



EAGLE
BIOSCIENCES

Complement Factor H-Related Protein 2 (FHR-2) ELISA Kit

Catalog Number:

HK3004-01 (1 x 96 wells)

HK3004-02 (2 x 96 wells)

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

v. 3.24

Eagle Biosciences, Inc.

20A Northwest Blvd., Suite 112, Nashua, NH 03063

Phone: 617-419-2019 Fax: 617-419-1110

www.EagleBio.com



INTENDED USE

The human FHR-2 ELISA kit is to be used for the in vitro quantitative determination of FHR-2 in serum and plasma samples. This kit is intended for laboratory research use only and is not for use in diagnostic or therapeutic procedures. The analysis should be performed by trained laboratory professionals.

INTRODUCTION

Complement Factor H-Related Protein 2 (FHR-2) is a significant member of the complement regulatory protein family, with a molecular weight of around 29 kDa. Synthesized mainly in the liver, FHR-2 is also produced by immune cells like monocytes, macrophages, and dendritic cells. Sharing structural homology with Complement Factor H (CFH), FHR-2 plays a role in regulating the complement cascade, although its exact biological function remains somewhat elusive and subject to debate. Research indicates that FHR-2 may inhibit the activity of the alternative pathway (AP) C3 convertase, which would theoretically prevent the formation of the membrane attack complex (MAC). Conversely, some studies suggest that FHR-2 might actually promote AP activation, highlighted by findings where a hybrid mutant protein comprising FHR-5 domains and full-length FHR-2 led to overactivation of the complement system.

FHR-2 has been linked to various diseases where abnormal complement activation is a contributing factor. Particularly, it is associated with atypical hemolytic uremic syndrome (aHUS) and age-related macular degeneration (AMD), where dysregulation in complement activation leads to detrimental tissue deposition. The involvement of FHR-2 in these pathologies, specifically in the alternative pathway, makes it a potential therapeutic target in managing diseases driven by complement dysregulation.

FHR-2 can be measured with our HK3004 assay, designed to precisely quantify this intriguing glycoprotein.

PRINCIPLE OF THE ASSAY

The human FHR-2 ELISA is a ready-to-use solid-phase enzyme-linked immunosorbent assay based on the sandwich principle with a working time of 1 hour and 15 minutes. The efficient format of a plate with twelve disposable 8-well strips allows free choice of batch size for the assay. Samples and standards are incubated in microtiter wells coated with antibodies recognizing human FHR-2. Peroxidase-conjugated antibody will bind to the captured FHR-2. Peroxidase-conjugate will react with the substrate, tetramethylbenzidine (TMB). The enzyme reaction is stopped by the addition of oxalic acid. The absorbance at 450 nm is measured with a spectrophotometer. A standard curve is obtained by plotting the absorbance (linear) versus the corresponding concentrations of the human FHR-2 standards (log). The human FHR-2 concentration of samples, which are run concurrently with the standards, can be determined from the standard curve.

KIT FEATURES

- Working time of 1 ¼ hours.
- Minimum concentration which can be measured is 0.88 ng/ml.
- Measurable concentration range of 0.88 to 10 ng/ml.
- Working volume of 100 µl/well.



CROSS REACTIVITY

Potential cross-reacting proteins detected in the Complement FHR-2 ELISA:

Cross reactant	Reactivity
FH (native)	Negative
FHR-1 (recombinant)	Negative
FHR-3 (recombinant)	Negative
FHR-4A (recombinant)	Negative
FHR-4B (recombinant)	Negative
FHR-5 (recombinant)	Negative

Table 1

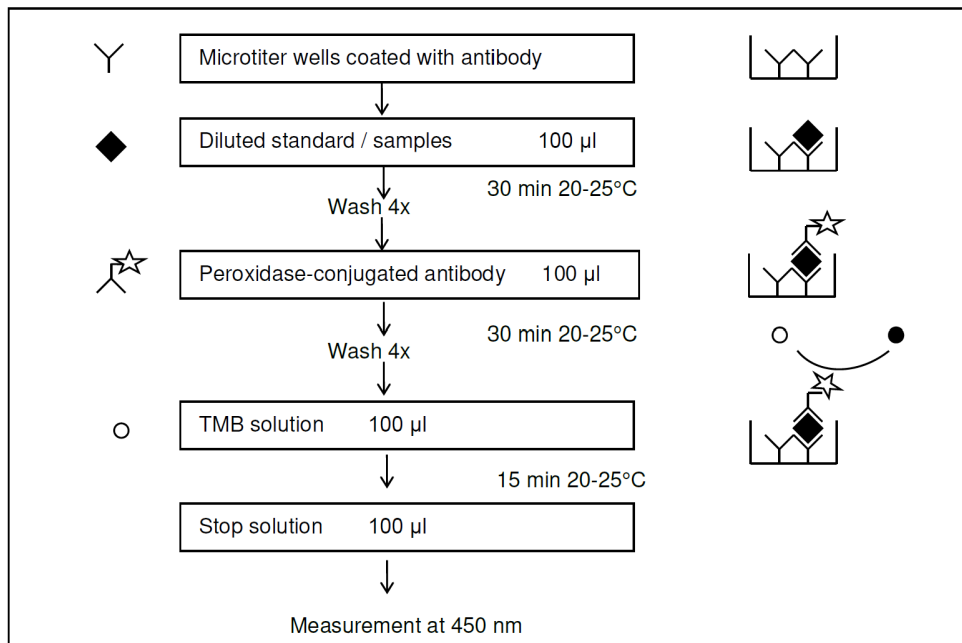
Potential cross-reacting species detected in the Complement FHR-2 ELISA:

Cross reactant	Reactivity
Rat	Negative
Mouse	Negative
Pig	Negative
Monkey (Cynomolgus)	Positive

Table 2

Cross-reactivity for other species or proteins/peptides has not been tested.

PROTOCOL OVERVIEW





- The human FHR-2 ELISA is a ready-to-use solid-phase enzyme-linked immunosorbent assay based on the sandwich principle with a working time of 1 hour and 15 minutes.
- The efficient format of a plate with twelve disposable 8-well strips allows free choice of batch size for the assay.
- Samples and standards are incubated in microtiter wells coated with antibodies recognizing human FHR-2.
- Peroxidase-conjugated antibody will bind to the captured FHR-2.
- Peroxidase-conjugate will react with the substrate, tetramethylbenzidine (TMB).
- The enzyme reaction is stopped by the addition of oxalic acid.
- The absorbance at 450 nm is measured with a spectrophotometer. A standard curve is obtained by plotting the absorbance (linear) versus the corresponding concentrations of the human FHR-2 standards (log).
- The human FHR-2 concentration of samples, which are run concurrently with the standards, can be determined from the standard curve.

KIT COMPONENTS AND STORAGE INSTRUCTIONS

Kit Component	Quantity HK3004-01	Quantity HK3004-02	Color Code
Wash buffer 20x	1 vial (60 ml)	1 vial (60 ml)	Colorless
Dilution Buffer 10x	1 vial (15 ml)	2 vials (30 ml)	Green
Standard	2 vials, lyophilized	4 vials, lyophilized	White
Peroxidase-conjugated antibody	1 vial, 1 ml lyophilized	2 vials, 1 ml lyophilized	Blue
TMB substrate	1 vial (11 ml)	1 vial (22 ml)	Brown
Stop solution	1 vial (22 ml)	1 vial (22 ml)	Red
12 Microtiter strips, pre-coated	1 plate	2 plates	
Certificate of Analysis	1	1	
Manual	1	1	
Data collection sheet	2	2	

Table 3

- Store the kit components at 2 - 8°C immediately upon receipt. Do not freeze.
- Ensure components are used before the expiration date indicated on the kit label.
- The standard and peroxidase-conjugated antibody in lyophilized form are stable until the expiration date indicated on the kit label, if stored at 2 - 8°C.
- The exact quantity of the standard is specified on the vial label and the Certificate of Analysis.
- The standard is intended for a single use; it cannot be stored after reconstitution.



- Ensure that the foil pouch containing the plate is vacuum-sealed and undamaged upon receipt. Any deviation may impact assay performance.
- Immediately return unused strips back into the foil pouch with the desiccant, sealing it fully along the entire edge of the zip-seal. Stored at 2 - 8°C, the quality is ensured for one month.

MATERIALS REQUIRED BUT NOT PROVIDED

- Calibrated micropipettes and disposable tips.
- Distilled or de-ionized water.
- Plate washer: automatic or manual.
- Polypropylene tubes.
- Calibrated ELISA plate reader capable of measuring absorbance at 450 nm.
- Centrifuge for 1 ml tubes.

WARNINGS AND PRECAUTIONS

- This product is intended for research purposes only and not for use in diagnostic or therapeutic procedures.
- Only qualified personnel trained in laboratory procedures should handle this kit.
- Under no circumstances should sodium azide be added to any component as a preservative.
- Refrain from using kit components beyond their expiration date.
- To ensure accuracy, do not interchange reagents from different kits or lots. Each kit and lot is calibrated as a complete unit; use only the reagents supplied by the manufacturer.
- The assay is specifically optimized for the stated standard range. Alterations to the standard range are not recommended.
- Exercise caution when opening vials as they are under vacuum.
- Avoid ingestion of any kit components.
- The kit reagents include 2-chloroacetamide, a preservative known for its harmful effects upon skin contact and toxicity if ingested. In the event of an accident or discomfort, immediate medical consultation is advised.
- Protect the TMB substrate from intense light exposure; it should remain colorless until utilized.
- The stop solution contains 2% oxalic acid, a substance that can irritate or burn the respiratory system, skin, and eyes. Avoid any direct contact, and in case of exposure, rinse thoroughly with water and seek medical attention.
- Deviations from the specified incubation times, temperatures, or pipetting volumes may result in inaccurate results.
- Once dispensed, avoid reusing microwells or returning reagents to their original bottles.
- Treat all biological samples as potentially hazardous or infectious and handle them under conditions that minimize the risk of disease transmission.
- Be aware that samples that are hemolyzed, hyperlipemic, heat-treated, or contaminated may yield inaccurate results.
- Utilize polypropylene tubes for the preparation of standards and samples, avoiding the use of polystyrene tubes or sample plates.



- The standard is derived from human sources and has been tested for various viruses with negative results. However, as no testing method can guarantee the complete absence of infectious agents, treat this reagent with the same precautions as you would any potentially infectious human serum or blood specimen. Follow established guidelines for preventing the transmission of blood-borne infections when handling materials in contact with this reagent.

SAMPLE PREPARATION

The recommended sample medium is EDTA plasma. All samples should be kept on ice whilst performing the assay.

Collection and handling

Serum and Plasma

Collect blood using normal aseptic techniques. Blood samples should be kept on ice. If serum is used, separate serum from blood after clotting at room temperature within one hour by centrifugation (1,500xg at 4°C for 15 min). Transfer the serum to a fresh polypropylene tube. If plasma is used, separate plasma from blood within 20 minutes after blood sampling by centrifugation (1,500xg at 4°C for 15 min). Transfer the plasma to a fresh polypropylene tube. Most reliable results are obtained if EDTA plasma is used.

Storage

Store samples below -20°C, preferably at -70°C in polypropylene tubes. Storage at -20°C can affect recovery of human FHR-2. Preferable use samples within 1-2 hours after thawing. Avoid multiple freeze-thaw cycles which may cause loss of human FHR-2 activity and give erroneous results. Do not use hemolyzed, hyperlipemic, heat-treated or contaminated samples. Before performing the assay, samples should be mixed gently. Prepare all samples (controls and test samples) prior to starting the assay procedure. Avoid foaming.

Dilution procedures

Serum or plasma samples

Due to expected high levels of FHR-2 it is advised to apply a 400 – 3200x dilution range for accurate measurement of serum or plasma samples with supplied dilution buffer in polypropylene tubes. This recommended dilution applies to healthy samples, the optimal dilution for diseased samples should be tested. Optimal dilution is dependent on sample quality and expected FHR-2 quantity and should be determined for the specific sample set/study.

Comment regarding recommended sample dilution

The mentioned dilution for samples is a minimum dilution and should be used as a guideline. The recovery of FHR-2 from an undiluted sample is not 100% and may vary from sample to sample. When testing less diluted samples it is advisable to run recovery experiments to determine the influence of the matrix on the detection of FHR-2.

Do not use polystyrene tubes or sample plates for preparation or dilution of the samples.



Guideline for dilution of samples

Please see Table 4 for recommended sample dilutions. Volumes are based on a total volume of at least 230 μ l of diluted sample, which is sufficient for one sample in duplicate in the ELISA. For dilution of samples we recommend to use at least 10 μ l of sample.

Dilution	Pre-Dilution	Amount of sample or pre-dilution required	Amount of dilution buffer required
100x	Not necessary	10 μ l (sample)	990 μ l
400x	Recommended: 100x (see nr. 1)	75 μ l (pre-dilution)	225 μ l
800x	Recommended: 100x (see nr. 1)	50 μ l (pre-dilution)	350 μ l
1200x	Recommended: 100x (see nr. 1)	50 μ l (pre-dilution)	550 μ l
1600x	Recommended: 100x (see nr. 1)	25 μ l (pre-dilution)	375 μ l
2000x	Recommended: 100x (see nr. 1)	25 μ l (pre-dilution)	475 μ l
2500x	Recommended: 100x (see nr. 1)	10 μ l (pre-dilution)	240 μ l
3200x	Recommended: 100x (see nr. 3)	10 μ l (pre-dilution)	310 μ l

Table 4

REAGENT PREPARATION

The standard should be kept on ice. Allow all remaining reagents to equilibrate to room temperature (20 – 25°C) prior to use.

Allow all reagents to equilibrate to room temperature (20 – 25°C) prior to use. Return to proper storage conditions immediately after use.

Wash buffer

Prepare wash buffer by mixing 60 ml of 20x wash buffer with 1140 ml of distilled or de-ionized water, which is sufficient for 2 x 96 tests. In case less volume is required, prepare the desired volume of wash buffer by diluting 1 part of the 20x wash buffer with 19 parts of distilled or deionized water.

Dilution buffer

Prepare dilution buffer by mixing 15 ml of the 10x dilution buffer with 145 ml of distilled or deionized water, which is sufficient for 1 x 96 tests. In the instance of 2 x 96 tests, prepare the dilution buffer by mixing 30 ml of the 10x dilution buffer with 270 ml of distilled or de-ionized water. In case less volume is required, prepare the desired volume of dilution buffer by diluting 1 part of the 10x dilution buffer with 9 parts of distilled or de-ionized water. Concentrated dilution buffer may contain crystals. In case the crystals do not disappear at room temperature within 1 hour, concentrated dilution buffer can be warmed up to 37°C. Do not shake the solution.



Standard solution

The standard is reconstituted by pipetting the amount of dilution buffer mentioned on the CoA in the standard vial. Use the standard vial as Tube 1 in Figure 2. Prepare each FHR-2 standard in polypropylene tubes by serial dilution of the reconstituted standard with dilution buffer as shown in Figure 2*. After reconstitution the standard cannot be stored for repeated use.

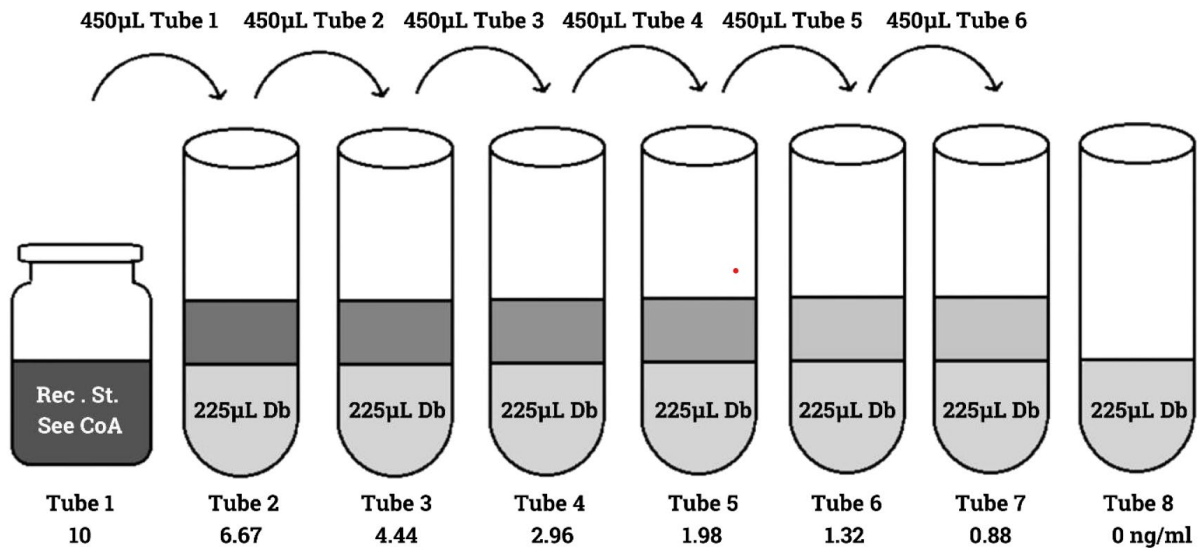


Figure 2

*) CoA: Certificate of Analysis, St: Standard, Db: Dilution buffer

Peroxidase-conjugated antibody

The peroxidase-conjugated antibody is reconstituted by pipetting 1 ml distilled or de-ionized water. Dilute the reconstituted 1 ml detection antibody with 11 ml dilution buffer, which is sufficient for 1 x 96 tests. In case less volume is required, prepare the desired volume of peroxidase-conjugated antibody by diluting 1 part of the reconstituted vial with 11 parts of dilution buffer.

ELISA PROTOCOL

Bring all reagents to room temperature (20 - 25°C) before use.

1. Determine the number of test wells required, put the necessary microwell strips into the supplied frame, and fill out the data collection sheet. Return the unused strips to the storage bag with desiccant, seal and store at 2 - 8°C.
2. Transfer 100 µl in duplicate of standard, samples, or controls into appropriate wells. Do not touch the side or bottom of the wells.
3. Cover the tray and tap the tray to eliminate any air bubbles. Be careful not to splash liquid onto the cover.
4. Incubate the strips or plate for 30 minutes at room temperature.



5. Wash the plates 4 times with wash buffer as follows*:
 - a. Carefully remove the cover, avoid splashing.
 - b. Empty the plate by inverting plate and shaking contents out over the sink, keep inverted and tap dry on a thick layer of tissues.
 - c. Add 200 μ l of 1x wash buffer to each well, wait 20 seconds, empty the plate as described in 5b.
 - d. Repeat the washing procedure 5b/5c three times.
 - e. Empty the plate and gently tap on thick layer of tissues.
6. Add 100 μ l of diluted peroxidase-conjugated antibody to each well using the same pipetting order as applied in step 2. Do not touch the side or bottom of the wells.
7. Cover the tray and incubate the tray for 30 minutes at room temperature.
8. Repeat the wash procedure described in step 5a-e.
9. Add 100 μ l of TMB substrate to each well, using the same pipetting order as applied in step 2. Do not touch the side or bottom of the wells.
10. Cover the tray and incubate the tray for 15 minutes at room temperature. It is advised to control the reaction on the plate regularly. In the case of strong development the TMB reaction can be stopped sooner. Avoid exposing the microwell strips to direct sunlight. Covering the plate with aluminum foil is recommended.
11. Stop the reaction by adding 100ul of the stop solution with the same sequence and timing as used in step 9. Mix solutions in the wells thoroughly by gently swirling the plate. Gently tap the tray to eliminate any air bubbles trapped in the wells.
12. Read the plate within 30 minutes after addition of stop solution at 450 nm using a plate reader, following the instructions provided by the instrument's manufacturer.

*) In case plate washer is used, please note: use of a plate washer can result in higher background and decrease in sensitivity. We advise validation of the plate washer with the manual procedure. Make sure the plate washer is used as specified for the manual method.

INTERPRETATION OF RESULTS

- Determine the average absorbance for each group of duplicate standards, controls, and samples.
- Discrepancies exceeding 15% from the mean absorbance value suggest potential inaccuracies, necessitating sample reanalysis.
- Ensure the mean absorbance of the zero standard does not surpass 0.3.
- Utilize specialized software to construct a standard curve, plotting mean absorbance values (Y-axis) against corresponding concentrations (X-axis) on a logarithmic scale.
- For diluted samples, adjust the concentration derived from the standard curve by the dilution factor.
- Samples yielding an average absorbance higher than that of the maximum standard concentration exceed the assay's scope and must be reanalyzed using a greater dilution factor.

TECHNICAL HINTS

- Technicians should be proficient and well-versed in ELISA assays and the specific test procedures before initiating the assay.



- For those unfamiliar with ELISA techniques, it is advisable to conduct a preliminary assay with a standard curve to ensure understanding and adherence to the protocol before proceeding with sample evaluations.
- Accurate and thorough washing is critical at all stages of the assay to prevent false positive or negative outcomes. Ensure complete removal of liquids from wells prior to adding wash buffer, adhere strictly to the specified volume for each washing cycle, and avoid allowing the wells to remain uncovered or dry for prolonged periods.
- A standard curve is essential for each assay run due to varying conditions; samples must be evaluated against a standard curve established on the same plate during that session.
- Do not interchange reagents from different kits or batches, including strips, and avoid combining remnants with contents from new vials.
- Prepare fresh dilutions of the standard, samples, biotinylated LPS, streptavidinperoxidase, and buffers each time the kit is utilized.
- Maintain cap-to-vial correspondence; caps are designed to fit their original vials and should not be swapped.
- Prevent cross-contamination by using new pipette tips for each addition across standards, samples, and reagents, and employ separate reservoirs for each reagent to ensure integrity.
- Dispose of all waste in accordance with the established laboratory safety protocols and regulations.

QUALITY CONTROL

The Certificate of Analysis accompanying this kit is specific to its lot number, intended to verify the results achieved in your facility. Please note that the absorbance values indicated on the Certificate serve merely as reference points; deviations in outcomes produced by your laboratory are to be expected. Designed to mitigate the impact of soluble receptors, binding proteins, and extraneous variables present in biological samples, this assay aims for precise measurement free from external interferences. Nonetheless, without exhaustive testing of all possible variables, the exclusion of interference cannot be entirely assured.

To ensure the highest efficacy of this kit, implementing good laboratory practices (GLP) is crucial.



Warranty Information

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

For further information about this kit, its application or the procedures in this kit, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 866-411-8023.