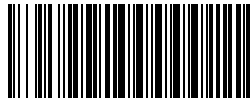


Certificate of Analysis

Product name	chemagic cfDNA 5k Kit	 cf - 01xxxxxx
Article No.	CMG-134	
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	10 mL	n/a	Brown suspension	Accord. spec.
Lysis Buffer 1	cf-01xxxxx	15 mL	n/a	Clear solution	Accord. spec.
Binding Buffer 2	cf-01xxxxx	2x 150 mL	5.3 – 5.9	Clear solution	Accord. spec.
Wash Buffer 3	cf-01xxxxx	200 mL	6.3 – 6.9	Clear solution	Accord. spec.
Wash Buffer 4	cf-01xxxxx	200 mL	n/a	Clear solution	Accord. spec.
Elution Buffer 5	cf-01xxxxx	10 mL	7.8 – 8.4	Clear solution	Accord. spec.
Poly(A) RNA	cf-01xxxxx	1 x 440 µL*	n/a	White powder	Accord. spec.
Poly(A) RNA Buffer	cf-01xxxxx	1 x 500 µL	n/a	Clear solution	Accord. spec.
Proteinase K	cf-01xxxxx	5 x 1.25 mL*	n/a	White powder	Accord. spec.

* = 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.