

# Cell-Based Tyrosine Kinase Assay Panel



CARNA BIOSCIENCES collaboration with



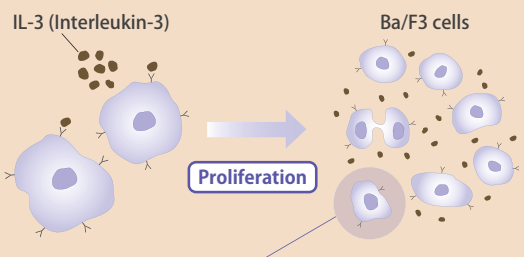
## Largest Commercially Available Panel of Tyrosine Kinase Cell-Based Assays

### Why ACD?



- Comparative cell-based analysis
- To discover direct inhibitory activity to targeted kinases
- Ready-to-run **80** Tyrosine Kinase (TK) Panel
- Time & cost saving solution for your in-house cellular assays

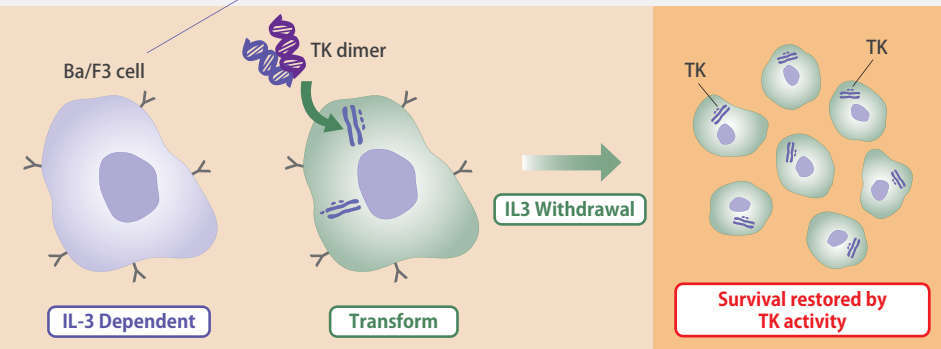
### Principle & Method of ACD Cell-Based Assays



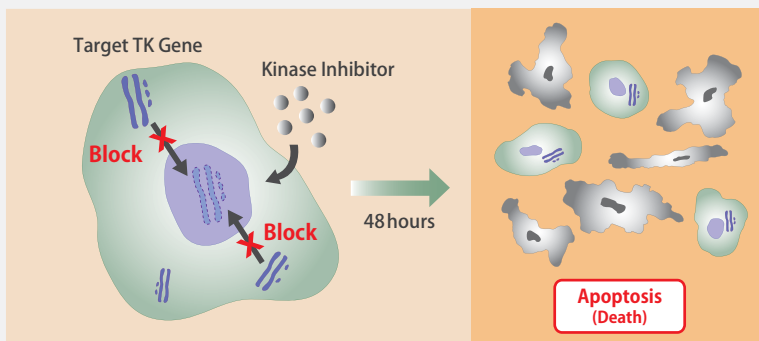
The assay principle builds upon the work of Daley & Baltimore (1988)\*.

In this system, IL3-dependent Ba/F3 cells are modified to express an activated recombinant kinase. Following removal of IL3, the modified cells are dependent on the activity of the recombinant kinase for survival and proliferation.

\* Daley and Baltimore; Proc. Natl. Acad. Sci. USA. 1988; 85(23):9312-6



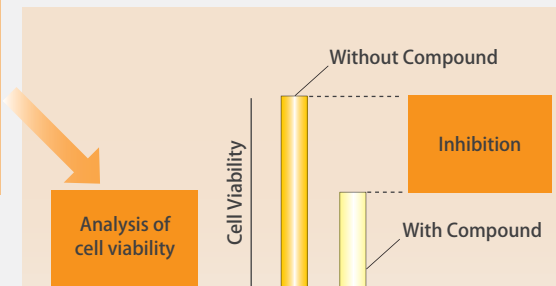
Ba/F3 cells are transformed by inducing target kinase dimerization via viral vectors. Activity of the transformed kinase overrides IL3 dependency for cellular proliferation and survival - modified cells no longer require IL3 for growth.



If the kinase inhibitor (compound) specifically blocks the activity of the recombinant kinase, the modified cells undergo programmed cell death (apoptosis).

### About ACD

Advanced Cellular Dynamics (San Diego, CA USA) is a leading provider of cell-based assay panel technologies and services to the life-sciences community. ACD develops and deploys families of cell-based screening assays, encompassing broad representations of important target gene families. Their assays are designed to simplify high-throughput screening and profiling of chemical entities in a physiologically relevant cellular environment.



Each assay is engineered to be dependent upon maintenance of the introduced kinase activity for survival. Inhibition of this activity results in a directly proportional decrease in cell viability.

Visit our website for more information:  
[www.carnabio.com](http://www.carnabio.com)

### Carna Biosciences, Inc.

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

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

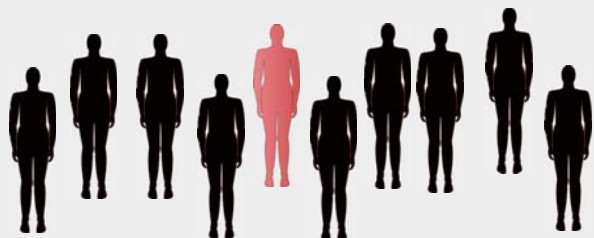
E-mail: [info@carnabio.com](mailto:info@carnabio.com)

# ACD's Cell-Based Tyrosine Kinase Assay Panel

Don't miss important biology using traditional assays.

## 1 EGFR and Lung Cancer

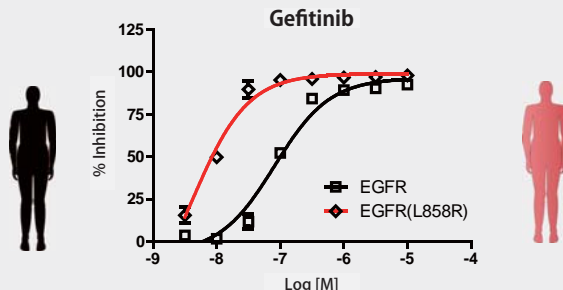
Gefitinib (Iressa™) was the first EGFR tyrosine kinase inhibitor for the treatment of Non-Small Cell Lung Cancer (NSCLC).



➤ Only 10% of the treated population responded.

## 2 Mutant EGF Receptors

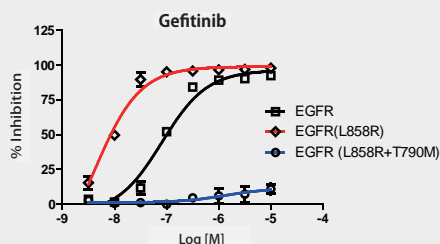
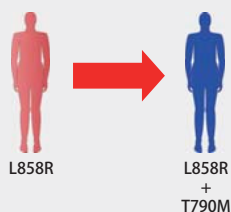
Responding NSCLC patients possess a mutant EGFR.



- Two "classical" mutations **L858R** and **Δ746-750**
- Mutant receptors are much more responsive to Gefitinib.

## 3 Gefitinib Resistance

Responsive patients become resistant over time.



- Resistance due to secondary "gatekeeper" mutation (T790M).
- Double mutant receptors (L858R + T790M) are much less responsive to Gefitinib.

## 4 Why Use Cell-based Assays?

Kinase biology can be complex.

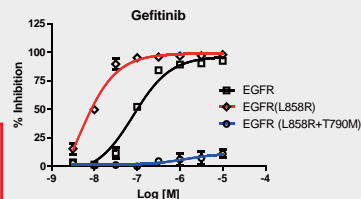
Biochemical Assay\*

ACD Cell-Based Assay

Kinase	K <sub>i</sub> , nM		K <sub>i</sub> /K <sub>i(ATP)</sub> , x 10 <sup>-3</sup>	
	Gefitinib	AEE788	Gefitinib	AEE788
WT	35.3 ± 0.4	5.3 ± 0.3	6.8	1
T790M	4.6 ± 0.1	27.6 ± 0.7	0.78	4.7
L858R	2.4 ± 0.1	1.1 ± 0.1	0.016	0.0074
L858R/T790M	10.9 ± 0.6	18.6 ± 0.5	1.3	2.2

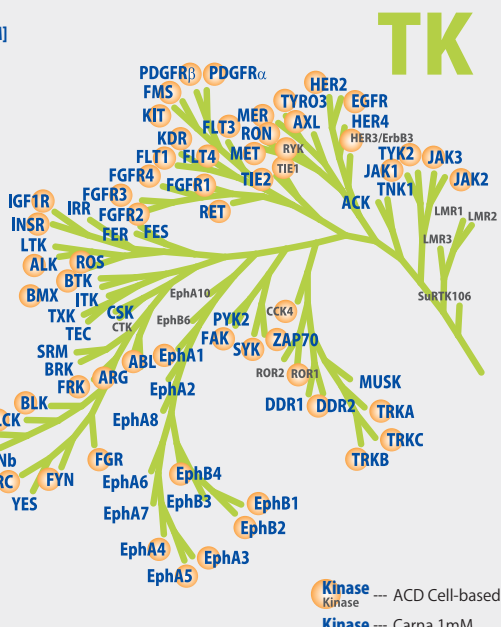
The ratio K<sub>i</sub>/K<sub>i(ATP)</sub> provides a relative estimate of inhibitor potency.

Notice that L858R/T790M affinity for Gefitinib is reduced < 5-fold relative to L858R. In tumors, the response differs by > 100-fold!



- Tumors bearing L858R/T790M respond poorly to Gefitinib. These differences can be missed when evaluation is performed using traditional biochemical assays, but are captured using ACD cell-based assays.

- ALK [L1196M]
- EGFR [Δ746-750]
- EGFR [Δ746-750+T790M]
- EGFR [L858R]
- EGFR [L858R+T790M]
- EGFR [L861Q]
- EGFR [L861Q+T790M]
- EGFR [T790M]
- FGFR1 [V561M]
- FGFR2 [K660E]
- FGFR2 [K660N]
- FGFR2 [N550K]
- FGFR2 [V565I]
- FGFR3 [K650M]
- FGFR4 [V550E]
- KIT [D816V]
- KIT [K642E]
- KIT [N822H]
- KIT [T670I]
- KIT [V654A]



ACD Cell-Based TK Assays Available for Screening Services

80 Total Kinases - Broad Coverage of the Tyrosine Kinome!

- |                       |               |             |             |
|-----------------------|---------------|-------------|-------------|
| ABL(BCR-ABL)          | EphB1         | FMS(CSF1R)  | MER(MERTK)  |
| ALK                   | EphB2         | FRK         | MET         |
| ALK [L1196M]          | EphB4         | FYN         | PDGFRa      |
| ARG(ABL2)             | FAK           | HCK         | PDGFRb      |
| AXL                   | FGFR1         | HER2(ERBB2) | RET         |
| BLK                   | FGFR1 [V561M] | HER3(ERBB3) | RON(MST1R)  |
| BMX                   | FGFR2         | IGF1R       | ROR1        |
| BTK                   | FGFR2 [K660E] | INSR        | ROS(ROS1)   |
| CCK4(PTK7)            | FGFR2 [K660N] | JAK1        | RYK         |
| DDR2                  | FGFR2 [N550K] | JAK2        | SRC         |
| EGFR                  | FGFR2 [V565I] | JAK3        | SYK         |
| EGFR [Δ746-750]       | FGFR3         | KDR         | TIE1        |
| EGFR [Δ746-750+T790M] | FGFR3 [K650M] | KIT         | TIE2        |
| EGFR [L858R]          | FGFR4         | KIT [D816V] | TRKA(NTRK1) |
| EGFR [L858R+T790M]    | FGFR4 [V550E] | KIT [K642E] | TRKB(NTRK2) |
| EGFR [L861Q]          | FGR           | KIT [N822H] | TRKC(NTRK3) |
| EGFR [L861Q+T790M]    | FLT1          | KIT [T670I] | TYK2        |
| EGFR [T790M]          | FLT3          | KIT [V654A] | TYRO3       |
| EphA1                 | FLT4          | LCK         | ZAP70       |
| EphA3                 |               | LYN         |             |
| EphA4                 |               |             |             |
| EphA5                 |               |             |             |

Updated: 2014/1/15