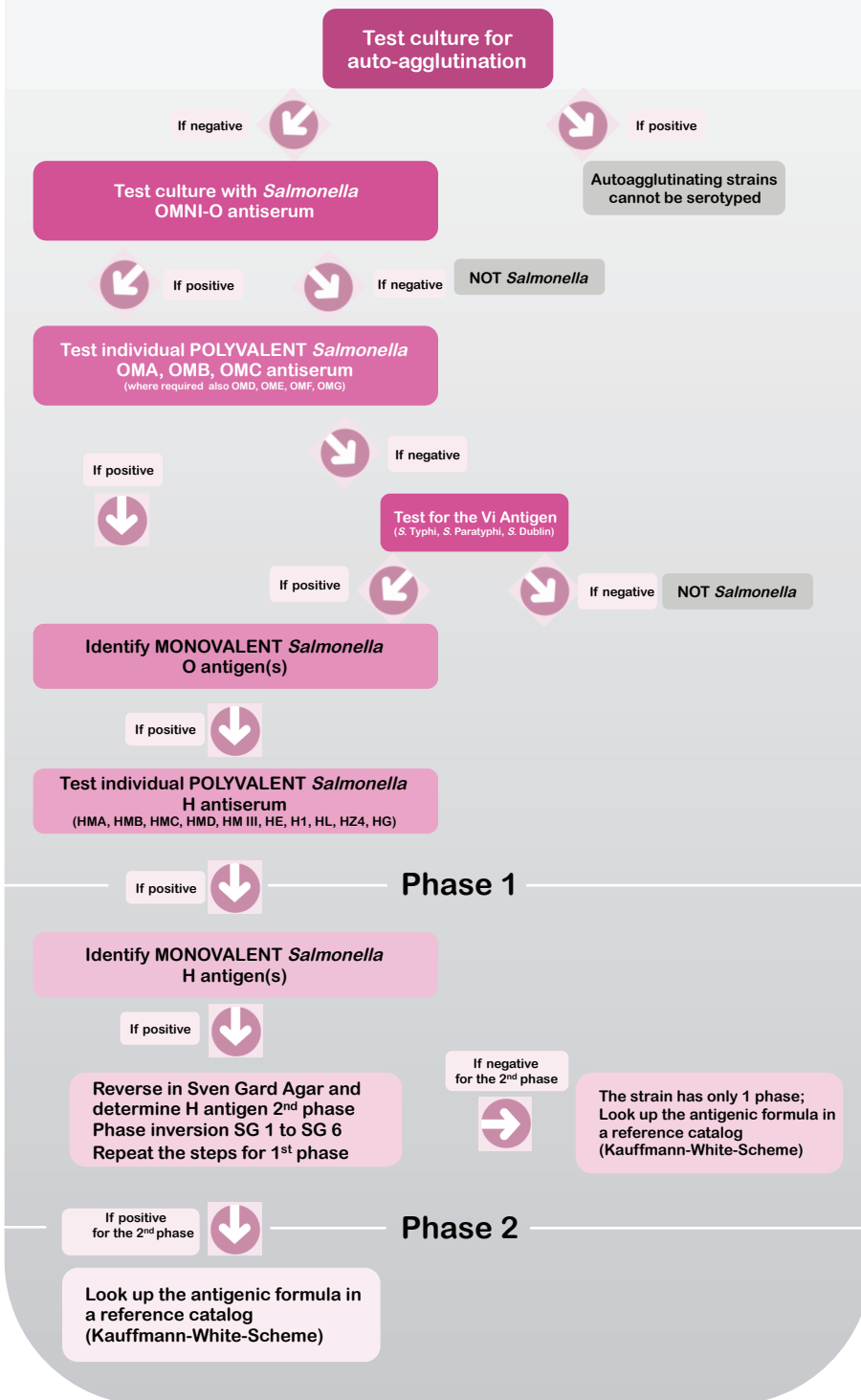


**No agglutination =  
NEGATIVE result**

**Agglutination =  
POSITIVE result**



## Salmonella Serotyping Flowchart



### 1. Test the culture for auto-agglutination

- Test the culture first in Physiological Water/Saline; strains that produce auto-agglutination cannot be serotyped.
- In addition, all strains should be tested with the *Salmonella* Omni-O antiserum which contains antigens A – 60 for the presumptive identification of O-agglutinable strains of *Salmonellae*.
- Agglutination of a strain of *Salmonella* with Omni-O antiserum indicates that the strain is O-agglutinable and can be serotyped with specific sera.

### 2. Test for the O antigens

- Begin by testing the isolate with polyvalent O antiserum. The majority (about 98%) of *Salmonella* encountered in warm-blooded animals possess an O antigen corresponding to the agglutinins contained in OMA, OMB and OMC sera.
- When agglutination occurs with one of these 3 groups, the isolate is positive for that group.
- The individual monovalent O antisera are used to identify the O antigen(s).
- Repeat the agglutination step by testing the isolate in each monovalent O antiserum present in the group.

**Example:** The polyvalent O antiserum OMA shows agglutination therefore the following monovalent O antisera must be tested: O:1,2; O:4,5; O:9, O:46; O:3,10,15; O:1,3,19

- When a strain does not agglutinate the OMA, OMB or OMC polyvalent sera, it is recommended to test this strain with Vi serum and the other polyvalent O sera.
- If a Vi positive reaction is observed, the bacterial suspension must then be heated to 100°C for 30 minutes, before repeating the test with polyvalent OMA, OMB and OMC sera and the corresponding monovalent sera to define the O antigen.

### 3. Test for the H Antigen – Phase 1

- Begin by testing the isolate with a polyvalent H antiserum (HMA – HG). When agglutination occurs with one of these groups, the isolate is positive for that group.
- The individual monovalent H antiserum is used to identify the H antigen.
- Repeat the agglutination step by testing the isolate in each monovalent H antiserum present in the group.

**Example:** The polyvalent H Antiserum HMA shows agglutination therefore the following monovalent H antisera must be tested: a; c; d; l; z10.

Once the first H antigen is identified, a phase inversion on the isolate must be performed to force the organism to repress its dominant H phase and grow in the second phase.

### 4. Phase Inversion – Sven Gard Method

Sven Gard medium is used during serotyping of *Salmonella* to demonstrate the inapparent H antigen phase of biphasic *Salmonella* (Sven Gard method). Sven Gard agar should be used with the following *Salmonella* antisera: SG 1 to SG 6.

#### Example:

H:1,2 antigens were identified in a culture → SG6 (1,2 + 1,5 + 1,7 + z6) serum is used for phase inversion.

H:r antigen is identified in a culture → SG4 (r + z) antiserum is used for phase inversion.

### 5. Test for the H Antigen – Phase 2

- A culture at the periphery of the invasion zone of the Sven Gard agar should be taken.
- Start testing again by using the H polyvalent antisera (HMA – HG). If there is no agglutination, this serotype contains only one phase.
- If one of these groups shows agglutination, define the specific H phase by using the relevant H monovalent antisera.

As the antigenic formula with O, H – phase 1 and H – phase 2 are identified, the serotype is now specified by referring to a reference catalog, such as the Kauffman-White scheme.

Serotype	O Antigens	H Antigens Phase 1	H Antigens Phase 2
Agona	1, 4, [5], 12	f, g, s	[1, 2]
Anatum	3, {10} {15} {15, 34}	e, h	1, 6
Bareilly	6, 7, 14	y	1, 5
Blockley	6, 8	k	1, 5
Bovis Morbificans	6, 8, 20	r, [i]	1, 5
Brandenburg	1, 4, 12	e, h	e, n, z15
Bredeney	1, 4, 12, 27	l, v	1, 7
Chester	1, 4, [5], 12	e, h	e, n, x
Derby	1, 4, [5], 12	f, g	[1, 2]
Dublin	1, 9, 12, [Vi]	g, p	-
Enteritidis	1, 9, 12	[f], g, m, [p]	[1, 7]
Gallinarium	1, 9, 12	-	-
Gloucester	1, 4, 12, 27	i	l, w
Hadar	6, 8	z10	e, n, x
Heidelberg	1, 4, [5], 12	R	1, 2
Indiana	1, 4, 12	z	1, 7
Infantis	6, 7, 14	R	1, 5
Javiana	1, 9, 12	l, z28	1, 5
Kentucky	8, 20	i	z6
Kottbus	6, 8	e, h	1, 5
Lagos	1, 4, [5], 12	i	1, 5
Lille	6, 7, 14	z38	-
Livingstone	6, 7, 14	d	l, w
Mbandaka	6, 7, 14	z10	e, n, z15
Meleagridis	3, {10} {15} {15, 34}	e, h	l, w
Montevideo	6, 7, 14	g, m, [p], s	[1, 2, 7]
Muenchen	6, 8	d	1, 2
Newport	6, 8, 20	e, h	1, 2
Orion	3, {10} {15} {15, 34}	y	1, 5
Paratyphi B	1, 4, [5], 12	b	1, 2
Saintpaul	1, 4, [5], 12	e, h	1, 2
Senftenberg	1, 3, 19	g, [s], t	-
Stanley	1, 4, [5], 12, [27]	d	1, 2
Thomson	6, 7, 14	k	1, 5
Typhimurium	1, 4, [5], 12	i	1, 2
Virchow	6, 7	r	1, 2
Weltervreden	3, {10} {15}	r	z6



Bio-Rad provides a complete solution for the detection and confirmation of *Salmonella* spp. and a large panel for the serological identification of a variety of *Salmonella* serotypes.

### Detect and Identify *Salmonella* in 4 Simple Steps



**Sample enrichment**

1



Standard Methods:  
EN ISO 6579/A1  
FDA BAM  
USDA FSIS MLG  
MFHPB-20

OR



OR



iQ-Check Prep, CFX96 Deep Well & and IQ-Check Kits

**Detection of *Salmonella***

2



***Salmonella* Latex  
Oxidase Test  
ONPG**

**Biochemical  
confirmation**

3



**Omni-O / Vi  
Polyvalent / Monovalent O  
Polyvalent / Monovalent H  
Phase inversion**

**Serological  
identification**

4

The product range includes the required enrichment media. Detection can be performed by either classical standard reference methods or by one of Bio-Rad's alternative validated rapid methods, iQ-Check® real-time PCR solution or RAPID Chromogenic media. Several easy-to-use confirmation and identification tests are also available.

# Detection and Identification of *Salmonella*

## Ordering Information



### Catalog # Description

#### Special *Salmonella* Antisera

- 356-0781 Total polyvalent *Salmonella* Omni-O antiserum (A60); 3 ml vial with 60 tests  
356-1261 *Salmonella* Kit/Typhi and Paratyphi fever; *Salmonella* polyvalent agglutinating antisera T, A, B, C et Vi; 5 dropper vials  
356-0951 *Salmonella* agglutinating antiserum Vi; 3 ml vial with 60 tests

#### *Salmonella* Monovalent O Antisera; 3 ml dropper bottle with 60 tests

- | Bio-Rad equivalent serum                                      | O group designated by Letters        |
|---|--------------------------------------|
| 355-9021 <i>Salmonella</i> monovalent antiserum O: 4, 5       | <i>Salmonella</i> O group B          |
| 355-9031 <i>Salmonella</i> monovalent antiserum O: 1, 2       | <i>Salmonella</i> O group A          |
| 355-9062 <i>Salmonella</i> monovalent antiserum O: 6, 7, 8    | <i>Salmonella</i> O group C1-C4      |
| 355-9081 <i>Salmonella</i> monovalent antiserum O: 7          | <i>Salmonella</i> O group C1, C4     |
| 355-9091 <i>Salmonella</i> monovalent antiserum O: 8          | <i>Salmonella</i> O group C2, C3     |
| 355-9101 <i>Salmonella</i> monovalent antiserum O: 9          | <i>Salmonella</i> O group D1         |
| 355-9112 <i>Salmonella</i> monovalent antiserum O: 3, 10, 15  | <i>Salmonella</i> O group E1, E2, E3 |
| 355-9127 <i>Salmonella</i> monovalent antiserum O: 15         | <i>Salmonella</i> O group            |
| 355-9131 <i>Salmonella</i> monovalent antiserum O: 1, 3, 19   | <i>Salmonella</i> O group E4         |
| 355-9141 <i>Salmonella</i> monovalent antiserum O: 11         | <i>Salmonella</i> O group F          |
| 355-9162 <i>Salmonella</i> monovalent antiserum O: 13, 22, 23 | <i>Salmonella</i> O group G1, G2     |
| 355-9171 <i>Salmonella</i> monovalent antiserum O: 6, 14, 24  | <i>Salmonella</i> O group H          |

#### *Salmonella* Monovalent H Antisera; 3 ml dropper bottle with 60 tests

- 356-1111 *Salmonella* monovalent H antiserum: 2  
356-1112 *Salmonella* monovalent H antiserum: 5  
356-1113 *Salmonella* monovalent H antiserum: 6  
356-1114 *Salmonella* monovalent H antiserum: 7  
356-0111 *Salmonella* monovalent H antiserum: a  
356-0121 *Salmonella* monovalent H antiserum: b  
356-0131 *Salmonella* monovalent H antiserum: c  
356-0141 *Salmonella* monovalent H antiserum: d  
356-1121 *Salmonella* monovalent H antiserum: g, m  
356-1122 *Salmonella* monovalent H antiserum: g, p  
356-1119 *Salmonella* monovalent H antiserum: h  
356-0161 *Salmonella* monovalent H antiserum: i  
356-0171 *Salmonella* monovalent H antiserum: k

#### Diluents and Physiological Water

- 355-4154 Distilled Sterile Water; 9 ml x 25 tubes  
355-4164 Physiological Sterile Water (0,85%); 9 ml x 25 tubes  
356-1723 Disposable Agglutination Cards; 3 x 15 cards

#### *Salmonella* Confirmation

- 356-4384 TSI (Triple Sugar Iron) Agar, 500 g  
356-4482 Nutrient Agar (2,1%) without NaCl, 500 g  
355-3834 Oxidase disks, 2 x 50 disks  
356-0781 Antiserum *Salmonella* Omnivalent; Omni-O (A-60), 3 ml dropper bottle  
355-3822 ONPG, 50 disks  
355-6710 *Salmonella* Latex kit, 75 tests  
355-6711 *Salmonella* Confirm Latex, 50 tests

#### RAPID *Salmonella* — Chromogenic medium

- 356-3961 90 mm x 20 dishes  
356-3963 90 mm x 120 dishes  
356-4705 500 g

#### Selective Enrichment

- 355-6140 MKTTn Broth (Muller-Kauffmann Tetrathionat Novobiocin), 10 ml x 50 tubes  
356-4714 MKTTn Broth (Muller-Kauffmann Tetrathionat Novobiocin), dehydrated base, 500 g  
355-6139 MSRV Medium (semi-solid Rappaport Vassiliadis), 200 ml x 6 bottles  
356-4325 MSRV Medium (semi-solid Rappaport Vassiliadis), dehydrated, 500 g  
355-5773 RVS Broth (Rappaport Vassiliadis Soy), 10 ml x 50 tubes  
356-4324 RVS Broth (Rappaport Vassiliadis Soy), dehydrated, 500 g

### Catalog # Description

#### *Salmonella* Polyvalent O Antisera; 3 ml dropper bottle with 60 tests

- 356-0801 *Salmonella* polyvalent O antiserum OMA Group A, B, D, E, L (1, 2, 12 + 4, 5, 12 + 9, 12 + 9, 46 + 3, 10 + 3, 15 + 1, 3, 19 + 21)  
356-0811 *Salmonella* polyvalent O antiserum OMB Group C, F, G, H (6, 7 + 6, 8 + 11 + 13, 22 + 13, 23 + 6, 14, 24 + 8, 20)  
356-0821 *Salmonella* polyvalent O antiserum OMC Group I, J, K, M, N, O, P (16 + 17 + 18 + 28 + 30 + 35 + 38)  
356-0831 *Salmonella* polyvalent O antiserum OMD Group Q, R, S, T, U, V, W (39 + 40 + 41 + 42 + 43 + 44 + 45)  
356-0841 *Salmonella* polyvalent O antiserum OME Group X, Y, Z, 51 to 53, 61 (47 + 48 + 50 + 51 + 52 + 53 + 61)  
356-0851 *Salmonella* polyvalent O antiserum OMF Group O:54 to O:59 54 + 55 + 56 + 57 + 58 + 59)

#### *Salmonella* Polyvalent H Antisera; 3 ml dropper bottle with 60 tests

- 356-0451 *Salmonella* polyvalent H antiserum HMA (a + c + d + i + z10 + z29)  
356-0461 *Salmonella* polyvalent H antiserum HMB (e, h + e, n, x + e, n, z15 + g)  
356-0471 *Salmonella* polyvalent H antiserum HMC (k + y + l + z4 + r)  
356-0481 *Salmonella* polyvalent H antiserum HMD (z35 + z36 + z38 + z39 + z41 + z42 + z44 + z60)  
356-0493 *Salmonella* polyvalent H antiserum HM III (z52 + z53 + z54 + z55 + z57 + z61)  
356-0391 *Salmonella* polyvalent HE antiserum (e, h + e, n, x + e, n, z15)  
356-0401 *Salmonella* polyvalent H1 antiserum (1, 2 + 1, 5 + 1, 6 + 1, 7 + z6)  
356-0411 *Salmonella* polyvalent HL antiserum (l, v + l, w + l, z13 + l, z28 + l, z40)  
356-0431 *Salmonella* polyvalent HZ4 antiserum (z4, z23 + z4, z24 + z4, z32)  
356-0441 *Salmonella* polyvalent HG antiserum (f, g + g, p + g, m, s + g, m + m, t)

#### *Salmonella* Monovalent H Antisera; 3 ml dropper bottle with 60 tests

- 356-1117 *Salmonella* monovalent H antiserum: m  
356-1118 *Salmonella* monovalent H antiserum: p  
356-1124 *Salmonella* monovalent H antiserum: q  
356-0201 *Salmonella* monovalent H antiserum: r  
356-0202 *Salmonella* monovalent H antiserum: s  
356-1115 *Salmonella* monovalent H antiserum: v  
356-1116 *Salmonella* monovalent H antiserum: w  
356-1123 *Salmonella* monovalent H antiserum: x  
356-0211 *Salmonella* monovalent H antiserum: y  
356-0221 *Salmonella* monovalent H antiserum: z  
356-0241 *Salmonella* monovalent H antiserum: z10  
356-1120 *Salmonella* monovalent H antiserum: z15

#### Phase Inversion/Sven Gard Method Antisera; 3 ml dropper bottle with 60 tests

- 356-1011 *Salmonella* phase inversion SG 1 antiserum (a + b + c + z10)  
356-1021 *Salmonella* phase inversion SG 2 antiserum (d + i + e, h)  
356-1031 *Salmonella* phase inversion SG 3 antiserum (k + y + l, v + l, w + l, z13 + z28)  
356-1041 *Salmonella* phase inversion SG 4 antiserum (r + z)  
356-1051 *Salmonella* phase inversion SG 5 antiserum (e, n, x + e, n, z15)  
356-1061 *Salmonella* phase inversion SG 6 antiserum (1, 2 + 1, 5 + 1, 6 + 1, 7 + z6)  
355-3430 Sven Gard, Semiagar; 25 ml x 25 tubes

#### Real-time PCR Detection tests

- 357-8123 iQ-Check™ *Salmonella* II kit, 96 reactions  
357-8142 iQ-Check S. Enteritidis kit, 96 reactions  
357-8141 iQ-Check S. Typhimurium kit, 96 reactions

#### Enrichment Supplements

- 356-4709 RAPID *Salmonella* capsules, 10 x conc., 100 capsules  
356-4710 RAPID *Salmonella* capsules, 1 x conc., 100 capsules  
356-4712 RAPID *Salmonella* supplement, 1 box qsp 100 analysis  
356-4610 Novobiocin supplement, 1 g

#### Non-Selective (Pre-)Enrichment

- 356-4684 Buffered Peptone Water, 500 g  
355-4179 Buffered Peptone Water, 225 ml x 6 bottles  
355-5795 Buffered Peptone Water, 3 l x 4 bags  
355-5790 Buffered Peptone Water, 5 l x 2 bags  
356-4426 mTSB (modified Tryptone-Soya)/Broth, 500 g, dehydrated base  
355-5426 mTSB (modified Tryptone-Soya)/Broth, 225 ml x 6 bottles



**Bio-Rad**  
Laboratories, Inc.

Life Science  
Group

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 03 710 53 00 **Brazil** 55 11 3065 7550  
**Canada** 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00  
**France** 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300  
**Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666  
**New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04  
**Singapore** 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026 674 55 05  
**Taiwan** 886 2 2578 7189 **Thailand** 1800 88 22 88 **United Kingdom** 020 8328 2000