

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 \_\_\_\_\_

Reagent / Réactif / Reagenz / Reagente / Reactivo	None / Aucun / Keine, Nessuno / Ninguno						RapID Yeast Plus Reagent A / Réactif A RapID Yeast Plus / RapIDYeast Plus Reagens A / RapID Yeast Plus Reagent A / Reactivo RapID Yeast Plus A								None / Aucun / Keine, Nessuno / Ninguno		RapID Yeast Plus Reagent B / Réactif B RapID Yeast Plus / RapID Yeast Plus Reagens B / RapID Yeast Plus Reagent B / Reactivo RapID Yeast Plus B			
Positive Reactions Réactions positives Positive Reaktionen Reazioni positive Reacciones positivas	Yellow Jaune Gelb Giallo Amarillo						Yellow Jaune Gelb Giallo Amarillo								Red or Dark red-orange Rouge ou rouge orangé foncé Rot oder dunkles Rotorange Rosso o rosso-arancione scuro Rojo o rojo-naranja oscuro		Purple, red, or dark pink Violacé, rouge ou rose foncé Purpur, rot oder dunkelrosa Porpora, rosso o rosa scuro Morado, rojo o rosa oscuro			
Cavity #, No. cavité / Kammer-Nr. / Cavità N. / N° de cavidad	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18		
Test Code Code du test Testcode Codice esame Código de prueba	GLU	MAL	SUC	TRE	RAF	LIP	NAGA	aGLU	BGLU	ONPG	aGAL	BFUC	PHS	PCHO	URE	PRO	HIST	LGY		
Value / Valeur / Wert Valore / Valor	1	2	4	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 / IDENTIFIZIERUNG / IDENTIFICAZIONE / IDENTIFICACIÓN \_\_\_\_\_



Microcode \_\_\_\_\_