


Certificate of Analysis

Product name	chemagic DNA Blood 4k Kit H24	 B4k - 01xxxxx
Article No.	CMG-1074	
Lot No.	B4k-01xxxxx	
Manufacturing date	yyyy-mm-dd	
Storage	At +2 to +25 °C (Proteinase K after dissolving: store at +2 to +8 °C)	
Stability	yyyy-mm-dd (expiry date of kit as supplied) (Proteinase K after dissolving: stable for 28 days at +2 to +8 °C)	

Kit Components

	Lot no.	Amount	pH	Appearance	Quality for DNA isolation
Magnetic Beads	B4k-01xxxxx	130 mL	n/a	Brown suspension	Accord. spec.
Lysis Buffer 1	B4k-01xxxxx	1.1 L	6.7 – 7.2	Clear solution, slightly yellow	Accord. spec.
Binding Buffer 2	B4k-01xxxxx	2 x 1.6 L	5.2 – 5.9	Clear solution	Accord. spec.
Wash Buffer 3	B4k-01xxxxx	1.6 L	5.0 – 5.6	Clear solution	Accord. spec.
Wash Buffer 4	B4k-01xxxxx	1.6 L	5.0 – 5.6	Clear solution	Accord. spec.
Wash Buffer 5	B4k-01xxxxx	1.6 L	n/a	Clear solution	Accord. spec.
Wash Buffer 6	B4k-01xxxxx	1.6 L	6.0 – 6.7	Clear solution	Accord. spec.
Elution Buffer 7	B4k-01xxxxx	200 mL	7.8 – 8.4	Clear solution	Accord. spec.
Proteinase K	B4k-01xxxxx	4 x 1.25 mL*	n/a	White powder	Accord. spec.
chemagic Deep Well Plate 24 XL	-	50	-	-	-
chemagic Elution Tube 13 mL	-	288	-	-	-
chemagic Tips XL	-	250	-	-	-

* = 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.