



The importance of assessing intrinsic target kinase activity

Carna Biosciences' long standing commitment to manufacture highly active protein kinases is based on our belief that drugs designed to inhibit disease related protein dysregulation should be evaluated in activity- based enzyme assays using the most active kinases. Nearly all kinases produced at Carna are expressed in insect cells, purified, and amino acid sequence-confirmed using peptide mass fingerprinting, however ...

One may ask whether it is the expressed Kinase (X) that catalyzes the phosphorylation reaction giving rise to the product, phosphorylated peptide substrate or ADP, in the presence of ATP. Alternatively, is it possibly another kinase derived from the insect cell host system, Kinase (Z)? Also, could ATPase activity be generating ADP in the assay?

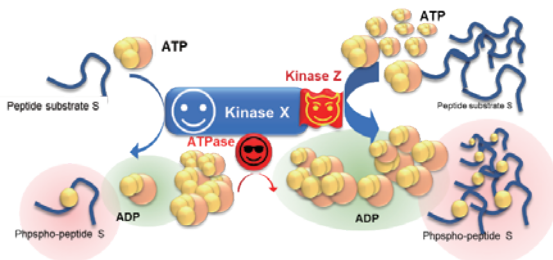


Fig1. If a small amount of Kinase Z exists in the preparation of Kinase X, is it possible to confirm that the measured substrate phosphorylation is the result of Kinase X? Additionally, can we establish that the ADP produced is from the target of interest as opposed to an ATPase activity?

Let's assume that the purity of Kinase X is 99% and that the remaining 1% is contributed from Kinase Z. What would happen if Kinase Z has a much higher affinity for ATP and peptide substrate S compared to Kinase X? In that case, the majority of the phosphorylation product may be generated from Kinase Z activity, as opposed to the target kinase, despite the fact that it only represents 1% of the kinase preparation. At the

same time, ATPase contamination could affect the detection of kinase activity of interest because of excess ADP production. Is there a way to determine whether the obtained phosphorylation activity is truly from the target kinase?

[Risks lie in production using the insect expression system and purification using a His tag]

When an E. coli expression system is utilized to produce a kinase, the result is often a product with no or minimal activity, making the insect cell expression system the more favorable alternative. Carna became aware of the fact that insect cells have their own unique, highly active kinases, which also possess the rather undesirable characteristic of low substrate selectivity. What's worse, Carna found that when a His tag is used for kinase purification, the insect cell-derived kinases tend to remain in the preparation, and become inseparable from the kinase of interest after completing the purification. In addition, there are cases where a tyrosine kinase has a strong serine-threonine kinase activity even if it has a purity of 95% or higher (Fig. 2).

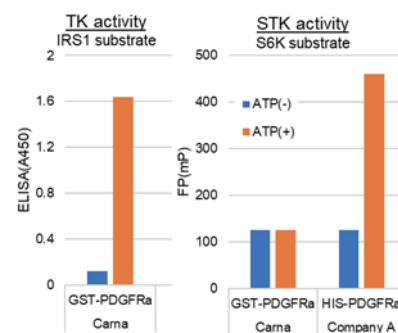


Fig 2. Comparison of 10 ng GST-PDGFRα (Carna) vs. 10ng HIS-PDGFRα (Company A). HIS-PDGFRα had the ability to phosphorylate Ser/Thr on S6K substrate, presumably from an insect cell-derived kinase activity. GST-PDGFRα showed Tyr phosphorylation of IRS1 peptide substrate, however no phosphorylation of the S6K



CARNA COFFEE BREAK - TECHNICAL NOTE - No.2

substrate was detected. Similar results were obtained using PKC ϵ pseudo-substrate as the substrate for PDGFR α .

Based on these observations, Carna has elected to use a GST tag as its first choice when producing kinases.

Insect cell-derived kinases have also demonstrated phosphorylation activity, particularly targeting the substrates shown in Table 1.

phosphorylation site	substrate
Tyr	Lyn-peptide
	Gastrin-peptide
Ser / Thr	PKC ϵ -pseudopeptide
	S6K-peptide
	SRPKtide
	MBP
	Casein

Table 1. Substrates susceptible to phosphorylation by insect cell-derived kinase

[The importance of using a specific substrate to the kinase of interest]

Additionally, phosphorylation by insect-derived kinases, when detected, gives rise to erroneous data for the target kinase, particularly when using non-specific substrates or when using certain antibodies to a substrate.

Let's illustrate one example here. Even when a protein kinase was prepared as a GST-kinase to eliminate contamination by insect cell-derived kinase(s), contaminant kinase activity was detected when using anti-phosphorylated threonine-histone H3 antibody with histone H3 as a substrate (Fig. 3).

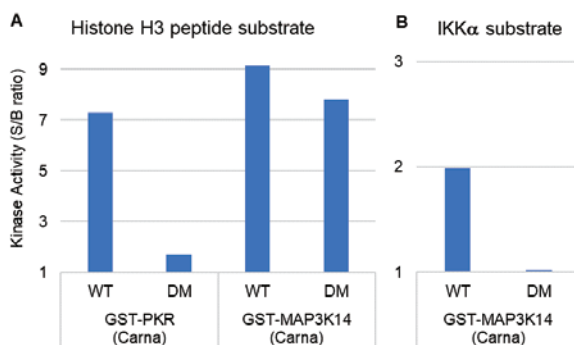


Fig 3. A) Kinase activity of PKR wild type and K296R (DM; dead mutant) on the left, and MAP3K14 wild type and K429A/K430A (DM) produced using GST tags (each kinase 10 nM) on the right. Detection was performed using Histone H3 peptide as a substrate and anti-phosphorylated H3 (Thr3) antibody.

B) Kinase activity of MAP3K14 wild type and K429A/K430A (DM) produced using GST tags (each kinase 12.5 nM) measured by use of IKK α as a substrate. The kinase activity of phosphorylated IKK α was detected by the phosphorylation of IKK α , a specific substrate of IKK α , in a cascade assay on the Mobility Shift Assay platform. GST-PKR was found to have phosphorylation activity to H3 peptide only in its WT form. Contrarily, MAP3K14 was found to have H3 phosphorylation activity even in its DM form, and therefore it was concluded that this activity was not from MAP3K14 itself. Currently, at Carna, IKK α , which is a specific substrate and not phosphorylated using DM, is utilized as a preferred substrate.

In summary, even if a kinase protein of interest is highly purified, the activity that one detects in experiments may be derived from a kinase other than that of the target kinase. Carna has been reflecting upon each lesson learned from a long production history with our products, and we strive to analyze activity via different approaches to make sure we validate that the activity is coming from the kinase of interest. This way, we can not only ensure that quality of our products, but we also protect the quality of our assay services that utilize these kinases.

Kinase List

Biotinykated Kinase List