

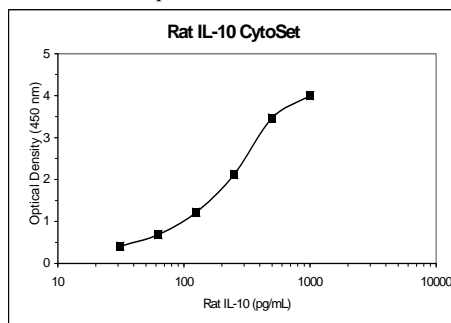
Catalog # CRC0103

Lot #: S030709

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

1. **Coating Antibody:** **Ms Anti-Rat IL-10 (0.125mg/0.125mL)**
 Part Number: ARC9104-D
 Lot Number: S042612
 Form: Liquid, 1 vial, contains 0.1% sodium azide
 Storage: Store at 2-8°C for 1 month. For longer periods, aliquot and store at ≤ -20°C.
 Recommended Dilution: Dilute to 1.25 µg/mL with Coating Buffer B (Cat. # CB01100, or see Recommended Buffers). For example, to make 10 mL (enough to coat 1 plate), add 12.5 µL coating antibody to 9.988 mL Coating Buffer B.
2. **Detection Antibody:** **Ms Anti-Rat IL-10 Biotin (0.025mg/0.125mL)**
 Part Number: ARC7109-D
 Lot Number: S042613
 Form: Liquid, 1 vial, contains 0.1% sodium azide
 Storage: Store at 2-8°C for 1 month. For longer periods, aliquot and store at ≤ -20°C.
 Recommended Dilution: Dilute to 0.125 µg/mL with Assay Buffer **supplemented with 5% calf serum** (Cat. # DS98200, or see Recommended Buffers). For example, to make enough for 1 plate, add 3.4 µL detection antibody to 5.49 mL Assay Buffer.
3. **Standard:** **Recombinant Rat IL-10**
 Part Number: SD048 (inquire regarding additional vials)
 Lot Number: R121916
 Form: Lyophilized, 3 vials (single use)
 Storage: Store at 2-8°C.
 Concentration of Reconstituted Standard: 20,000 pg/mL.
 Reconstitution: Reconstitute in Assay Buffer **supplemented with 50% calf serum** (Cat. # DS98200 or see Recommended Buffers) according to instructions on vial label. Allow standard to rehydrate for approximately 10 minutes before dilutions. If the standard stock is not being used immediately, please aliquot into polypropylene tubes and freeze at -80°C. *Do not store at room temperature or at 4°C for any extended time or subject to more than one freeze-thaw cycle.*
 Recommended Starting Standard Curve: Dilutions of the standard should be made in Assay Buffer **supplemented with 50% calf serum**. Dilute standard stock to 1000 pg/mL (a 1:20 dilution) followed by six 1:2 serial dilutions using at least 300 µL of buffer. Mix thoroughly between dilutions. Avoid foaming. To an empty tube add 300 µL of buffer and label as zero standard.
4. **Streptavidin-HRP:** **0.025 mg/0.125 mL**
 Part Number: SNN4004Y
 Lot Number: R102526C
 Form: Liquid, 1 vial, contains animal serum and 50% glycerol in phosphate buffered saline with 0.05% thymol as a preservative.
 Storage: Store concentrate at 2-8°C for 1 month. For longer periods, aliquot and store at ≤ -20°C. Diluted streptavidin-HRP should not be stored; discard remaining solution after use.
 Recommended Dilution: Dilute to 0.2 µg/mL. For example, to make enough for 1 plate, add 10 µL of streptavidin-HRP to 9.990 mL of Assay Buffer (Cat. # DS98200 or see Recommended Buffers).

Following the recommended assay procedure using the **Invitrogen CytoSet™ Buffer Set** (Cat. # CNB0011), the representative standard curve was generated.



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

Intended Use and Materials Provided

The CytoSet™ for Rat IL-10 contains components required to construct an enzyme-linked immunoassay for the specific and quantitative measurement of IL-10. Sufficient quantities of all reagents are provided to yield 10 plates of 96 wells if the recommended assay procedure and recommended storage and handling of materials are followed as specified on this insert. The materials provided are **FOR RESEARCH USE ONLY**.

Recommended Buffers and Solutions

The Invitrogen CytoSet™ Buffer Set (Cat. # CNB0011) containing Coating Buffers A and B, Assay Buffer, Substrate Solution (TMB), Stop Solution, and Wash Buffer is recommended.

- Coating Buffer A:** *Coating Buffer A (Cat. # CB07100) from Invitrogen is recommended. Alternate buffer choice listed below.*
8.0 g NaCl, 1.13 g Na₂HPO₄, 0.2 g KH₂PO₄, 0.2 g KCl; q.s. to 1.0 L with distilled H₂O, pH to 7.4.
- Coating Buffer B:** *Coating Buffer B (Cat. # CB01100) from Invitrogen is recommended. Alternate buffer choice listed below.*
4.3 g NaHCO₃, 5.3 g Na₂CO₃; q.s. to 1.0 L with distilled H₂O, pH to 9.4.
- Assay Buffer:** *Assay Buffer (Cat. # DS98200) from Invitrogen is recommended. Alternate buffer choice listed below.*
8.0 g NaCl, 1.13 g Na₂HPO₄, 0.2 g KH₂PO₄, 0.2 g KCl, 5.0 g bovine serum albumin (fraction V), 1 mL Tween 20; q.s. to 1.0 L with distilled H₂O, pH to 7.4.
- Wash Buffer:** *Wash Buffer (Cat. # WB01) from Invitrogen is recommended. Alternate buffer choice listed below.*
9.0 g NaCl, 1 mL Tween 20; q.s. to 1.0 L with distilled H₂O, pH to 7.4.
- Substrate Solution:** *TMB (Cat. # SB01) from Invitrogen is recommended. Alternate solution choice listed below.*
Tetramethylbenzidine (TMB) and Hydrogen Peroxide.
- Stop Solution:** *Stop Solution (Cat. # SS01100) from Invitrogen is recommended. Alternate solution choice listed below.*
1.8 N H₂SO₄.

Assay Optimization

CytoSets™ from Invitrogen are designed to be very flexible for your experiments. Consequently, the assay procedure contains only recommendations. The assay procedure has been optimized for use with tissue culture samples. However, serum and plasma samples may be used but may require that certain assay parameters be modified. Investigators are advised to determine optimal buffer formulations, concentrations and incubation times for individual applications.

Recommended Assay Procedure

- Prepare coating solution by diluting the coating antibody. See “coating antibody” section for the recommended coating antibody dilution.
- Coat plates with 100 µL per well of the coating solution. Cover plates and incubate overnight (12-18 hr.) at 4°C.
- Aspirate wells and wash 1 time with > 400 µL of Wash Buffer per well. Following wash, invert and tap on absorbent paper to remove excess liquid.
- Block plate with 300 µL per well of Assay Buffer for 1 hour at room temperature.
- Aspirate, invert, and tap on absorbent paper to remove excess liquid.
- Prepare standards and sample dilutions in Assay Buffer (or in a diluent that most closely matches the matrix of your sample).
- Pipette 100 µL of standards (in duplicate), samples and controls into designated wells.
- Immediately following step 7, add 50 µL of the working detection antibody into each well. For recommended dilutions, see “detection antibody” section. Gently tap the plate on the side 10 times to mix. **Cover plate and incubate for 2 hours at room temperature.**
- Aspirate and wash 5 times using the method in step 3.
- Add 100 µL of the working streptavidin-HRP solution into each well. For recommended dilutions, see “streptavidin-HRP” section. **Cover plate and incubate for 30 minutes at room temperature.**
- Aspirate and wash 5 times using the method in step 3.
- Add 100 µL of the TMB substrate to each well. **Incubate plate without a plate cover for 30 minutes in the dark at room temperature.**
- Add 100 µL of Stop Solution to each well.
- Measure absorbance at 450 nm (reference absorbance: 650 nm) within 30 minutes of adding Stop Solution. Calculate results using a log-log or 4-parameter curve fit.

Additional Materials Required

- 96 well NUNC MaxiSorp microplates; NUNC Cat. # 434797.
- Pipettes; plate covers or plate sealers and timer.
- Microplate reader with a detector that can measure absorbance at 450 nm.
- 1 L graduated cylinder; plate washer or wash bottle.
- Polypropylene tubes for standards and sample dilutions, if needed.

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