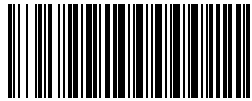


## Certificate of Analysis

Product name	chemagic cfDNA 5k Kit H24	 cf - 01xxxxxx
Article No.	CMG-1304	
Lot No.	cf-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	cf-01xxxxx	50 mL	n/a	Brown suspension	Accord. spec.
Lysis Buffer 1	cf-01xxxxx	70 mL	n/a	Clear solution	Accord. spec.
Binding Buffer 2	cf-01xxxxx	2.0 L	n/a	Clear solution	Accord. spec.
Wash Buffer 3	cf-01xxxxx	800 mL	n/a	Clear solution	Accord. spec.
Wash Buffer 4	cf-01xxxxx	800 mL	n/a	Clear solution	Accord. spec.
Elution Buffer 5	cf-01xxxxx	100 mL	7.8 – 8.4	Clear solution	Accord. spec.
Poly(A) RNA	cf-01xxxxx	2 x 440 µL*	n/a	White powder	Accord. spec.
Poly(A) RNA Buffer	cf-01xxxxx	2 x 500 µL	n/a	Clear solution	Accord. spec.
Proteinase K	cf-01xxxxx	32 x 1.25 mL*	n/a	White powder	Accord. spec.
chemagic Tips XL	-	250	-	-	-
chemagic Deep Well Plate 24 XL	-	12	-	-	-
chemagic Elution Tube 4 mL	-	250	-	-	-
chemagic Elution Tube 13 mL	-	1008	-	-	-

\* = 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).