


Certificate of Analysis

Product name	chemagic RNA Tissue 10mg Kit H96	 T i R N A - 0 1 x x x x x
Article No.	CMG-1212	
Lot No.	TiRNA-01xxxxx	
Manufacturing date	yyyy-mm-dd	
Storage	At +2 to +25 °C (except DNase I at +2 to +8 °C) (DNase I after dissolving: store at -15 to -20 °C)	
Stability	yyyy-mm-dd (expiry date of kit as supplied)	

Kit Components

	Lot no.	Amount	pH	Appearance	Quality for RNA isolation
Magnetic Beads	TiRNA-01xxxxx	85 mL	n/a	Brown suspension	Accord. spec.
Lysis Buffer 1	TiRNA-01xxxxx	640 mL	n/a	Clear solution	Accord. spec.
Binding Buffer 2	TiRNA-01xxxxx	1.9 L	5.2 – 6.1	Clear solution	Accord. spec.
Wash Buffer 3	TiRNA-01xxxxx	2.2 L	4.4 – 5.2	Clear solution	Accord. spec.
Wash Buffer 4	TiRNA-01xxxxx	1.6 L	n/a	Clear solution	Accord. spec.
DNase I Buffer 5	TiRNA-01xxxxx	65 mL	7.8 – 8.1	Clear solution	Accord. spec.
Elution Buffer 6	TiRNA-01xxxxx	220 mL	7.8 – 8.4	Clear solution	Accord. spec.
DNase I	TiRNA-01xxxxx	30.000 U *	n/a	White powder	Accord. spec.
chemagic Tips 96 Racks	-	10	-	-	-
chemagic Deep Well Plate 2 mL	-	72	-	-	-

* = 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.