

AQT0661 – JAK3 (JH1) Assay Validation

PhosphoSens[®]-Red Assay Format

Outline for this Study

PhosphoSens-Red Assay Validation



Enzyme Source, Construct, and Lot Information:

Carna JAK3(JH1) (Cat/Lot #, 08-046/19CBS-0798B) amino acids 795-1124, N-term His tag

Experiments to be run:

PhosphoSens-Kinetic Assay

PhosphoSens-Red TRF End Point S:B Determination

Has this Sensor Peptide been fully validated in the PhosphoSens-Kinetic Assay Format?

No

Red-Shift TRF S:B Determination

Reaction Conditions and Set Up

Reaction Conditions:

54 mM HEPES, pH 7.5

1 mM ATP

1.2 mM DTT

0.012% Brij-35

1% glycerol

0.2 mg/ml BSA

0.55 mM EGTA

10 mM MgCl₂

15 μM AQT0661

2 nM JAK3(JH1)

Notes:

Enzyme Dilution Buffer (EDB): 20 mM HEPES, pH 7.5, 0.01% Brij-35, 5% Glycerol, 0.5 mM EGTA, 1 mM DTT, 1 mg/ml Bovine Serum Albumin.

Reaction Set Up:

2 or 2.5 μL

10x Sensor Peptide

14 or 17.5 μL

Reaction Mix with ATP & DTT

4 or 5 μL

1x EDB or Kinase dilutions (5x in EDB)

20 or 25 μL

Final reaction volume

Reactions were run at 30°C for 240 minutes in either Corning, low volume 384-well, white flat round bottom polystyrene NBS microplates (Cat. #3824) at 25 μL final well volume or in PerkinElmer, ProxiPlate-384 Plus, white shallow well microplates (Cat. #6008280) at 20 μL final well volume after sealing using optically-clear adhesive film (TopSealA-Plus plate seal, PerkinElmer [Cat. #6050185]) in a Biotek Synergy Neo 2 microplate reader with excitation (360 nm) and emission (485 nm) wavelengths.

Red-Shift Reaction Set Up:

Following completion of the kinase reaction, remove plate from the reader and remove plate seal and add Europium to a final concentration of 5 mM.

4 or 5 μL

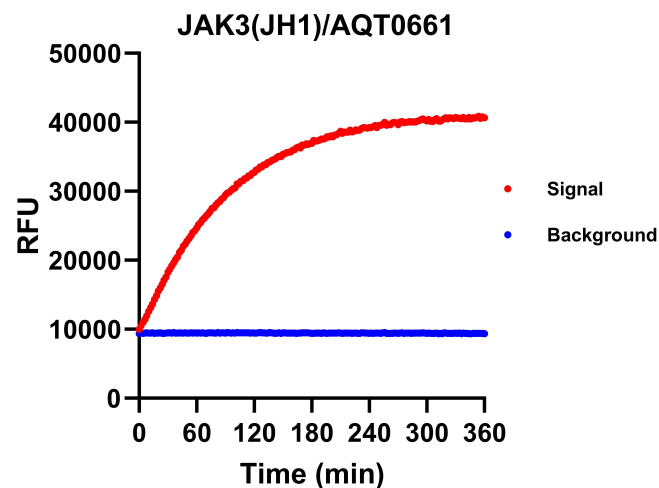
30 mM Europium (6x)

Incubate for 5 minutes at room temperature (RT), return to plate reader without seal and read in TRF mode with excitation (360 nm) and emission (620 nm) wavelengths.

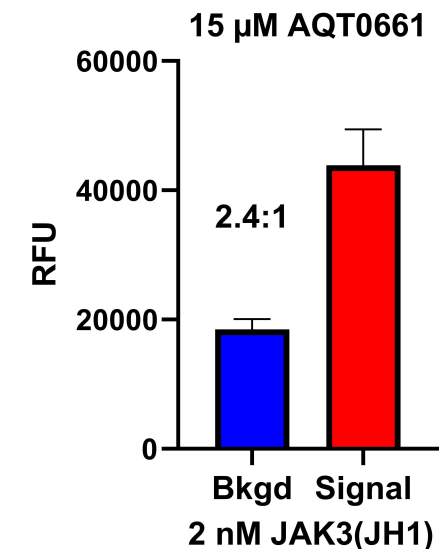
Red-Shift TRF S:B Determination

Progress and S:B Curves

PhosphoSens-Kinetic Progress Curves



PhosphoSens-Red S:B Chart



Using the reactions conditions previously described, the TRF S:B for this assay is 2.4