



RapID™ ANAII

Report Form

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/Morphology, Souche Gram/Morphologie, Gram-Färbung/Morphologie, Colorazione di Gram/Morfologia, Tinción de Gram/Morfologia _____

Reagent / Réactif / Reagenz / Reagente / Reactivo	None / Aucun / Keine / Nessuno / Ninguno										RapID ANA II Reagent / Réactif RapID ANA II / RapID ANA II Reagens / RapID ANA II Reagent / Reactivo RapID ANA II						RapID Spot Indole		
Positive Reactions Réactions positives Positive Reaktionen Reazioni positive Reacciones positivas	Red or purple Rouge ou violacé Rot oder purpur Rosso o porpora Rojo o púrpura		Medium or bright yellow jaune moyen ou vif mittel-oder hellgelb Giallo di intensità media o brillante Amarillo medio o brillante										Purple, violet, red, or dark pink violacé, violet, rouge ou rose soutenu purpur, violett, rot oder dunkelrosa Porpora, violetto, rosso o rosa scuro Púrpura, violeta, rojo o rosa oscuro						Blue or blue-green bleu ou bleu-vert blau oder blaugrün Blu o blu-verde Azul o verde azulado
Cavity # / No. cavité / Kammer-Nr. / Cavità N. / N° de cavidad	1	2	3	4	5	6	7	8	9	10	3	4	5	6	7	8	9	10	
Test Code Code du test Testcode Codice esame Código de prueba	URE	BLTS	aARA	ONPG	aGLU	BGLU	aGAL	aFUC	NAG	PO4	LGY	GLY	PRO	PAL	ARG	SER	PYR	IND	
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 _____