

Reference # / No. de référence / Referenz-Nr. / Riferimento N. / N° de referencia _____

Date / Date / Datum / Data / Fecha _____

Tech / Tech / Techn. / Tech / Tec _____

Source / Source / Quelle / Origine / Origen _____

Gram Stain / Souche Gram / Gram Färbung / Colorazione di Gram / Tinción de Gram _____

Lancefield Group / Groupe Lancefield / Lancefield-Gruppe / Gruppo di Lancefield / Grupo de Lancefield _____

Reagent / Réactif / Reagenz / Reagente / Reactivo	None / Aucun / Keine / Nessuno / Ninguno										RapID STR Reagent / Réactif RapID STR RapID STR Reagens / Reagente RapID STR / Reactivo RapID STR				None
Positive Reactions Réactions positives Positive Reaktionen Reazioni positive Reacciones positivas	Red or Dark Orange Rouge ou orange foncé Rot oder dunkelorange Rosso o arancione scuro Rojo o naranja oscuro	Black Noir Schwarz Nero Negro	Yellow or Yellow-orange Jaune ou Jaune orangé Gelb oder gelb-orange Giallo o giallo-arancio Amarillo o amarillo-naranja		Yellow, yellow-orange or orange Jaune, jaune orangé ou orange Gelb, gelb-orange oder orange Giallo, giallo-arancio o arancio Amarillo, amarillo-naranja o naranja		Yellow Jaune Gelb Giallo Amarillo				Light Purple or purple Violacé léger ou franc Helles Violett oder Violett Porpora chiaro o Porpora Morado claro o morado		Very Dark Purple Violacé très foncé Sehr dunkles Violett Porpora molto scuro Morado muy oscuro		Beta Beta Beta Beta Beta
Cavity # / No. cavité / Kammer-Nr. / Cavità N. / N° de cavidad	1	2	3	4	5	6	7	8	9	10	7	8	9	10	
Test Code Code du test Testcode Codice esame Código de prueba	ARG	ESC	MNL	SBL	RAF	INU	GAL	GLU	NAG	PO4	TYR	HPR	LYS	PYR	HEM
Value / Valeur / Wert Valore / Valor	1	2	4	1	2	4	1	2	4	1	2	4	1	2	4
Result / Résultat Ergebnis / Risultato Resultado															
Value Total Total des valeurs Gesamtwert Valore totale Valor total															

IDENTIFICATION / IDENTIFICATION / IDENTIFICAZIONE / IDENTIFICACIÓN _____

