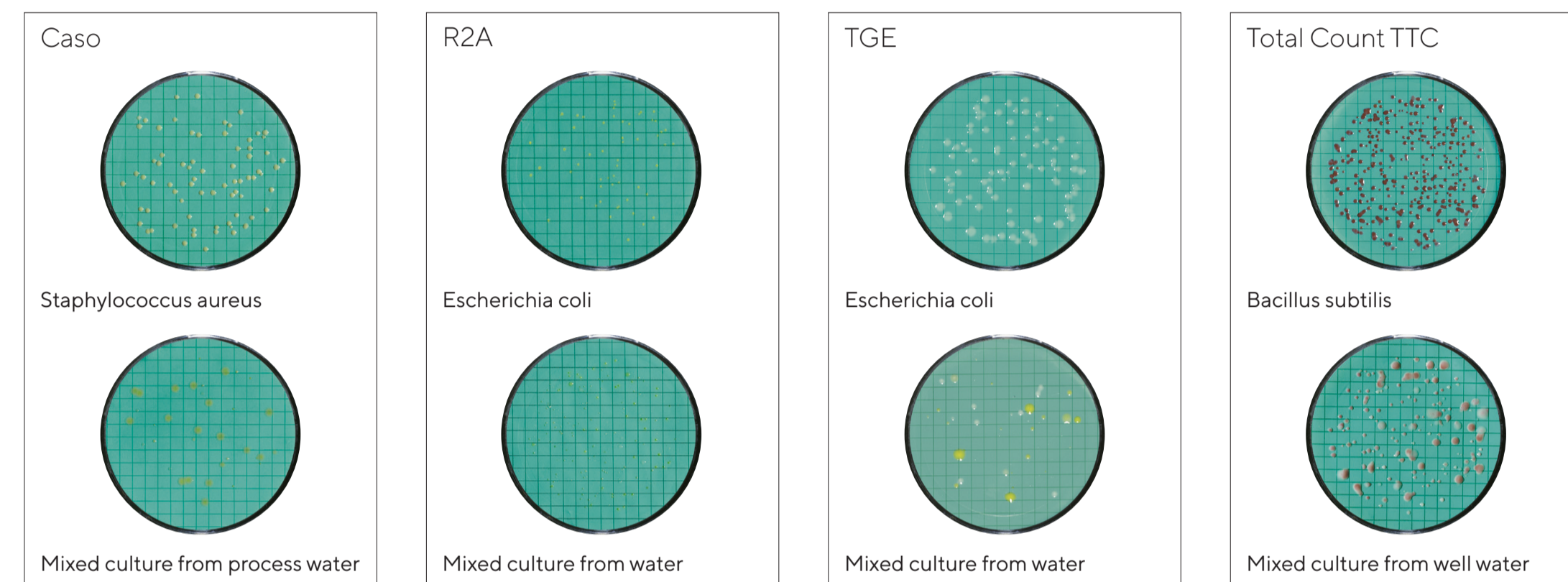
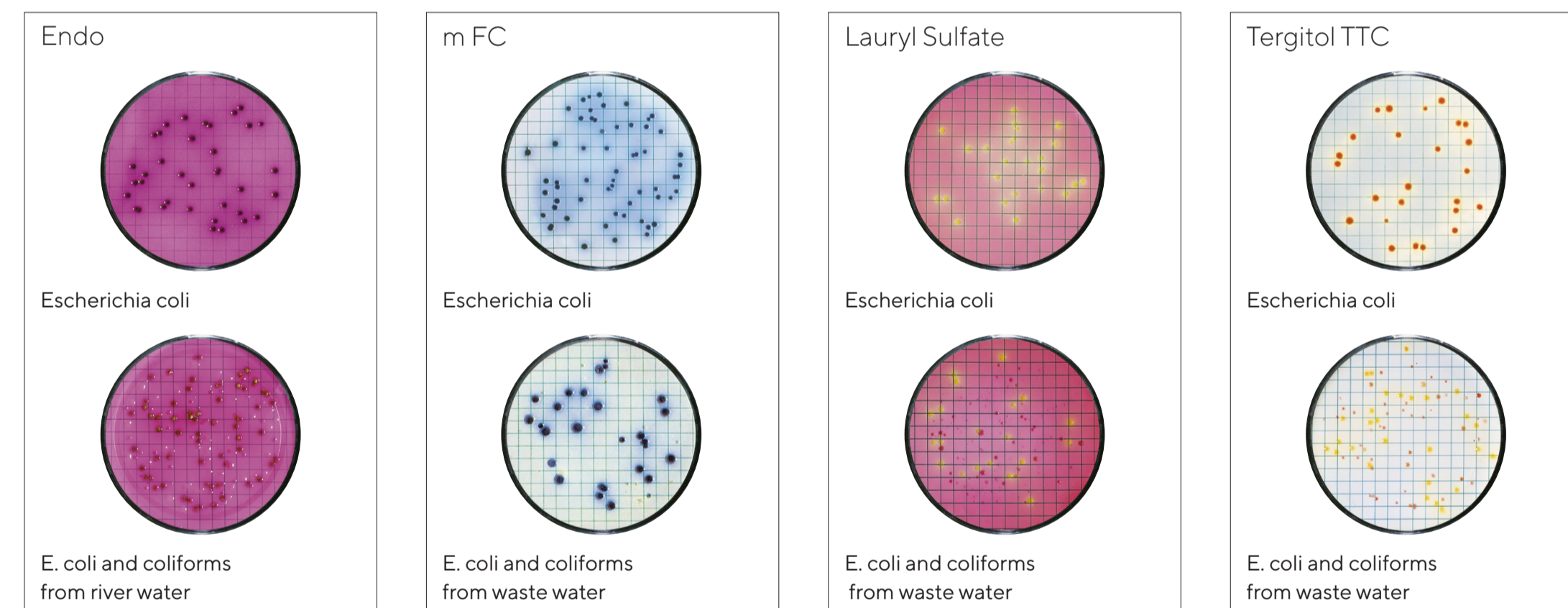


Biosart® 100 Nutrient Media

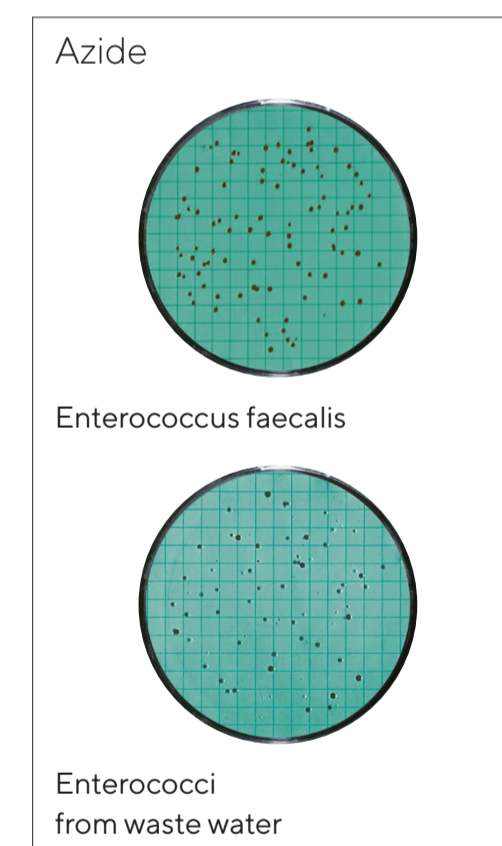
Total colony count



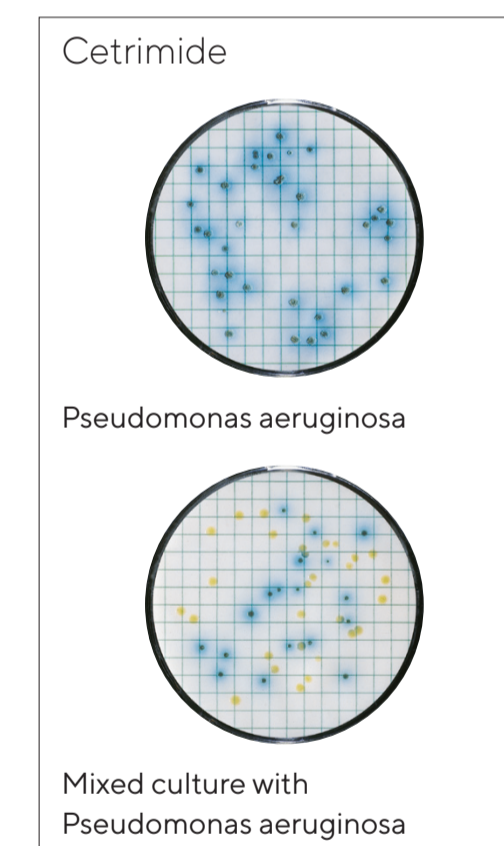
E. coli and coliforms, Enterobacteria



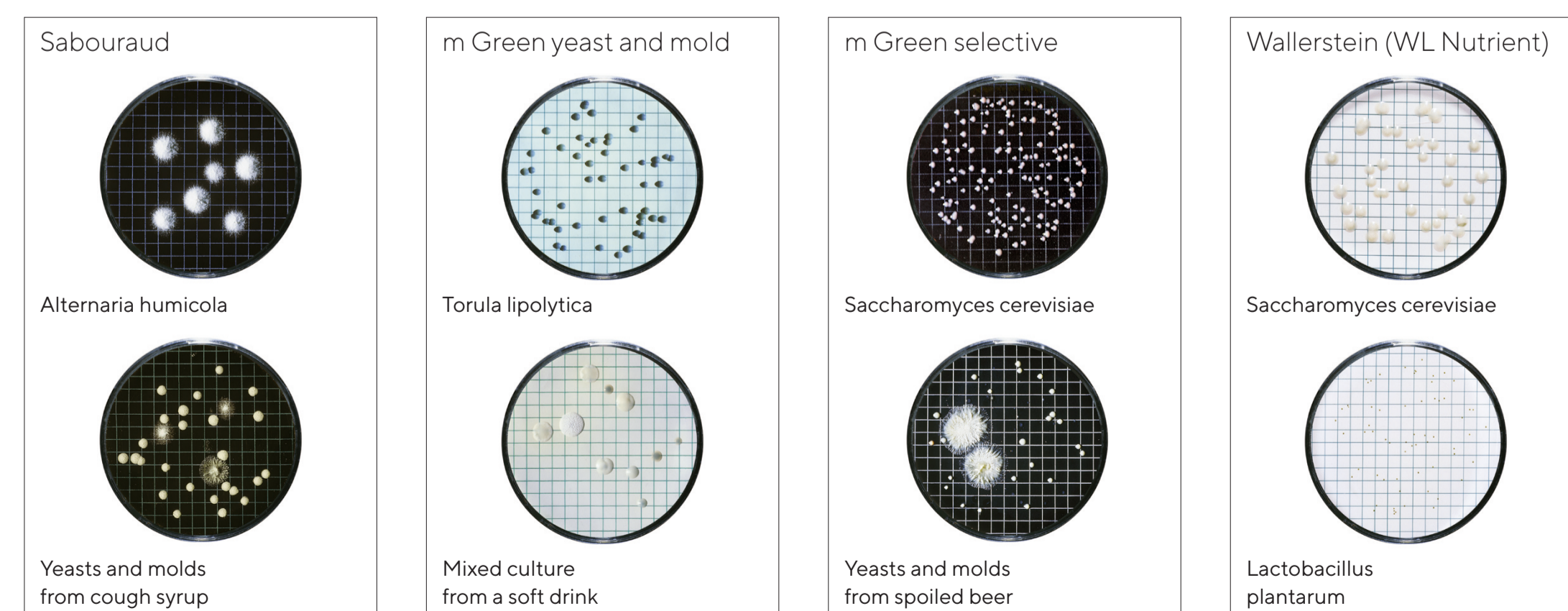
Other faecal bacteria



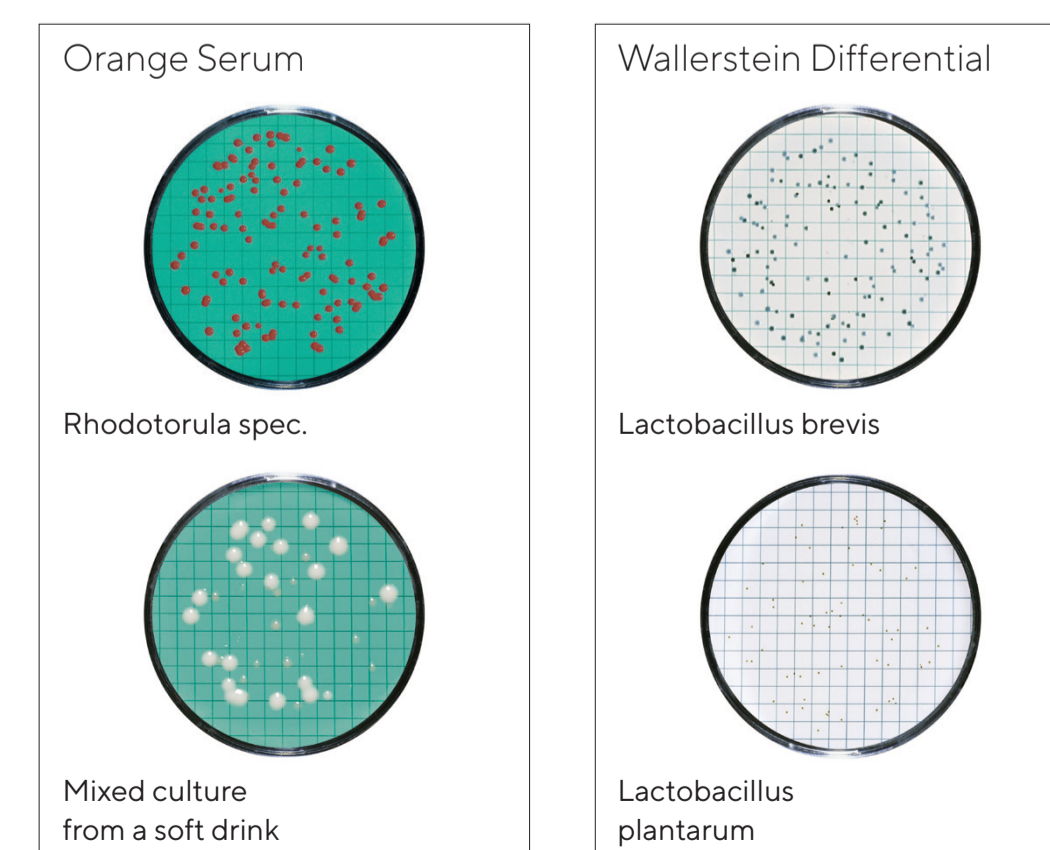
Non-faecal, pathogenic bacteria



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	Lactobacilli and Pediococci and
	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 Lauryl Sulphate, 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
	Enterococci, Enterococcus faecalis	Azide KF Strep
	Pseudomonas aeruginosa	Cetrimide
	Total colony count	Caso, R2A
	Yeasts and molds, Candida albicans	Sabouraud
Soft drinks, concentrates	Acid-tolerant microorganisms, Lactic-acid bacteria	Orange Serum, Wallerstein Differential
	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 Lauryl Sulphate, Tergitol TTC
	Enterococci, Enterococcus faecalis	Azide KF Strep
Water (general quality), mineral water, natural water, waste water	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)
	Acid-tolerant microorganisms, Lactic-acid bacteria	Orange Serum, Wallerstein Differential
Wine	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

- 1 | Pour the sample
- 2 | Apply vacuum and filter the sample
- 3 | Add the Biosart® 100 Nutrient Media
- 4 | Close the outlet
- 5 | Remove the funnel
- 6 | Incubate the petri dish

Remarks
The pictures show typical appearance of the mentioned microorganisms. In particular cases, color and shape of the colonies could vary from the expected habits. Further tests may be necessary to validate the result.

Sartorius Stedim Biotech shall not be liable for consequential and/or incidental damage sustained by any customer from the use of its products.

Biosart® 100 Nutrient Media are subject to continuous product improvement as part of our product development program to align our products with changing application requirements. For current specifications and lot release criteria please visit our homepage under: www.sartorius-stedim.com/BiosartMediaSearch.