

distributed in the US/Canada by:

EAGLE BIOSCIENCES, INC.

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

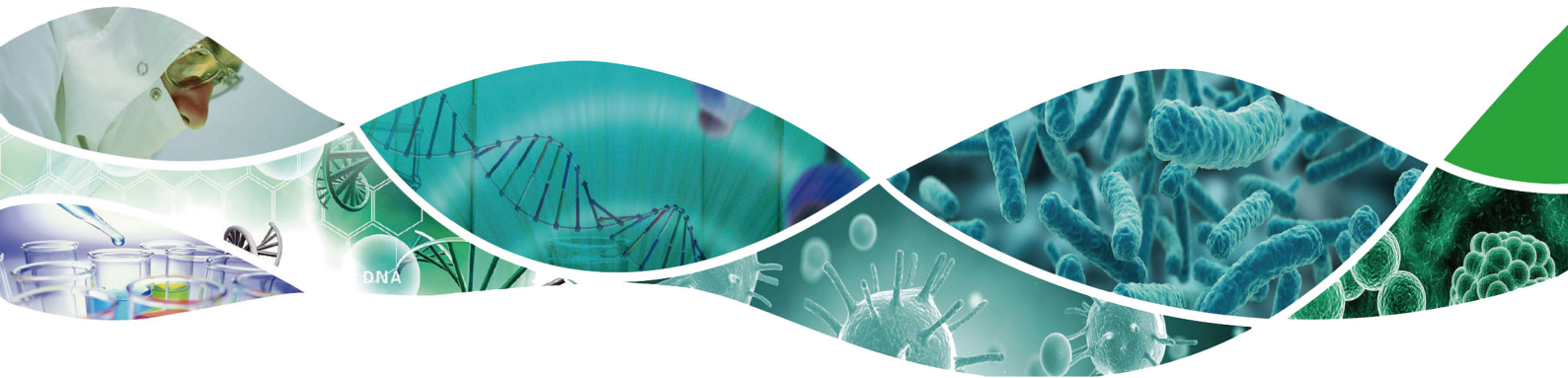
www.EagleBio.com • info@eaglebio.com



EAGLE
BIOSCIENCES

PRODUCT CATALOGUE

SSNP-3000A Nucleic Acid Extraction System



FEATURES

- 10.1" touch screen, user-friendly software
- Easy to use, fully automatic operation
- Magnetic bead method nucleic acid extraction kit results high purity and yield efficiency.
- High throughput: 64 samples simultaneously
- High efficiency: 15-45 minutes result
- Avoid biological contamination: UV light to disinfect, HEPA filter, Independent air filter



SSNP-3000A (64 channels)

SPECIFICATION

Model	SSNP-3000A
Sample Capacity	20μL -1000μL
Sample Throughput	1-64
Efficiency of Magnetic Bead Recycle	≥98%
Plate Type	96-well plate
Size (Length×Width×Height)	435×320×575mm
Weight	28kg
Reagent Type	Reagent for magnetic bead
Oscillating Mixing	Multi-mode and multi-speed for adjustment
Auxiliary Reagent	Viral Nucleic Acid Extraction Kit (Magnetic Bead Method) (Cat.No.SDK60104)
Program Management	New build, edit, delete mode programs, free and flexibly edit extraction system

Heating Temperature

Cracking heating temperature: room temperature to 120℃
Elution heating temperature: room temperature to 120℃

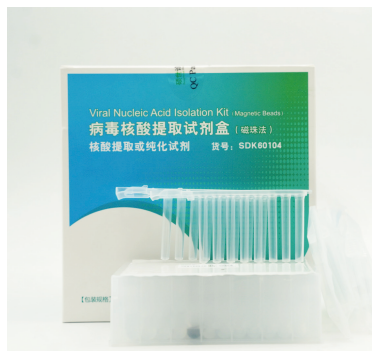
Contamination Control

The experimental cabin is equipped with an external HEPA filter independent air path, in which the biological filter cotton can absorb the nucleic acid aerosol.

EASY TO OPERATE

- **Step1:** Power on the instrument
- **Step2:** Add the samples in the 96-Well plate
- **Step3:** Place the 96-Well plate in the instrument
- **Step4:** Select and start the program
- **Step5:** Extract automatically and get high-purity nucleic acid

AUXILIARY REAGENT



Viral Nucleic Acid Extraction Kit (Magnetic Bead Method)

- Sample volume: 200μL
- Format: 96-Well plate
- Storage temperature: 4-30℃
- Transportation temperature: 4-30℃

SAMPLE TYPE

- Serum, plasma, virus culture fluid, cerebrospinal fluid, urine, swab wash, fecal supernatant

PACKAGING

CATALOG	PRODUCT	PACKAGE
SDK60104	Viral Nucleic Acid Extraction Kit (Magnetic Bead Method)	32 Tests/Kit 48 Tests/Kit 96 Tests/Kit



JIANGSU BIOPERFECTUS TECHNOLOGIES CO.,LTD

OFFICE: 2nd floor, Block 10, No.188 Xinjun Ring Road, Pujiang High-Tech Park, Caohejing Development Area, Shanghai, China.

HQ: 3rd & 4th floors of Building A(G19), 4th floor of Building F(G14), Ground floor of Building G20, Shuaiyu Village, Fuye village, Sixiang town, Taizhou National Medical Hi tech Development Zone, 225300 Taizhou, Jiangsu P.R. China.

EMAIL: info@s-sbio.com

WEBSITE: www.s-sbio.com

AUTHORIZED DISTRIBUTOR

Kindly follow us "bioperfectus"

