


Certificate of Analysis

Product name	chemagic DNA Saliva 4k Kit H24	 Sa l 4 k - 0 1 x x x x x
Article No.	CMG-1081	
Lot No.	Sal4k-01xxxxx	
Manufacturing date	yyyy-mm-dd	
Storage	At +2 to +25 °C (Proteinase K after dissolving: store at +2 to +8 °C)	
Stability	yyyy-mm-dd (expiry date of kit as supplied) (Proteinase K after dissolving: stable for 28 days at +2 to +8 °C)	

Kit Components

	Lot no.	Amount	pH	Appearance	Quality for DNA isolation
Magnetic Beads	Sal4k-01xxxxx	130 mL	n/a	Brown suspension	Accord. spec.
Binding Buffer 2	Sal4k-01xxxxx	2.2 L	5.2 – 5.9	Clear solution	Accord. spec.
Wash Buffer 3	Sal4k-01xxxxx	1.6 L	5.0 – 5.6	Clear solution	Accord. spec.
Wash Buffer 4	Sal4k-01xxxxx	1.6 L	5.0 – 5.6	Clear solution	Accord. spec.
Wash Buffer 5	Sal4k-01xxxxx	2 x 1.5 L	n/a	Clear solution	Accord. spec.
Wash Buffer 6	Sal4k-01xxxxx	2 x 1.5 L	6.0 – 6.7	Clear solution	Accord. spec.
Elution Buffer 7	Sal4k-01xxxxx	200 mL	7.8 – 8.4	Clear solution	Accord. spec.
chemagic Tips XL	-	250	-	-	-
chemagic Deep Well Plate 24 XL	-	50	-	-	-
chemagic Elution Tube 13 mL	-	288	-	-	-

* = after dissolving the lyophilisate

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.