

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 \_\_\_\_\_

Oxidase / Oxydase / Oxidase / Ossidasi / Oxidasa \_\_\_\_\_

Growth Media / Milieu de croissance / Nährböden / Terreno di crescita / Medios de crecimiento \_\_\_\_\_

Reagent/ Réactif/ Reagenz/ Reagente/ Reactivo	None / Aucun / Keine / Nessuno / Ninguno						IDS RapID SS/u Reagent / Réactif RapID SS/u / RapID SS/u Reagens /RapID SS/u Reagent / Reactivo RapID SS/u	Spot indole				
Positive Reactions Réactions positives Positive Reaktionen Reazioni positive Reacciones positivas	Medium or bright yellow Jaune moyen ou vif Mittleres oder starkes Gelb Giallo di intensità media o brillante Amarillo medio o brillante						Dark red, red, red-orange or orange Rouge foncé, rouge, rouge orange ou orange Dunkelrot, Rot, Rot-Orange oder orange Rosso scuro, rosso, rosso-arancio o arancio Rojo oscuro, rojo, rojo-naranja o naranja	Purple, violet, red, or dark pink Violacé, violet, rouge ou rose soutenu Purpur, Violett, Rot oder Dunkelrosa Porpora, violetto, rosso o rosa scuro Morado, violeta, rojo o rosa oscuro	Black, brown, or tan Noir, marron ou légère coloration Schwarz, Braun oder dunkel getönt Nero, marrone o beige Negro, marrón o tostado			
Cavity # / No. cavité / Kammer-Nr. / Cavità N. / N° de cavidad	1	2	3	4	5	6	7	8	9	10	7	
Test Code / Code du test / Testcode / Codice esame Código de prueba	GMS	ONPG	G1	G2	G3	PHS	URE	A1	A2	A3	IND	OXI
Value / Valeur / Wert Valore / Valor	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 \_\_\_\_\_

