




Certificate of Analysis

Product name	chemagic Viral DNA/RNA 300 Kit H96	 V300 - 01xxxxx
Article No.	CMG-1033-S	
Lot No.	V300-01xxxxx	
Manufacturing date	yyyy-mm-dd	
Storage	At +2 to +25 °C (Proteinase K and Poly(A) RNA after dissolving: store at +2 to +8 °C)	
Stability	yyyy-mm-dd (expiry date of kit as supplied) (Proteinase K and Poly(A) RNA after dissolving: stable for 28 days at +2 to +8 °C)	

Kit Components

	Lot no.	Amount	pH	Appearance	Quality for DNA isolation
Magnetic Beads	V300-01XXXXX	150 mL	n/a	Brown suspension	Accord. spec.
Lysis Buffer 1	V300-01XXXXX	320 mL	n/a	Clear solution	Accord. spec.
Binding Buffer 2	V300-01XXXXX	1.3 L	5.2 – 6.1	Clear solution	Accord. spec.
Wash Buffer 3	V300-01XXXXX	700 mL	4.8 – 5.6	Clear solution	Accord. spec.
Wash Buffer 4	V300-01XXXXX	700 mL	n/a	Clear solution	Accord. spec.
Wash Buffer 5	V300-01XXXXX	700 mL	6.0 – 6.7	Clear solution	Accord. spec.
Elution Buffer 6	V300-01XXXXX	200 mL	7.8 – 8.4	Clear solution	Accord. spec.
Proteinase K	V300-01XXXXX	11 mL*	n/a	White powder	Accord. spec.
Poly(A) RNA	V300-01XXXXX	10 x 440 µL*	n/a	White powder	Accord. spec.
Poly(A) RNA Buffer	V300-01XXXXX	10 x 500 µL	n/a	Clear solution	Accord. spec.
chemagic Deep Well Plate 2 mL	-	50	-		
chemagic Low Well Plates	-	10	-		
chemagic Tips 96 Racks	-	10	-		

* = 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.