

Protocol report

Blood_RNA_1mL_Flex

2013.02.21 08:37:04+02:00

1/7

General info

Protocol information

Protocol name

Blood_RNA_1mL_Flex

Modified by

admin

Kit name

MagJET Whole Blood RNA Purification Kit

Description

MagJET Whole Blood RNA Kit Protocol for RNA purification from 1mL of blood using KingFisher Flex instrument.

Protocol report

Blood_RNA_1mL_Flex

2013.02.21 08:37:04+02:00

Sample layout

| | 1 | 2 | 3 | 4 | 5 | 6 |
|---|---|---|---|---|---|---|
| A | | | | | | |
| B | | | | | | |
| C | | | | | | |
| D | | | | | | |








Reagent info

| Sample | | KingFisher 24 DW plate | | |
|----------------------|------------------|---------------------------|---------|--|
| Name | Well volume [µl] | Total reagent volume [µl] | Type | |
| Blood Sample | 1000 | - | Sample | |
| Lysis Buffer | 1000 | - | Reagent | |
| Proteinase K | 80 | - | Reagent | |
| Wash 1 | | KingFisher 24 DW plate | | |
| Name | Well volume [µl] | Total reagent volume [µl] | Type | |
| Wash Buffer 1 | 2800 | - | Reagent | |
| Wash 2 | | KingFisher 24 DW plate | | |
| Name | Well volume [µl] | Total reagent volume [µl] | Type | |
| Wash Buffer 2 | 2800 | - | Reagent | |
| Wash 3 | | KingFisher 24 DW plate | | |
| Name | Well volume [µl] | Total reagent volume [µl] | Type | |
| Wash Buffer 2 | 2800 | - | Reagent | |
| Elution | | KingFisher 24 DW plate | | |
| Name | Well volume [µl] | Total reagent volume [µl] | Type | |
| Water, nuclease free | 200 | - | Reagent | |
| Tip plate | | KingFisher 24 DW plate | | |
| Name | Well volume [µl] | Total reagent volume [µl] | Type | |
| - | - | - | - | |
| Prewash 1 | | KingFisher 24 DW plate | | |
| Name | Well volume [µl] | Total reagent volume [µl] | Type | |
| Wash Buffer 1 | 2800 | - | Reagent | |
| DNase I | | KingFisher 24 DW plate | | |
| Name | Well volume [µl] | Total reagent volume [µl] | Type | |
| DNase I solution | 200 | - | Reagent | |

Dispensed reagents

| Sample | | KingFisher 24 DW plate | | |
|----------------|------------|------------------------|---------------------------|--|
| Name | Step | Well volume [µl] | Total reagent volume [µl] | |
| Magnetic Beads | Dispense 1 | 60 | - | |
| Isopropanol | Dispense 1 | 2000 | - | |
| DNase I | | KingFisher 24 DW plate | | |
| Name | Step | Well volume [µl] | Total reagent volume [µl] | |


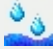




Steps data

| | | | |
|---|-------------------|------------------------|--|
|  | Tip | 24 DW tip comb | |
|  | Pick-Up | Tip plate | |
|  | Lysis | Sample | |
| | Beginning of step | Precollect | No |
| | | Release beads | No |
| | Mixing / heating: | Mixing time, speed | 00:15:00, Medium |
| | | Heating during mixing | No |
| | End of step | Postmix | No |
| | | Collect beads | No |
|  | Dispense 1 | Sample | |
| | | Message | Add Magnetic Beads 60 µl and Isopropanol 2000 µl |
| | Reagent(s) | Dispensing volume [µl] | 2060 |
| | | Name | Magnetic Beads |
| | | Volume [µl] | 60 |
| | | Name | Isopropanol |
| | | Volume [µl] | 2000 |
|  | Bind | Sample | |
| | Beginning of step | Precollect | No |
| | | Release beads | No |
| | Mixing / heating: | Mixing time, speed | 00:05:00, Bottom mix |
| | | Heating during mixing | No |
| | End of step | Postmix | No |
| | | Collect count | 5 |
| | | Collect time [s] | 3 |
|  | Pre_WS1 | Prewash 1 | |
| | Beginning of step | Precollect | No |
| | | Release beads | Yes |
| | Mixing / heating: | Shake 1 time, speed | 00:00:05, Bottom mix |
| | | Shake 2 time, speed | 00:01:00, Fast |
| | | Heating during mixing | No |
| | End of step | Postmix | No |
| | | Collect count | 3 |
| | | Collect time [s] | 1 |
|  | Dry | Prewash 1 | |
| | | Dry time | 00:02:00 |
| | | Tip position | Outside well / tube |

Protocol report

Blood_RNA_1mL_Flex





2013.02.21 08:37:04+02:00

| | | | |
|---|-------------------|--------------------------|------------------------------|
|  | DNase I | DNase I | |
| | Beginning of step | Precollect | No |
| | | Release beads | Yes |
| | Mixing / heating: | Shake 1 time, speed | 00:00:15, Bottom mix |
| | | Shake 2 time, speed | 00:15:00, Medium |
| | | Heating temperature [°C] | 37 |
| | End of step | Preheat | Yes |
| | | Postmix | No |
| | | Collect beads | No |
|  | Dispense 2 | DNase I | |
| | | Message | Add 1.4ml Rebinding Solution |
| | | Dispensing volume [µl] | 1400 |
| | Reagent(s) | Name | Rebinding Buffer |
| | | Volume [µl] | 1400 |
|  | Rebinding | DNase I | |
| | Beginning of step | Precollect | No |
| | | Release beads | Yes |
| | Mixing / heating: | Mixing time, speed | 00:05:00, Fast |
| | | Heating during mixing | No |
| | End of step | Postmix | No |
| | Collect count | 3 | |
| | Collect time [s] | 1 | |
|  | Wash 1 | Wash 1 | |
| | Beginning of step | Precollect | No |
| | | Release time, speed | 00:00:05, Bottom mix |
| | Mixing / heating: | Mixing time, speed | 00:01:00, Fast |
| | | Heating during mixing | No |
| | End of step | Postmix | No |
| | Collect count | 3 | |
| | Collect time [s] | 1 | |
|  | Wash 2 | Wash 2 | |
| | Beginning of step | Precollect | No |
| | | Release time, speed | 00:00:05, Bottom mix |
| | Mixing / heating: | Mixing time, speed | 00:01:00, Fast |
| | | Heating during mixing | No |
| | End of step | Postmix | No |
| | Collect count | 3 | |
| | Collect time [s] | 1 | |
|  | Wash 3 | Wash 3 | |
| | Beginning of step | Precollect | No |
| | | Release time, speed | 00:00:05, Bottom mix |
| | Mixing / heating: | Mixing time, speed | 00:01:00, Fast |
| | | Heating during mixing | No |
| | End of step | Postmix | No |
| | Collect count | 3 | |
| | Collect time [s] | 2 | |

Protocol report

Blood_RNA_1mL_Flex

2013.02.21 08:37:04+02:00

| | | | |
|---|-------------------|--------------------------|---------------------------|
|  | Dry 2 | Wash 3 | |
| | | Dry time | 00:05:00 |
| | | Tip position | Above well / tube surface |
|  | Elution | Elution | |
| | Beginning of step | Precollect | No |
| | | Release time, speed | 00:00:05, Fast |
| | Mixing / heating: | Mixing time, speed | 00:05:00, Fast |
| | | Heating temperature [°C] | 60 |
| | | Preheat | Yes |
| | End of step | Postmix | No |
| | | Collect count | 5 |
| | | Collect time [s] | 30 |
|  | ReleaseBeads | Wash 3 | |
| | | Release time, speed | 00:00:01, Fast |
|  | Leave | Tip plate | |

Lot info

No lot numbers have been defined.