




Certificate of Analysis

Product name	chemagic DNA Blood 400 Kit H96	 B400 - 01xxxxx
Article No.	CMG-1091	
Lot No.	B400-01xxxxx	
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	B400-01XXXXX	160 mL	n/a	Brown suspension	Accord. spec.
Lysis Buffer 1	B400-01XXXXX	750 mL	6.7 – 7.2	Clear solution	Accord. spec.
Binding Buffer 2	B400-01XXXXX	1.6 L	5.2 – 5.9	Clear solution	Accord. spec.
Wash Buffer 3	B400-01XXXXX	1.4 L	5.0 – 5.6	Clear solution	Accord. spec.
Wash Buffer 4	B400-01XXXXX	1.4 L	5.0 – 5.6	Clear solution	Accord. spec.
Wash Buffer 5	B400-01XXXXX	1.4 L	n/a	Clear solution	Accord. spec.
Wash Buffer 6	B400-01XXXXX	1.4 L	6.0 – 6.7	Clear solution	Accord. spec.
Elution Buffer 7	B400-01XXXXX	2 x 160 mL	7.8 – 8.4	Clear solution	Accord. spec.
Proteinase K	B400-01XXXXX	9 x 1.25 mL*	n/a	White powder	Accord. spec.
chemagic Tips 96 Racks	-	10	-	-	-
chemagic Deep Well Plate 2 mL	-	62	-	-	-
chemagic Low Well Plates	-	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.