

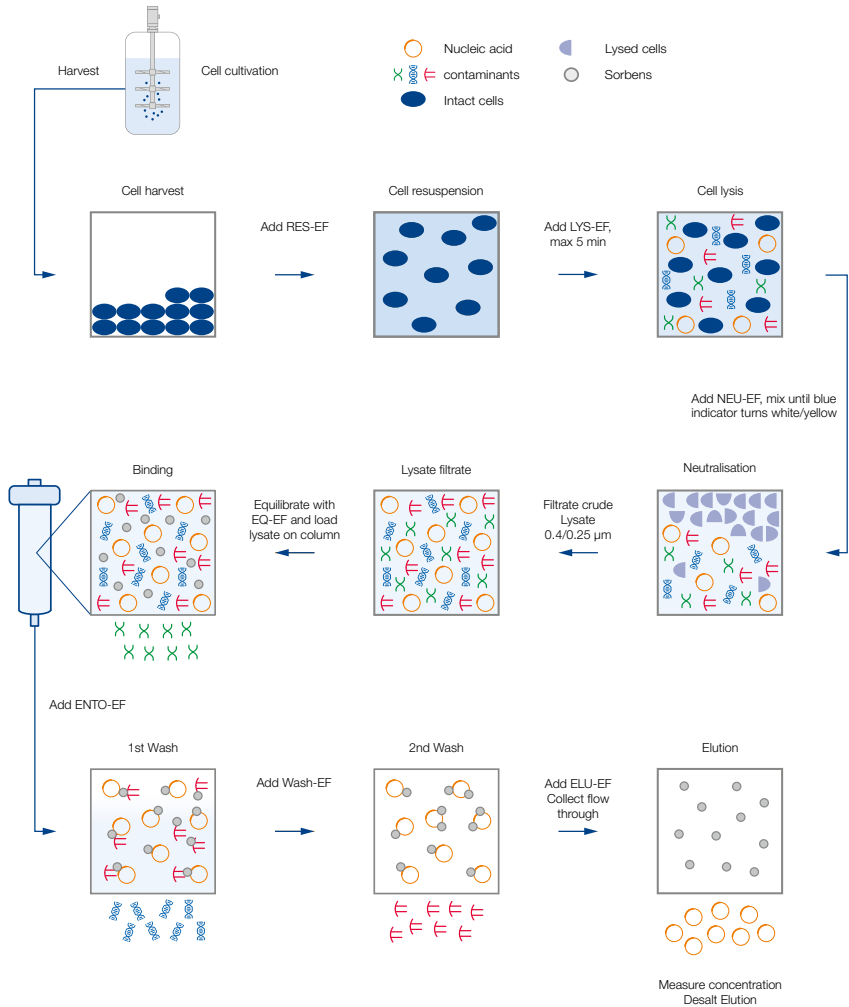
NucleoBond® RS purification workflow

Protocol at a glance (Rev. 01)

	RS 10	RS 50	RS 100	RS 200
1–3 Cell cultivation and harvest	4 °C, 15 min 4,500–6,000 × g			
	max. recom: ODV: 3,000 CDW: 8 g typ. Yield: ~10 mg	max. recom: ODV: 15,000 CDW: 40 g typ. Yield: ~50 mg	max. recom: ODV: 30,000 CDW: 80 g typ. Yield: ~100 mg	max. recom: ODV: 54,000 CDW: 145 g typ. Yield: ~190 mg
4–5 Cell lysis <i>(Important: Check Buffer LYS-EF for precipitated SDS)</i>	60 mL Buffer RES-EF	300 mL Buffer RES-EF	600 mL Buffer RES-EF	1080 mL Buffer RES-EF
	60 mL Buffer LYS-EF	300 mL Buffer LYS-EF	600 mL Buffer LYS-EF	1080 mL Buffer LYS-EF
RT, max 5 min				
6 Neutralization	60 mL Buffer NEU-EF	300 mL Buffer NEU-EF	600 mL Buffer NEU-EF	1080 mL Buffer NEU-EF
	Mix thoroughly until colorless Incubate 15 min on ice!			
<i>Maximum recommended pump speed</i>	5 mL/min	10 mL/min	10 mL/min	10 mL/min
7 Equilibration of the column	30 mL EQ-EF	120 mL EQ-EF	250 mL EQ-EF	400 mL EQ-EF
8 Clarification of the lysate	Clarify and filtrate the crude lysate to 0.45/0.22 µm			
9 Binding	Load cleared lysate on NucleoBond® RS column			
10 1st Wash	30 mL Buffer ENTO-EF	150 mL Buffer ENTO-EF	300 mL Buffer ENTO-EF	550 mL Buffer ENTO-EF
11 2nd Wash	30 mL Buffer WASH-EF	150 mL Buffer WASH-EF	300 mL Buffer WASH-EF	550 mL Buffer WASH-EF
13 Elution	30 mL Buffer ELU-EF	120 mL Buffer ELU-EF	250 mL Buffer ELU-EF	400 mL Buffer ELU-EF
14 Precipitation	21 mL Isopropanol	84 mL Isopropanol	175 mL Isopropanol	280 mL Isopropanol
	Mix thoroughly 4,500–15,000 × g 4 °C, 15 min			
15 Washing and drying	5 mL 70 % EtOH 4,500–15,000 × g RT, 5 min	15 mL 70 % EtOH 4,500–15,000 × g RT, 5 min	25 mL 70 % EtOH 4,500–15,000 × g RT, 5 min	50 mL 70 % EtOH 4,500–15,000 × g RT, 5 min
	Approx. RT, 10–15 min	Approx. RT, 30–60 min	Approx. RT, 45–90 min	Approx. RT, 75–120 min
16 Reconstitution	Appropriate volume of TE-EF or H ₂ O-EF.			
	It is recommended to choose the resuspension volume according to the requirements of the downstream application. (for var., @ 1 mg/mL)			
	8–10 mL	40–50 mL	80–100 mL	170–190 mL

NucleoBond® RS purification workflow

Protocol at a glance (Rev. 01)



This product distributed by

Takara Bio USA, Inc.

Order now
1.800.662.2566
ordersUS@takarabio.com

Talk to a scientist
1.800.662.2566
technical_support@takarabio.com



takarabio.com

For research use only. Not for resale. Unless otherwise indicated, all trademarks are the property of Takara Bio Inc. ©2023

