

# Quick Guide for *Salmonella* Serotyping

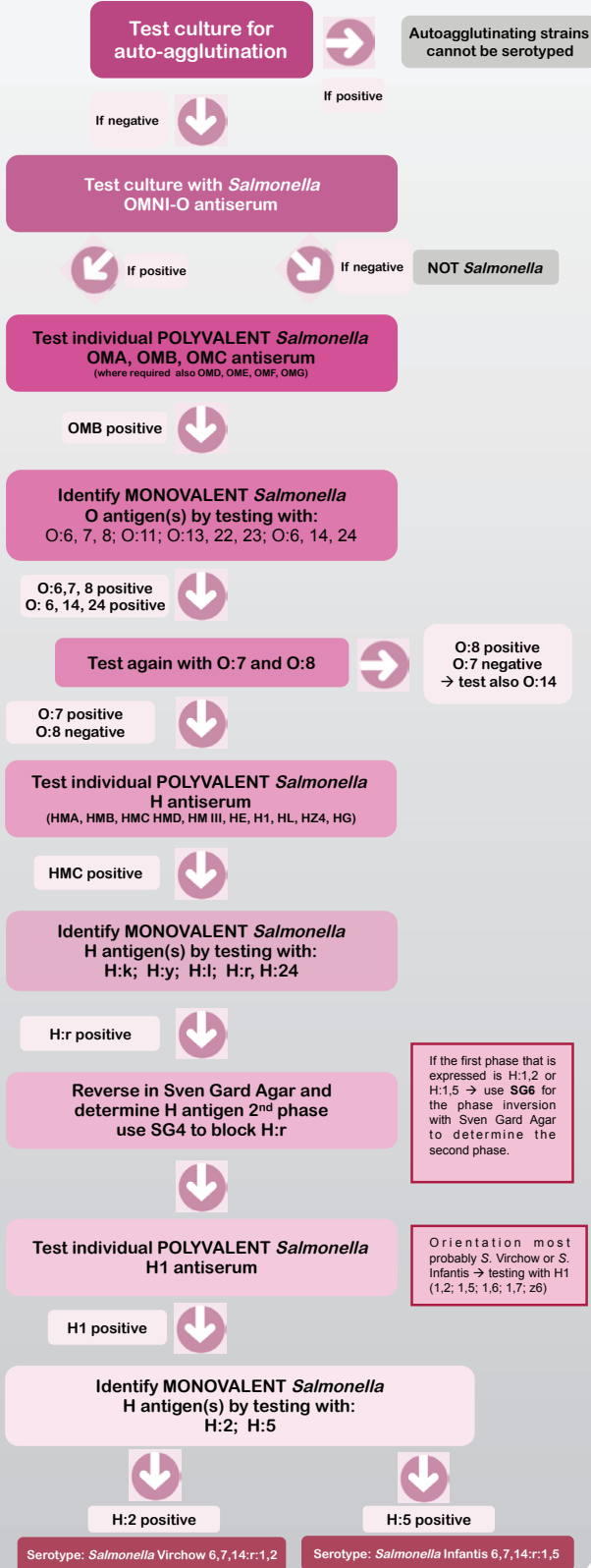
FOOD SCIENCE

Serotype	O Antigens	H Antigen Phase 1	H Antigen 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
Weltevreden	3, {10} {15}	r	z6

**Table 1:** Examples of some typical *Salmonella* serotypes isolated in food products.

# Quick Guide for *Salmonella* Serotyping Food Science Division

## Serotyping Flowchart – Example I



## Serotyping Flowchart – Example II

