

## QS S Assist **KINASE**\_TR-FRET Kit

### **Description**

**KINASE** TR-FRET kit is designed for use in pharmacological assays for **KINASE** based on Time-Resolved Fluorescence Resonance Energy Transfer (TR-FRET). The kit includes assay buffer, human protein kinase, ATP/ULight™ labeled substrate peptide/Metal and a protocol to perform 384 well plate assay.

### **Components (800 dpt x 1 set)**

| Materials               | Volume    | Storage |
|-------------------------|-----------|---------|
| 10 x Assay Buffer       | 5 mL x 1  | -80°C   |
| <b>KINASE</b> *         | 60 µL x 1 | -80°C   |
| 5 x ATP/Substrate/Metal | 1 mL x 1  | -80°C   |

\* Concentration of the kinase; 200 µg/mL

Please avoid repeated freeze-thaw cycles.

### **Reagent Preparation (per 400 dpt)**

Bring all reagents (except kinases) to room temperature before use.

### **Materials provided**

#### **Assay Buffer**

Thaw 10 x Assay Buffer and take 2 mL. Dilute with 18 mL of distilled water. Adjusted Assay Buffer is able to keep room temperature before use. Please do not carry over this buffer on the next day, because the buffer component DTT is unstable.

#### **ATP/Substrate/Metal Solution**

Thaw 5 x ATP/Substrate/Metal component and dilute 0.45 mL into 1.8 mL of Assay Buffer (total volume: 2.25 mL). Bring the solution to room temperature until use.

#### **Enzyme Solution**

Thaw **KINASE** and dilute it appropriate-fold with Assay Buffer. Please keep it on ice before use.

## **Materials required**

### **Compound Solution**

Prepare a hundred times concentrated compound stock solution with DMSO. Dilute the solution 25 times with Assay Buffer. For the vehicle control, prepare 4% DMSO-Assay Buffer solution.

### **Detection Mixture**

Prepare the Detection Mixture containing appropriate concentrations of Tris-HCl (pH7.5), Tween20, EDTA, and Eu-labeled Antibody in distilled water. The Detection Mixture is able to keep room temperature before use with light shielding.

Minimal 24 mL of the Detection Mixture is needed for 400 dpt (60 µL/well).

The final concentrations of Tris-HCl (pH7.5), Tween20, EDTA, and Eu-labeled Antibody in the detection mixture is 15 mM, 0.01 %, 20 mM, and 0.53 nM, respectively in the protocol.

### **Example of Reaction Mixture**

| Sample | Compound Solution<br>(µL) | Vehicle<br>(µL) | ATP/Substrate/Metal<br>Solution<br>(µL) | Enzyme<br>Solution<br>(µL) | Assay Buffer<br>(µL) |
|--------|---------------------------|-----------------|-----------------------------------------|----------------------------|----------------------|
| A      | —                         | 5               | 5                                       | —                          | 10                   |
| B      | —                         | 5               | 5                                       | 10                         | —                    |
| C      | 5                         | —               | 5                                       | 10                         | —                    |

Where A equals negative control, B equals positive control and C equals test sample.

Calculate the percent inhibition of compound as follows;

$$\text{Inhibition (\%)} = (1 - (C - A) / (B - A)) \times 100$$

### **Final Concentrations of Components in Reaction Mixture**

15 mM Tris-HCl (pH7.5), 0.01 % Tween 20, 2 mM DTT

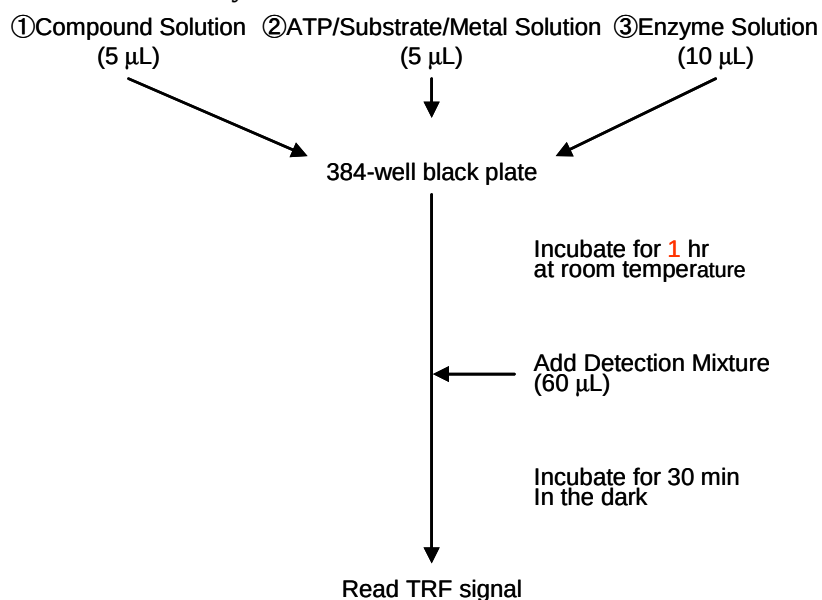
50 nM ULight™ labeled substrate peptide, 5 µM ATP, 5 mM Metal

### ASSAY PROCEDURE:

All procedures are performed at room temperature.

1. Add 5  $\mu$ L of Vehicle (4% DMSO) to wells of “A” and “B” and Compound Solution to wells of , “C” of a 384-well assay plate.
2. Add 5  $\mu$ L of ATP/Substrate/Metal Solution to each well.
3. Add 10  $\mu$ L of Assay Buffer to wells of “A” and Enzyme Solution to wells of “B” and “C” to start kinase reaction. Cover the plate and incubate for **1** hour at room temperature.
4. Add 60  $\mu$ L of Detection Mixture to each well. Incubate for 30 minutes at room temperature with light shielding.
5. Measure the TR-FRET signal with a plate reader (excitation 360 nm, emission 665 nm).

### Illustration of Assay Procedures:



### The settings for the instrument (2104EnVision D, PerkinElmer)

| Parameter          | Setting          |
|--------------------|------------------|
| Light source       | Laser            |
| Top mirror         | LANCE/DELFIABias |
| Excitation filter  | TRF 320nm        |
| Emission filter    | APC 665nm        |
| Emission filter 2  | Europium 615nm   |
| Measurement height | 7 mm             |
| Cycle              | 16600            |
| Delay              | 50 $\mu$ s       |
| Number of flashes  | 40               |
| Window time        | 400 $\mu$ s      |

### Assay result example

The inhibitory effect of **Reference compound** on **KINASE** evaluated with **KINASE** TR-FRET kit is shown below.

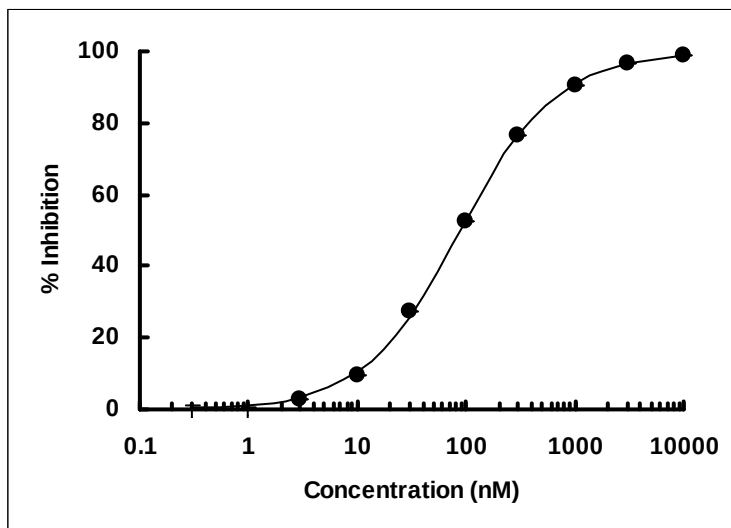


Figure.1. **Reference compound** inhibition curve.