

# Preincubation Kinase Profiling Service

Available for  
**165**  
Kinases\*



## Preincubation Assay is now available from Carna!

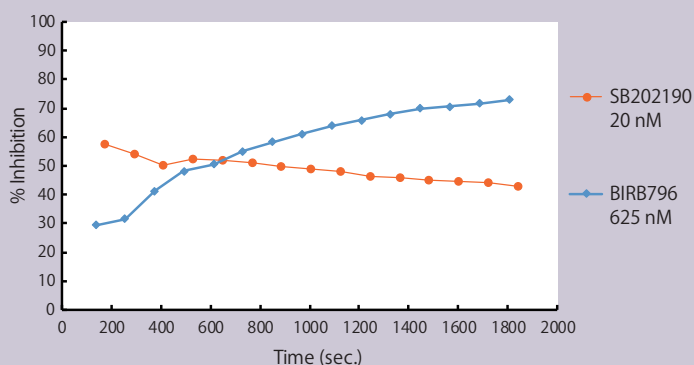
Certain classes of kinase inhibitors exhibit slow binding kinetics as shown in Figure 1.

For such compounds, preincubation (Figure 2) is necessary to assess their maximal effect (Figure 3 and 4).

We are pleased to announce a new kinase profiling service that incorporates preincubation with the test compound(s). A thirty (30) minute preincubation at room temperature is typically sufficient to capture such slow binding compounds, after which their activity is evaluated using our standard Mobility Shift Assay.

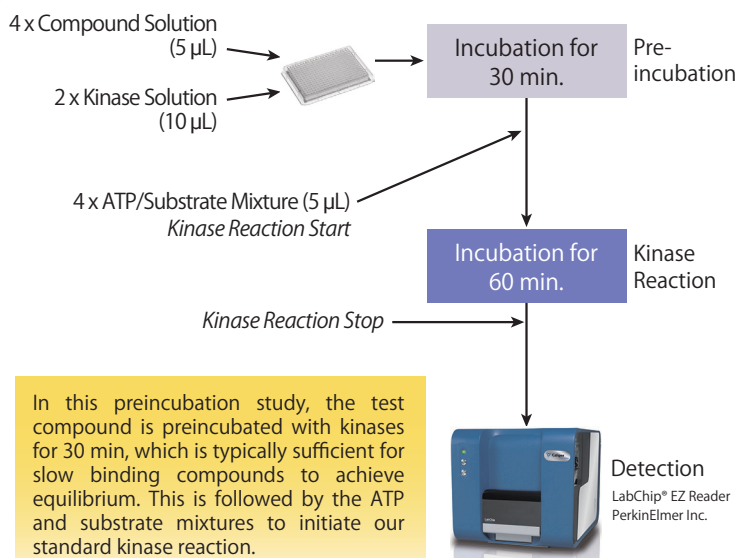
This service is specifically designed to measure compounds with slow association kinetics, and the corresponding studies are conducted once per month. Please inquire for more detailed information.

\* As of June 30, 2014



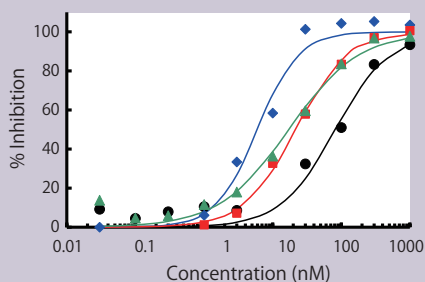
**Fig. 1 Time course of p38 $\alpha$  inhibitors**

Structurally unrelated SB202190 and BIRB796 are potent inhibitors of p38 $\alpha$ . BIRB796 interacts with p38 $\alpha$  in a manner different from SB202190, and its binding induces a slow conformational change that locks the protein into an inactive conformation. Therefore, the potency of BIRB796 increased with the period of incubation. The assay was performed at an ATP concentration of 1 mM.



In this preincubation study, the test compound is preincubated with kinases for 30 min, which is typically sufficient for slow binding compounds to achieve equilibrium. This is followed by the ATP and substrate mixtures to initiate our standard kinase reaction.

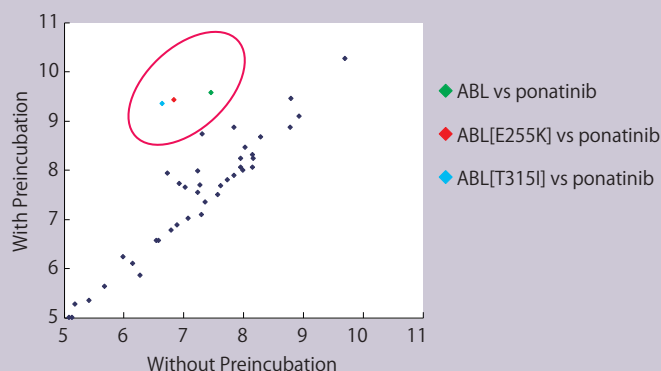
**Fig. 2 Preincubation Study Methodology**



|          | Preincubation |        |
|----------|---------------|--------|
|          | +             | -      |
| BIRB796  | 9.5 (◆)       | 82 (●) |
| SB202190 | 19 (▲)        | 22 (■) |

**Fig. 3 IC<sub>50</sub> Comparison with or without Preincubation for p38 $\alpha$  Inhibitors**

Preincubation did not affect the IC<sub>50</sub> of SB202190, however, a significant change was observed for the activity of BIRB796. The assay was performed at an ATP concentration of 1 mM.



**Fig. 4 IC<sub>50</sub> Comparison with or without Preincubation for Clinical Inhibitors**

Several clinical kinase inhibitors and their corresponding target kinases were evaluated in a preincubation study; the resulting pIC<sub>50</sub> values are compared above. In most cases, good correlations were observed. However, Ponatinib showed enhanced inhibition of native and mutant isoforms of the ABL kinase following preincubation. All assays were performed at an ATP concentration of 1 mM.

## Carna Biosciences, Inc.

BMA 3F, 1-5-5 Minatojima-Minamimachi, Chuo-ku, Kobe 650-0047 JAPAN

TEL: 078-302-7091 / FAX: 078-302-7086

E-mail: info@carnabio.com

