

# Activity- and Affinity- based assay of FMS tyrosine kinase: the impact of kinase phosphorylation state on the activities of various kinase inhibitory compounds.

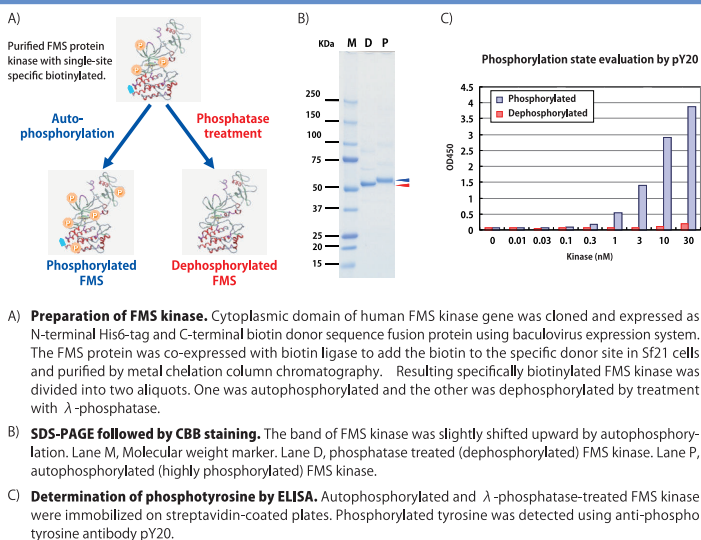
Daisuke Kitagawa, Masaki Gouda, Masanori Nakamura, Kensaku Akita, and Koichi Yokota

Carna Biosciences, Inc., BMA 3F 1-5-5 Minatojima-Minami, Chuo, Kobe 650-0047 Japan

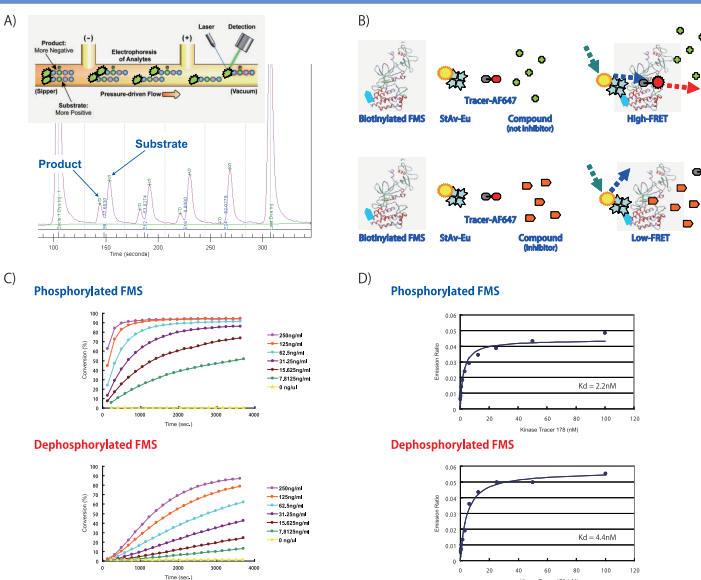
## 1. Introduction

It is well-known that the sensitivity of certain kinase inhibitors is influenced by the degree at which the kinase is phosphorylated (phosphorylation state). Although it is clear that such inhibitory compounds can discriminate between phosphorylated and nonphosphorylated protein, the biological consequences of this phosphorylation sensitivity are poorly defined. To further understand the relationship between selectivity, efficacy and safety of kinase inhibitors, we investigated effect of kinase protein phosphorylation state on various kinase inhibitors. We generated FMS (also known as CSF1R) tyrosine kinase protein to yield two differently phosphorylated batches; one was highly phosphorylated by its autophosphorylation and the other was dephosphorylated by phosphatase treatment. These two forms were then subjected to evaluation using nine kinase inhibitors in both activity-based and affinity-based assays.

## 2. Preparation of FMS protein

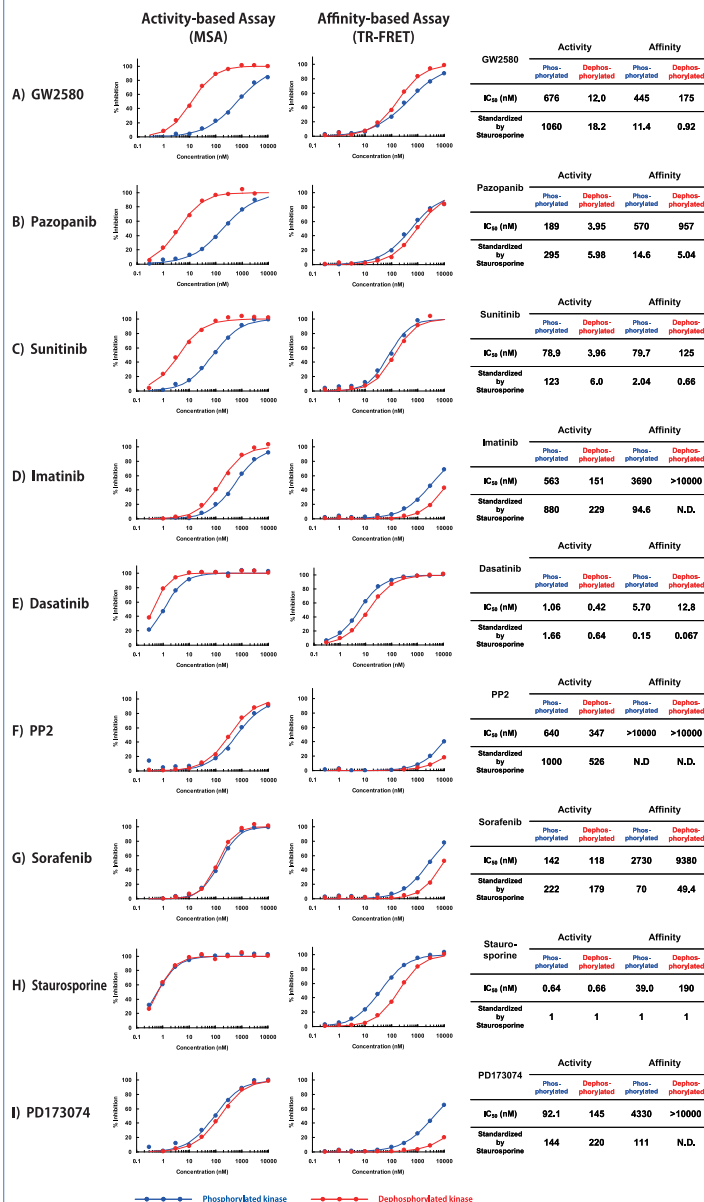


## 3. Assay platform



- A) Activity-based assay:** FMS kinase activity was determined by ability to phosphorylate a peptide substrate. The peak ratio of phosphosubstrate to substrate (P/(P+S)) was measured by off-chip mobility shift assay (MSA) using LabChip<sup>®</sup>3000 (Caliper Life Science).
- B) Affinity-based assay:** Europium-labeled streptavidin (StAv-Eu, FRET donor) is coupled to the FMS kinase through an interaction with its C-terminal biotin. Alexa-Fluor 647-conjugated compound (kinase tracer 178, FRET acceptor) is used as tracer. Binding of the tracer to FMS kinase results in increased TR-FRET signal between the fluorophores and the tracer.
- C) Kinase activity of autophosphorylated and dephosphorylated FMS.** The activity of the autophosphorylated FMS was about 5-times stronger than that of dephosphorylated FMS by continuous monitoring the production of phospho-peptides. The assay was performed at the ATP concentration of 1mM.
- D) TR-FRET signal of autophosphorylated and dephosphorylated FMS.** Both types of FMS kinase showed the similar TR-FRET signal.

## 4. Results



A) ~ I) Potency of inhibitors toward two types of FMS tyrosine kinase was assessed by MSA and TR-FRET competition assay. IC<sub>50</sub> values were calculated from concentration vs. %inhibition curves by fitting to a four-parameter logistic curve. Each point is an average value of at least two independent experiments.

Some compounds (GW2580, pazopanib, and sunitinib) showed different potency between phosphorylated and dephosphorylated FMS in MSA. TR-FRET assay was insensitive to the difference of phosphorylation state of FMS.

## 5. Conclusions

We have successfully generated single-site specific biotinylated FMS kinase using the baculovirus expression system, that can be easily adapted for use in affinity-based assay systems.

Dephosphorylation of FMS resulted in an inactive form that possessed approximately one fifth activity of phosphorylated FMS.

Surprisingly, GW2580, pazopanib, and sunitinib showed strong inhibitory activity toward the dephosphorylated FMS compared to phosphorylated FMS, however, such a difference was not observed in TR-FRET.

Further study on the effects of phosphorylation state on potency using SPR methods will be in due course.