

Using FRET-based Measurements of Protein Complexes to Determine Stoichiometry with the Job Plot

Francesca Mattioli¹, Yajie Gu^{1,2}, and Karolin Luger¹

¹HHMI, University of Colorado Boulder, Dept. Chem. & Biochem., Boulder, CO, USA; ²Colorado State University, Dept. Biochem. & Mol. Biol., Fort Collins, CO, USA

- Determination of stoichiometry of protein-protein interactions is based on observed FRET maximum
- The CLARIOstar was used in a 384-well plate format to measure FRET and relevant controls for up to 4 interactions
- Determination of stoichiometry is assisted by a MARS data analysis template

Introduction

Measurement of stoichiometry is important for understanding biochemical reactions since these reactions are dependent on the interaction of at least 2 cellular components. Measuring these interactions has been limited to approaches that are useful for observing large molecular weight changes such as size-exclusion chromatography. Furthermore, these approaches require a large amount of purified cellular components.

With these limitations in mind we sought to provide a platform where protein-protein interactions could be observed. Herein we describe an adaptation of the Job plot¹. Our approach uses a microplate reader, which was used to read FRET and relevant acceptor and donor control intensities. Using a microplate reader enables the use of small sample volumes (40 µl) in 384-well plates. Furthermore, 4 separate interactions can be studied on an individual plate.

Assay Principle

A key factor of the Job plot, also called the continuous variation method, is that the total concentration of the 2 molecules is held constant². It is the ratio of the 2 molecules that is changed. An observable parameter that is proportional to complex formation is employed, in this case FRET between the labelled proteins. Figure 1 depicts the expected assay results for a 1:1 protein-protein interaction.

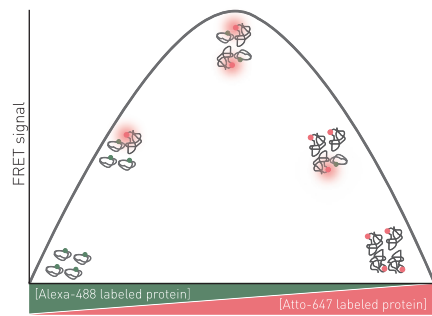


Fig. 1: Assay Principle for FRET Job plot

In the conditions where 100% donor labelled protein are present there is no FRET signal as the ratio shifts to larger amounts of acceptor the FRET signal increases until maximum FRET is obtained. Maximum FRET will indicate the stoichiometry of the interaction. Continued changes in the ratio toward greater acceptor will be associated with a decrease in FRET until zero FRET is again seen with 100% acceptor labelled protein.

Materials & Methods

- Black 384-well microplates (Corning)
- BMG LABTECH CLARIOstar
- For a complete list of reagents and procedure please refer to Mattioli et al.¹

Experimental Procedure

Stock solutions with a 1 µM concentration were prepared for each protein, both labeled and unlabeled. Further dilution of these stocks were created such that diluted stocks have 2X concentration needed for the well reactions. Final well reactions are created by combining 20 µl each of the appropriate diluted stocks for proteins 1 and 2. This can be done according to the scheme depicted in Figure 2.

	1		2		3		4		5		6		
C	B	A	B	A	B	A	B	B	A	B	A		
D	S1A	S1A	S1B	S1B	S1C	S1C							Buffer control Protein 1 Protein 2 Protein 1/Protein 2 ratio 1:0 11:1 5:1 3:1 2:1 1.4:1 1:1 1:1.4 1:2 1:3 1:5 1:11
E	S2A	S2A	S2B	S2B	S2C	S2C							
F	S3A	S3A	S3B	S3B	S3C	S3C							
G	S4A	S4A	S4B	S4B	S4C	S4C							
H	S5A	S5A	S5B	S5B	S5C	S5C							
I	S6A	S6A	S6B	S6B	S6C	S6C							
J	S7A	S7A	S7B	S7B	S7C	S7C							
K	S8A	S8A	S8B	S8B	S8C	S8C							
L	S9A	S9A	S9B	S9B	S9C	S9C							
M	S10A	S10A	S10B	S10B	S10C	S10C							
N	S11A	S11A	S11B	S11B	S11C	S11C							
O	S12A	S12A	S12B	S12B	S12C	S12C							
P	S13A	S13A	S13B	S13B	S13C	S13C							
	donor only control		acceptor only control		FRET experiment								

Fig. 2: 384-well plate preparation layout



Instrument settings

Because of the multichromatic nature of the test performed appropriate setting of the gain for each chromatic was an important consideration. To measure the acceptor fluorescence gain was set using a sample with highest acceptor dye and no donor, for example well P3. Similarly, gain for donor fluorescence measurement was set on well with highest donor dye and no acceptor dye, such as well D1. For FRET measurement plate is first read and the well with max FRET signal is found. This well is used to perform gain adjustment.

Optic settings	Fluorescence, multichromatic, endpoint	
	Chromatic 1: Alexa 488 preset	
	LVF Ex	488-14
	Dichroic	Auto: 507.5
	LVF Em	535-30
	Chromatic 2: Atto 647 preset	
	LVF Ex	625-30
	Dichroic	Auto: 647.5
	LVF Em	680-40
	Chromatic 3: Alexa 488/Atto 647 FRET	
	LVF Ex	488-15
	Dichroic	Auto: 577.8
LVF Em	680-40	
Gain	Adjusted as described	
General settings	Number of flashes	50
	Settling time	0.1 s

Results & Discussion

As proof of principle we first looked at the interaction between histone binding protein and histone H3-H4. In the layout described in figure 2 histone binding protein is protein 1 and H3-H4 is protein 2. Figure 3 show the expected 1:1 binding interaction.

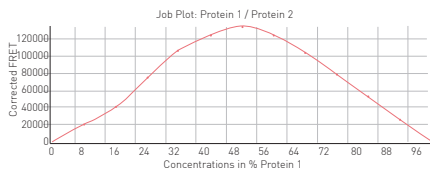


Fig. 3: Job plot for interaction between H3-H4 and histone binding protein

Figure 4 shows the results of 2 additional protein interaction tests. As you can see one of the interactions also exhibits a 1:1 stoichiometry while the other is an example of a 2:1 interaction stoichiometry.

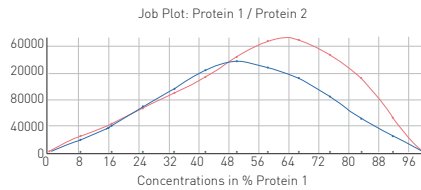


Fig. 4: Job plot depicting 1:1 and 2:1 protein-protein interaction stoichiometry

Conclusion

The Job plot to assess protein-protein interaction stoichiometry has been successfully adapted to a microplate reader-based system. This enables both miniaturization to save on protein components and improved throughput, up to 4 protein pairs can be studied in one 384-well plate.

References

1. Mattioli, F. et al. FRET-based Stoichiometry Measurements of Protein Complexes *in vitro*. *Bio. Protoc.* (2018) 8: e2713. DOI: 10.21769/BioProtoc.2713
2. Huang, C.Y. Determination of Binding Stoichiometry by the Continuous Variation Method: The Job Plot. *Methods Enzymol.* (1982) 87: 509-525

