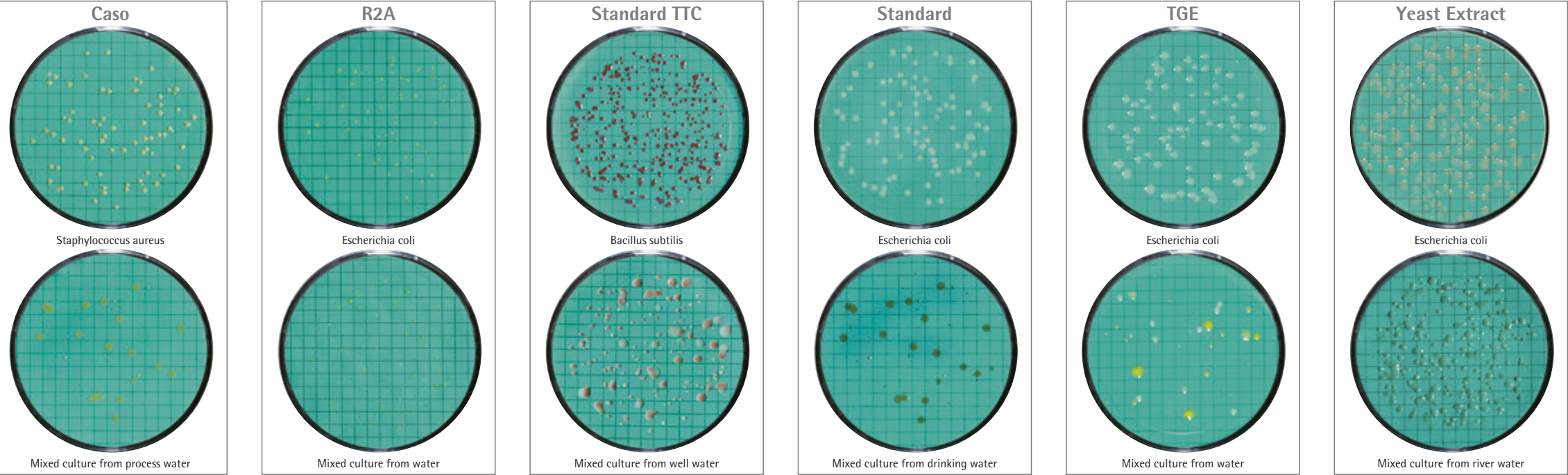


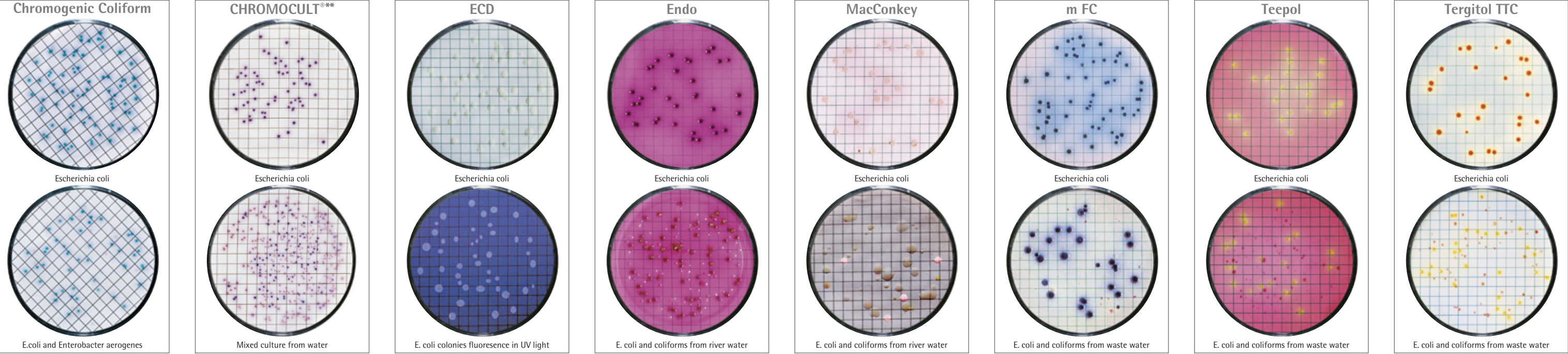
Nutrient Pad Sets



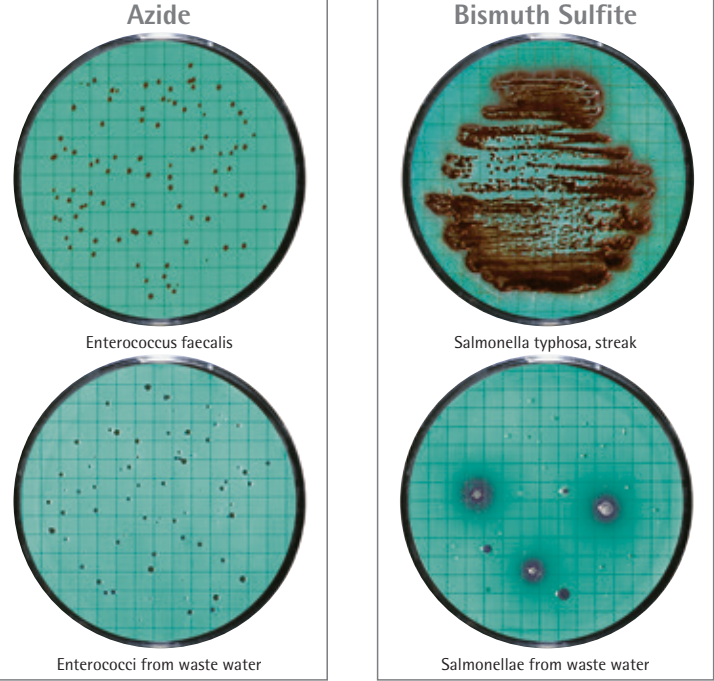
Total colony count



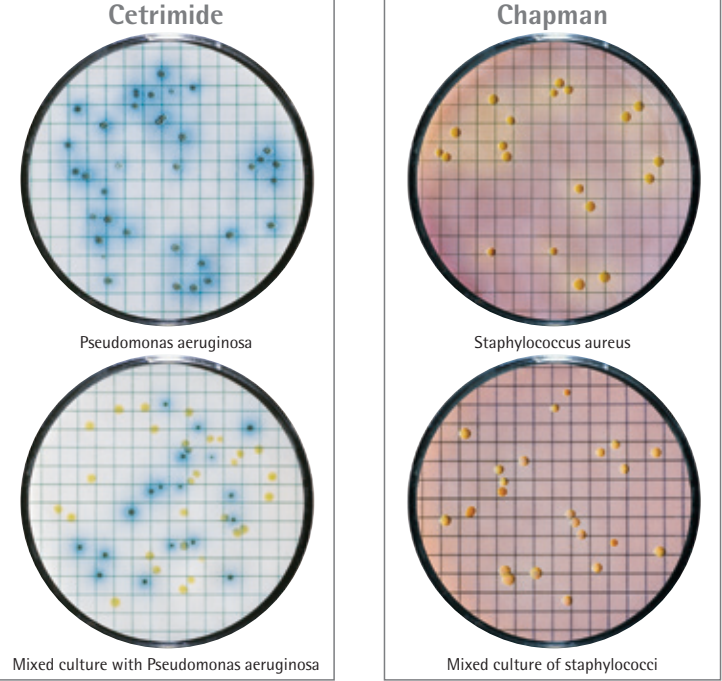
E. coli and coliforms, Enterobacteria



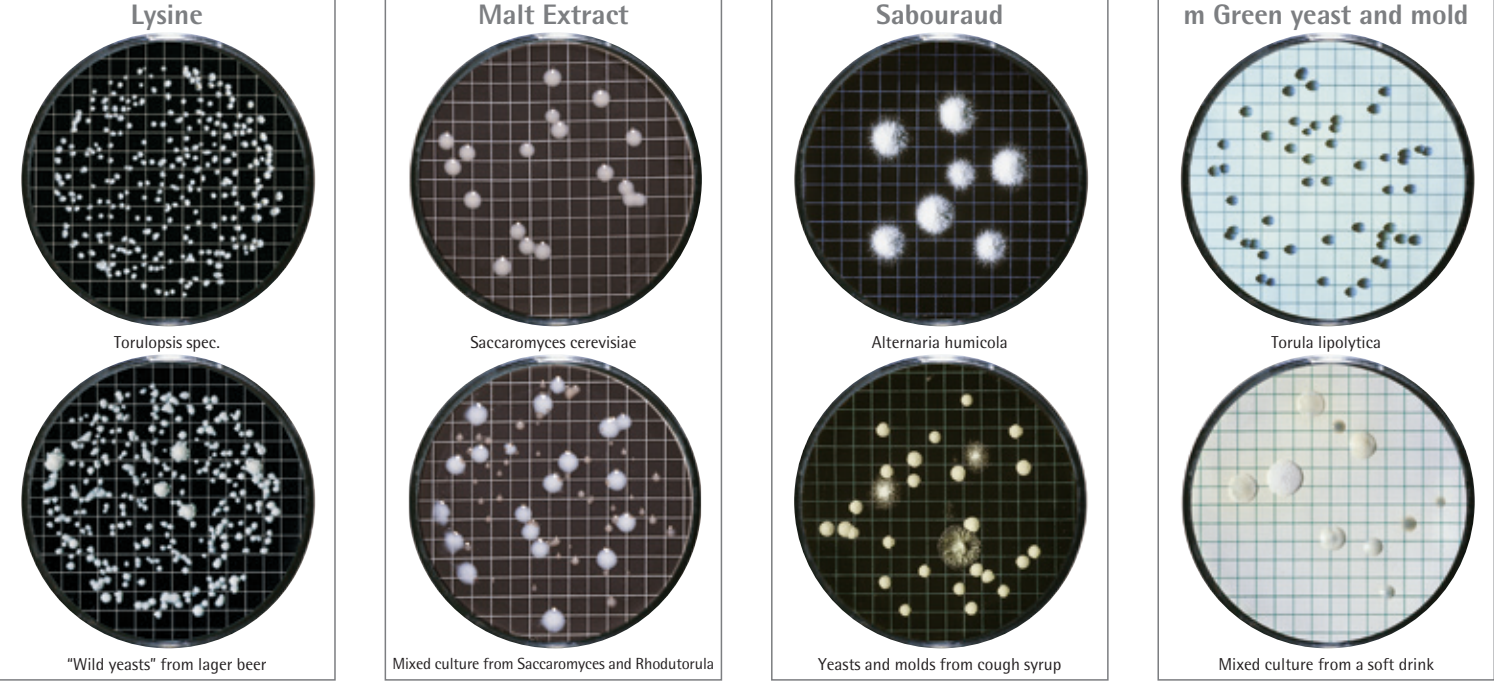
Other faecal bacteria



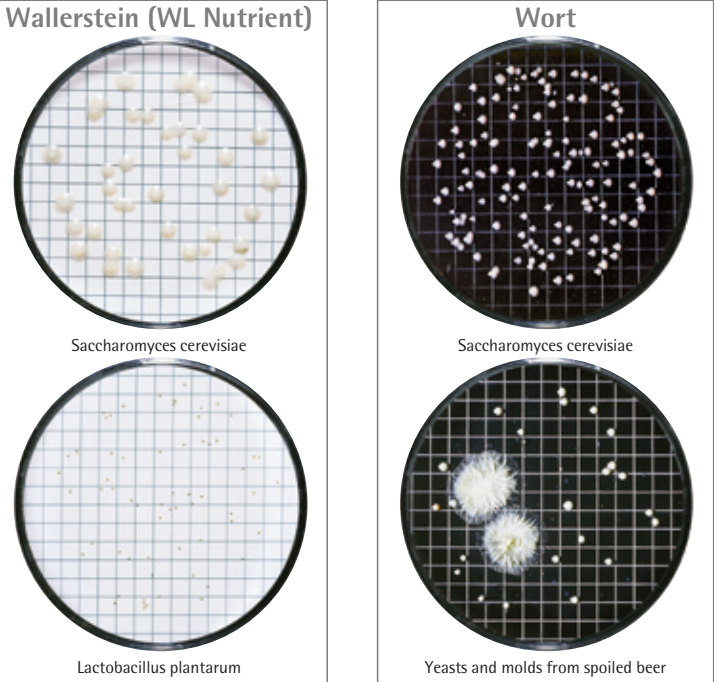
Non-faecal, pathogenic bacteria



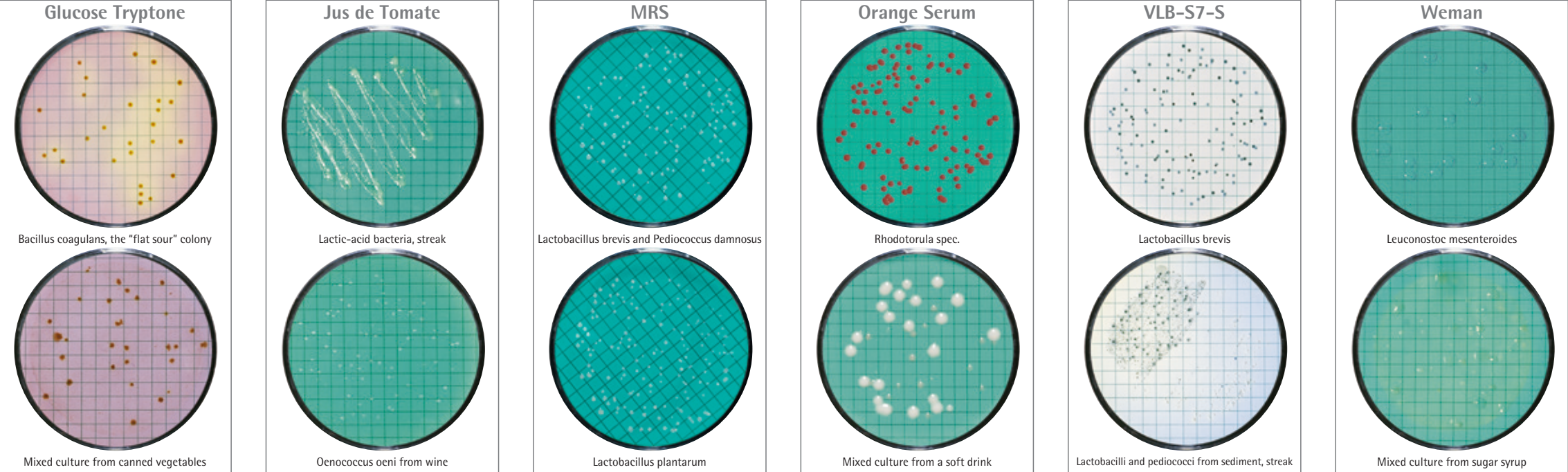
Yeasts and molds



Yeasts and molds



Product Spoiling Microorganisms



Typical Application Examples

Product	Detection and enumeration of...	Nutrient Pad type
Beer	Lactobacilli and Pediococi and other beer spoiling organisms	VLB-S7-S
	Total colony count	Standard, Standard TTC,
	Wild yeasts	Lysine
	Yeasts and molds	Malt Extract*, Wallerstein Nutrient, Wort
Diary products	Lactobacilli	MRS
Foods	Acid-tolerant microorganisms	Orange Serum
	Enterobacteria, E. coli and coliforms	CHROMOCULT***, ECD, Endo, (MacConkey), m FC, Teepol   Lauryl Sulphate, Tergitol TTC
	Enterococci, Enterococcus faecalis	Azide   KF Strep
	Lactobacilli	MRS
	Pseudomonas aeruginosa	Cetrimide
	Salmonellae	Bismuth Sulfite
	Staphylococci, Staphylococcus aureus	Chapman
	Thermophilic spore formers and mesophilic bacteria	Glucose Tryptone
	Total colony count	Caso, Standard, Standard TTC, TGE   Tryptone Glucose Extract
	Yeasts and molds	Malt Extract, Wort
Food and Beverages	Lactobacilli	MRS
Fruit juice	Enterobacteria, E. coli and coliforms	Endo, (MacConkey) Tergitol TTC*
	Oenococcus and other product spoiling organisms	Jus de Tomate   Tomato Juice, Orange Serum
	Lactobacilli	MRS
	Yeasts and molds	Malt Extract, m Green yeast and mold   Schaufus Pottinger, Wallerstein Nutrient, Wort
Milk	E. coli and coliforms	Endo
	Enterococci, Enterococcus faecalis	Azide   KF Strep
	Salmonellae	Bismuth Sulfite
Pharmaceuticals, WFI, raw materials, and cosmetics	Enterobacteria, E. coli	MacConkey
	Enterococci, Enterococcus faecalis	Azide   KF Strep
	Pseudomonas aeruginosa	Cetrimide (cosmetics only)
	Salmonellae	Bismuth Sulfite
	Staphylococci, Staphylococcus aureus	Chapman
	Total colony count	Caso, R2A
Soft drinks, concentrates	Yeasts and molds, Candida albicans	Sabouraud
	Acid-tolerant microorganisms, Lactic-acid bacteria	Orange Serum, VLB-S-7-S
	Enterobacteria, E. coli and coliforms	Endo, MacConkey
	Lactobacilli	MRS
	Mesophilic slime-forming bacteria, Leuconostoc	Weman
	Total colony count	Standard*, Standard TTC*, TGE   Tryptone Glucose Extract
Sugar, sugar products	Yeasts and molds	Malt Extract, m Green yeast and mold   Schaufus Pottinger, Wallerstein Nutrient, Wort
	E. coli and coliforms	Endo
	Mesophilic slime-forming bacteria, Leuconostoc	Weman
	Thermophilic spore formers and mesophilic bacteria	Glucose Tryptone
Water (general quality), mineral water, natural water, waste water	Yeasts and molds	Malt Extract*, Schaufus Pottinger   m Green yeast and mold, Wort*
	Acid-tolerant microorganisms, Lactic-acid bacteria	Orange Serum
	Enterobacteria, E. coli and coliforms	Chromogenic Coliform (according to new DIN EN ISO 9308-1), CHROMOCULT***, ECD, Endo, (MacConkey), m FC, Teepol   Lauryl Sulphate, Tergitol TTC
	Enterococci, Enterococcus faecalis	Azide   KF Strep
	Pseudomonas aeruginosa	Cetrimide
	Salmonellae	Bismuth Sulfite
	Staphylococci, Staphylococcus aureus	Chapman
	Total colony count	Caso, R2A, Standard, Standard TTC, TGE   Tryptone Glucose Extract, Yeast Extract
	Yeasts and molds, Candida albicans	Sabouraud
	Acetobacter	Orange Serum, Wort (both wetted with 5-8% ethanol)
Wine	Acid-tolerant microorganisms, Lactic-acid bacteria	Orange Serum
	Lactobacilli	MRS
	Oenococcus and other wine spoiling microorgan.	Jus de Tomate   Tomato Juice
	Yeasts and molds	Malt Extract, m Green yeast and mold   Schaufus Pottinger, Wallerstein Nutrient, Wort

\* These NPS types are suitable for the determination of the mentioned microorganisms, although the media are not explicit declared in references. The description of the typical results or any pictures show typical appearance of the mentioned microorganisms. In particular cases, color and shape of the colonies could vary from the expected habitus. 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. Nutrient Pad Sets (NPS) 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/NPSearch](http://www.sartorius-stedim.com/NPSearch). \*\* Trade mark owner and manufacturer is Merck KGaA