



RapID™ NF Plus

Report Form

Reference # / No. de référence / Referenz-Nr. _____

Date / Date / Datum _____

Tech / Tech / Techn. _____

Source / Source / Quelle _____

MacConkey / MacConkey / MacConkey _____

Pigment / Pigment / Pigment _____

Reagent / Réactif / Reagenz	None, Aucun, Keine										RapID NF Plus Reagent / Réactif RapID NF Plus / RapID NF Plus Reagens					RapID Spot Indole	Nitrate A	
	Red Rouge Rot	Yellow, gold, or yellow-orange / Jaune, or ou jaune orangé / Gelb, goldgelb oder Gelborange	Yellow, gold, yellow-orange, or light orange / Jaune, doré, jaune-orange ou orange clair / Gelb, Goldgelb, Gelborange oder Hellorange	Yellow Jaune Gelb					Red Rouge Rot	Yellow Jaune Gelb	Purple, violet, red, dark orange or dark pink / Violacé, violet, rouge, orange soutenu ou rose soutenu / Purpur, Violet, Rot, Dunkelorange oder Dunkelrosa	Brown or Black / Marron ou noir Braun oder Schwarz				Red or orange / Rouge ou orange / Rot oder Orange		
Cavity #, No. cavité / Kammer-Nr.	1	2	3	4	5	6	7	8	9	10	4	5	6	7	8	9	10	
Test Code / Code du test / Testcode	ADH	TRD	EST	PHS	NAG	aGLU	BGLU	ONPG	URE	GLU	PRO	PYR	GGT	TRY	BANA	IND	NO3	OXI
Value / Valeur / Wert	1	2	4	1	2	4	1	2	4	1	2	4	1	2	4	1	2	4
Result / Résultat Ergebnis																		
Value Total Total des valeurs Gesamtwert																		

 GLU (color / couleur / Farbe)

IDENTIFICATION / IDENTIFICATION / IDENTIFIZIERUNG _____



Microcode _____

REMEL Inc

800-255-6730

Printed in USA 9/03

Riferimento N. / N° de referencia _____

Data / Fecha _____

Tech / Tec _____

Origine / Origen _____

MacConkey / MacConkey _____

Pigmento / Pigmento _____

Reagente Reactivo	Nessuno / Ninguno										RapID NF PlusReagente / Reactivo RapID NF Plus				RapID Spot Indole	Nitrate A		
	Rosso / Rojo	Giallo, oro o giallo- arancio / Amarillo, dorado o amarillo-naranja	Giallo, oro, giallo- arancio o arancio chiaro / Amarillo, dorado amarillo- naranja o naranja claro	Giallo / Amarillo				Rosso / Rojo	Giallo / Amarillo	Porpora, violetto, rosso, arancio scuro o rosa scuro / Morado, violeta, rojo, naranja oscuro o rosa oscuro				Marrone o nero / Marrón o negro	Rosso o arancione / Rojo o naranja			
Cavità N. N° de cavidad	1	2	3	4	5	6	7	8	9	10	4	5	6	7	8	9	10	
Codice esame Código de prueba	ADH	TRD	EST	PHS	NAG	aGLU	BGLU	ONPG	URE	GLU	PRO	PYR	GGT	TRY	BANA	IND	NO3	OXI
Valore Valor	1	2	4	1	2	4	1	2	4	1	2	4	1	2	4	1	2	4
Risultato Resultado																		
Valore totale Valor total																		

GLU (colore / color)

Identificazione / Identificación _____

