

ab316915 GFP SimpleStep ELISA® Kit – Extracellular

For the quantitative measurement of GFP in human and mouse cell culture supernatant.
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit: www.abcam.com/ab316915

Storage and Stability: Store kit at 2-8°C immediately upon receipt. Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Standard Preparation and Reagent preparation sections.

Materials Supplied

Item	Quantity	Storage Condition
GFP Capture Antibody 10X	600 µL	+4°C
GFP Detector Antibody 10X	600 µL	+4°C
GFP Lyophilized Recombinant Protein	2 Vials	+4°C
Antibody Diluent CPR2	6 mL	+4°C
Cell Extraction Enhancer Solution 50X	1 mL	+4°C
Sample Diluent NS	50 mL	+4°C
Wash Buffer PT 10X	20 mL	+4°C
TMB Development Solution	12 mL	+4°C
Stop Solution	12 mL	+4°C
SimpleStep Pre-Coated 96-Well Microplate	96 wells	+4°C
Plate Seal	1	+4°C

Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

Microplate reader capable of measuring absorbance at 450 or 600 nm.
Deionized water.
Multi- and single-channel pipettes.
Tubes for standard dilution.
Plate shaker for all incubation steps.
Optional: Phenylmethylsulfonyl Fluoride (PMSF) (or other protease inhibitors).

Reagent Preparation

Equilibrate all reagents to room temperature (18-25°C) prior to use. The kit contains enough reagents for 96 wells. The sample volumes below are sufficient for 48 wells (6 x 8-well strips); adjust volumes as needed for the number of strips in your experiment.

Prepare only as much reagent as is needed on the day of the experiment. Capture and Detector Antibodies have only been tested for stability in the provided 10X formulations.

The provided Cell Extraction Enhancer Solution 50X may precipitate when stored at + 4°C. To dissolve, warm briefly at + 37°C and mix gently. The Cell Extraction Enhancer Solution 50X can be stored at room temperature to avoid precipitation.

Sample Diluent NS + 1X Enhancer: Prepare Sample Diluent NS + 1X Enhancer by combining Sample Diluent NS and 50X Cell Extraction Enhancer Solution. To make 10 mL Sample Diluent NS + 1X Enhancer, combine 9.8 mL Sample Diluent NS and 200 µL Cell Extraction Enhancer Solution 50X. Mix thoroughly and gently.

1X Wash Buffer PT: Prepare 1X Wash Buffer PT by diluting Wash Buffer PT 10X with deionized water. To make 50 mL 1X Wash Buffer PT combine 5 mL Wash Buffer PT 10X with 45 mL deionized water. Mix thoroughly and gently.

Antibody Cocktail: Prepare Antibody Cocktail by diluting the capture and detector antibodies in Antibody Diluent CPR2. To make 3 mL of the Antibody Cocktail combine 300 µL 10X Capture Antibody and 300 µL 10X Detector Antibody with 2.4 mL Antibody Diluent CPR2. Mix thoroughly and gently.

Standard Preparation

Always prepare a fresh set of standards for every use. Discard working standard dilutions after use as they do not store well. The following section describes the preparation of a standard curve for duplicate measurements (recommended).

1. Reconstitute the GFP standard sample by adding the volume of **Sample Diluent NS + 1X Enhancer** indicated on the protein vial label. Hold at room temperature for 10 minutes. Mix thoroughly and gently. This is the 9,600 pg/mL **Stock Standard** Solution.
2. Label eight tubes, Standards 1–8.
3. Add 375 µL of **NS + 1X Enhancer** into tube number 1 and 150 µL of **Sample Diluent NS + 1X Enhancer** into numbers 2-8.
4. Use the **Stock Standard** to prepare the following dilution series. Standard #8 contains no protein and is the Blank control:

Standard #	Dilution Sample	Volume to Dilute (µL)	Volume of Diluent (µL)	Starting Conc. (pg/mL)	Final Conc. (pg/mL)
1	Stock Standard	25	375	9,600	600
2	Standard#1	150	150	600	300
3	Standard#2	150	150	300	150
4	Standard#3	150	150	150	75
5	Standard#4	150	150	75	37.5
6	Standard#5	150	150	37.5	18.75
7	Standard#6	150	150	18.75	9.38
8	Blank Control	0	150	0	0

Sample Preparation

Typical Sample Dynamic Range	
Sample Type	Range
HeLa Cell Supernatant	≤50%
Neuro 2A Cell Supernatant	≤50%
Mouse Lung Tissue Supernatant	≤50%

Note: This kit is compatible with samples containing 0.1% SDS. If sample contains higher concentration of SDS, dilute initially the sample with Sample Diluent NS without Enhancer to lower SDS concentration to 0.1%. Following dilutions can be done in **Sample Diluent NS + 1X Enhancer**.

Cell Culture Supernatants Centrifuge cell culture media at 2,000 x g for 10 minutes to remove debris. Collect supernatants. Dilute samples at least 1:2 into **Sample Diluent NS + 1X Enhancer** and assay. Store un-diluted samples at -20°C or below. Avoid repeated freeze-thaw cycles.

Plate Preparation

The 96 well plate strips included with this kit are supplied ready to use. It is not necessary to rinse the plate prior to adding reagents.

Unused plate strips should be immediately returned to the foil pouch containing the desiccant pack, resealed and stored at 4°C.

For each assay performed, a minimum of two wells must be used as the zero control.

For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates).

Differences in well absorbance or “edge effects” have not been observed with this assay.

Assay Procedure

Equilibrate all materials and prepared reagents to room temperature prior to use.

We recommend that you assay all standards, controls and samples in duplicate.

1. Prepare all reagents, working standards, and samples as directed in the previous sections.
2. Remove excess microplate strips from the plate frame, return them to the foil pouch containing the desiccant pack, resealed and return to 4°C storage.
3. Add 50 µL of all sample or standard to appropriate wells.
4. Add 50 µL of the Antibody Cocktail to each well.
5. Seal the plate and incubate for 1 hour at room temperature on a plate shaker set to 400 rpm.
6. Wash each well with 3 x 350 µL 1X Wash Buffer PT. Wash by aspirating or decanting from wells then dispensing 350 µL 1X Wash Buffer PT into each well. Wash Buffer PT should remain in wells for at least 10 seconds. Complete removal of liquid at each step is essential for good performance. After the last wash invert the plate and tap gently against clean paper towels to remove excess liquid.
7. Add 100 µL of TMB Development Solution to each well and incubate for 10 minutes in the dark on a plate shaker set to 400 rpm.

Given variability in laboratory environmental conditions, optimal incubation time may vary between 5 and 20 minutes.

Note: The addition of Stop Solution will change the color from blue to yellow and enhance the signal intensity about 3X. To avoid signal saturation, proceed to the next step before the high concentration of the standard reaches a blue color of O.D.600 equal to 1.0.

8. Add 100 µL of Stop Solution to each well. Shake plate on a plate shaker for 1 minute to mix. Record the OD at 450 nm. This is an endpoint reading.
9. Alternative to 7 – 8: Instead of the endpoint reading at 450 nm, record the development of TMB Substrate kinetically. Immediately after addition of TMB Development Solution begin recording the blue color development with elapsed time in the microplate reader prepared with the following settings:

Mode	Kinetic
Wavelength:	600 nm
Time:	up to 20 min
Interval:	20 sec - 1 min
Shaking:	Shake between readings

Note that an endpoint reading can also be recorded at the completion of the kinetic read by adding 100 µL Stop Solution to each well and recording the OD at 450 nm.

Download our ELISA guide for technical hints, results, calculation, and troubleshooting tips:

www.abcam.com/protocols/the-complete-elisa-guide

For technical support contact information, visit: www.abcam.com/contactus

ab316915 – GFP SimpleStep ELISA® Kit

Additional information

ASSAY SPECIFICITY

This kit is designed for the quantification of GFP.

The standard protein in this kit is full length enhanced GFP (eGFP).

This kit recognizes the GFP, eGFP, as well as the SDS-denatured eGFP.

500 pg/mL of recombinant GFP was tested for reactivity. 92% reactivity was observed.

600 pg/mL of recombinant SDS-denatured EGFP was tested for reactivity. 89% reactivity was observed.

Spiked protein experiments were used to validate human and mouse cell and tissue culture supernatant sample types.

50% HeLa, Neuro 2A, and mouse lung cell culture supernatants were measured in duplicate. All values were below the detectable range of the assay.

Serum, plasma (citrate), plasma (EDTA), plasma (heparin), urine, milk, saliva, CSF, cell extract, and tissue extract samples have not been tested with this kit.

For the measurement of GFP in cell extract and tissue extract sample types, use GFP ELISA kit ab316916.

CROSS REACTIVITY

6,000 pg/mL of recombinant YFP and 6,000 pg/mL of recombinant mCherry were tested for cross reactivity. No cross reactivity was observed.

INTERFERENCE

6,000 pg/mL of recombinant YFP and 6,000 pg/mL of recombinant mCherry were tested for interference with 600 pg/mL of recombinant GFP. No interference was observed.

CALCULATION

- Calculate the average absorbance value for the blank control (zero) standards. Subtract the average blank control standard absorbance value from all other absorbance values.
- Create a standard curve by plotting the average blank control subtracted absorbance value for each standard concentration (y-axis) against the target protein concentration (x-axis) of the standard. Use graphing software to draw the best smooth curve through these points to construct the standard curve.
 Δ Note: Most microplate reader software or graphing software will plot these values and fit a curve to the data. A four parameter curve fit (4PL) is often the best choice; however, other algorithms (e.g. linear, semi-log, log/log, 4 parameter logistic) can also be tested to determine if it provides a better curve fit to the standard values.
- Determine the concentration of the target protein in the sample by interpolating the blank control subtracted absorbance values against the standard curve. Multiply the resulting

value by the appropriate sample dilution factor, if used, to obtain the concentration of target protein in the sample.

- Samples generating absorbance values greater than that of the highest standard should be further diluted and reanalyzed. Similarly, samples which measure at absorbance values less than that of the lowest standard should be retested in a less dilute form.

TYPICAL DATA

Typical standard curve – data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

Standard Curve Measurements			
Concentration (pg/mL)	O.D 450 nm		Mean O.D
	1	2	
0	0.040	0.040	0.040
9.38	0.095	0.097	0.096
18.75	0.156	0.155	0.156
37.5	0.239	0.225	0.232
75	0.407	0.400	0.404
150	0.713	0.704	0.708
300	1.586	1.536	1.561
600	2.851	2.750	2.800

Table 1. Example of species GFP standard curve in **Sample Diluent NS + 1X Enhancer**. The GFP standard curve was prepared as described in the Standard Preparation section. The table shows raw data values.

TYPICAL SAMPLE VALUES

Sensitivity:

The calculated minimal detectable dose (MDD) is 1.8 pg/mL. The MDD was determined by calculating the mean of zero standard replicates (n=16) and adding 2 standard deviations then extrapolating the corresponding concentration.

Recovery

Three concentrations of eGFP were spiked in duplicate to the indicated biological matrix to evaluate signal recovery in the working range of the assay.

Sample Type	Average % Recovery	Range (%)
50% HeLa Cell Culture Supernatant*	98	94 - 104
50% Neuro 2A Cell Culture Supernatant*	101	91 - 108
50% Mouse Lung Tissue Culture Supernatant**	86	79 - 93

*Media is DMEM containing 10% fetal bovine serum

** Media is RPMI 1640 containing 10% fetal bovine serum

Linearity of Dilution

Linearity of dilution is determined based on interpolated values from the standard curve. Linearity of dilution defines a sample concentration interval in which interpolated target concentrations are directly proportional to sample dilution.

Recombinant eGFP was spiked into the following biological samples in a 2-fold dilution series. Sample dilutions are made in **Sample Diluent NS + 1X Enhancer**.

Dilution Factor	Interpolated value	50% HeLa Cell Supernatant	50% Neuro 2A Cell Supernatant	50% Mouse Lung Tissue Supernatant
Undiluted	pg/mL	237.0	279.8	287.0
	% Expected value	100%	100%	100%
2	pg/mL	111.2	164.3	146.1
	% Expected value	94%	117%	102%
4	pg/mL	57.6	82.7	69.9
	% Expected value	97%	118%	97%
8	pg/mL	33.0	39.6	37.7
	% Expected value	111%	113%	105%
16	pg/mL	16.8	16.6	15.5
	% Expected value	113%	95%	86%

Precision

Mean coefficient of variations of interpolated values of GFP from two concentrations of HeLa cell culture supernatant within the working range of the assay.

	Intra-assay	Inter-assay
N=	8	3
CV (%)	8.8	7.4

Download our ELISA guide for technical hints, results, calculation, and troubleshooting tips:

www.abcam.com/protocols/the-complete-elisa-guide

For technical support contact information, visit: www.abcam.com/contactus

Copyright © 2024 Abcam. All Rights Reserved. All information / detail is correct at time of going to print.
Version 1 | 2024-03-05