

Selected PHOTOPROBE® Biotin references:

- Saito M, et al. 2013. *Leptospira idonii* sp. nov., isolated from environmental water. Int J Syst Evol Microbiol. 63:2457-62. Doi: 10.1099/ijs.0.047233-0
- Arai R, et al. 2012. Diversity of *Melissococcus plutonius* from honeybee larvae in Japan and experimental reproduction of European foulbrood with cultured atypical isolates. PLoS ONE. 7(3):e33708. doi: 10.1371/journal.pone.0033708.
- Stearns NA et al. 2012. The inhibition of anti-DNA binding to DNA by nucleic acid binding polymers. PLoS ONE. 7(7):e40862. doi: 10.1371/journal.pone.0040862
- Boothby TC and Wolniak SM. 2011. Masked mRNA is stored with aggregated nuclear speckles and its asymmetric redistribution requires a homolog of mago nashi. BMC Cell Biol. 12:45.
- Ohkuni K and Kitagawa K. 2011. Endogenous transcription at the centromere facilitates centromere activity in budding yeast. Cur. Biol. 21:1695-1703. Doi: 10.1016/j.cub.2011.08.056
- Tsutsui KM et al. 2011. Nuclear protein LEDGF/p75 recognizes supercoiled DNA by a novel DNA-binding domain. Nucleic Acids Res. 39(12):5067-81.
- Fuxman Bass JI et al. 2010. Extracellular DNA: a major proinflammatory component of *Pseudomonas aeruginosa* biofilms. J Immunol. 184:6386-95.
- Matsunaga F, et al. 2010. Localized melting of duplex DNA by Cdc6/Orc1 at the DNA replication origin in the hyperthermophilic archaeon *Pyrococcus furiosus*. Extremophiles. 14:21-31.
- Tomida J, et al. 2010. *Streptococcus fryi* sp. Nov., a novel species with Lancefield group M antigens. FEMS Microbiol. Lett. 314(1) doi:10.1111/j.1574-6968.2010.02148.x
- Evrin C, et al. 2009. A double-hexameric MCM2-7 complex is loaded onto origin DNA during licensing of eukaryotic DNA replication. Proc. Natl. Acad. Sci. USA. 106:20240-5.
- Ford CT, et al. 2009. Liver X receptor β : maintenance of epidermal expression in intrinsic and extrinsic skin aging. AGE. 31:365-72.
- Hagi T, and Hoshino T. 2009. Screening and characterization of potential probiotic lactic acid bacteria from cultured common carp intestine. Biosci. Biotechnol. Biochem. 73(7):1479-83.

Nakamura N et al. 2009. Characterization of the interaction between serum mannan-binding protein and nucleic acid ligands. *J. Leukocyte Biol.* 86:1-13.

Singh M, et al. 2009. Chromosomal localization of 18S and 5S rDNA using FISH in the genus *Tor* (Pisces, Cyprinidae). *Genetica.* 137:245-52.

Wang K, Mahalingam G, Imai Y, Pesnicak L, Margolis TT, Straus SE, Cohen JI. 2009. Cell type specific accumulation of the major latency-associated transcript (LAT) of herpes simplex virus type 2 in LAT transgenic mice. *Virology.* 386:79-87.

Yang J-N et al. 2009. Study on HBV vertical transmission via the *in vitro* fertilization (IVF) technique. *J Reproduct. Contracept.* 20(2):73-9.

Bass JIF et al. 2008. Characterization of bacterial DNA binding to human neutrophil surface. *Lab. Invest.* 88:926-37.

Fossati-Jimack L et al. 2008. C1q deficiency promotes the production of transgenic-derived IgM and IgG3 autoantibodies in anti-DNA knock-in transgenic mice. *Mol. Immunol.* 45:787-95.

Hagi T and Hoshino T. 2008. Screening and characterization of potential probiotic lactic acid bacteria from cultured common carp intestine. *Biosci. Biotechnol. Biochem.* 73(7):1479-83.

Rameshkumar N et al. 2008. *Vibrio porteresiae* sp. nov., a diazotrophic bacterium isolated from a mangrove-associated wild rice (*Porteresia coarctata* Tateoka). *Int. J. Syst. Evol. Microbiol.* 58:1608-15.

Voynova E et al. 2008. An antibody-based construct carrying DNA-mimotope and targeting CR1 (CD35) selectively suppresses human autoreactive B-lymphocytes. *Immunol. Letters.* 116:168-73.

Yamahira K, et al. 2008. *Acinetobacter* sp strain Ths, a novel psychrotolerant and alkalitolerant bacterium that utilizes hydrocarbon. *Extremophiles.* 12:729-34.

Huhtamella S et al. 2007. RNA-based sandwich hybridization method for detection of lactic acid bacteria in brewery samples. *J. Microbiol. Methods.* 68:543-53.

Jiang C, Foley J, Clayton N, Kissling G, Jokinen M, Herbert R and Diaz M. 2007. Abrogation of lupus nephritis in activation-induced deaminase-deficient MRL/lpr mice. *J Immunol.* 178:7422-31.

Kitamura T et al. 2007. *Helicobacter cinaedi* cellulitis and bacteremia in immunocompetent hosts after orthopedic surgery. *J. Clin Microbiol.* 45(1):31-8.

Nishiwaki H et al. 2007. Insecticidal bacteria isolated from predatory larvae of the antlion species *Myrmeleon bore* (Neuroptera:Myrmeleontidae). J. Invertebrate Pathol.. 96:80-88.

Hamilton KJ et al. 2006. The binding of sera of patients with SLE to bacterial and mammalian DNA. Clin. Immunol. 118:209-18.

Kawasaki Y, Kim H-D, Kojima A, Seki T, Sugino A. 2006. Reconstitution of *Saccharomyces cerevisiae* prereplicative complex assembly *in vitro*. Genes to Cells. 11:745-56.

Manderson AP, Carlucci F, Lachmann PJ, Lazarus RA, Restenstein RJ, Cook HT, Walport MJ, Botto M. 2006. The *in vivo* expression of actin/salt-resistant hyperactive DNase I inhibits the development of anti-ssDNA and anti-histone autoantibodies in a murine model of systemic lupus erythematosus. Arthritis Res. and Ther. 8:(3)R68.

Maukonen JN, Matto J, Satokari R, Soderlund H, Mattila-Sandholm T, Saarela M. 2006. PCR DGGE and RT-PCR DGGE show diversity and short-term temporal stability in the *Clostridium coccooides-Eubacterium rectale* group in the human intestinal microbiota. FEMS Microbiol Ecol. 58:517-28.

Maukonen J, Satokari R, Matto J, Soderlund H, Mattila-Sandholm T, Saarela M. 2006. Prevalence and temporal stability of selected clostridial groups in irritable bowel syndrome in relation to predominant faecal bacteria. J Med Microbiol. 55:625-33.

Nandivada H et al. 2006. Reactive polymer coatings that “click”. Agnew. Chem. Int. Ed. 45:3360-63.

Shirey KA, Jung Y-Y, Maeder GS, Carlin JM. 2006. Upregulation of EFN- γ receptor expression by proinflammatory cytokines influences IDO activation in epithelial cells. J Interferon and Cytokine Res. 26:53-62.

Will C, Schewe C, Schluns K, Petersen I. 2006. HPV typing and CGH analysis for the differentiation of primary and metastatic squamous cell carcinomas of the aerodigestive tract. Cellular Oncol. 28(3):97-105.

Hirota K, Nodasaka Y, Orikasa Y, Okuyama H, Yumoto I. 2005. *Shewanella pneumatophori* sp. nov., an eicosapentaenoic acid-producing marine bacterium isolated from the intestines of Pacific mackerel (*Pneumatophorus japonicus*). Int J of Systematic Evol Microbiol. 55:2355-9.

Kandasamy K, Joseph K, Subramaniam K, Raymond JR, Tholanikunnel BG. 2005. Translational control of β_2 -adrenergic receptor mRNA by T-cell-restricted

intracellular antigen-related protein. J Biol Chem. 280(3):1931-43.

Katsuta A, Adachi K, Matsuda S, Shizuri Y, Kasai H. 2005. *Ferrimonas marina* sp. nov. Int J of Systematic Evol Microbiol. 55:1851-5.

Korinth D, Pacyna-Gengelbach M, Deutchmann N, Hattenberger S, Bockmuhl U, Dietel M, Schroeder H-G, Donhuijsen K, Petersen I. 2005. Chromosomal imbalances in wood dust-related adenocarcinomas of the inner nose and their associations with pathological parameters. J Pathol. 207:207-15.

Satokari RM, Kataja K, Soderlund H. 2005. Multiplexed quantification of bacterial 16S rRNA by solution hybridization with oligonucleotides probes and affinity capture. Microbial Ecol. 50:120-27.

Tan J, Tannock GA. 2005. Role of viral load in the pathogenesis of chicken anemia virus. J Gen Vir. 86:1327-33.

Yumoto I, Hirota K, Nodasaka Y, Nakajima K. 2005. *Oceanobacillus oncorhynchi* sp. nov., a halotolerant obligate alkaliphile isolated from the skin of a rainbow trout (*Oncorhynchus mykiss*), and emended description of the genus *Oceanobacillus*. Int J of Systematic Evol Microbiol. 55:1521-4.

Gruber HE, Leslie K, Ingram J, Norton HJ, Hanley EN. 2004. Cell-based tissue engineering for the intervertebral disc: in vitro studies of human disc cell gene expression and matrix production within selected cell carriers. Spine J. 4:44-55.

Kawakami M, Kimura T, Kishimoto Y, Tatekawa T, Baba Y, Nishizaki T, Matsuzaki N, Taniguchi Y, Yoshihara S, Idegame K, Shirakata T, Nishida S, Masuda T, Hosen N, Tsuboi A, Oji Y, Oka Y, Ogawa H, Sonoda Y, Sugiyama H, Kawase I, Soma T. 2004. Preferential expression of the vasoactive intestinal peptide (VIP) receptor VPAC1 in human cord blood-derived CD34⁺CD38⁻ cells: possible role of VIP as a growth-promoting factor for hematopoietic stem/progenitor cells. Leukemia. 18:912-21.

Korinth D, Donhuijsen K, Bockmuhl U, Petersen I. 2004. PHOTOPROBE® biotin: An alternative method for labeling archival DNA for comparative genomic hybridization. Cellular Oncology. 26:329-34.

Krzywinski J, Nusskern DR, Kern MK, Besansky NJ. 2004. Isolation and characterization of Y chromosome sequences from the African malaria mosquito *Anopheles gambiae*. Genetics. 166:1291-1302.

Lighthall GK, Hamilton BP, Hamlyn JM. 2004. Identification of salt-sensitive genes in the kidneys of Dahl rats. J Hypertension. 22:1487-94.

- Matsuno Y, Adachi T, Koh K-B, Fukata H, Sugimura A, Sakurai K, Shibayama T, Iguchi T, Komiyama M, Mori C. 2004. Effect of neonatal exposure to diethylstilbestrol on testicular gene expression in adult mouse: comprehensive analysis with cDNA subtraction method. *Int J Androl.* 27:115-22.
- Rigby RJ, Rozzo SJ, Boyle JJ, Lewis M, Kotzin BL, Vyse TJ. 2004. New loci from New Zealand black and New Zealand white on chromosomes 4 and 12 contribute to lupus-like disease in the context of BALB/c¹. *J Immunol.* 172:4609-17.
- Sengupta TK, Bandyopadhyay S, Fernandes DJ, Spicer EK. 2004. Identification of nucleolin as an AU-rich element binding protein involved in bcl-2 mRNA stabilization. *J Biol Chem.* 279(12):10855-63.
- Shepherd BR, Chen HYS, Smith CM, Gruionu G, Williams SK, Hoying JB. 2004. Rapid perfusion and network remodeling in a microvascular construct after implantation. *Arterioscler Thromb Vasc Biol.* 24:898-904.
- Tan MGK, Ooi LLPJ, Aw SE, Hui KM. 2004. Cloning and identification of hepatocellular carcinoma down-regulated mitochondrial carrier protein, a novel liver-specific uncoupling protein. *J Biol Chem.* 279(43):45235-44.
- Gruber HE, Ingram JA, Leslie K, Norton HJ, Hanley EN. 2003. Cell shape and gene expression in human intervertebral disc cells: in vitro tissue engineering studies. *Biotechnic and Histochem.* 78(2):109-17.
- Imai KS. 2003. Isolation and characterization of β -catenin downstream genes in early embryos of the ascidian *Ciona savignyi*. *Differentiation.* 71:346-60.