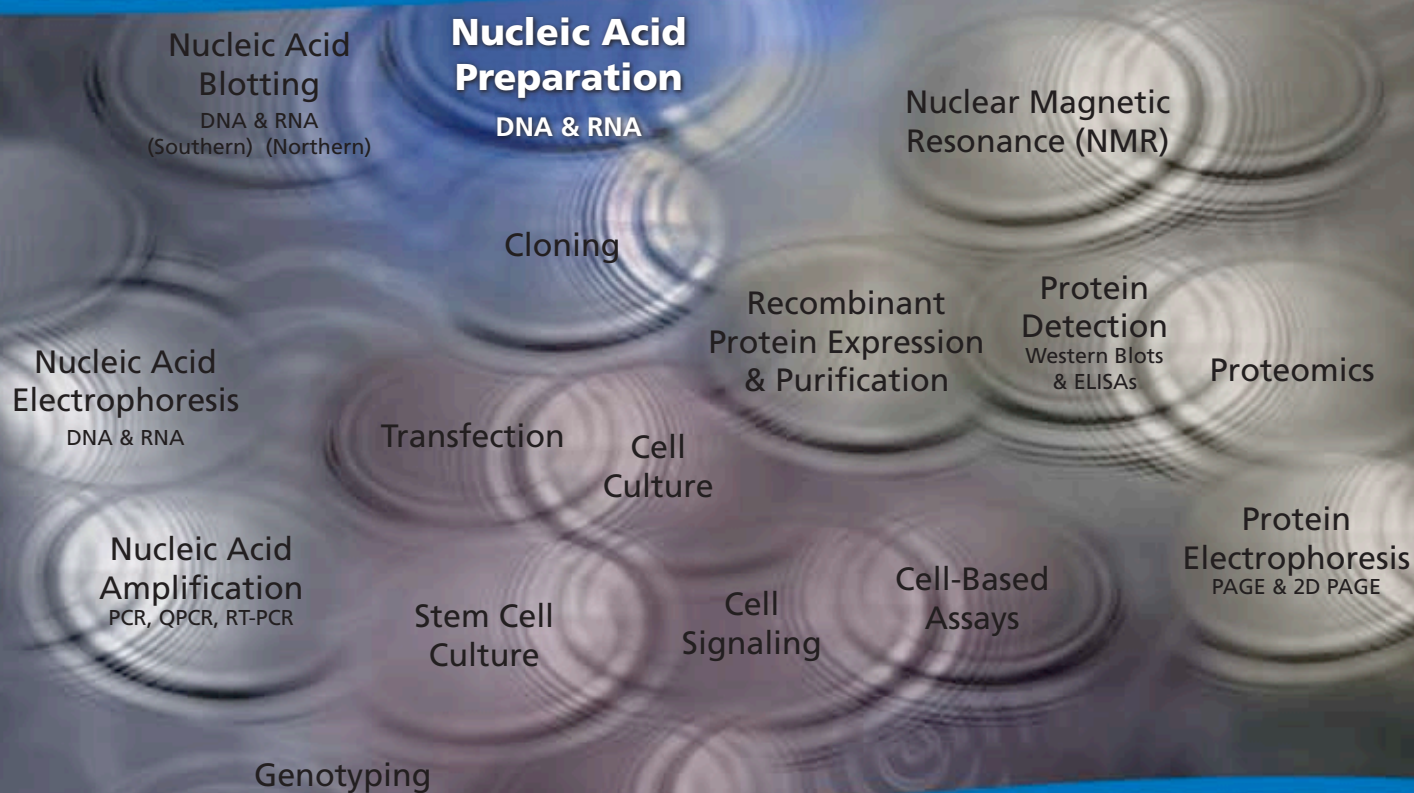


Life Science Reagents for Nucleic Acid Purification from Sigma-Aldrich®



We Ripple Through Everything You Do!

- Extensive reagent offering to support your Nucleic Acid Purification based research needs
- Full selection of Purification kits including mammalian, plant, blood, and bacterial kits
- Quality and availability assured
- Market leading Technical Service
- Leading Web support tools:

DNA & RNA Purification — sigma-aldrich.com/dnapurification
Plasmid Partner — sigma-aldrich.com/plasmidpartner
Extract-N-Amp™ Kits — sigma-aldrich.com/xna

Life Science Reagents for Nucleic Acid Purification from Sigma-Aldrich®

Sigma offers a full line of purification kits and reagents!

• Plasmid Kits

- For use with large and small scale plasmid, endotoxin-free plasmid, and BAC DNA
- Technology uses advanced silica bind and elute with vacuum and spin format
- For use with transfections, sequencing, restriction digests, cloning, and gene therapy

• Genomic DNA Kits

- Kits for isolating mammalian, plant, blood, and bacterial DNA.
- Useful with applications such as PCR, restriction digests, cloning, southern blots, and sequencing

• RNA Kits

- Kits for the purification of total RNA and mRNA
- Mammalian kits may be used with multiple tissue types including brain, heart, kidney and more
- Isolated RNA is suitable for northern blotting, RT-PCR and other common applications

To view our full line of purification products, visit sigma-aldrich.com/dnapurification.

Cat. No.	Description
Acids & Bases	
H1758	Hydrochloric acid Molecular Biology Grade, 36.5-38.0 %
438081	Phosphoric acid ACS reagent, ≥85 wt. % in H ₂ O
S2770	Sodium hydroxide solution 1.0 N, cell culture tested
Agarose	
A9539	Agarose for molecular biology
A4718	Agarose High resolution, DNase, RNase, NICKase, none detected
A5093	Agarose Low EEO, for molecular biology
A6013	Agarose Type I, Low EEO
A0169	Agarose Type I-A Low EEO
A6877	Agarose Type II, Medium EEO
A2576	Agarose Type IX-A, Ultra-low Gelling Temperature
A0701	Agarose Type VII-A, Low Gelling Temperature
A7431	Agarose Wide range/Standard 3:1, DNase, RNase, NICKase, none detected
Buffers	
A2706	Ammonium acetate solution for molecular biology, 7.5 M
A4418	Ammonium sulfate for molecular biology, ≥99.0%
B6768	Boric acid for molecular biology, ~99%
M2933	MES hydrate Biotechnology Performance Certified, plant cell culture tested, ≥99.5%
M3058	MES sodium salt Biotechnology Performance Certified
M1317	MES solution Biotechnology Performance Certified, 1 M, cell culture tested
P5493	Phosphate buffered saline 10× Concentrate, Biotechnology Performance Certified, cell culture tested
P5368	Phosphate buffered saline pH 7.4, Biotechnology Performance Certified
P1851	PIPES Biotechnology Performance Certified, cell culture tested, ≥99%
P5833	Potassium carbonate SigmaUltra, ≥99.0%
P9791	Potassium phosphate monobasic for molecular biology, ≥98%
S2889	Sodium acetate anhydrous for molecular biology
S3014	Sodium chloride for molecular biology, ≥98%
S6316	Sodium chloride solution 5 M, Biotechnology Performance Certified, cell culture tested
T9650	Tris Acetate-EDTA buffer 10× Concentrate, DNase and RNase, none detected
T4415	Tris-Borate-EDTA buffer 10× Concentrate
T6066	Trizma® base Biotechnology Performance Certified, meets EP, USP testing specifications, cell culture tested, ≥99.9%
T5941	Trizma hydrochloride for molecular biology, Biotechnology Performance Certified, cell culture tested, ≥99.0%
Detergents	
P5413	Poly(ethylene glycol) for molecular biology, average mol. wt. 8,000
L4390	Sodium dodecyl sulfate for molecular biology, ≥98.5%
T8787	Triton® X-100 for molecular biology
P9416	TWEEN® 20 for molecular biology, viscous liquid

Cat. No.	Description
General Reagents	
M3148	2-Mercaptoethanol for molecular biology, ≥98%
B4287	Albumin from bovine serum for molecular biology, powder
B2518	Albumin, Acetylated from bovine serum lyophilized powder, for molecular biology
A9434	Ammonium chloride for molecular biology, ≥99.5%
A3678	Ammonium persulfate for electrophoresis, ≥98%
C3306	Calcium chloride dihydrate for molecular biology, ~99%
C3032	Cesium chloride for molecular biology, ≥99%
C2432	Chloroform ≥99%, contains amylenes as stabilizer
C0549	Chloroform:Isoamyl alcohol 24:1
D2438	Dimethyl sulfoxide Biotechnology Performance Certified, sterile-filtered, hybridoma tested, meets EP, USP testing specifications
D0632	DL-Dithiothreitol ≥98% (TLC), ≥99% (titration)
D9779	DL-Dithiothreitol for molecular biology, ≥98% (TLC), ≥99% (titration)
M9546	D-Mannitol SigmaUltra™
E5134	EDTA disodium salt dihydrate for molecular biology, ≥99%
E7889	EDTA disodium salt solution for molecular biology, 0.5 M
E3889	EGTA for molecular biology, ≥97.0%
E7023	Ethanol 200 proof (absolute), for molecular biology
E1510	Ethidium bromide solution 10 mg/mL, for molecular biology, aqueous solution
G7663	Gelatin blocking buffer for molecular biology, powder blend
G5516	Glycerol for molecular biology, ≥99%
G3272	Guanidine hydrochloride for molecular biology, ≥99%
G9277	Guanidine thiocyanate for molecular biology, ≥99%
L7026	Lithium chloride solution 8 M, for molecular biology, ≥99%
M1028	Magnesium chloride solution for molecular biology, 1.00 M ±0.01 M
P1037	Phenol for molecular biology
P2069	Phenol:Chloroform:Isoamyl Alcohol 25:24:1 Saturated with 10 mM Tris, pH 8.0, 1 mM EDTA. for molecular biology
P1190	Potassium acetate for molecular biology, ≥99%
383112	Sodium iodide ACS reagent, ≥99.5 %
S0389	Sucrose for molecular biology, ≥99.5%
T9281	TEMED for electrophoresis, ~99%
U5378	Urea for molecular biology, ≥98%
W4502	Water Molecular Biology Reagent
534056	Xylenes histological grade
T9424	TRI Reagent® For processing tissues, cells cultured in monolayer or cell pellets
R0901	RNAlater®
P3803	Phenol:Chloroform:Isoamyl Alcohol 25:24:1 Saturated with 10 mM Tris, pH 8.0, 1 mM EDTA for molecular biology
R2020	RNaseZAP®
P1944	Phenol:Chloroform 5:1 for molecular biology
T3809	TRI Reagent BD, For processing whole blood, plasma, or serum

Cat. No.	Description
Kits	
56500	GenElute™ Agarose Spin Columns
56501	GenElute Minus EtBr Spin Column
NA2110	GenElute Bacterial Genomic DNA Kit
PFM250	GenElute Five-Minute Plasmid Miniprep Kits
NA0410	GenElute HP Endotoxin-Free Plasmid Maxiprep Kit
NA0600	GenElute HP Endotoxin-Free Plasmid Megaprep Kit
NA0310	GenElute HP Plasmid Maxiprep Kit
NA0500	GenElute HP Plasmid Megaprep Kit
NA0200	GenElute HP Plasmid Midiprep Kit
NA0160	GenElute HP Plasmid Miniprep Kit
NA0800	GenElute HP Select Plasmid Gigaprep Kit
G1N70	GenElute Mammalian Genomic DNA Miniprep Kit
G1N350	GenElute Mammalian Genomic DNA Miniprep Kit
RTN70	GenElute Mammalian Total RNA Miniprep Kit
NA0100	PhasePrep BAC DNA Kit

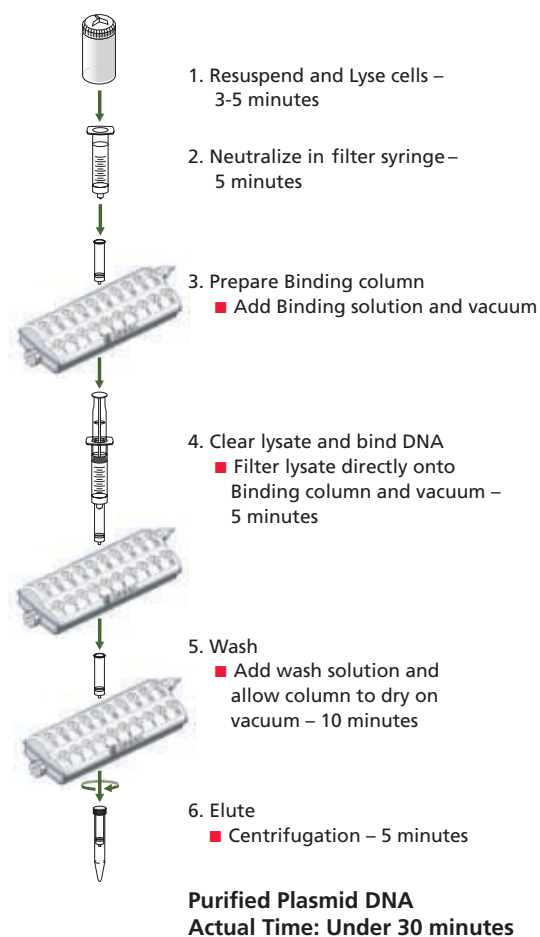


GenElute™ HP Plasmid Midiprep Kit
(Cat. No. NA0200)

Plasmid Purification

More Plasmid DNA in Less Time

HP Midi or Maxiprep



GenElute™ HP Plasmid Purification Kits offer the fastest, most efficient, and highest yielding plasmid purification in the market.

Our kits feature an optimized silica-based membrane technology that eliminates the need for alcohol precipitation. They are also compatible with both vacuum and spin purification formats for maximum flexibility.

Features & Benefits

- Saves time and provides increased yield through an optimized silica-based membrane technology
- Purified DNA is suitable for all downstream applications including Sequencing, PCR, QPCR, Transfection
- Compatible with vacuum or spin format
- No phenol/chloroform extraction or alcohol precipitation required

Performance Information

Kit Type	Prep Time w/vacuum manifold	Yield (µg)
GenElute Five-Minute Plasmid Miniprep Kit	5 minutes	5
GenElute HP Plasmid Miniprep	25 minutes	25
GenElute HP Plasmid Midiprep Kit	30 minutes	350
GenElute HP Plasmid Maxiprep Kit	45 minutes	1200
GenElute HP Endotoxin-Free Plasmid Maxiprep Kit	45 minutes	1200

Learn more at sigma-aldrich.com/dnapurification.

World Headquarters

3050 Spruce St., St. Louis, MO 63103
(314) 771-5765
sigma-aldrich.com

Order/Customer Service (800) 325-3010 • Fax (800) 325-5052

Technical Service (800) 325-5832 • sigma-aldrich.com/techservice

Development/Bulk Manufacturing Inquiries **SAFC**™ (800) 244-1173

*Accelerating Customers'
Success through Leadership
in Life Science, High
Technology and Service*

©2008 Sigma-Aldrich Co. All rights reserved. SIGMA, SAFC, SAFC, SIGMA-ALDRICH, ALDRICH, FLUKA, and SUPELCO are trademarks belonging to Sigma-Aldrich Co. and its affiliate Sigma-Aldrich Biotechnology, L.P. Sigma brand products are sold through Sigma-Aldrich, Inc. Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.

Trizma, REDTaq, Restorase, Sigma and Sigma-Aldrich are registered trademarks of Sigma-Aldrich Biotechnology LP and Sigma-Aldrich Co. eAMV, JumpStart, AccuTag, ReadyMix, GenElute and SigmaUltra are trademarks of Sigma-Aldrich Biotechnology LP and Sigma-Aldrich Co. SYBR is a registered trademark of Molecular Probes, Inc. KlenTaq is a registered trademark of Wayne Barnes. Triton is a registered trademark of Union Carbide Corp. TWEEN is a registered trademark of Uniqema, a business unit of ICI Americas, Inc. TRI Reagent are registered trademarks of Molecular Research Center, Inc. RNAlater and RNaseZAP are registered trademarks of Ambion, Inc.

KSY
04653
0078