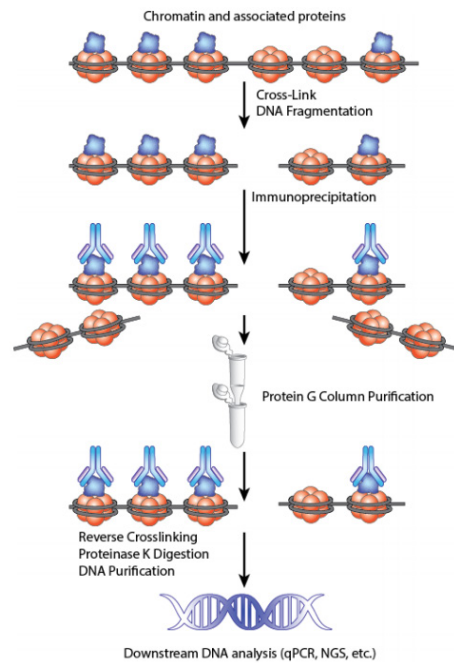


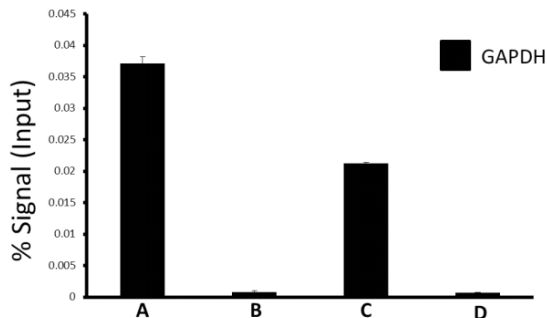
Go-ChIP-Grade™ Protein G Enzymatic Kit

BioLegend is proud to introduce the Go-ChIP-Grade™ Protein G Enzymatic Kit that provides a quicker, easier, and more efficient way of performing Chromatin Immunoprecipitation (ChIP) assays. The kit offers all of the major components required from chromatin sample preparation to immunoprecipitation (IP), protein-DNA complex purification, reverse-crosslinking, and DNA purification. The kit is based on the solid state technology from Chromatrap® that allows high-throughput screening from small cell numbers and low chromatin concentrations. The columns used in the kit contain discs of an inert, porous polymer to which Protein G has been covalently bound to maximize the capture efficiency of the target chromatin/antibody complex.

Principle of the ChIP assay



Anti-RNA Polymerase II Antibody



ChIP was performed using Go-ChIP-Grade™ Protein G Enzymatic Kit by loading 3 µg of crosslinked chromatin samples from HeLa cells with either A) 1:300 dilution of Go-ChIP-Grade™ Purified anti-RNA Polymerase II Antibody (Clone 8WG16), B) equal amount of Purified Mouse IgG2a Isotype Control Antibody, or C) competitor's ChIP-grade Purified anti-RNA Polymerase II Antibody and D) equal amount of matched Isotype Control Antibody as recommended by the manufacturer. The enriched DNA was purified and quantified by real-time qPCR using primers targeting human GAPDH gene region. The amount of IP'd DNA in each sample is represented as signal relative to the 5% of total amount of input chromatin.

Advantages of the Kit

- Compatible for use with ChIP-qPCR, ChIP-on-chip, and ChIP-seq.
- Suitable for ChIP assays from as little as 1 µg, to up to 50 µg of chromatin.
- ChIP assay can be completed in 1-2 days.
- The polymer disc is inert, reducing nonspecific binding.
- Provides reduced IP incubation times.
- Elution chemistry is optimized for high quality and quantity of IP'd DNA. Sensitive and provides selective enrichment of low chromatin loading.

Learn more at: biolegend.com/chip

Come and Meet Us!

January 27-30, 2018
Midwinter Conference of Immunologists
Pacific Grove, CA

February 14-18, 2018
Translational Research Cancer Centers Consortium
Seven Springs, PA

March 14-17, 2018
WIRM 2018
Davos, Switzerland

March 21-26, 2018
ThymOz International Workshop on T Lymphocytes
Heron Island, Australia

Talkin' Immunology Podcast

In our latest podcast, we speak with a special guest, Dr. Vijay Kuchroo, covering cancer research, innate lymphoid cells, and the integration of immune cells into multiple systems.



Listen to the podcast: biolegend.com/podcast

Neuroscience Spotlight

New Tau Antibodies

Tau is a family of neuronal proteins that bind to microtubules and stabilize their formation and maintenance. In the human brain, Tau proteins have 6 isoforms produced by alternative splicing of a single gene called MAPT. Research interest in tau proteins began to grow when tangled forms of these proteins were found to make up the paired helical filaments in brains of Alzheimer's disease (AD) patients. BioLegend has several new purified and conjugated formats of antibodies against Tau, its modifications, and isoforms.

| Description | Size | Cat. No. |
|-------------------------------------|----------------|-----------------|
| Biotin anti-Tau, 1-100 Antibody | 25 µg 100 µg | 816611 816612 |
| Purified anti-Tau, 1-223 Antibody | 25 µg 100 µg | 851301 851302 |
| Purified anti-Tau, 359-373 Antibody | 25 µg 100 µg | 851401 851402 |
| Purified anti-Tau, 368-441 Antibody | 25 µg 100 µg | 851201 851202 |
| Purified anti-Tau, 419-433 Antibody | 25 µg 100 µg | 851001 851002 |
| Purified anti-Tau, 425-441 Antibody | 25 µg 100 µg | 851101 851102 |

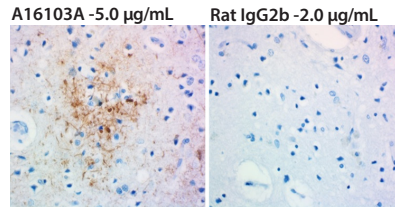
Learn more at: biolegend.com/tau

New Human LEGENDplex™ Kits

Hematopoietic stem cells (HSCs) give rise to myeloid, erythroid, and lymphoid lineages of cells, which further differentiate into specialized blood cells with the help of cytokines, growth factors, and chemokines. Dysregulation of these lineage-specific markers can lead to a hematological condition. HSCs develop in the bone marrow, where bone metabolism is a complex equilibrium between bone formation (by osteoblasts) and resorption (by osteoclasts). This is regulated through numerous hormones, steroids, growth factors, and cytokines. Disruptions to the homeostasis between formation and resorption results in diseases such as osteoporosis and Paget's disease. As research reveals the importance of these cytokines, our selection of LEGENDplex™ panels expands to match researchers' needs and analyze multiple cytokines simultaneously in stem cell and bone homeostasis studies.

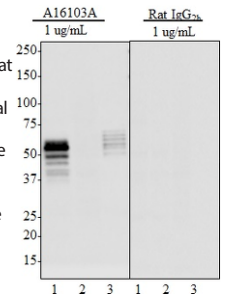
Learn more about LEGENDplex™ and watch a full protocol video at: biolegend.com/legendplex

Purified anti-Tau, 1-223 Antibody

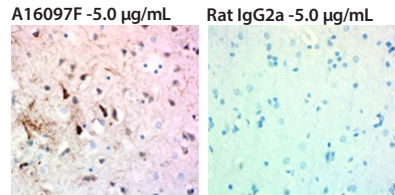


IHC staining of anti-Tau, 1-223 antibody (clone A16103A) and rat IgG2b isotype control on formalin-fixed paraffin-embedded Alzheimer's disease brain tissue. Following antigen retrieval using Sodium Citrate, the tissues were incubated with anti-Tau or rat IgG2b at 0.2 µg/mL for 1 hour at room temperature. Biotinylated anti-rat IgG, HRP Streptavidin, and DAB substrate were used as the detection system. Slides were counterstained with hematoxylin.

Western blots of anti-Tau, 1-223 antibody (clone A16103A) and rat IgG2b isotype control. Lane 1: 20 µg of normal human brain lysate; Lane 2: 20 µg of mouse brain lysate; Lane 3: 50 ng of recombinant human Tau ladder. The blots were incubated overnight at 4°C with 1 µg/mL of anti-Tau or rat IgG2b, followed by incubation with horseradish peroxidase labeled goat anti-rat IgG secondary antibody. Enhanced chemiluminescence was used as the detection system.

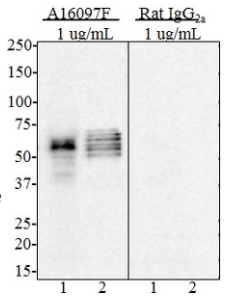


Purified anti-Tau, 368-441 Antibody



IHC staining of anti-Tau, 368-441 antibody (clone A161097F) and rat IgG2a isotype control on formalin-fixed paraffin-embedded Alzheimer's disease brain tissue. Following antigen retrieval using Sodium Citrate, the tissues were incubated with 5.0 µg/mL of anti-Tau or rat IgG2a overnight at 4°C. Biotinylated anti-rat IgG, HRP Streptavidin, and DAB substrate were used as the detection system. Slides were counterstained with hematoxylin.

Western blots of anti-Tau, 368-441 antibody (clone A161097F) and rat IgG2a isotype control. Lane 1: 20 µg of normal human brain lysate; Lane 2: 50 ng of recombinant human Tau ladder. The blots were incubated overnight at 4°C with 1 µg/mL of anti-Tau or rat IgG2a, followed by incubation with horseradish peroxidase labeled goat anti-rat IgG secondary antibody. Enhanced chemiluminescence was used as the detection system.



Human Hematopoietic Stem Cell Panel

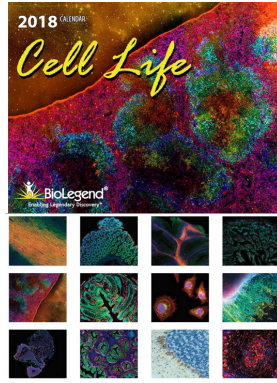
| Target | Bead ID | Hematopoietic Stem Cell Panel | HSC Lymphoid Panel | HSC Erythroid Panel | HSC Myeloid Panel |
|---------------|--------------------------|-------------------------------|--------------------|---------------------|-------------------|
| IL-6 | A4 | ✓ | ✓ | ✓ | |
| M-CSF | B7 | ✓ | | | ✓ |
| IL-15 | B6 | ✓ | ✓ | | |
| CXCL12 (SDF1) | B5 | ✓ | ✓ | ✓ | ✓ |
| LIF | B4 | ✓ | ✓ | | |
| SCF | B3 | ✓ | ✓ | ✓ | ✓ |
| IL-11 | B2 | ✓ | ✓ | ✓ | |
| IL-34 | A10 | ✓ | | | ✓ |
| IL-3 | A8 | ✓ | ✓ | ✓ | ✓ |
| GM-CSF | A6 | ✓ | | ✓ | ✓ |
| FLT3L | A5 | ✓ | ✓ | | ✓ |
| IL-7 | B9 | ✓ | ✓ | | |
| Options | Kit with Filter Plates | 740610 | 740612 | 740614 | 740616 |
| Options | Kit with V-bottom Plates | 740611 | 740613 | 740615 | 740617 |

Human Bone Metabolism Panel

| Target | Bead ID | Bone Metabolism | Bone Resorption | Bone Formation |
|---------|--------------------------|-----------------|-----------------|----------------|
| OPG | A4 | ✓ | | ✓ |
| OPN | A5 | ✓ | ✓ | |
| ALPL | A7 | ✓ | | ✓ |
| ACP5 | A8 | ✓ | ✓ | |
| Leptin | A10 | ✓ | | ✓ |
| RANKL | B2 | ✓ | ✓ | |
| TNF-α | B3 | ✓ | ✓ | |
| IL-6 | B4 | ✓ | ✓ | |
| PTH | B5 | ✓ | ✓ | |
| IL-1β | B6 | ✓ | ✓ | |
| BMP-2 | B7 | ✓ | | ✓ |
| DKK-1 | B9 | ✓ | ✓ | |
| Options | Kit with Filter Plates | 740655 | 740674 | 740672 |
| Options | Kit with V-bottom Plates | 740656 | 740675 | 740673 |

2018 Cell Life Calendar

Plan out your 2018 with our new Cell Life Calendar. Each image is generated using BioLegend microscopy reagents and features extraordinary and unique staining.



Request your calendar at:
biolegend.com/literature/brochures

New Year Promotion

LAB TIMER PROMOTION

Get a complimentary Lab Timer when using promo code **TIME18** on orders over \$300 USD. For European orders, value must be over €250 EURO, £250 GBP, or 300 CHF.



Valid for direct end users in USA, Canada, UK, Ireland, Germany, Austria, Switzerland, France, Belgium, Netherlands, and Luxembourg. Promotion and discount codes must be referenced when placing your order. Order must total \$300+ after discounts, before shipping charges. Not valid for distributors or re-sellers. One gift per order, valid until 3/31/18 or while supplies last. Void where prohibited.

View more promotions at: biolegend.com/promotions

BioLegend | New Products Released

Human Flow Cytometry Antibodies

| Description | Clone | Format | Applications |
|-------------------------------------|-------------------------------------|--|--------------|
| BTLA | MIH26 | Purified | FC |
| CD3/CD4/CD8/CD16/CD19/CD45/CD56 | SK7/5.1H11/B73.1/2D1/SK3/SJ25C1/SK1 | Cocktail | FC |
| CD4 | A161A1 | Biotin, Brilliant Violet 421™ | FC |
| Human CD4 Treg Surface ID Panel | A019D5, BC96, UCHT1, A161A1 | | FC |
| Human Activated CD8 T Cell ID Panel | UCHT1, RPA-T8, BC96, UCHL1 | | FC |
| CD9 | H19a | APC | FC |
| CD11a | H1111 | PE/Dazzle™ 594, PerCP/Cy5.5 | FC |
| CD14 | 63D3 | Alexa Fluor® 488, Alexa Fluor® 647, Brilliant Violet 605™ | FC |
| CD18 | m24 | APC/Fire™ 750, PE, PerCP/Cy5.5 | FC |
| CD22 | HIB22 | Alexa Fluor® 700, APC/Fire™ 750 | FC |
| CD33 | P67.6 | Alexa Fluor® 647, Brilliant Violet 711™ | FC |
| CD36L1 | m1B9 | APC, PerCP/Cy5.5 | FC |
| CD42b | HIP1 | PE/Dazzle™ 594 | FC |
| CD44 | BJ18 | Pacific Blue™ | FC |
| CD45 | 2D1 | Alexa Fluor® 488, Biotin | FC |
| CD45 | OX-1 | Pacific Blue™, PE/Dazzle™ 594 | FC |
| CD47 | CC2C6 | Brilliant Violet 421™, Brilliant Violet 605™, Pacific Blue™ | FC |
| CD74 | Pin.1 | PE/Cy7, PerCP/Cy5.5 | ICFC |
| CD85a (ILT5) | MKT5.1 | APC | FC |
| CD86 | BU63 | APC, FITC, PE | FC |
| CD99 | 3B2/TA8 | PE/Cy7, PerCP/Cy5.5 | FC |
| CD107b (LAMP-2) | H4B4 | Alexa Fluor® 647, Brilliant Violet 421™ | ICFC |
| CD117 (c-kit) | 104D2 | Brilliant Violet 650™ | FC |
| CD122 | TU27 | PE/Dazzle™ 594 | FC |
| CD155 (PVR) | SK1L4 | Alexa Fluor® 647, Biotin | FC |
| CD169 | 7-239 | Brilliant Violet 605™ | FC |
| CD172a | 15-414 | FITC | FC |
| CD195 (CCR5) | J418F1 | Brilliant Violet 510™ | FC |
| CD209 (DC-SIGN) | 9E9A8 | APC/Fire™ 750 | FC |
| CD223 | 11C3C65 | Alexa Fluor® 488, Biotin, Brilliant Violet 605™ | FC |
| CD226 (DNAM-1) | 11A8 | Brilliant Violet 605™, Brilliant Violet 785™, PE/Dazzle™ 594 | FC |
| CD258 (LIGHT) | T5-39 | PE/Cy7 | FC |
| CD279 | NAT105 | Alexa Fluor® 647, Biotin, Brilliant Violet 421™, Brilliant Violet 510™ | FC |
| CD282 | W15145C | Purified | FC |
| CD360 | 17A12 | PE/Cy7 | FC |
| FcεR1α | AER-37 (CRA-1) | Alexa Fluor® 488 | FC |
| Folate Receptor β | 94b/FOLR2 | APC, PE | FC |
| Galectin-9 | 9M1-3 | Alexa Fluor® 647 | ICFC, FC |
| GL7 | GL7 | Biotin | FC |
| Granzyme A | CB9 | APC, PE/Cy7 | ICFC |
| HLA-A2 | BB7.2 | Brilliant Violet 421™, Brilliant Violet 785™ | FC |
| HLA-DR | LN3 | PerCP/Cy5.5 | FC |
| HLA-E | 3D12 | Brilliant Violet 421™ | FC |

Human Flow Cytometry Antibodies

| Description | Clone | Format | Applications |
|--------------------|------------|-----------------------|--------------|
| Ig light chain λ | 1-155-2 | Purified | FC |
| IgG Fc | M1310G05 | Biotin, FITC | FC |
| IL-22 | 2G12A41 | FITC | ICFC |
| LAP | TW4-2F8 | Alexa Fluor® 647 | FC |
| Mac-2 (Galectin-3) | M3/38 | APC, PE/Cy7 | FC |
| MERTK | 590H11G1E3 | APC/Fire™ 750, Biotin | FC |
| NTAL | W15102A | PE | ICFC |
| SSEA-3 | MC-631 | FITC | FC |
| TCR Vβ8 | JR2 (JR.2) | APC | FC |
| T-bet | 4B10 | Brilliant Violet 785™ | ICFC |
| Vβ13.1 | H131 | PE/Cy7 | FC |

Mouse Flow Cytometry Antibodies

| Specificity | Clone | Format | Applications |
|-------------------------|------------|---|--------------|
| β2-microglobulin | A16041A | APC, PE | FC |
| CD1d (CD1.1, Ly-38) | 1B1 | PE/Cy7 | FC |
| CD14 | Sa14-2 | Alexa Fluor® 647, Brilliant Violet 421™, PE/Dazzle™ 594 | FC |
| CD20 | SA275A11 | Alexa Fluor® 700 | FC |
| CD31 | MEC13.3 | APC/Fire™ 750 | FC |
| CD34 | HM34 | PE/Cy7 | FC |
| CD125 | DIH37 | APC | FC |
| CD140a | APA5 | Brilliant Violet 605™ | FC |
| CD146 | ME-9F1 | Alexa Fluor® 647, Biotin | FC |
| CD184 | QA16A08 | Ultra-LEAF™ | FC, Block |
| CD186 | SA051D1 | Alexa Fluor® 647, Biotin, Brilliant Violet 711™, PE/Cy7, PE/Dazzle™ 594 | FC |
| CD215 | 6B4C88 | PE, Purified | FC |
| CD355 | 11-5/CRTAM | Biotin, PE/Cy7, PerCP/Cy5.5 | FC |
| CX3CR1 | QA16A03 | Ultra-LEAF™ | FC, Block |
| Folate Receptor β | 10/FR2 | APC, PE | FC |
| FR4 (Folate Receptor 4) | 12A5 | APC/Fire™ 750, PE/Dazzle™ 594, PerCP/Cy5.5 | FC |
| GL7 | GL7 | Biotin | FC |
| GM-CSF | MP1-22E9 | APC, PE/Cy7 | ICFC |
| H-2Kb | AF6-88.5 | PE/Cy7 | FC |
| I-Ab | AF6-120.1 | APC/Fire™ 750 | FC |
| IgG2a | RMG2a-62 | Alexa Fluor® 488, APC/Fire™ 750 | FC |
| LAP (TGF-β1) | TW7-16B4 | FITC | FC |
| Mac-2 (Galectin-3) | M3/38 | APC, PE/Cy7 | FC |
| P2X7R | 1F11 | PE/Cy7 | FC |
| Perforin | S16009A | APC, PE | ICFC |
| Perforin | S16009B | APC, PE, Purified | ICFC |
| SSEA-3 | MC-631 | FITC | FC |
| STAT3 Phospho (Ser727) | A16089B | PE | ICFC |
| TNF-α | TN3-19.12 | APC | ICFC |
| VISTA | MIH63 | Biotin | FC |
| TCR Va2 | B20.1 | PE/Cy7 | FC |
| ZAP-70 | A16043B | Alexa Fluor® 647 | ICFC |

Rat Flow Cytometry Antibodies

| Description | Clone | Format | Applications |
|-------------|-----------|--------|--------------|
| SiglecE | M1304A01 | FITC | FC |
| TNF-α | TN3-19.12 | APC | ICFC |

BioLegend | New Products Released

LEGENDplex™ Products

| Description | Applications |
|---|--------------|
| Buffer Set H | Multiplex |
| Human Hematopoietic Stem Cell Panel (12-plex) with Filter Plate | Multiplex |
| Human Hematopoietic Stem Cell Panel (12-plex) with V-bottom Plate | Multiplex |
| Human HSC Lymphoid Panel (9-plex) with Filter Plate | Multiplex |
| Human HSC Lymphoid Panel (9-plex) with V-bottom Plate | Multiplex |
| Human HSC Erythroid Panel (6-plex) with Filter Plate | Multiplex |
| Human HSC Erythroid Panel (6-plex) with V-bottom Plate | Multiplex |
| Human HSC Myeloid Panel (7-plex) with Filter Plate | Multiplex |
| Human HSC Myeloid Panel (7-plex) with V-bottom Plate | Multiplex |
| Human Bone Metabolism Panel 1 (12-plex) with Filter Plate | Multiplex |
| Human Bone Metabolism Panel 1 (12-plex) with V-bottom Plate | Multiplex |
| Human Bone Resorption Panel (8-plex) with Filter Plate | Multiplex |
| Human Bone Resorption Panel (8-plex) with V-bottom Plate | Multiplex |
| Human Bone Formation Panel (4-plex) with Filter Plate | Multiplex |
| Human Bone Formation Panel (4-plex) with V-bottom Plate | Multiplex |

Human Recombinant Proteins

| Description | Format | Applications |
|------------------|-------------|--------------|
| ALCAM/CD166 | Recombinant | BA |
| CD6 | Recombinant | BA |
| ICAM-2 | Recombinant | BA |
| Serpin E2 | Recombinant | BA |
| TIGIT-Fc Chimera | Recombinant | BA |

Mouse Recombinant Proteins

| Description | Format | Applications |
|--------------------------|-------------|--------------|
| BMPRI1A/ALK-3-Fc Chimera | Recombinant | BA |
| PDGFRβ | Recombinant | BA |
| TRANCE | Recombinant | BA |

Microscopy Products

| Specificity | Clone | Format | Applications |
|-----------------|----------|------------------------------------|--------------|
| ATF7 | W16152A | Alexa Fluor® 594 | IF |
| Cadherin 11 | 16G5 | Alexa Fluor® 647 | IHC-P |
| CD107b (LAMP-2) | H4B4 | Alexa Fluor® 594 | IHC-P |
| c-Myc | 9E10 | Alexa Fluor® 594 | IF |
| Cytokeratin 7 | W16155A | Alexa Fluor® 594 | IF |
| Cytokeratin 17 | W16131A | Alexa Fluor® 594 | IF |
| Galectin-9 | 9M1-3 | Alexa Fluor® 594 | IHC-P |
| IκB-α | 3D6C02 | Alexa Fluor® 647 | IF |
| Keratin 5 | Poly9059 | Purified | IHC-P |
| Ki-67 | 11F6 | Alexa Fluor® 594 | IHC-F |
| Rat IgG1 | MRG1-58 | Alexa Fluor® 594 | IHC-F |
| Rat IgG2b | MRG2b-85 | Alexa Fluor® 488, Alexa Fluor® 647 | IHC-F |
| Lypd8 | SA025A4 | Alexa Fluor® 594, Alexa Fluor® 647 | IHC-F |
| Lypd8 | SA024B4 | Alexa Fluor® 594, Alexa Fluor® 647 | IHC-F |
| MAP2 | SMI 52 | Alexa Fluor® 594 | IHC-P |
| NR1D2 | W15136A | Alexa Fluor® 594 | IF |
| PARP | 5A5 | Alexa Fluor® 647 | IF, IHC-P |

Alexa Fluor® and Pacific Blue™ are registered trademarks of Life Technologies Corporation. Brilliant Violet™ is a trademark of Sirigen Group Ltd. Chromatrap™ Technology: This patented format is unique and has been granted in the UK (Patent No. GB2482209), the US (Patent No. 9523681), China (Patent No. ZL 2011 8 0067254.X) Japan (Patent No. JP 6088434) and Australia (Patent No. AU 2011340263).

Go-CHIP-Grade™ Products

| Description | Clone | Format | Applications |
|-------------------------|-----------|-------------------------|---------------------|
| Bcl-6 | IG191E/A8 | Go-CHIP-Grade™ Purified | ChIP, WB, IHC-P, IP |
| Ikaros | 16B5C71 | Go-CHIP-Grade™ Purified | ChIP, WB |
| IRF3 | 12A4A35 | Go-CHIP-Grade™ Purified | ChIP, WB, IF |
| IRF9 | 5A3A39 | Go-CHIP-Grade™ Purified | ChIP, WB, IF |
| NFATc1 | 7A6 | Go-CHIP-Grade™ Purified | ChIP, WB, ICC, ICFC |
| Spt16 (FACT140 complex) | 8D2 | Go-CHIP-Grade™ Purified | ChIP, ICC, WB |
| STAT2 | W15088A | Go-CHIP-Grade™ Purified | ChIP, WB |
| Protein G Enzymatic Kit | | Go-CHIP-Grade™ Purified | ChIP |

Isotype Controls

| Specificity | Clone | Format | Applications |
|----------------------------------|---------|----------------|--------------|
| Syrian Hamster IgG, Isotype Ctrl | SHG-1 | PE/Dazzle™ 594 | FC, ICFC |
| Human IgG1, Isotype Ctrl | QA16A12 | APC, FITC, PE | FC |
| Human IgG2, Isotype Ctrl | QA16A13 | FITC, PE | FC |
| Human IgG4, Isotype ctrl | QA16A15 | APC, FITC, PE | FC |
| Mouse IgG2b, κ Isotype Ctrl | 27-35 | APC/Cy7, FITC | FC |
| Mouse IgM, κ Isotype Ctrl | MM-30 | PE/Cy7 | FC |

Human ELISA Products

| Description | Clone | Format | Applications |
|-------------|----------|----------|-----------------|
| CCL20 | L2412F11 | Purified | ELISA Capture |
| CCL20 | L2418G3 | Biotin | ELISA Detection |

Neuroscience Products

| Specificity | Clone | Format | Applications |
|-----------------------------|-----------|--|---------------------------|
| α-Synuclein Phospho (Ser12) | P-syn/81A | Alexa Fluor® 594 | IHC-P, IHC-F, WB |
| β-Amyloid, 17-24 | 4G8 | Alexa Fluor® 488, Alexa Fluor® 594, Alexa Fluor® 647 | IHC-P |
| Connexin 43 | 2F5 | Purified | WB, IHC-P |
| DJ-1 | A16125E | Purified | WB, Direct ELISA, IHC-P |
| SNAP-25 | SMI 81 | Biotin | WB, IHC-P, EM, FC, IF, IP |
| Tau, 1-100 | 43D | Alexa Fluor® 488, Alexa Fluor® 647, Biotin | WB, IHC-P, IHC, WB |
| Tau, 1-223 | A16103A | Purified | WB, Direct ELISA, IHC-P |
| Tau, 359-373 | A16097B | Purified | WB, Direct ELISA, IHC-P |
| Tau, 368-441 | A16097F | Purified | WB, Direct ELISA, IHC-P |
| Tau, 419-433 | A16097D | Purified | WB, Direct ELISA, IHC-P |
| Tau, 425-441 | A16097E | Purified | WB, Direct ELISA, IHC-P |
| Tubulin β 3 (TUBB3) | TUJ1 | Biotin | WB, IHC-P, FC, IF |

Western Blot Antibodies

| Specificity | Clone | Format | Applications |
|--|---------|----------------------------|--------------|
| <i>E. coli</i> RNA Polymerase Antibody Sampler Kit | | | WB |
| β-actin | W16197A | Purified | WB |
| BTK Phospho (Tyr223)/ITK Phospho (Tyr180) | A16128C | Purified | WB, ICFC |
| His Tag | J095G46 | Alexa Fluor® 647, Purified | WB, IF |
| RUNX1 | W16118A | Purified | WB |
| SIRT2 | W16205A | Purified | WB |
| Stx2A | 1C6A6G9 | Purified | WB |
| Tata Binding Protein | 1TBP18 | Purified | WB |



Chromatrap

BioLegend products are manufactured in an ISO 13485:2003 certified facility following GMP compliant procedures, ensuring the highest quality and stability.

9727 Pacific Heights Blvd. | San Diego, CA 92121

1.877.246.5343 (877-BIOLEGEND)

International: 1.858.768.5800

biolegend.com

02-0001-58