

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

Date / Date / Datum: _____

Tech / Tech / Techn: _____

Source / Source / Quelle: _____

Reagent / Réactif / Reagenz	None / Aucun / Keine											STAPH PLUS REAGENT / RÉACTIF RapID STAPH PLUS / RapID STAPH PLUS REAGENS					Nitrate A&B	
	Purple, blue, bluish-gray or greenish-gray / Violet, bleu, gris-bleu ou gris-vert / Purpurrot, Blau, Blaugrau oder Grüngrau			Yellow or yellow-orange / Jaune ou jaune-orangé / Gelb oder Gelborange			Yellow or pale yellow / Jaune ou jaune pale / Gelb oder Blassgelb		Any yellow or light yellow / Tous les tons de jaune ou jaune clair / Jegliches Gelb oder Hellgelb			Red or red-orange / Rouge ou rouge orange / Rot oder Rot-orange		Purple or dark red / Violet ou rouge foncé / Purpurrot oder Dunkelrot				Purple, red, light red or dark pink / Violet, rouge, rouge clair ou rose foncé / Purpur, Rot, Hellrot oder Dunkelrosa
Cavity # / No. cavité / Kammer-Nr.	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18
Test Code Code du test Testcode	ADH	ODC	LIP	SUC	MANO	PO4	αGLU	βGLU	ONPG	GUR	NAGA	UREA	PYR	ARG	ALA	LEU	LGLY	NIT
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 (Microcode)																		

IDENTIFICATION / IDENTIFICATION / IDENTIFIZIERUNG



Riferimento N. / N° de referencia: _____

Data / Fecha: _____

Tech / Tec: _____

Origine / Origen: _____

Reagente Reactivo	Nessuno / Ninguno												RapID STAPH PLUS REAGENT / REACTIVO RapID STAPH PLUS				Nitrate A & B	
Reazioni positive Reacciones positivas	Porpora, blu, grigio-blu o grigio-verde / Morado, azul, gris azulado o gris verdoso		Giallo o giallo arancione / Amarillo o amarillo-naranja			Giallo o giallo chiaro / Amarillo o amarillo pálido	Qualsiasi giallo o giallo chiaro / Cualquier amarillo o amarillo claro					Rosso o rosso arancione/ Rojo o rojo- naranja	Porpora o rosso scuro / Morado o rojo oscuro	Porpora, rosso, rosso chiaro o rosa scuro / Morado, rojo, rojo claro o rosa oscuro			Rosso ciliegia o rosa scuro / Rojo cereza o rosa oscuro	
Cavità N. / N° de cavidad	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18
Codice esame Código de prueba	ADH	ODC	LIP	SUC	MANO	PO4	αGLU	βGLU	ONPG	GUR	NAGA	UREA	PYR	ARG	ALA	LEU	LGLY	NIT
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 (Microcode)																		

IDENTIFICAZIONE / IDENTIFICACIÓN

