

Comprehensive Protein Kinase Profiling Panel for Selective Inhibitor Screening

provided by Carna Biosciences "The Kinase company"

Masaki Gouda*, Hiroshi Ohmoto, Etsuko Ishibushi, Tetsuya Myojin, Kensaku Akita, Yoshimasa Inoue, Ryoko Nakai, Hiroshi Ishiguro and Koichi Yokota, Carna Biosciences, Inc., Japan.

Abstract

Kinase selectivity profiling has become an invaluable tool to assess the potential of kinase inhibitors in the drug discovery process. To support drug discovery programs at biotechnology and pharmaceutical companies, a Japanese biotechnology company, Carna provides 255 active protein kinases and profiling services against over 200 kinases under its optimized assay conditions. In this presentation, we describe a new kinase profiling panel using an off-chip mobility shift assays leveraged on the Caliper LifeSciences' LabChip® 3000 (LC3000) systems. Peptide substrates were identified for approximately 200 kinases. The peak-separation conditions of the fluorescence-labeled substrates and their phosphorylated products were optimized. Carna's active kinases phosphorylated their peptide substrates on dose- and time-dependent manners. The Km value for ATP was determined for each kinase. Consequently more than 150 kinase assays are available in the profiling panel on the LC3000. These kinases belong to TK, TKL, STE, CK1, AGC, CMGC, and CAMK families in the kinome. Therefore, our kinase profiling panel is useful for optimizing the lead compounds.

Procedure Off-Chip mobility shift assay for profiling

Characterization of recombinant kinase (Fig. 1).

Optimization of peak separation (Fig. 2).

Km determination (Fig. 3 and Fig. 4).

Relationship between % conversion and IC50 (Fig. 5).

IC50 determination (Fig. 6 and Specificity profiles for clinical kinase inhibitors).

Fig. 1 Characterization of recombinant kinase

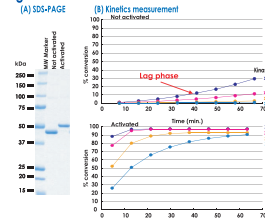
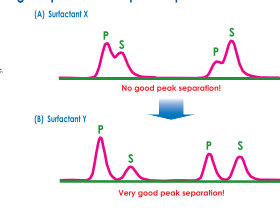


Fig. 2 Optimization of peak separation



Full-length GST-tagged kinase protein was produced with baculovirus-infected insect cells. Kinases were purified with several steps chromatography (without any contaminated kinase derived from host cell). Characterization of recombinant protein kinase by SDS-PAGE (A) and kinetic measurement of non-activated kinase (B, top) and activated kinase after activation by auto-phosphorylation or upstream kinases (B, bottom).

Peak separation is most importance for analysis on the mobility shift assay. Figures show peak separation condition used (A) surfactant X (B) surfactant Y in running buffer. Peaks show ratio of substrate to phosphorylated product from same wells. We examined that peak separation was deemed acceptable for use, conditions for validation were optimized.

Fig. 3 Km determination

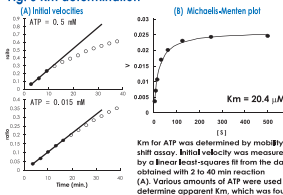


Fig. 4 Distribution of Km values for ATP

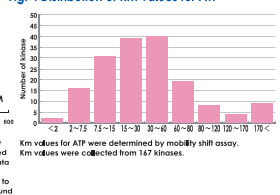


Fig. 5 Relationship between % conversion and IC50

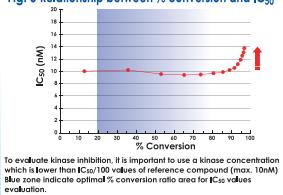
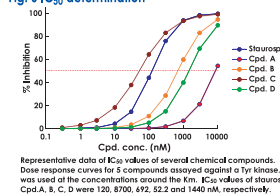


Fig. 6 IC50 determination



Protocol

ATP/Substrate solution (5 μL) + Compound solution (5 μL) + Kinase solution (10 μL) → 384 well plate

Incubate at room temperature 1 or 5 hr

Add stop buffer (40 μL)

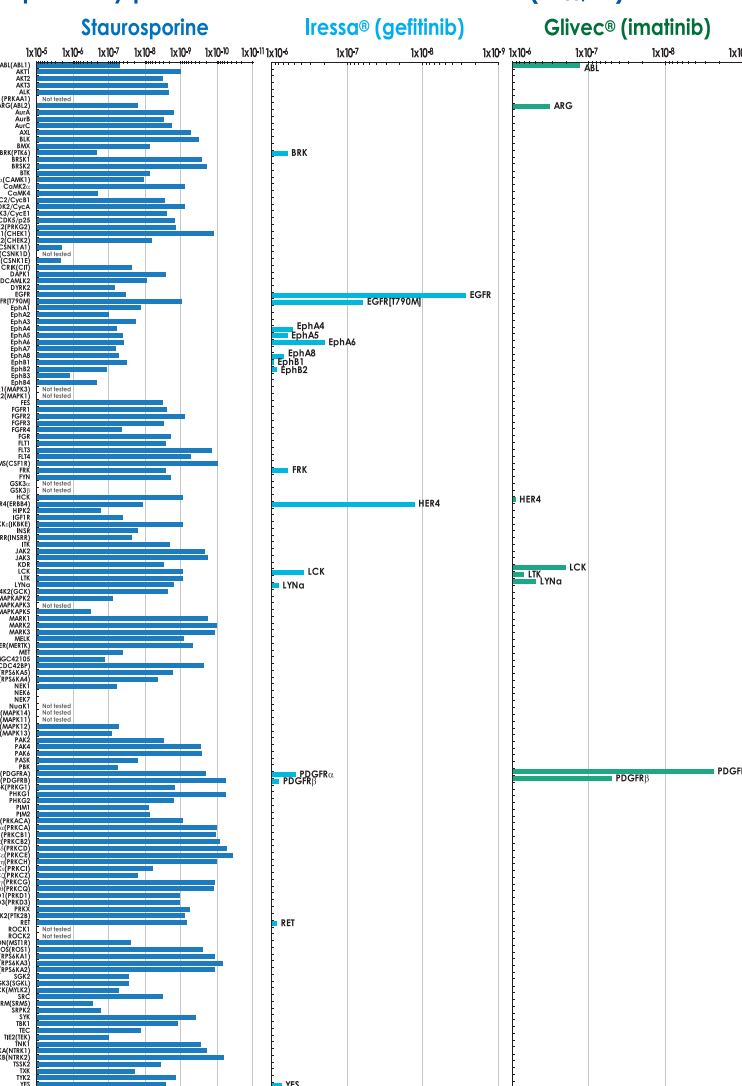
Read Signal

-Final assay condition-
Hepes buffer (pH 7.5)
5 mM MgCl2
2 mM DTT
0.01 % Surfactant (lipid activator, Ca/CaM, cGMP etc.)
ATP (nearest to ATP Km app.)
Kinase protein (max. 10nM)
Compound (Staurosporine, Iressa®, Glivec®)

M: Marker Peaks
S: Substrate Peaks
P: Product Peaks

Example of peak separation result showing varying amounts of substrate conversion to product.
Clear separation and detection of "Product" and "Substrate" peaks allow accurate quantification of inhibition.

Specificity profiles for clinical kinase inhibitors (IC50, M)



Conclusion

Carna Biosciences has developed assays for kinase profiling panel using mobility shift assay, LabChip® microfluidic platform, which is applicable to various kinases.

We have obtained the high reproducibility of kinase selectivity profiling by Iressa® and Glivec® which can prove the accuracy of new assay platform.

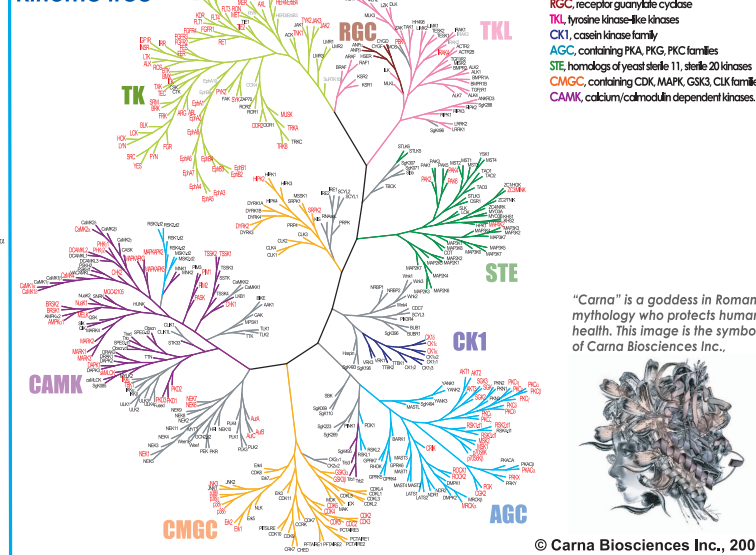
Profiling panel with 167 kinases (71 TKs and 96 STKs) has been developed by using Mobility Shift Assay!

Kinases represented in this presentation are shown in kinome tree, as well as in Kinase List.

Kinase List

| Tyr kinase | | | | | | | | | | Ser/Thr kinase | | | | | | | | | |
|-------------|------|------------|----------|--------|-------|-----------|-------|---------|------------|----------------|----------|------------|------------|------------|-------------|----------|-----------|--|--|
| Cytoplasmic | | | Receptor | | | AGC group | | | CAMK group | | | CMGC group | | | Other group | | CK1 group | | |
| ABL | BLK | BTK | ALK | EPHA3 | FGFR3 | ROS | AKT1 | PKCβ2 | CRIK | CaMK1α | MG42105 | PIM2 | CDC2/CycB1 | Erk1 | AurA | CK1α | | | |
| ABL3133 | IRK | BRK | ITK | LTK | EPHA4 | FGFR4 | TIE2 | AKT2 | PKCγ | RSK1 | CaMK1b | Nvak | PKD1 | CDC2/CycA | Erk2 | CK1β | | | |
| ARG | FGR | TEC | AXL | EPHA5 | IGF1R | TRKA | AKT3 | PKCδ | RSK2 | CaMK2α | PASK | PKD2 | CDC3/CycE1 | p38α | AurC | CK1ε | | | |
| TNK1 | FRK | TKX | MER | EPHA6 | INSR | TRKB | MRCKα | PKCε | RSK3 | CaMK4 | DAPK1 | CHK2 | CDC4/CycD3 | p38β | IKKα | | | | |
| PYK2 | FYN | TYRO3 | EPHA7 | IEI | FLT1 | FLT3 | MSK1 | PKCζ | SGK | AMPKα1 | DCAMK2 | CHK1 | CDK5/p25 | p38γ | TBK1 | MAPKAPK2 | | | |
| FES | HCK | DDR2 | EPHA8 | AET | FLT4 | FLT4 | MSK2 | PKCθ | SGK2 | BRSK1 | MAPKAPK2 | TSSK1 | CDK6/CycD3 | p38δ | NEK1 | MAPKAPK3 | | | |
| FER | LCK | EGFR | EPHA11 | RON | KDR | | CGK2 | PKCη | SGK3 | BRSK2 | MAPKAPK3 | TSSK2 | DYRK2 | CDK6/CycD3 | NEK6 | PAK2 | | | |
| JAK2 | LYNα | EGFRIT/TPM | EPHB2 | FLT3 | | | PKG | PKCι | | CHK1 | MAPKAPK5 | | HIPK2 | | NEK7 | PAK4 | | | |
| JAK3 | LYNβ | | MUSK | EPHB3 | FMS | | PKAα | p70S6K | | MARK1 | sMLCK | | SRPK2 | | PBK | PAK4 | | | |
| TYK2 | SRC | HER4 | EPHA4 | PDGFRα | | | PRKX | p70S6Kβ | | MARK2 | PHKG1 | | GSK3α | | BRK4 | PAK6 | | | |
| SRM | SYK | EPHA1 | FGFR1 | PDGFRβ | | | PKCα | ROCK1 | | MARK3 | PHKG2 | | GSK3β | | | | | | |
| YES | BMX | EPHA2 | FGFR2 | RET | | | PKCβ1 | ROCK2 | | MELK | PIM1 | | JNK3 | | | | | | |

Kinome tree



"Carna" is a goddess in Roman mythology who protects human health. This image is the symbol of Carna Biosciences Inc.,

151 kinases (42 Tyr kinases, 1 drug-resistant EGFR(T790M) and 88 Ser/Thr kinases) for specificity profiling were picked up from kinase list. Profiling was conducted at 10 concentrations (10, 3, 1, 0.3, 0.1, 0.03, 0.01, 0.003, 0.001 and 0.0003 μM) in duplicate for each compound.

Please contact us; info@carnabio.com Visit our web site; www.carnabio.com