

Amplex® Red Uric Acid/Uricase Assay Kit

Catalog no. A22181

Table 1 Contents and storage

Material	A22181	Concentration	Storage*
Amplex® Red reagent (MW = 257, Component A)	2 vials of 0.26 mg	—	<ul style="list-style-type: none"> • -20°C • Desiccate • Protect from light
Dimethylsulfoxide (DMSO), anhydrous (Component B)	500 µL	—	
Horseradish peroxidase (Component C)	20 U**	—	
Hydrogen peroxyde (H ₂ O ₂) (MW = 34, Component D)	500 µL	~3%†	
5X Reaction Buffer (Tris-HCl, pH 7.5; Component E)	10 mL	0.5 M	
Uricase (Component F)	10 U‡	—	
Uric acid (MW = 168, Component G)	3 vials of 1.5 mL	5 mM in 1X Reaction Buffer	

*These storage conditions are appropriate when storing the entire kit upon receipt. For optimal storage of each component, see vial labels. When stored as directed this product is stable for 6 months. **One unit is defined as the amount of enzyme that will form 1.0 mg purpurogallin from pyrogallol in 20 seconds at pH 6.0 and 20°C. †The actual concentration is indicated on the component label. ‡One unit is define as the amount of enzyme that will convert 1.0 µmole of uric acid to allantoin per minute at pH 8.5 and 25°C

Approximate fluorescence excitation/emission maxima of reaction product: 571/585 in nm.

Introduction

The Amplex® Red Uric Acid/Uricase Assay Kit provides an ultrasensitive method for detecting uric acid or for monitoring uricase activity. In the assay, uricase catalyzes the conversion of uric acid to allantoin, hydrogen peroxide (H₂O₂) and carbon dioxide. The H₂O₂ then, in the presence of horseradish peroxidase (HRP), reacts stoichiometrically with Amplex® Red reagent to generate the red-fluorescent oxidation product, resorufin.¹ Resorufin has absorption and fluorescence emission maxima of approximately 571 nm and 585 nm, respectively (Figure 1), and because the extinction coefficient is high (54,000 cm⁻¹M⁻¹), the assay can be performed either fluorometrically or spectrophotometrically.

Serum uric acid is the end product of purine metabolism in the body tissues and is cleared through the kidneys by glomerular filtration. Most animals can metabolize uric acid to more readily excreted products, but humans lack the necessary enzyme, urate oxidase (uricase), as a result of the presence of two “nonsense mutations” in the human gene for uricase.² Increased uric acid levels may result from leukemia, polycythemia, ingestion of foods high in nucleoproteins (e.g. liver and kidney) or impaired renal function. Gout results from the deposit of uric acid in body joints.³ Previous literature reports colormetric detection limits at 3.6 µM,⁴ whereas the Amplex® Red Uric Acid/Uricase Assay Kit can be used to detect as little as 100 nM of uric acid in a purified system (Figure 2). The kit can also be used to detect a little as 0.2 mU/mL uricase in a purified system (Figure 3).

For Research Use Only. Not for use in diagnostic procedures.

Figure 1 Normalized absorption and fluorescence emission spectra of resorufin, the product of the Amplex® Red reagent.

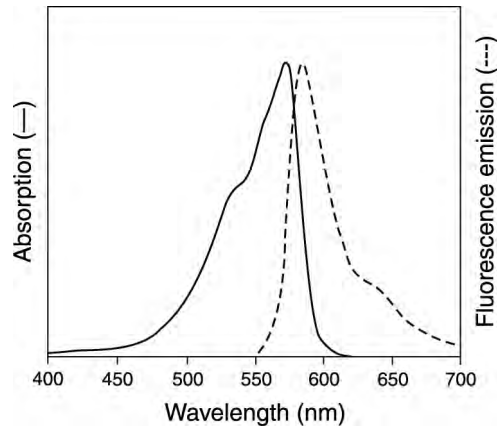


Figure 2 Detection of uric acid using the Amplex® Red reagent-based assay. Each reaction contained 50 μ M Amplex® Red reagent, 0.2 U/mL HRP, 0.2 U/mL uricase, and the indicated amount of uric acid in 1X Reaction Buffer. After 30 minutes incubation at 37°C, fluorescence was measured in a fluorescence microplate reader using excitation at 530 ± 12.5 nm and fluorescence detection at 590 ± 17.5 nm. A background fluorescence of 26 fluorescence units was subtracted from each data point.

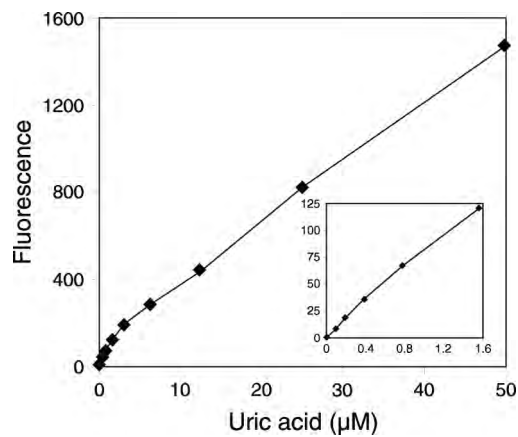
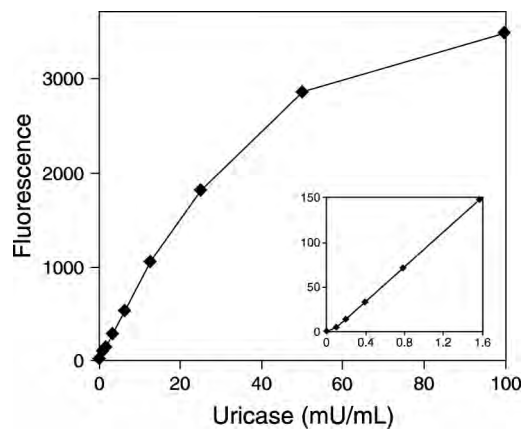


Figure 3 Detection of uric acid using the Amplex® Red reagent-based assay. Each reaction contained 50 μ M Amplex® Red reagent, 0.2 U/mL HRP, 1 mM uric acid, and the indicated amount of uricase in 1X Reaction Buffer. After 30 minutes incubation at 37°C, fluorescence was measured in a fluorescence microplate reader using excitation at 530 ± 12.5 nm and fluorescence detection at 590 ± 17.5 nm. A background fluorescence of 26 fluorescence units was subtracted from each data point.



Before Starting

Storage and Handling Upon receipt, the kit should be stored frozen at -20°C , protected from light. Stored properly, the kit components should remain stable for at least six months. Allow reagents to warm to room temperature before opening vials. The Amplex[®] Red reagent is somewhat air sensitive. Once a vial of Amplex[®] Red reagent is opened, the reagent should be used promptly. PROTECT THE Amplex[®] RED REAGENT FROM LIGHT.

Preparing Stock Solutions

- 1.1 Prepare a 10 mM stock solution of Amplex[®] Red reagent: Allow one vial of Amplex[®] Red reagent (Component A) and DMSO (Component B) to warm to room temperature. Just prior to use, dissolve the contents of the vial of Amplex[®] Red reagent (0.26 mg) in 100 μL of DMSO. Each vial of Amplex[®] Red reagent is sufficient for approximately 200 assays, with a final reaction volume of 100 μL per assay. This stock solution should be stored frozen at -20°C , protected from light.
- 1.2 Prepare a 1X working solution of Reaction Buffer by adding 4 mL of 5X Reaction Buffer stock solution (Component E) to 16 mL of deionized water (dH_2O). This 20 mL volume of 1X Reaction Buffer is sufficient for approximately 100 assays of 100 μL each with a 10 mL excess for making stock solutions and dilutions.
- 1.3 Prepare a 100 U/mL stock solution of horseradish peroxidase (HRP) by dissolving the contents of the vial of HRP (Component C) in 200 μL of 1X Reaction Buffer. After use, the remaining solution should be divided into small aliquots and stored frozen at -20°C .
- 1.4 Prepare a 20 mM H_2O_2 working solution by diluting the $\sim 3\%$ H_2O_2 stock solution (Component D) into the appropriate volume of dH_2O . The actual H_2O_2 concentration is indicated on the component label. For instance, a 20 mM H_2O_2 working solution can be prepared from a 3.0% H_2O_2 stock solution by diluting 23 μL of 3.0% H_2O_2 into 977 μL of dH_2O . Please note that although the $\sim 3\%$ H_2O_2 stock solution has been stabilized to slow degradation, the 20 mM H_2O_2 working solution will be less stable and should be used promptly.
- 1.5 Prepare a 100 U/mL uricase stock solution by dissolving the contents of the vial of uricase (Component F) in 100 μL of dH_2O . After use the remaining solution should be stored at -20°C .

Experimentals Protocols

The following procedure is designed for use with a fluorescence or absorbance multiwell plate scanner. For use with a standard fluorometer or spectrophotometer, volumes must be increased accordingly.

Please note that resorufin, the product of the Amplex[®] Red reaction, is unstable in the presence of thiols such as dithiothreitol (DTT) and 2-mercaptoethanol. For this reason, the final DTT or 2-mercaptoethanol concentration in the reaction should be no higher than 10 μ M.

The absorption and fluorescence of resorufin are pH-dependent. Below the pK_a (~6.0), the absorption maximum shifts to ~480 nm and the fluorescence quantum yield is markedly lower. In addition, the Amplex[®] Red reagent is unstable at high pH (>8.5). For these reasons, the reactions should be performed at pH 7–8. We recommend using the included Reaction Buffer (pH 7.5) for optimal performance of the Amplex[®] Red reagent.

Uric Acid Assay The following protocol describes the assay of uric acid in a total volume of 100 μ L per microplate well. The volumes recommended here are sufficient for ~100 assays. The kit provides sufficient material for ~400 assays.

- 2.1 Prepare a uric acid standard curve by diluting the appropriate amount of 5 mM uric acid (Component G) into 1X Reaction Buffer to produce uric acid concentrations of 0 to 100 μ M. Use 1X Reaction Buffer without uric acid as a negative control. A volume of 50 μ L will be used for each reaction. Note that the final concentration will be twofold lower in the final reaction.
- 2.2 If desired, prepare a positive control by diluting the 20 mM H₂O₂ working solution (prepared in Step 1.4) to 10 μ M in 1X Reaction Buffer.
- 2.3 Dilute the uric acid-containing samples in 1X Reaction Buffer. A volume of 50 μ L will be used for each reaction.
- 2.4 Pipet 50 μ L of the diluted samples, standards and controls into separate wells of a microplate.
- 2.5 Prepare a working solution of 100 μ M Amplex[®] Red reagent containing 0.4 U/mL HRP and 0.4 U/mL uricase according to Table 2.

Table 2 Amplex[®] Red reagent working solution for uricase assay

Reaction Components	Amount needed (for ~100 assays)
Amplex [®] Red reagent stock solution (prepared in Step 1.1)	50 μ L
HRP stock solution (prepared in Step 1.3)	20 μ L
Uricase stock solution (prepared in Step 1.5)	20 μ L
1X Reaction Buffer	4.91 mL
Total volume	5 mL

Note that the final concentration of each component will be two-fold lower in the final reaction.

- 2.6 Begin the reactions by adding 50 μL of the Amplex[®] Red reagent/HRP/uricase working solution to each microplate well containing the samples and controls.
- 2.7 Incubate the reactions for 30 minutes or longer at 37°C, protected from light. Because the assay is continuous (not terminated), fluorescence or absorbance may be measured at multiple time points to follow the kinetics of the reactions.
- 2.8 Measure the fluorescence or absorbance in a microplate reader using excitation in the range of 530–560 nm and emission detection at ~590 nm or absorbance at ~560 nm (see Figure 1).
- 2.9 For each point, correct for background fluorescence or absorbance by subtracting the value derived from the no-uric acid control.

Uric Acid Assay The following protocol describes the assay of uricase activity in a total volume of 100 μL per microplate well. The volumes recommended here are sufficient for ~100 assays. The kit provides sufficient material for ~400 assays.

- 3.1 Prepare a uricase standard curve by diluting the 100 U/mL uricase stock solution (prepared in step 1.5) in 1X Reaction Buffer to produce uricase concentrations of 0 to 100 mU/mL. Use 1X Reaction Buffer without uricase as a negative control. A volume of 50 μL will be used for each reaction. Note that the uricase concentration will be twofold lower in the final reaction volume.
- 3.2 Dilute the uricase-containing samples in 1X Reaction Buffer. A volume of 50 μL will be used for each reaction.
- 3.3 If desired, prepare a positive control by diluting the 20 mM H₂O₂ work solution (prepared in step 1.4) to 10 μM in 1X Reaction Buffer.
- 3.4 Pipet 50 μL of the diluted uricase-containing standards, controls and samples into separate wells of a microplate.
- 3.5 Prepare a working solution of 100 μM Amplex[®] Red reagent containing 0.4 U/mL HRP and 1.0 mM uric acid according to Table 3.

Table 3 Amplex[®] Red reagent working solution for uric acid assays

Reaction Components	Amount needed (for ~100 assays)
Amplex [®] Red reagent stock solution (prepared in Step 1.1)	50 μL
HRP stock solution (prepared in Step 1.3)	20 μL
5 mM uric acid stock solution (Component G)	1 mL
1X Reaction Buffer	3.93 mL
Total volume	5 mL

Note that the final concentration of each component will be two-fold lower in the final reaction.

- 3.6 Begin the reaction by adding 50 μL of the Amplex[®] Red reagent/HRP/uric acid working solution to each microplate well containing the samples and controls.
- 3.7 Incubate the reaction for 30 minutes or longer at 37°C, protected from light. Because the assay is continuous (not terminated), fluorescence or absorbance may be measured at multiple time points to follow the kinetics of the reactions.

- 3.8 Measure the fluorescence or absorbance in a microplate reader using excitation in the range of 530–560 nm and emission detection at ~590 nm (see Figure 1).
- 3.9 For each point, correct for background fluorescence or absorbance by subtracting the value derived from the no-uricase control.

References

1. J Immunol Methods 202, 133 (1997); 2. J Mol Evol 34, 78 (1992); 3. Anal Chem 71, 1928 (1999); 4. Clin Biochem 27, 93 (1994).

Product List

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

Cat. no.	Product Name	Unit Size
A12222	Amplex® Red reagent (10-acetyl-3,7-dihydroxyphenoxazine).....	5 mg
A22177	Amplex® Red reagent*packaged for high-throughput screening*	10 x 10 mg
A22181	Amplex® Red Uric Acid/Uricase Assay Kit *400 assays*.....	1 kit
A36006	Amplex® UltraRed reagent.....	5 x 1 mg

Purchaser Notification

Corporate Headquarters

5791 Van Allen Way
Carlsbad, CA 92008
USA
Phone: +1 760 603 7200
Fax: +1 760 602 6500
Email: techsupport@lifetech.com

European Headquarters

Inchinnan Business Park
3 Fountain Drive
Paisley PA4 9RF
UK
Phone: +44 141 814 6100
Toll-Free Phone: 0800 269 210
Toll-Free Tech: 0800 838 380
Fax: +44 141 814 6260
Tech Fax: +44 141 814 6117
Email: euroinfo@invitrogen.com
Email Tech: eurotech@invitrogen.com

Japanese Headquarters

LOOP-X Bldg. 6F
3-9-15, Kaigan
Minato-ku, Tokyo 108-0022
Japan
Phone: +81 3 5730 6509
Fax: +81 3 5730 6519
Email: jpinfo@invitrogen.com

Additional international offices are listed at
www.lifetechnologies.com

These high-quality reagents and materials must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Read the Safety Data Sheet provided for each product; other regulatory considerations may apply.

Obtaining Support

For the latest services and support information for all locations, go to www.lifetechnologies.com.

At the website, you can:

- Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support (techsupport@lifetech.com)
- Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents
- Obtain information about customer training
- Download software updates and patches

SDS

Safety Data Sheets (SDSs) are available at www.lifetechnologies.com/sds.

Certificate of Analysis

The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to www.lifetechnologies.com/support and search for the Certificate of Analysis by product lot number, which is printed on the product packaging (tube, pouch, or box).

Limited Product Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

Disclaimer

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

Important Licensing Information

This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliate(s) or their respective owners.

©2014 Life Technologies Corporation. All rights reserved.

