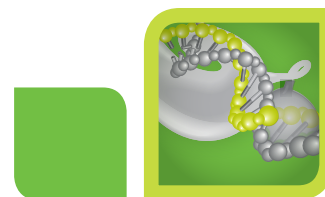


PrimePCR™ Publications List



PrimePCR Assays and Panels

Bulletin 6734



Top Ten PrimePCR Publications

Baker BA et al. (2014).

Ontology analysis of global gene expression differences of human bone marrow stromal cells cultured on 3D scaffolds or 2D films.
Biomaterials 35, 6,716–6,726.



Miotke L et al. (2014).

High sensitivity detection and quantitation of DNA copy number and single nucleotide variants with single color droplet digital PCR.
Anal Chem 86, 2,618–2,624.



Oestreich KJ et al. (2014).

Bcl-6 directly represses the gene program of the glycolysis pathway.
Nat Immunol 15, 957–964.



Poloni A et al. (2015).

Biosafety evidence for human dedifferentiated adipocytes.
J Cell Physiol 230, 1,525–1,533.



Rivas DA et al. (2014).

Diminished skeletal muscle microRNA expression with aging is associated with attenuated muscle plasticity and inhibition of IGF-1 signaling.
FASEB J 28, 4,133–4,147.



Sahler J et al. (2014).

Microparticles engineered to highly express peroxisome proliferator-activated receptor- γ decreased inflammatory mediator production and increased adhesion of recipient monocytes.
PLoS One 9, e113189.



SEQC/MAQC-III Consortium (2014).

A comprehensive assessment of RNA-seq accuracy, reproducibility and information content by the Sequencing Quality Control Consortium.
Nat Biotechnol 32, 903–914.



Supernat A et al. (2013).

Epithelial-mesenchymal transition and cancer stem cells in endometrial cancer.
Anticancer Res 33, 5,461–5,469.



Van Deun J et al. (2014).

The impact of disparate isolation methods for extracellular vesicles on downstream RNA profiling.
J Extracell Vesicles 3, 10.3402.



Zhang X et al. (2014).

Flavonoid apigenin inhibits lipopolysaccharide-induced inflammatory response through multiple mechanisms in macrophages.
PLoS One 9, e107072.





PrimePCR Publications by Date

2015

Arkatkar T et al. (2015).

Murine microRNA-214 regulates intracellular adhesion molecule (ICAM1) gene expression in genital *Chlamydia muridarum* infection. Immunology 145, 534–542.

**Baldrige MG et al. (2015).**

Very low-dose (femtomolar) 2,3,7,8-tetrachlorodibenzo-p-dioxin (TCDD) disrupts steroidogenic enzyme mRNAs and steroid secretion by human luteinizing granulosa cells. Reprod Toxicol 52, 57–61.

**Bol GM et al. (2015).**

Targeting DDX3 with a small molecule inhibitor for lung cancer therapy. EMBO Mol Med 7, 648–669.

**Bouch S et al. (2015).**

Neonatal exposure to mild hyperoxia causes persistent increases in oxidative stress and immune cells in the lungs of mice without altering lung structure. Am J Physiol Lung Cell Mol Physiol, Jul 2 [Epub ahead of print]. Accessed August 18, 2015.

**Croci S et al. (2015).**

Interleukin-15 is required for immunosurveillance and immunoprevention of HER2/neu-driven mammary carcinogenesis. Breast Cancer Res 17, 70.

**Ding Z et al. (2015).**

Cross-talk between LOX-1 and PCSK9 in vascular tissues. Cardiovasc Res 107, 556–567.

**Ding Z et al. (2015).**

Hemodynamic shear stress via ROS modulates PCSK9 expression in human vascular endothelial and smooth muscle cells and along the mouse aorta. Antioxid Redox Signal 22, 760–761.

**Fu J et al. (2015).**

STING agonist formulated cancer vaccines can cure established tumors resistant to PD-1 blockade. Sci Transl Med 7, 283ra52.

**Gibellini L et al. (2015).**

Different origin of adipogenic stem cells influences the response to antiretroviral drugs. Exp Cell Res, Jul 31 [Epub ahead of print]. Accessed August 18, 2015.

**Gosau M et al. (2015).**

Evaluation of implant-materials as cell carriers for dental stem cells under in vitro conditions. Int J Impl Dent 1, 2.

**Hui CC et al. (2015).**

Thymic stromal lymphopoietin (TSLP) secretion from human nasal epithelium is a function of TSLP genotype. Mucosal Immunol 8, 993–999.

**Jones AV et al. (2015).**

Evaluation of methods to detect CALR mutations in myeloproliferative neoplasms. Leuk Res 39, 82–87.

**Kozicky LK et al. (2015).**

Intravenous immunoglobulin skews macrophages to an anti-inflammatory, IL-10-producing activation state. J Leukoc Biol, Jul 27 [Epub ahead of print]. Accessed August 18, 2015.



**Merna N et al. (2015).**

Differential $\beta 3$ integrin expression regulates the response of human lung and cardiac fibroblasts to extracellular matrix and its components.
Tissue Eng Part A 21, 2,195–2,205.

**Molina SA et al. (2015).**

Junctional abnormalities in human airway epithelial cells expressing delF508 CFTR.
Am J Physiol Lung Cell Mol Physiol, Jun 26 [Epub ahead of print]. Accessed August 18, 2015.

**Opata MM et al. (2015).**

Early effector cells survive the contraction phase in malaria infection and generate both central and effector memory T cells.
J Immunol 194, 5,346–5,354.

**Qu T et al. (2015).**

Complete pulpodentin complex regeneration by modulating the stiffness of biomimetic matrix.
Acta Biomater 16, 60–70.

**Sung J et al. (2015).**

Garlic (*Allium sativum*) stimulates lipopolysaccharide-induced tumor necrosis factor-alpha production from J774A.1 murine macrophages.
Phytother Res 29, 288–294.

**Suzuki A et al. (2015).**

WNT/ β -catenin signaling regulates multiple steps of myogenesis by regulating step-specific targets.
Mol Cell Biol 35, 1,763–1,776.

**Uchino Y et al. (2015).**

Alteration of galectin-3 in tears of patients with dry eye disease.
Am J Ophthalmol 159, 1,027–1,035.

**Wenzel UA et al. (2015).**

CD103+CD11b+ dendritic cells induce Th17 T cells in Muc2-deficient mice with extensively spread colitis.
PLoS One 10, e0130750.

**Xie G et al. (2015).**

A West Nile virus NS4B-P38G mutant strain induces cell intrinsic innate cytokine responses in human monocytic and macrophage cells.
Vaccine 33, 869–878.

**Zhang JD et al. (2015).**

Pathway reporter genes define molecular phenotypes of human cells.
BMC Genomics 16, 342.

**2014****Galoian K et al. (2014).**

Lost miRNA surveillance of Notch, IGFR pathway — road to sarcomagenesis.
Tumour Biol 35, 483–492.

**Kroll CM et al. (2014).**

Enhanced resistance of CXCR3 deficient mice to ocular HSV-1 infection is due to control of replication in the brain ependyma.
J Neuroimmunol 276, 219–223.

**Landi A et al. (2014).**

Genome-wide shRNA screening identifies host factors involved in early endocytic events for HIV-1-induced CD4 down-regulation.
Retrovirology 11, 118.



**Leigh NJ et al. (2014).**

Expression of resolvin D1 biosynthetic pathways in salivary epithelium.
J Dent Res 93, 300–305.

**Lewarchik CM et al. (2014).**

The ryanodine receptor is expressed in human pancreatic acinar cells and contributes to acinar cell injury.
Am J Physiol Gastrointest Liver Physiol 307, G574–G581.

**Li S et al. (2014).**

Multi-platform assessment of transcriptome profiling using RNA-seq in the ABRF next-generation sequencing study.
Nat Biotechnol 32, 915–925.

**Petrie MA et al. (2014).**

A minimal dose of electrically induced muscle activity regulates distinct gene signaling pathways in humans with spinal cord injury.
PLoS One 9, e115791.

**Ramon S et al. (2014).**

Lipoxin A₄ modulates adaptive immunity by decreasing memory B-cell responses via an ALX/FPR2-dependent mechanism.
Eur J Immunol 44, 357–369.

**Schacter JL et al. (2014).**

Estrogen regulation of anti-apoptotic Bcl-2 family member Mcl-1 expression in breast cancer cells.
PLoS One 9, e100364.

**Woodward AM and Argüeso P (2014).**

Expression analysis of the transmembrane mucin MUC20 in human corneal and conjunctival epithelia.
Invest Ophthalmol Vis Sci 55, 6,132–6,138.

**Yang AX et al. (2014).**

Molecular characterization of antigen-peptide pulsed dendritic cells: Immature dendritic cells develop a distinct molecular profile when pulsed with antigen peptide.
PLoS One 9, e86306.

**2013****Minor EA et al. (2013).**

AMD-associated variants at the chromosome 10q26 locus and the stability of ARMS2 transcripts.
Invest Ophthalmol Vis Sci 54, 5,913–5,919.



Visit bio-rad.com/web/PrimePCRpublished for a wide range of videos, technical reports, published findings, and other resources describing Bio-Rad's PrimePCR Assays and Panels.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 **Belgium** 03 710 53 00 **Brazil** 55 11 3065 7550
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 49 89 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026674 55 05
Taiwan 886 2 2578 7189 **Thailand** 1800 88 22 88 **United Kingdom** 020 8328 2000