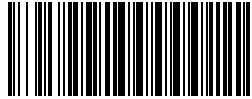




Certificate of Analysis

Product name	chemagic cfNA 1.5k Kit H96	 cf - 01xxxxxx
Article No.	CMG-1396	
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	95 mL	n/a	Brown suspension	Accord. spec.
Lysis Buffer 1	cf-01xxxxx	80 mL	n/a	Clear solution	Accord. spec.
Binding Buffer 2	cf-01xxxxx	2.5 L	n/a	Clear solution	Accord. spec.
Wash Buffer 3	cf-01xxxxx	1.5 L	n/a	Clear solution	Accord. spec.
Wash Buffer 4	cf-01xxxxx	1.5 L	n/a	Clear solution	Accord. spec.
Deionized Water	cf-01xxxxx	1.0 L	6.0 – 6.7	Clear solution	Accord. spec.
Elution Buffer 5	cf-01xxxxx	100 mL	7.8 – 8.4	Clear solution	Accord. spec.
Poly(A) RNA	cf-01xxxxx	10 x 440 µL*	n/a	White powder	Accord. spec.
Poly(A) RNA Buffer	cf-01xxxxx	10 x 500 µL	n/a	Clear solution	Accord. spec.
Proteinase K	cf-01xxxxx	38 x 2.5 mL*	n/a	White powder	Accord. spec.
chemagic Tips 96 Racks	-	10	-	-	-
chemagic Deep Well Plate 2 mL	-	70	-	-	-

* = 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.