



Monkey IL-17 Antibody Bead Kit

INFORMATION SHEET

Catalog #: LPC0171 Description: Monkey IL-17 Lot:* A10671

*Note: A letter at the end of the lot number signifies an additional packaging of this same lot.

Intended Use

This reagent set comprises the analyte specific components for the measurement of monkey IL-17 in serum, plasma, buffered solution or tissue culture supernatant. The assay may be run alone or in combination with other Antibody Bead Kits from Invitrogen. Buffer reagents needed to complete the reaction are sold separately under Catalog #LHB0001. These reagents are intended for use in the **Luminex® 100™ or 200™ System** only. **This kit is configured for research use only and is not to be used in diagnostic procedures.**

Reagents Provided

1. Antibody Bead Concentrate (10x):

Catalog #: LM208 Description: Ms x Mky IL-17 Lot: A10672 Size: 0.25 mL-100 tests

Bead Region: 25

Form: 0.25 mL 10x bead concentrate solution in storage buffer. Contains 7.5 mM sodium azide as preservative.

Storage: Store at 2 to 8°C until the expiration date indicated on the kit.

2. Biotinylated Antibody Concentrate (10x):

Catalog #: BN208 Description: Ms x Mky IL-17 biotin Lot: A10673 Size: 1 mL-100 tests

Form: 1 mL of a 10x stock of Biotinylated Antibody Concentrate in Biotin Diluent. Contains 15 mM sodium azide as preservative. Concentration of antibody is matched to this lot of beads. Do not mix lots of Coated Beads and Detection Antibody.

Storage: Store at 2 to 8°C until the expiration date indicated on the kit.

3. Monkey Cytokine II Five-Plex Standard (IL-1 β , IL-6, IL-12, IL-17, and G-CSF) (2 vials).

Catalog #: SM206 Description: Rec. Mky Cytokine II Five-Plex Std Lot: A10638 Size: Single Use

Form: Lyophilized. The proteins in this standard have been calibrated against the masses of highly purified recombinant proteins, and the respective Luminex kits.

Storage: Store at 2 to 8°C. Use within 1 hour after reconstitution. Discard immediately after use.

Concentration of Reconstituted Standards**:

IL-1 β (13,750 pg/mL)

IL-6 (7,580 pg/mL)

IL-12 (62,000 pg/mL)

IL-17 (7,360 pg/mL)

G-CSF (27,560 pg/mL)

****Important note: The concentrations of reconstituted standards are lot-specific. Please verify all concentration values entered in data analysis software.**

Reconstitution: Reconstitute with 1.0 mL Assay Diluent when measuring IL-17 in serum or plasma samples. For other sample types, such as tissue culture supernatants, reconstitute the standard in 1.0 mL of a solution consisting of 50% Assay Diluent + 50% of sample matrix. Allow standard to rehydrate for approximately 10 minutes before further dilution.

Recommended Starting Concentration for Standard Curve: Upon reconstitution, the starting concentration of standard is the value cited above. Make serial 1:3 dilutions in Assay Diluent (serum/plasma samples) or other appropriate matrix. Use 100 μ L per assay. If establishing a Multiplex Assay, this same standard can be used to measure the other related cytokines cited above in a Multiplex Assay format. See the Product Insert included in the Buffer Reagent Kit for further information.

This product is for research use only. Not for use in diagnostic procedures.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PILPC0171 (Rev 03/09)

(Page 1 of 2) DCC-09-0372

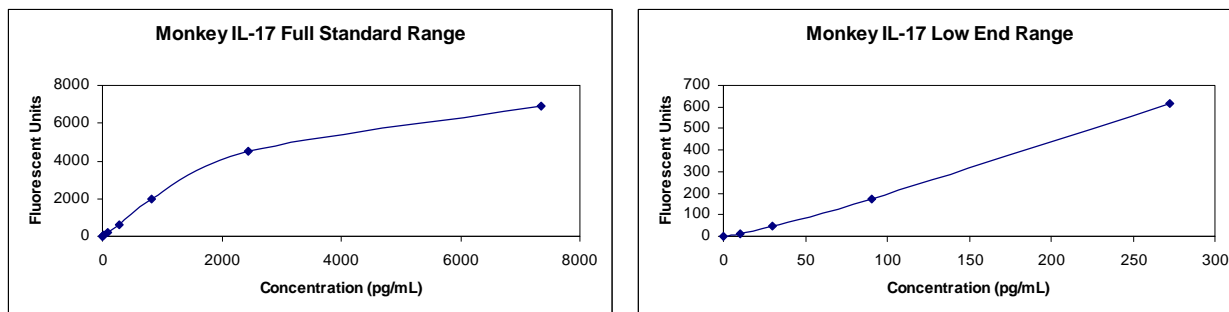
Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.

Monkey IL-17 Antibody Bead Kit

INFORMATION SHEET

Performance Characteristics

Analytical Sensitivity: The minimum detectable dose of monkey IL-17 is 5 pg/mL. This was determined by adding two standard deviations to the mean FI obtained when the zero standard was assayed 32 times.



Typical Standard Curve

Specificity: Buffered solutions of panels of substances at 5 to 10 ng/mL were assayed with the Invitrogen Monkey IL-17 Antibody Bead Kit. The following substances were tested and all were found to have no cross-reactivity: monkey IL-1 β , IL-2, IL-4, IL-6, IL-10, IL-12, G-CSF, IFN- γ , TNF- α , MCP-1, MIP-1 α , MIP-1 β , IL-8, RANTES; human adiponectin, BDNF, GDNF, GM-CSF, insulin, leptin, IFN- α , IFN- γ , IL-1 α , IL-1 β , IL-3, IL-4, IL-5, IL-6, IL-7, IL-8, IL-9, IL-10, IL-11, IL-12, IL-13, IL-15, IL-16, eotaxin, IP-10, MCP-1, MIG, MIP-1 α , MIP-1 β , RANTES, TNF- α , TGF- β 1, Gro- α , MCP-2, MCP-3, PDGF-BB, EGF, FGF basic, G-CSF, HGF, VEGF, DR5, IL-1RA, IL-2R, IL-4R, IL-6R, TNF-RI, TNF-RII; mouse FGF basic, GM-CSF, IFN- γ , IL-1 α , IL-1 β , IL-2, IL-4, IL-5, IL-6, IL-10, IL-12p40/p70, IL-13, IL-17, IP-10, KC, MCP-1, MIG, MIP-1 α , TNF- α , and VEGF; rat GM-CSF, IFN- γ , IL-1 α , IL-1 β , IL-2, IL-4, IL-6, IL-10, IL-12 and TNF- α .

Precision:

	Intra-assay (n=16)	Inter-assay (n=32)
Mean (pg/mL)	3437.6	3208.2
SD	222.2	308.5
%CV	6.46	9.62

Linearity: Monkey serum and tissue culture medium containing 10% fetal calf serum were spiked with monkey IL-17 and serially diluted in *Assay Diluent* and a solution consisting of 50% *Assay Diluent* and 50% tissue culture medium containing 10% fetal calf serum, respectively, over the range of the assay. Linear regression analysis of samples versus the expected concentration yielded a correlation coefficient of 0.99 and 0.98 for serum and tissue culture, respectively.

Recovery:

Cynomolgus monkey serum averaged 73.5%.

Cynomolgus citrate plasma averaged 73.6%; Heparinized plasma averaged 81.3%; EDTA plasma averaged 66.4%.

Rhesus monkey serum averaged 85.8%.

Rhesus citrate plasma averaged 90.1%; Heparin plasma averaged 76.7%; EDTA plasma averaged 84.9%.

Tissue culture medium containing 10% fetal calf serum averaged 105.7%.

By purchasing this Kit, which contains fluorescently labeled microsphere beads authorized by Luminex® Corporation ("Luminex®"), you, the customer, acquire the right under Luminex's patent rights to use this Kit or any portion of this Kit, including without limitation the microsphere beads contained herein, only with Luminex's laser based fluorescent analytical test instrumentation marketed under the name Luminex® 100™ or 200™. This product is covered by one or more of the following U.S. patents: 6,046,807.

This product is for research use only. Not for use in diagnostic procedures.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PILPC0171 (Rev 03/09)

(Page 2 of 2) DCC-09-0372

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.