



DNA from Practically Anything

Citations Isolating Bacterial DNA with the Wizard® Genomic DNA Purification Kit

The Wizard® Genomic DNA Purification Kit is a gentle, scalable solution-based method for the isolation of high molecular weight genomic DNA (>50kb) from a variety of starting materials. Includes protocols for the isolation of gDNA from Gram-positive and Gram-negative bacteria as well as fungi and yeast. This is a partial list of citations describing the Wizard® Genomic DNA Purification Kit use with bacterial sources. The system has been cited for isolation of gDNA from other bacteria including ***Bordetella*, *Campylobacter*, *Enterobacter*, *Klebsiella*, *Lactobacillus*, *Leptospira*, *Mycobacterium*, *Desulfovibrio*, *Escherichia*, *Flavobacterium*, *Methanococcus*, *Microbacterium*, *Mycoplasma*, *Paracoccus*, *Prevotella*, *Proteus*, *Rhodopseudomonas*, *Rickettsia*, *Salmonella*, *Serratia*, *Shigella*, *Sphingomonas*, *Treponema* and *Vibrio***. See the protocol online at: www.promega.com/tbs/tm050/tm050.html

Acinetobacter

Zarrilli, R. *et al.* (2004) Molecular epidemiology of sequential outbreaks of *Acinetobacter baumannii* in an intensive care unit shows the emergence of carbapenem resistance. *J. Clin. Microbiol.* **42**, 946–53.

Aeromonas

Yu, H.B. *et al.* (2004) A type III secretion system is required for *Aeromonas hydrophila* AH-1 pathogenesis. *Infect. Immun.* **72**, 1248–56.

Agrobacterium

Weller, S.A., Stead, D.E. and Young, J.P.W. (2004) Acquisition of an *Agrobacterium* Ri plasmid and pathogenicity of other α -*proteobacteria* in cucumber and tomato crops affected by root mat. *Appl. Environ. Microbiol.* **70**, 2779–85.

Bacillus

van Schaik, W. *et al.* (2004) The alternative sigma factor σ^B of *Bacillus cereus*: Response to stress and role in heat adaptation. *J. Bacteriol.* **186**, 316–25.

Bordetella

Brockmeier, S.L. *et al.* (2002) Role of the dermonecrotic toxin of *Bordetella bronchiseptica* in the pathogenesis of respiratory disease in swine. *Infect. Immun.* **70**, 481–90.

Brucella

Baek, S.-H. *et al.* (2004) Denitrification genes regulate *Brucella* virulence in mice. *J. Bacteriol.* **186**, 6025–31.

Burkholderia

Reckseidler-Zenteno, S.L., DeVinney, R. and Woods, D.E. (2005) The capsular polysaccharide of *Burkholderia pseudomallei* contributes to survival in serum by reducing complement factor C3b deposition. *Infect. Immun.* **73**, 1106–15.

Candidatus Glomeribacter gigasporarum

Jargeat, P. *et al.* (2004) Isolation, free-living capacities, and genome structure of "Candidatus Glomeribacter gigasporarum," the endo cellular bacterium of the mycorrhizal fungus *Gigaspora margarita*. *J. Bacteriol.* **186**, 6876–84.

Chlamydomphila

Slepenkin, A., de la Maza, L.M. and Peterson, E.M. (2005) Interaction between components of the type III secretion system of the *Chlamydiaceae*. *J. Bacteriol.* **187**, 473–9.

Clostridium

Zhou, Y. and Singh, B.R. (2004) Cloning, high-level expression, single-step purification, and binding activity of His₆-tagged recombinant type B botulinum neurotoxin heavy chain transmembrane and binding domain. *Prot. Exp. Purif.* **34**, 8–16.

Desulfatibacillum

Cravo-Lareau, C. *et al.* (2004) *Desulfatibacillum aliphaticivorans* gen. Nov., sp. nov., an n-alkane- and n-alkene-degrading, sulfate-reducing bacterium. *Int. J. Syst. Evol. Microbiol.* **54**, 77–83.

Enterococcus

Arbeloa, A. *et al.* (2004) Role of class A penicillin-binding proteins in PBP5-mediated β -lactam resistance in *Enterococcus faecalis*. *J. Bacteriol.* **186**, 1221–8.

Haemophilus

Saeed-Kothe, A., Yang, W. and Mills, S.D. (2004) Use of the riboflavin synthase gene (*ribC*) as a model for development of an essential gene disruption and complementation system for *Haemophilus influenzae*. *Appl. Environ. Microbiol.* **70**, 4136–43.

Helicobacter

Moore, J.M. and Salama, N.R. (2005) Mutational analysis of metronidazole resistance in *Helicobacter pylori*. *Antimicrob. Agents Chemother.* **49**, 1236–7.

Legionella

De Buck, E. *et al.* (2004) A putative twin-arginine translocation pathway in *Legionella pneumophila*. *Biochem. Biophys. Res. Comm.* **317**, 654–61.

Listeria

Rodríguez-Lázaro, D. *et al.* (2004) Quantitative detection of *Listeria monocytogenes* and *Listeria innocua* by real-time PCR: Assessment of *hly*, *iap*, and *lin02483* targets and AmpliFluor technology. *Appl. Environ. Microbiol.* **70**, 1366–77.

Myxococcus

Pérez-Marín, M. *et al.* (2004) The N terminus of *Myxococcus xanthus* CarA repressor is an autonomously folding domain that mediates physical and functional interactions with both operator DNA and antirepressor protein. *J. Biol. Chem.* **279**, 33093–103.

Neisseria

Lynn, F. *et al.* (2005) Genetic typing of the porin protein of *Neisseria gonorrhoeae* from clinical noncultured samples for strain characterization and identification of mixed gonococcal infections. *J. Clin. Microbiol.* **43**, 368–75.

Pseudomonas

Lian, W. *et al.* (2004) Ultrasensitive detection of biomolecules with fluorescent dye-doped nanoparticles. *Anal. Biochem.* **334**, 135–44.

Rhodobacter

Morehouse, K.A., Goodfellow, I.G. and Sockett, R.E. (2005) A chimeric N-terminal *Escherichia coli*-C-terminal *Rhodobacter sphaeroides* FliG rotor protein supports bidirectional *E. coli* flagellar rotation and chemotaxis. *J. Bacteriol.* **187**, 1695–701.

Spiroplasma

Boutaread, A. *et al.* (2004) Disruption of a gene predicted to encode a solute binding protein of an ABC transporter reduces transmission of *Spiroplasma citri* by the leafhopper *Circulifer haematoceps*. *Appl. Environ. Microbiol.* **70**, 3960–7.

Staphylococcus

Muthaiyan, A., Jayaswal, R.K. and Wilkinson, B.J. (2004) Intact *mutS* in laboratory-derived and clinical glycopeptide-intermediate *Staphylococcus aureus* strains. *Antimicrob. Agents Chemother.* **48**, 623–5.

Streptococcus

Brown, J.S. *et al.* (2004) A locus contained within a variable region of pneumococcal pathogenicity island 1 contributes to virulence in mice. *Infect. Immun.* **72**, 1587–93.

Yersinia

Flashner, Y. *et al.* (2004) Generation of *Yersinia pestis* attenuated strains by signature-tagged mutagenesis in search of novel vaccine candidates. *Infect. Immun.* **72**, 908–15.

Product	Size	Cat. #
Wizard® Genomic DNA Purification Kit	100 isolations × 300µl	A1120
	500 isolations × 300µl	A1125
	100 isolations × 10ml	A1620

For Laboratory Use.

Note: Quantity "Isolations" for the Wizard® Genomic DNA Purification Kit refers to isolations from the indicated quantity of whole human blood. Cat.# A1120 will process ~80ml of bacterial cell culture; Cat. #A1125 will process ~400ml of bacterial cell culture; and Cat. #A1620 will process ~1,600ml of bacterial cell culture.