



Human IL-12 p70 Antibody Bead Kit

INFORMATION SHEET

Catalog #: LHC9121 Description: Human IL-12 p70 Lot:* 489604

*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 human IL-12 p70 in serum, plasma 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 #: LM086 Description: Ms x Hu IL-12 p70 Lot: 390178 Size: 0.25 mL-100 tests

Bead Region: 42

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

Storage: **Light-sensitive material.** Store at 2 to 8°C in the dark, until the expiration date indicated on the kit.

2. Biotinylated Antibody Concentrate (10x):

Catalog #: BN086 Description: Ms x Hu IL-12 p70 biotin Lot: 390185 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. Hu Cytokine II Standard (2 vials):

Catalog #: SM0073 Description: Rec. Hu Cytokine II Std. Lot: 481485 Size: Single use

Form: Lyophilized. The proteins in this standard have been calibrated against the mass of highly purified recombinant protein, with the respective Invitrogen ELISA kit, and NIBSC calibration standard (if available). See the Product Insert included in the Buffer Reagent Kit for further information.

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

Concentration of Reconstituted Standard**:

IL-12 (22,290 pg/mL) IL-13 (21,810 pg/mL) IL-15 (20,950 pg/mL)
IL-17 (15,000 pg/mL) IFN- α (13,150 pg/mL) IL-16 (28,160 pg/mL)

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

One nanogram of Invitrogen recombinant human IL-12 equals 10 IU of WHO reference preparation 95/544 (NIBSC, Hertfordshire, UK, EN6 3QG).

NOTE: Human IL-12 has been recalibrated to NIBSC reference material and the above value for IL-12 reflects this recalibration. Correlation of results using the previous standard (Cat. # SM0073, Lot # R102104) with results obtained with this standard may be obtained by using a conversion factor.

[results obtained with previous standard] x 1.5 = [results obtained with this standard]

Reconstitution: Reconstitute with 1 mL Assay Diluent when measuring IL-12 p70 in serum or plasma samples. For other sample types, such as tissue culture supernatants, reconstitute the standard in 1 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

PILHC9121 (Rev 10/08)

(Page 1 of 2) DCC-08-1089/1288

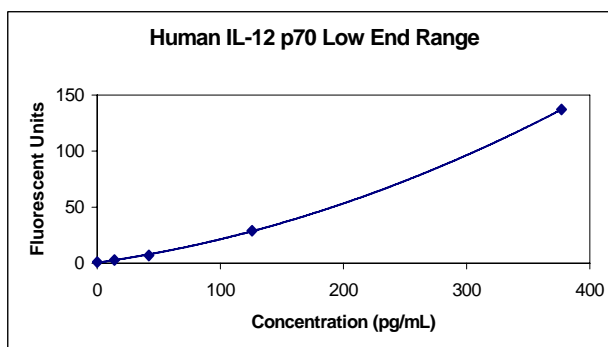
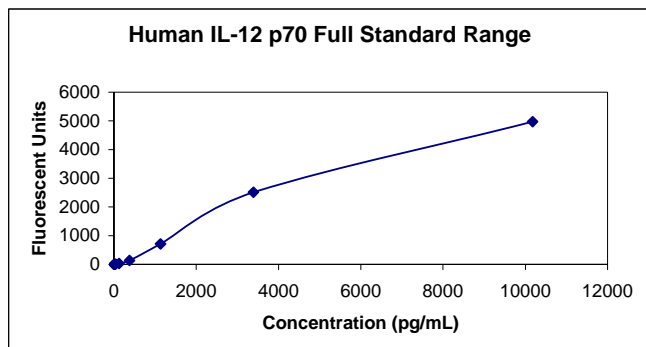
Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated,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

Human IL-12 p70 Antibody Bead Kit

INFORMATION SHEET

Performance Characteristics

Analytical Sensitivity: The minimum detectable dose of Hu IL-12 p70 is <10 pg/mL. This was determined by adding two standard deviations to the mean FI obtained when the zero standard was assayed 30 times.



Typical Standard Curve

Specificity: Buffered solutions of a panel of substances at 10 or 50 ng/mL were assayed with the Invitrogen Human IL-12 p70 Antibody Bead Kit. The following substances were tested and all were found to have no cross-reactivity: human IL-1 β , IL-2, IL-3, IL-4, IL-5, IL-6, IL-8, IL-10, IL-13, IL-15, IL-17, G-CSF, GM-CSF, IFN- γ , TNF- α , VEGF, FGF basic, HGF, EGF, IP-10, MIG, MCP-1, MIP-1 α , MIP-1 β , EOTAXIN, RANTES, IL-12 p40; mouse IL-12; rat IL-12.

Precision:

	Intra-assay (n=16)	Inter-assay (n=32)
Mean (pg/mL)	880.1	876.3
SD	41.6	35.2
%CV	4.7	4.0

Linearity: Human serum and tissue culture medium containing 10% fetal calf serum were spiked with human IL-12 p70 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 for both serum and tissue culture.

Recovery:

Human serum averaged 97% (range: 95% to 99.2%).

Human heparin plasma averaged 100% (range: 90% to 108.4%); EDTA plasma averaged 91.5% (range: 85.7% to 97.4%);

Citrate plasma is not recommended.

Tissue culture medium containing 10% fetal calf serum averaged 106.6% (range: 100% to 110.5%).

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

PILHC9121 (Rev 10/08)

(Page 2 of 2) DCC-08-1089/1288

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.