

**PRODUCT INSERT**

**CYTO-IC MONOCLONAL ANTIBODIES**

**RAT anti-MOUSE IL-12**

Product Code	Clone	Form	Volume	Antibody*	Excitation (nm)	Peak Emission (nm)	Matching Isotype Controls
RM9300	C17.8	Purified	1.0 ml	200 µg	N/A	N/A	Mouse IgG2a Purified Code MG2a00
RM9315	C17.8	Biotin	1.0 ml	100 µg	N/A	N/A	Mouse IgG2a Biotin Code MG2a15
RM9315-3	C17.8	Biotin	3.0 ml	300 µg			

**PRODUCT DESCRIPTION**

Rat monoclonal antibody to mouse IL-12

**Clone:** C17.8

**Isotype:** Rat IgG2a

**Lot No:** See label      **Expiration:** See label

**Immunogen:** CHO-expressed recombinant mouse IL-12 p70

**Preservatives:** 0.1% *sodium azide*. Sodium azide is an extremely toxic and dangerous compound particularly when combined with acids or metals. Solutions containing sodium azide should be disposed of properly.

**Stabilizer:** For conjugated products only, a highly purified grade of BSA has been added as a stabilizing protein.

**STORAGE AND HANDLING**

Store reagents at 2-8°C. Light exposure should be avoided when using fluorochrome-conjugated antibodies. Use dim light during handling, incubation with cells and prior to analysis. It is recommended that investigators dilute only the quantity of antibody to be used within one week.

**PRODUCT CHARACTERIZATION**




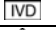



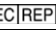
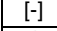
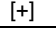



**Antigen Specificity:** The C17.8 monoclonal antibody (mAb) reacts with the p40 subunit of IL-12 p70, a heterodimeric cytokine<sup>1</sup>. C17.8 is a neutralizing mAb<sup>1</sup> that can also be used in ELISA and intracellular immunofluorescent staining for flow cytometry.

**REFERENCES**

1. Wysocka., M., M. Kubin, L. Q. Vieira, L. Ozmen, G. Garotta, P. Scott, and G. Trinchieri. 1995. Interleukin-12 is required for interferon-γ production and lethality in lipopolysaccharide-induced shock in mice. *Eur. J. Immunol.* 25: 672-676

\* The amount of antibody is determined by measuring the optical density using a spectrophotometer. The antibody titer is verified by immunofluorescent staining and flow cytometric analysis.

**Explanation of symbols**

Symbol	Description	Symbol	Description
	Catalogue Number		Batch code
	Research Use Only		<i>In vitro</i> diagnostic medical device
	Use by		Temperature limitation
	Manufacturer		European Community authorised representative
	Without, does not contain		With, contains
	Protect from light		Consult accompanying documents
	Directs the user to consult instructions for use (IFU), accompanying the product.		

**For research use only. CAUTION: Not for human or animal therapeutic or diagnostic use**

[www.invitrogen.com](http://www.invitrogen.com)

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

PI: L11309

(Rev 02/10) DCC-10-0338

**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](http://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.

**PRODUCT INSERT  
MONOCLONAL ANTIBODIES PAIR FOR ELISA TESTING**

**RAT anti-MOUSE IL-12**

Product Code	Clone	Form	Volume	Antibody*	Excitation (nm)	Peak Emission (nm)	Matching Isotype Controls	
RM9200	C15.6	Purified	1.0 ml	200 µg	N/A	N/A	Rat IgG-1 Purified	Code R100
RM9315	C17.8	Biotin	1.0 ml	100 µg	N/A	N/A	Mouse IgG-2a Biotin	Code MG2a15
RM9315-3	C17.8	Biotin	3.0 ml	300 µg				

**PRODUCT DESCRIPTION**

**Capture Antibody:** Rat monoclonal antibody to mouse IL-12

**Clone:** C15.6

**Isotype:** Rat IgG1

**Detection Antibody:** Rat monoclonal antibody to mouse IL-12

**Clone:** C17.8

**Isotype:** IgG2a

**Lot No.:** See label      **Expiration:** See label

**Buffer:** Phosphate buffered saline (PBS)

**Preservatives:** 0.1% *sodium azide*. Sodium azide is an extremely toxic and dangerous compound particularly when combined with acids or metals. Solutions containing sodium azide should be disposed of properly.

**Stabilizer:** For conjugated products only, a highly purified grade of BSA has been added as a stabilizing protein to bring the final protein concentration to 4-5 mg/ml after conjugation.

**STORAGE AND HANDLING**

Store reagents at 2-8°C. Light exposure should be avoided when using fluorochrome-conjugated antibodies. Use dim light during handling, incubation with cells and prior to analysis. It is recommended that investigators dilute only the quantity of antibody to be used within one week.

**PRODUCT QUALITY CONTROL**

**Use In Elisa:**

Clone C15.6 can be used as a capture antibody in a sandwich ELISA when paired with Caltag biotinylated IL-12 detection antibody clone C17.8 (RM9315). The antibody can be used in the range of 0.25 to 2.0 µg/ml. IL-12 capture antibody is known to have a substantial background. A special application booklet for ELISA procedures is available on request (L20025).

**Elisa Testing:**

To assure consistent activity and low background in sandwich ELISA assays, each lot of Caltag IL-12 capture antibody is tested with approved lots of Caltag biotinylated IL-12 detection antibody (RM9315) with and without murine IL-12. Quality control data is available on request.








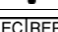
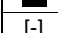
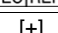

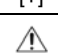
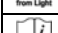
For conjugated products only, a highly purified grade of BSA has been added as a stabilizing protein to bring the final protein concentration to 4-5 mg/ml after conjugation.

**REFERENCES**

1. Wysocka, M., M. Kubin, L. Q. Vieira, L. Ozmen, G. Garotta, P. Scott, and G. Trinchieri. 1995. Interleukin-12 is required for interferon-γ production and lethality in lipopolysaccharide-induced shock in mice. *Eur. J. Immunol.* 25: 672-676

\* The amount of antibody is determined by measuring the optical density using a spectrophotometer. The antibody titer is verified by immunofluorescent staining and flow cytometric analysis.

**Explanation of symbols**

Symbol	Description	Symbol	Description
	Catalogue Number		Batch code
	Research Use Only		<i>In vitro</i> diagnostic medical device
	Use by		Temperature limitation
	Manufacturer		European Community authorised representative
	Without, does not contain		With, contains
	Protect from light		Consult accompanying documents
	Directs the user to consult instructions for use (IFU), accompanying the product.		

**For research use only. CAUTION: Not for human or animal therapeutic or diagnostic use**

[www.invitrogen.com](http://www.invitrogen.com)

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

PI: L11309

(Rev 02/10) DCC-10-0338

**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](http://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.