

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

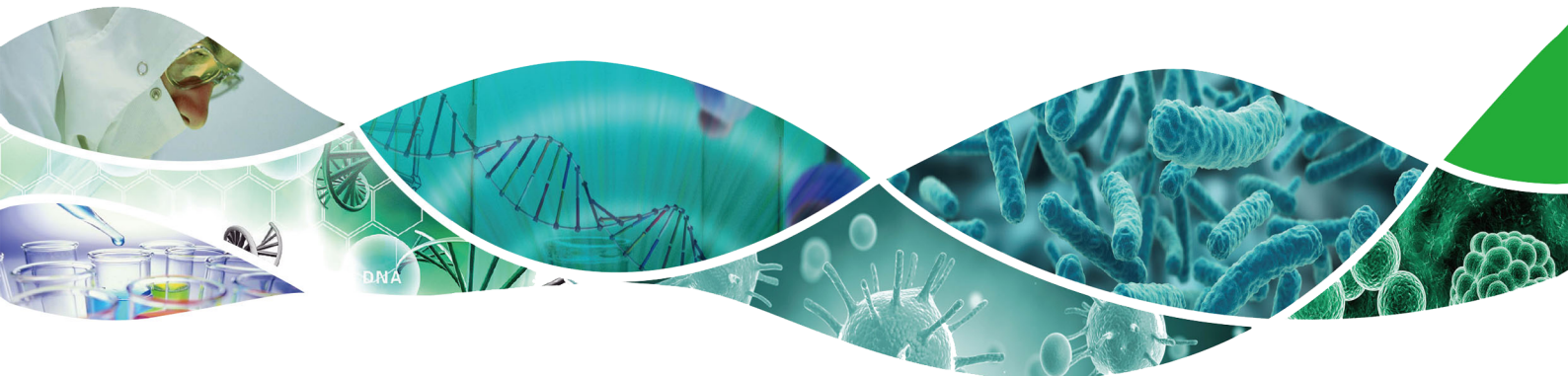
www.EagleBio.com • info@eaglebio.com



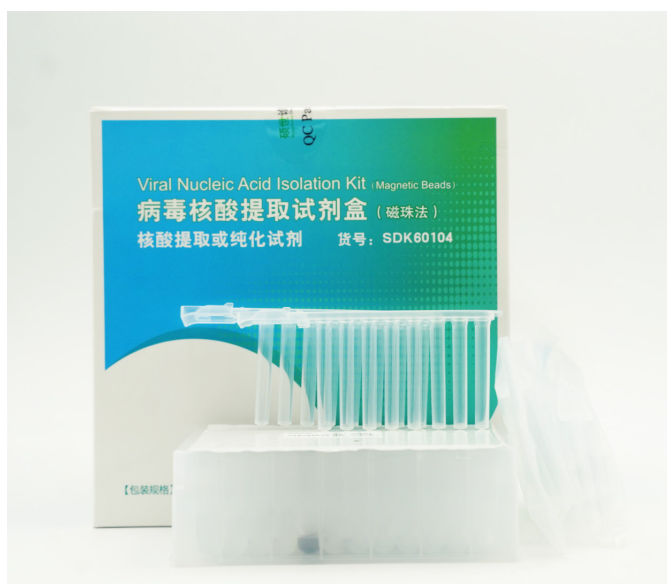
EAGLE
BIOSCIENCES

PRODUCT CATALOGUE

Viral Nucleic Acid Extraction Kit (Magnetic Bead Method)



FEATURES



- **Fast and high-quality nucleic acid**
- **Format: 96-Deep-Well plate**
- **Time per run (1-96 samples): 30-45 mins**
- **Application: Real-time PCR**
RT-PCR
PCR
Gene Sequencing
- **Storage Temperature: 4~30°C**
- **Transportation Temperature: 4~30°C**

SAMPLE TYPE

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

COMPONENT (Per Plate)

COLUMN	SOLUTION NAME	COMPONENTS	VOLUME
Column 1/7	Lysis binding solution	Guanidinium thiocyanate	650 µL
Column 2/8	Washing liquid 1	Guanidinium thiocyanate	500 µL
Column 3/9	Washing liquid 1	Guanidinium thiocyanate	500 µL
Column 4/10	Washing liquid 2	Ethanol	600 µL
Column 5/11	Magnetic bead solution	Magnetic beads	200 µL
Column 6/12	Eluent	RNase-Free water	80 µL

INSTRUCTION FOR USE

- **Step1:** Power on the instrument and self-check
- **Step2:** Add the samples in the 96-Deep-Well plate
- **Step3:** Place the 96-Deep-Well plate in the instrument
- **Step4:** Select a program
- **Step5:** Start the program and extract automatically
- **Step6:** Obtain high purity nucleic acid

PACKAGING

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

APPLICABLE MODELS

NUCLEIC ACID EXTRACTION SYSTEMS



SSNP-2000B(32 channels)

SSNP-3000A(64 channels)

SSNP-9600A(96 channels)

Performing fully automatic extraction of nucleic acid with Viral Nucleic Acid Extraction Kit (Magnetic Bead Method), above user-friendly instruments could simplify process and increase throughput.



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"

