

Amplex® ELISA Development Kit for Rabbit IgG with Amplex® UltraRed Reagent

Catalog no. A33852

Table 1. Contents and storage information.

Material	Amount	Storage*	Stability
Amplex® UltraRed reagent (Component A)	5 vials, each containing 180 µg	<ul style="list-style-type: none"> • ≤-20°C • Desiccate • Protect from light 	When stored as directed, the kit components are stable for at least 6 months.
Dimethyl sulfoxide (DMSO), anhydrous (Component B)	1.75 mL		
10X Phosphate-buffered saline (PBS) pH 7.2 (Component C)	200 mL		
Goat anti-rabbit IgG (H+L), horseradish peroxidase conjugate (Component D)	2 vials, each containing 100 µg		
Amplex® stop reagent (Component E)	20 mg		
Hydrogen peroxide (H ₂ O ₂), stabilized ~3% solution (Component F)	500 µL		
0.1 M sodium bicarbonate buffer, pH ~9.3 (Component G)	50 mL		
Bovine serum albumin (BSA) (Component H)	1.2 g		
Tween® 20 (Component I)	1.5 mL		
Nunc-Immuno™ MaxiSorp™ U96 plate (Component J)	5 each		
*The kit can be stored under the conditions listed. For optimal storage conditions of individual components, refer to the labels on the vials.			
Number of assays: Sufficient material is supplied for 500 reactions in 96-well microplates at 100 µL per well, based on the protocol below..			
Approximate fluorescence excitation and emission maxima: 568/581 nm for the reaction product.			

Introduction

The Amplex® ELISA Development Kit for Rabbit IgG provides a comprehensive set of components for creating a fluorescence-based ELISA using a rabbit primary antibody. The assay is based on Amplex® UltraRed reagent, a fluorogenic substrate for horseradish peroxidase (HRP) that reacts with hydrogen peroxide (H₂O₂) in a 1:1 stoichiometric ratio to produce Amplex® UltroRed product, a brightly fluorescent and strongly absorbing reaction product (excitation/emission maxima ~568/581 nm) (Figure 1). Because the Amplex® UltroRed product has long-wavelength spectra, there is little interference from the blue or

green autofluorescence found in most biological samples. With a high extinction coefficient, good quantum efficiency, and resistance to autooxidation, the fluorescence-based Amplex® UltraRed reagent delivers better sensitivity and a broader assay range than colorimetric reagents. In a sandwich ELISA format using C-reactive protein, it is possible to routinely detect 1 pg/ml of antigen (Figure 2). Using TMB (3,3',5,5'-Tetramethylbenzidine, a common colorimetric reagent) in the same sandwich ELISA format, the assay was 25-fold less sensitive.

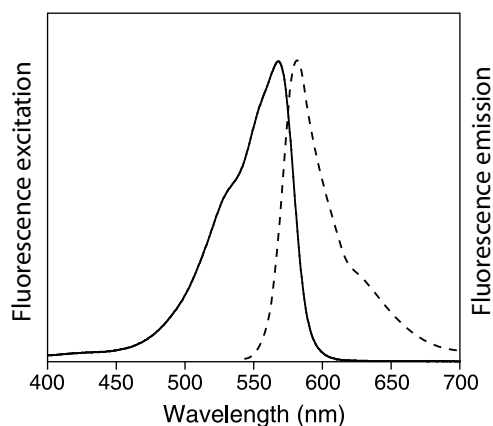


Figure 1. Normalized absorption and fluorescence emission spectra for the Amplex® UltraRed product.

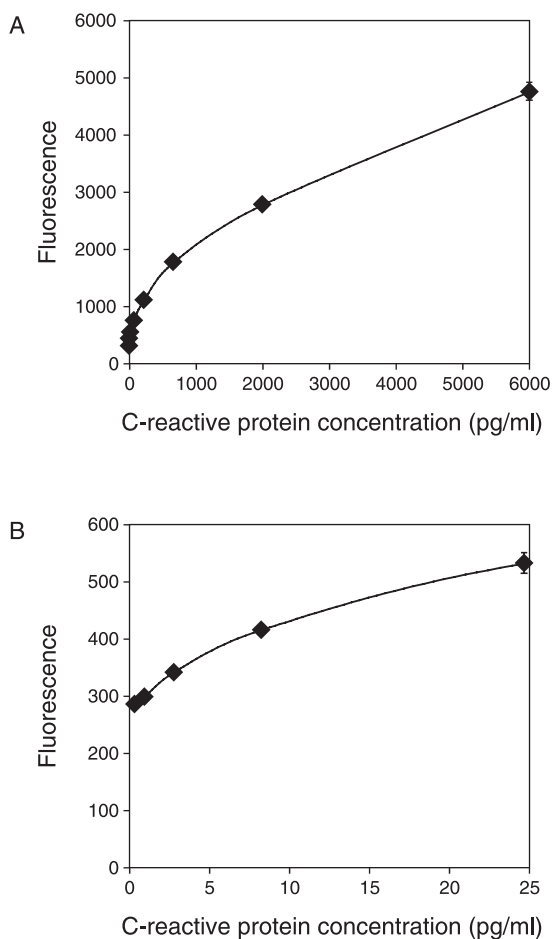


Figure 2. Detection range of C-reactive protein (CRP) using the Amplex® ELISA Development Kit for Rabbit IgG. The sandwich ELISA was carried out as described in the protocol using a mouse anti-CRP capture antibody, C-reactive protein in a concentration range from 6,000 pg/mL to 0.10 pg/mL, and a rabbit polyclonal anti-CRP primary antibody (100 µL per well of a 50 ng/mL solution). The Z' factor¹ analysis of the data obtained gives a lower limit of detection for CRP in this system of 1 pg/mL or 0.1 pg/well (based on a well volume of 100 µL in the sandwich ELISA).

Before you Begin

Materials Required but Not Provided

- Rabbit-derived antibody against target antigen
- Capture antibody against target antigen (e.g., mouse)
- Antigen (for generating a standard curve; see protocol step 2.5)
- Distilled and deionized water
- Single-channel and multichannel pipettes (1 μ L to 1 mL range)
- Fluorescence microplate reader capable of excitation/emission settings of 530 nm/590 nm

Caution

No data are currently available addressing the mutagenicity or toxicity of the Amplex® UltraRed reagent (Component A).

DMSO (Component B) is hazardous; avoid contact with skin and eyes and do not swallow. Handle reagents containing DMSO using equipment and practices appropriate for the hazards posed by such materials.

Amplex® stop reagent (Component E) is irritating to eyes, respiratory system, and skin, and may be harmful if swallowed. Avoid prolonged or repeated exposure. If eye or skin contact occurs, wash affected area with water for 15 minutes and seek medical advice. If inhaled, move individual to fresh air and seek medical advice. If swallowed, seek medical advice.

Preparing Stock Solutions

Allow components warm to room temperature before opening the vials and preparing various stock solutions.

- 1.1 Prepare 500 mL of 1X PBS** by adding 50 mL of 10X PBS (Component C) to 450 mL distilled and deionized water. You will use this stock of 1X PBS in preparing other buffers as well as in the final Amplex® UltraRed reaction.
- 1.2 Prepare 300 mL of 1X PBST** by adding 300 μ L of Tween® 20 (Component I) to 300 mL of 1X PBS. Shake well to mix. This solution is sufficient for 100 assays using the protocol below. Storage at 4°C is not required, but will not harm this solution.
- 1.3 Prepare 100 mL of 1X PBS-BSA** by adding 1 g of BSA (Component H) to 100 mL of 1X PBS. Dissolve completely. Store at 4°C when not in use.
- 1.4 Prepare a 200 μ g/mL stock solution of the goat anti-rabbit IgG HRP conjugate** (Component D) by adding 0.5 mL of PBS-BSA directly to the vial. Store this stock solution at 4°C after adding thimerosal to a final concentration of 0.02%.
- 1.5 Prepare a ~10 mM stock solution of Amplex® UltraRed** by adding 60 μ L of DMSO (Component B) to one vial of Amplex® UltraRed reagent (Component A). Vortex well to dissolve. Protect from light and moisture.
- 1.6 Prepare 10 mL of Amplex® stop solution** by resuspending the dried Amplex® stop reagent (Component E) in 1 mL of 1 M NaOH, and once resuspended, adding it to 9 mL of 1X PBS. This solution is stable for one month at 4°C when protected from light. If this solution begins to turn amber, it is no longer good and should be discarded.

Experimental Protocols

The following procedure describes a typical sandwich ELISA, and is designed for use with a fluorescence microplate reader. The procedure has been optimized for use with 96-well microplates and reaction volumes of 100 μ L per assay.

You may replace the protocol with any standard sandwich ELISA protocol at your discretion; however, you must follow steps 2.8–2.14 of this protocol as described.

- 2.1** Using 0.1 M sodium bicarbonate (Component G), prepare 10 mL of a 10 μ g/mL solution of the desired capture antibody and aliquot 100 μ L of this solution into each microplate well. You can use any capture antibody but it must **not** cross-react with the goat anti-rabbit IgG or the rabbit anti-target antibodies used for detection.

Incubate at room temperature for at least four hours, or overnight at 4°C (preferred).

Note: Overnight deposition of capture antibody provides optimal detection. If desired, you can incubate the plate at room temperature for 8 hours, and then block with 1X PBS-BSA overnight at 4°C.

- 2.2** Discard or shake contents of the plate into the sink, and wash the wells three times with 200 μ L of 1X PBST (prepared in step 1.2).
- 2.3** Add 200 μ L of 1X PBS-BSA (prepared in step 1.3) to each well of the microplate, and incubate at room temperature for at least four hours, or overnight at 4°C (preferred).
- 2.4** Shake plate contents into the sink, and wash the wells three times with 200 μ L of 1X PBST.
- 2.5** Add 100 μ L of 0.1X PBS-BSA (prepared by diluting the 1X PBS-BSA ten-fold in 1X PBS) to each well of the microplate. Add antigen to the first well of each row and serially dilute across the plate to achieve the desired range of concentrations. Leave the last well of the row as a no-antigen control. Incubate the plate at room temperature for one hour.

Example: Prepare 150 μ L of 60 ng/mL antigen in the first well and serially dilute 50 μ L into each successive well in that row to make a three-fold dilution series ranging from 6 ng to 0.1 pg antigen.

- 2.6** Following incubation, shake plate contents into the sink and wash the wells three times with 200 μ L of 1X PBST.
- 2.7** Prepare 10 mL of 50 ng/mL secondary capture antibody in 0.1X PBS-BSA and add 100 μ L to each well of the microplate. Incubate at room temperature for 30 minutes.
- 2.8** Shake plate contents into the sink, and wash the wells three or more times with 200 μ L of 1X PBST.
- 2.9** Prepare 10 mL of 50 ng/mL goat anti-rabbit IgG HRP by adding 2.5 μ L of the stock goat anti-rabbit IgG HRP solution (prepared in step 1.4) to 10 mL of 0.1X PBS-BSA. Add 100 μ L of this solution to each well of the microplate and incubate at room temperature for 30 minutes.
- 2.10** Shake plate contents into the sink, and wash the wells three times with 200 μ L of 1X PBST. You may adjust the stringency of the assay by washing more or fewer times with PBST, or by incubating or agitating PBST in the wells for a time during the wash steps. **Protect the plate from light at all times from this point onward.**

Note: Antibodies exposed to UV light can produce trace amounts of singlet oxygen, which can interfere with detection of Amplex® UltraRed reagent. Protecting the plate from light after the final wash provides optimal sensitivity for the assay.

- 2.11** Prepare 10 mL of reaction mixture by adding 50 µL of the 10 mM Amplex® UltraRed stock solution (prepared in step 1.5) and 22.7 µL of 3% H₂O₂ (Component F) to 10 mL of 1X PBS. If necessary, adjust the volume for the actual hydrogen peroxide concentration (check the label on Component F for actual H₂O₂ concentration). Protect the reaction mixture from light and use within 4 hours of preparation.
- 2.12** Using a multichannel pipette, add 100 µL of the reaction mixture to each assay well.
- Note:** Adding the reaction mixture (containing Amplex® UltraRed reagent) to the wells initiates the reaction.
- 2.13** Incubate the plate at room temperature, protected from light, until the fluorescence measurement is taken. For most reactions, a 30 minute incubation is sufficient. The plate can also be read continuously for up to an hour.
- If desired, you may add 20 µL of Amplex® stop solution (prepared in step 1.6) to each assay well. This will arrest the reaction, providing a stable signal that may be read for at least two hours if the plate is protected from light and kept at room temperature.
- 2.14** Measure the fluorescence in a microplate reader using filters for 530 nm (excitation) and 590 nm (emission)

Reference

1. J Biomol Screen 4, 67 (1999).

Product List

Current prices may be obtained from our website or from our Customer Service Department.

Cat. no.	Product Name	Unit Size
A33852	Amplex® ELISA Development Kit for Rabbit IgG *with Amplex® UltraRed reagent* *500 assays*	1 kit
Related Products		
A33851	Amplex® ELISA Development Kit for Mouse IgG *with Amplex® UltraRed reagent* *500 assays*	1 kit
A33855	Amplex® Red/UltraRed stop reagent *500 tests*	set of 5 vials
A36006	Amplex® UltraRed reagent.....	5 × 1 mg

Contact Information

Molecular Probes, Inc.

29851 Willow Creek Road
Eugene, OR 97402
Phone: (541) 465-8300
Fax: (541) 335-0504

Customer Service:

6:00 am to 4:30 pm (Pacific Time)
Phone: (541) 335-0338
Fax: (541) 335-0305
probesorder@invitrogen.com

Toll-Free Ordering for USA:

Order Phone: (800) 438-2209
Order Fax: (800) 438-0228

Technical Service:

8:00 am to 4:00 pm (Pacific Time)
Phone: (541) 335-0353
Toll-Free (800) 438-2209
Fax: (541) 335-0238
probetech@invitrogen.com

Invitrogen European Headquarters

Invitrogen, Ltd.
3 Fountain Drive
Inchinnan Business Park
Paisley PA4 9RF, UK
Phone: +44 (0) 141 814 6100
Fax: +44 (0) 141 814 6260
Email: euroinfo@invitrogen.com
Technical Services: eurotech@invitrogen.com

**For country-specific contact information,
visit www.invitrogen.com.**

Further information on Molecular Probes products, including product bibliographies, is available from your local distributor or directly from Molecular Probes. Customers in Europe, Africa and the Middle East should contact our office in Paisley, United Kingdom. All others should contact our Technical Service Department in Eugene, Oregon.

Molecular Probes products are high-quality reagents and materials intended for research purposes only. These products must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Please read the Material Safety Data Sheet provided for each product; other regulatory considerations may apply.

Limited Use Label License No. 223: Labeling and Detection Technology

The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) to not transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Molecular Probes, Inc., Business Development, 29851 Willow Creek Road, Eugene, OR 97402, Tel: (541) 465-8300. Fax: (541) 335-0354.

Several Molecular Probes products and product applications are covered by U.S. and foreign patents and patents pending. All names containing the designation ® are registered with the U.S. Patent and Trademark Office.

Copyright 2009, Molecular Probes, Inc. All rights reserved. This information is subject to change without notice.