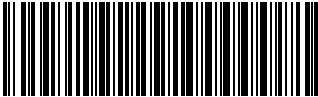




## Certificate of Analysis

Product name	chemagic Viral NA /gDNA 200 Kit H96	 V/B200-01xxxxx
Article No.	CMG-1049	
Lot No.	V/B200-01xxxxx	
Manufacturing date	yyyy-mm-dd	
Storage	At +2 to +25 °C (Proteinase K and Poly(A) after dissolving: store at +2 to +8 °C)	
Stability	yyyy-mm-dd (expiry date of kit as supplied) (Proteinase K and Poly(A) after dissolving: stable for 28 days at +2 to +8 °C)	

### Kit Components

	Lot no.	Amount	pH	Appearance	Quality for DNA isolation
Magnetic Beads	V/B200-01xxxxx	150 mL	n/a	Brown suspension	Accord. spec.
Lysis Buffer 1 (blood)	V/B200-01xxxxx	700 mL	6.9 – 7.4	Clear solution	Accord. spec.
Lysis Buffer 1 (plasma)	V/B200-01xxxxx	700 mL	n/a	Clear solution	Accord. spec.
Binding Buffer 2 (blood)	V/B200-01xxxxx	1.2 L	5.2 – 5.9	Clear solution	Accord. spec.
Binding Buffer 2 (plasma)	V/B200-01xxxxx	1.2 L	5.2 – 6.1	Clear solution	Accord. spec.
Wash Buffer 3	V/B200-01xxxxx	750 mL	5.0 – 5.6	Clear solution	Accord. spec.
Wash Buffer 4	V/B200-01xxxxx	750 mL	5.0 – 5.6	Clear solution	Accord. spec.
Wash Buffer 5	V/B200-01xxxxx	750 mL	n/a	Clear solution	Accord. spec.
Elution Buffer 6	V/B200-01xxxxx	220 mL	7.8 – 8.4	Clear solution	Accord. spec.
Proteinase K	V/B200-01xxxxx	9 x 1.25 mL*	n/a	White powder	Accord. spec.
Poly(A) RNA	V/B200-01xxxxx	10 x 440 µL*	n/a	White powder	Accord. spec.
Poly(A) RNA Buffer	V/B200-01xxxxx	10 x 500 µL	n/a	Clear solution	Accord. spec.
chemagic Deep Well Plate 2 mL	-	62	-	-	-
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](mailto:support.chemagen@revvity.com).