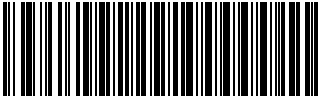




Certificate of Analysis

Product name	chemagic DNA Saliva 600 Kit H96	 Sa l 600 - 01 x x x x x
Article No.	CMG-1037	
Lot No.	Sal600-01xxxxx	
Manufacturing date	yyyy-mm-dd	
Storage	At +2 to +25 °C	
Stability	yyyy-mm-dd (expiry date of kit as supplied)	

Kit Components

	Lot no.	Amount	pH	Appearance	Quality for DNA isolation
Magnetic Beads	Sal600-xxxxxxx	160 mL	n/a	Brown suspension	Accord. spec.
Binding Buffer 2	Sal600-xxxxxxx	1.5 L	5.2 – 5.9	Clear solution	Accord. spec.
Wash Buffer 3	Sal600-xxxxxxx	1.4 L	5.0 – 5.6	Clear solution	Accord. spec.
Wash Buffer 4	Sal600-xxxxxxx	1.4 L	5.0 – 5.6	Clear solution	Accord. spec.
Wash Buffer 5	Sal600-xxxxxxx	1.4 L	n/a	Clear solution	Accord. spec.
Wash Buffer 6	Sal600-xxxxxxx	530 mL	6.0 – 6.7	Clear solution	Accord. spec.
Elution Buffer 7	Sal600-xxxxxxx	320 mL	7.8 – 8.4	Clear solution	Accord. spec.
chemagic Deep Well Plate 2 mL	-	62	-	-	-
chemagic Low Well Plates	-	10	-	-	-
chemagic Tips 96 Racks	-	10	-	-	-

This Certificate of Analysis has been produced electronically and is valid without a signature. For questions regarding this Certificate of Analysis please contact support.chemagen@revvity.com.