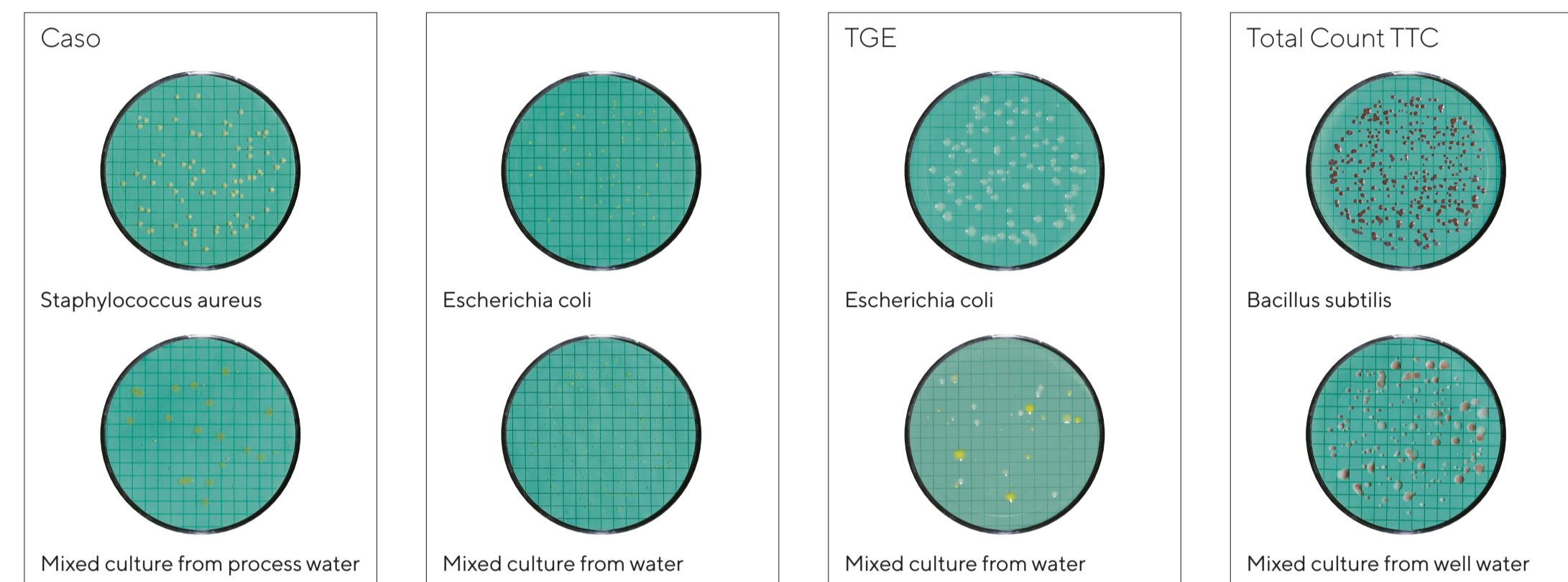
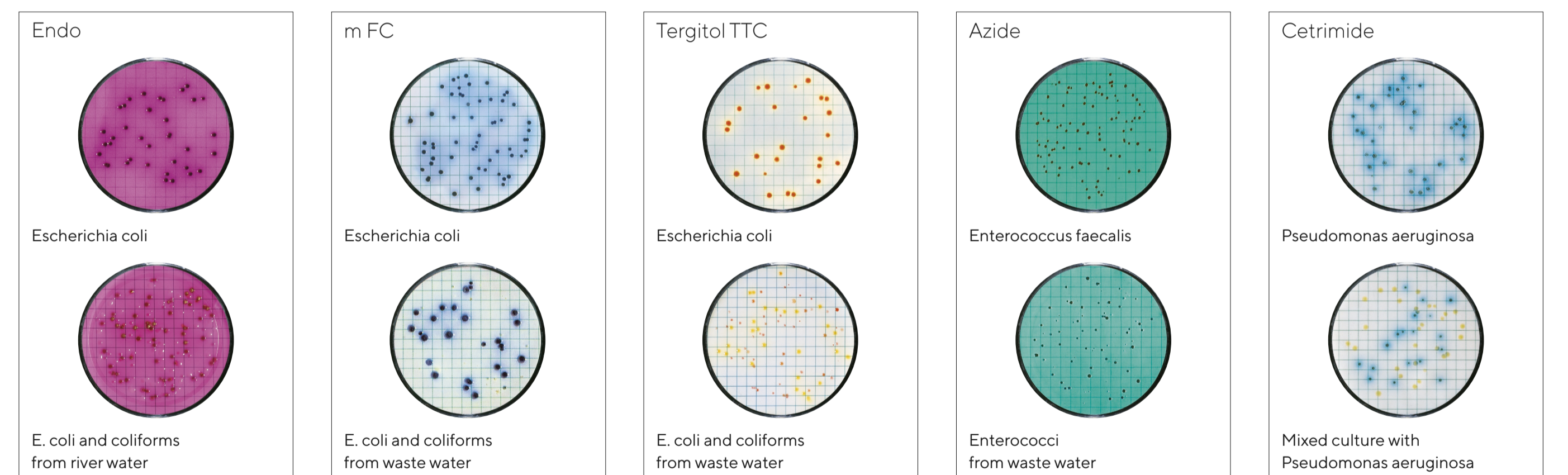


## Biosart® 100 Nutrient Media

### Total colony count



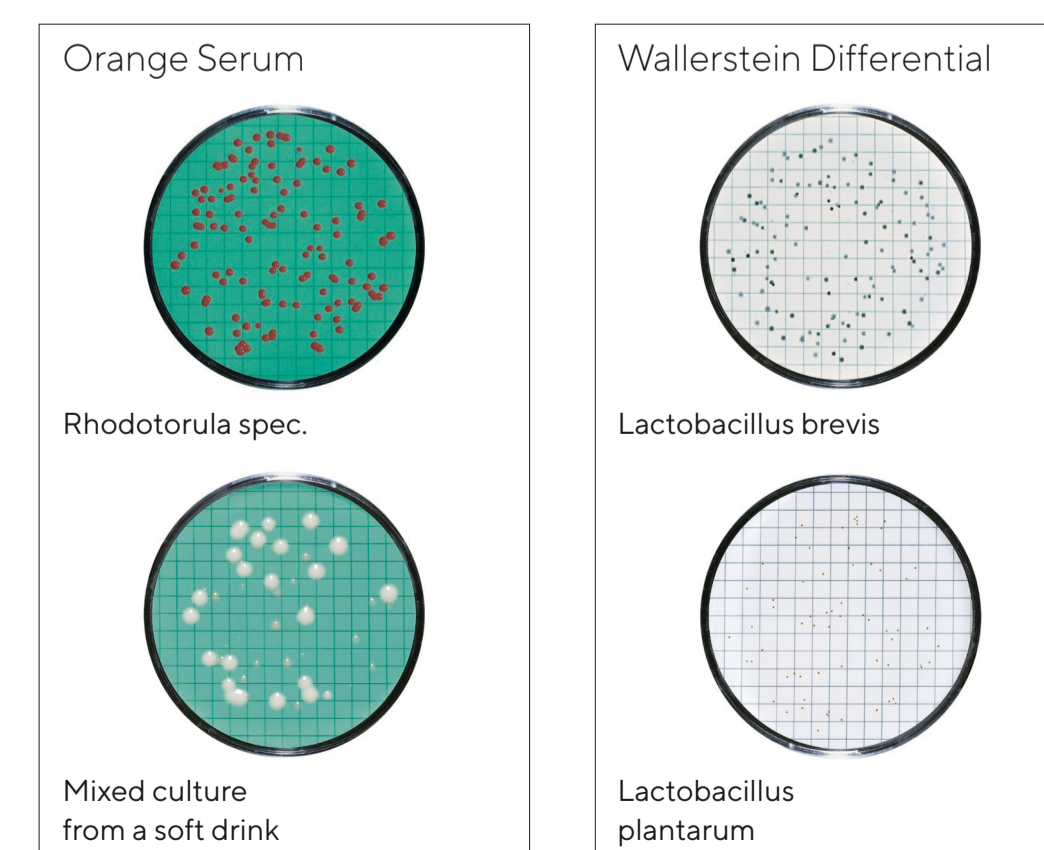
### E. coli and coliforms, Enterobacteria



### Yeasts and molds



### Product-spoiling microorganisms



### Typical Application Examples

Product	Detection and enumeration of ...	Biosart® 100 Nutrient Media Type
Beer	Lactobacilli and Pediococci other beer spoiling organisms	Wallerstein Differential
	Total colony count	Total Count TTC
	Yeasts and molds	Wallerstein Nutrient, Wort
Foods	Acid-tolerant microorganisms	Orange Serum
	Enterobacteria, E. coli and coliforms	Endo, m FC, Teepol   Tergitol TTC
	Enterococci, Enterococcus faecalis	Azide   KF Strep
	Pseudomonas aeruginosa	Cetrimide
	Total colony count	Caso, TGE   Tryptone Glucose Extract
Fruit juice	Yeasts and molds	Wort
	Enterobacteria, E. coli and coliforms	Endo, Tergitol TTC*
	Oenococcus and other product spoiling organisms	Orange Serum, Wallerstein Differential
Milk	Yeasts and molds	m Green yeast and mold   Schaufus Pottinger, Wallerstein Nutrient
	E. coli and coliforms	Endo
	Enterococci, Enterococcus faecalis	Azide   KF Strep
Pharmaceuticals, WFI, raw materials and cosmetics	Enterococci, Enterococcus faecalis	Azide   KF Strep
	Pseudomonas aeruginosa	Cetrimide
	Total colony count	Caso, R2A
	Yeasts and molds, Candida albicans	Sabouraud
	Acid-tolerant microorganisms, Lactic-acid bacteria	Orange Serum, Wallerstein Differential
Soft drinks, concentrates	Enterobacteria, E. coli and coliforms	Endo
	Total colony count	TGE   Tryptone Glucose Extract, Total Count TTC
	Yeasts and molds	m Green yeast and mold   Schaufus Pottinger, m Green yeast and mold selective, Wallerstein Nutrient, Wort
	E. coli and coliforms	Endo
Sugar, sugar products	Total colony count	Total Count TTC
	Yeasts and molds	m Green yeast and mold   Schaufus Pottinger, m Green yeast and mold selective, Wort
	Acid-tolerant microorganisms, Lactic-acid bacteria	Orange Serum
	Enterobacteria, E. coli and coliforms	Endo, m FC, Teepol   Tergitol TTC
Water (general quality), mineral water, natural water, waste water	Enterococci, Enterococcus faecalis	Azide   KF Strep
	Pseudomonas aeruginosa	Cetrimide
	Total colony count	Caso, R2A, TGE   Tryptone Glucose Extract
	Yeasts and molds, Candida albicans	Sabouraud
	Acetobacter	Orange Serum (by adding 5-8% ethanol)
Wine	Acid-tolerant microorganisms, Lactic-acid bacteria	Orange Serum, Wallerstein Differential
	Yeasts and molds	m Green yeast and mold   Schaufus Pottinger, Wallerstein Nutrient, Wort

\* These Biosart® 100 Media types are suitable for the determination of the mentioned microorganisms, although the media are not explicit declared in references.

### Easy work flow – reliable results

