

QuickScout™ Selectivity Profiling Service Application Form

Date: Enter mm/dd/yy

Customer Information

Customer Name	Dr. John Carna
Institution/Company	
Department	
Street Address	
City	
State/Zip	
Country	
Tel Number	
Fax Number	
Email Address	

Enter your information

Study Information

Assay Type	<input type="radio"/> % Inhibition at fixed concentration <input checked="" type="radio"/> IC50 determination	Choose the type of study
Physical State	<input checked="" type="radio"/> Solution <input type="radio"/> Solid	Choose the physical state of compound. Solution must be in DMSO
Unit of Concentration	<input checked="" type="radio"/> μ M <input type="radio"/> μ g/mL	Choose the unit of Concentration
Residual materials	<input type="radio"/> Dispose after completion of study. <input checked="" type="radio"/> Return to the customer (at customer's cost)	Choose how you would like us to handle the residual materials

Additional Information

For special request, please write here.

Notice

Detailed information about each kinase and assay condition can be checked on the latest Kinase Profiling Book.

http://www.carnabio.com/output/pdf/ProfilingProfilingBook_en.pdf

Sample Preparation

The volume of samples required for testing varies with the number and types of assays selected.

Please supply samples in 100% DMSO at 100X the highest test concentration.

Volume required : ≤ 100 target kinases = 500 μ L , > 100 target kinases = 1000 μ L

If your samples do not meet our requirements (i.e., 10 mM DMSO solution, solid form), please consult us. Also, please provide us the Safety Data Sheet (SDS, or cautions for sample handling), disposal instruction, and any special solubilization instructions if available.

Samples submitted as solids must be shipped pre-weighed with the accurate molecular weight, quantity shipped and purity recorded on this Application Form. Please provide sample(s) in a vessel deep enough to accommodate 5 mM* DMSO solution. An additional charge may apply if a large number of samples are supplied as powders. We do not weigh customer samples. Please consult us prior to shipment.

(*in case of >50 μ M test concentration, please contact us in advance)

Ship to: CARNA BIOSCIENCES, INC.□
 Attention: Yusuke Kawase
 BMA 3F 1-5-5 Minatojima-Minamimachi, Chuo-ku, Kobe 650-0047 Japan
 TEL: +81 78-302-7091 / FAX: +81 78-302-7086
 E-mail: info@carnabio.com

Storage

Unless otherwise directed by the client, samples are solubilized and diluted with dimethylsulfoxide (DMSO) to achieve 100-fold higher concentration than the final test concentrations. This Solution is stored at -10 to -30 Celsius. This solution is further diluted into 4% DMSO to achieve a 4-fold higher concentration than the final test concentrations. The addition of reagents used in the assay further dilutes the sample solution by 4-fold, so that the final DMSO concentration in the assay is 1%. If you have determined that your samples are insoluble in 100% DMSO, 4% DMSO, or require special handling, please consult us in prior to shipping.

Disposal

Unless otherwise indicated on the Application Form by the client, residuals will be disposed after three (3) months from the completion of study. Residuals can be returned to the client upon an request at the client's expense.

Please feel free to contact us if you have any questions or comments.

Compound and Assay Information

ATTENTION: Please do not DRAG & DROP (or PASTE) when completing cells. To change a cell entry, please delete the information then re-enter.

	Name	Lot #	Prep. Date	Conc. (mM)	Volume (μL)	Purity (%)	Storage Temp.	Start Concentration (μM)						
								Conc1						
1														
2														
3														
4														
5														
6														
7														
8														
9														
10														
11														
12														
13														
14														
15														
16														
17														
18														
19														
20														
21														
22														
23														
24														
25														
26														
27														
28														
29														
30														
31														
32														
33														
34														
35														
36														
37														
38														
39														
40														
41														
42														
43														
44														
45														
46														
47														
48														
49														
50														

Enter information on one (1) compound per line.
The colored cells must be filled in.

Enter if multiple concentrations.
For IC50 study, the top concentration must be
entered in "Conc 1" column.

For more than fifty (50) compounds,
please copy this sheet .

If you are going to choose a large number of kinases, please use the Kinase List function on the right!

Target Kinase Selection

ATP Conc.: K_m = around K_m , 1mM = 1mM ATP

Checked or red colored kinase is selected one.

QuickScout Panel

The total number of assys you have chosen

52 assays selected

☐ Hide All Unselected Kinases.

- TK Panel
- STK Panel
- Cell Cycle Panel
- MAPK Cascade Panel

- Km assay all ☐
- 1mM assay all ☐
- Cascade assay all ☐

[Click here for the pre-selected series.](#)

Check the box of kinase you would like to request.

Choose the ATP concentration of your request

~~Tyrosine Kinases~~

Km			1mM			Km			1mM			Km			1mM		
ABL	<input checked="" type="checkbox"/>	<input type="checkbox"/>	ABL [E255K]	<input checked="" type="checkbox"/>	<input type="checkbox"/>	ABL [T315I]	<input checked="" type="checkbox"/>	<input type="checkbox"/>	ACK	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
ALK	<input type="checkbox"/>	<input type="checkbox"/>	ALK [C1156Y]	<input type="checkbox"/>	<input type="checkbox"/>	ALK [F1174L]	<input type="checkbox"/>	<input type="checkbox"/>	ALK [G1202R]	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
ALK [G1269A]	<input type="checkbox"/>	<input type="checkbox"/>	ALK [L1152insT] (ALK [T1151_L1152insT])	<input type="checkbox"/>	<input type="checkbox"/>	ALK [L1196M]	<input type="checkbox"/>	<input type="checkbox"/>	ALK [R1275Q]	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
EML4-ALK	<input type="checkbox"/>	<input type="checkbox"/>	NPM1-ALK	<input type="checkbox"/>	<input type="checkbox"/>	ARG	<input type="checkbox"/>	<input type="checkbox"/>	AXL	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
BLK	<input type="checkbox"/>	<input type="checkbox"/>	BMX	<input type="checkbox"/>	<input type="checkbox"/>	BRK	<input type="checkbox"/>	<input type="checkbox"/>	BTK	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
BTK [C481S]	<input type="checkbox"/>	<input type="checkbox"/>	CSK	<input checked="" type="checkbox"/>	<input type="checkbox"/>	DDR1	<input type="checkbox"/>	<input type="checkbox"/>	DDR2	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
EGFR	<input checked="" type="checkbox"/>	<input type="checkbox"/>	EGFR [d746-750]	<input type="checkbox"/>	<input type="checkbox"/>	EGFR [d746-750/T790M]	<input type="checkbox"/>	<input type="checkbox"/>	EGFR [L858R]	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
EGFR [L861Q]	<input type="checkbox"/>	<input type="checkbox"/>	EGFR [T790M]	<input type="checkbox"/>	<input type="checkbox"/>	EGFR [T790M/L858R]	<input type="checkbox"/>	<input type="checkbox"/>	EPHA1	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
EPHA2	<input checked="" type="checkbox"/>	<input type="checkbox"/>	EPHA3	<input type="checkbox"/>	<input type="checkbox"/>	EPHA4	<input type="checkbox"/>	<input type="checkbox"/>	EPHA5	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
EPHA6	<input type="checkbox"/>	<input type="checkbox"/>	EPHA7	<input type="checkbox"/>	<input type="checkbox"/>	EPHA8	<input type="checkbox"/>	<input type="checkbox"/>	EPHB1	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
EPHB2	<input type="checkbox"/>	<input type="checkbox"/>	EPHB3	<input type="checkbox"/>	<input type="checkbox"/>	EPHB4	<input checked="" type="checkbox"/>	<input type="checkbox"/>	FAK	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
FER	<input type="checkbox"/>	<input type="checkbox"/>	FES	<input type="checkbox"/>	<input type="checkbox"/>	FGFR1	<input checked="" type="checkbox"/>	<input type="checkbox"/>	FGFR1 [V561M]	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
FGFR2	<input type="checkbox"/>	<input type="checkbox"/>	FGFR3	<input type="checkbox"/>	<input type="checkbox"/>	FGFR3 [K650E]	<input type="checkbox"/>	<input type="checkbox"/>	FGFR3 [K650M]	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
FGFR4	<input type="checkbox"/>	<input type="checkbox"/>	FGFR4 [N535K] ※			FGFR4 [V550E]	<input type="checkbox"/>	<input type="checkbox"/>	FGFR4 [V550L]	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
FGR	<input type="checkbox"/>	<input type="checkbox"/>	FLT1	<input type="checkbox"/>	<input type="checkbox"/>	FLT3	<input checked="" type="checkbox"/>	<input type="checkbox"/>	FLT4	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
FMS	<input type="checkbox"/>	<input type="checkbox"/>	FRK	<input type="checkbox"/>	<input type="checkbox"/>	FYN [isoform a]	<input type="checkbox"/>	<input type="checkbox"/>	FYN [isoform b]	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
HCK	<input type="checkbox"/>	<input type="checkbox"/>	HER2	<input type="checkbox"/>	<input type="checkbox"/>	HER4	<input type="checkbox"/>	<input type="checkbox"/>	IGF1R	<input checked="" type="checkbox"/>	<input type="checkbox"/>		<input checked="" type="checkbox"/>	<input type="checkbox"/>		<input checked="" type="checkbox"/>	
INSR	<input type="checkbox"/>	<input type="checkbox"/>	IRR	<input type="checkbox"/>	<input type="checkbox"/>	ITK	<input checked="" type="checkbox"/>	<input type="checkbox"/>	JAK1	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
JAK2	<input type="checkbox"/>	<input type="checkbox"/>	JAK3	<input checked="" type="checkbox"/>	<input type="checkbox"/>	KDR	<input checked="" type="checkbox"/>	<input type="checkbox"/>	KIT	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
KIT [D816E]	<input type="checkbox"/>	<input type="checkbox"/>	KIT [D816V]	<input type="checkbox"/>	<input type="checkbox"/>	KIT [D816Y]	<input type="checkbox"/>	<input type="checkbox"/>	KIT [T670I]	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
KIT [V560G]	<input type="checkbox"/>	<input type="checkbox"/>	KIT [V654A]	<input type="checkbox"/>	<input type="checkbox"/>	LCK	<input checked="" type="checkbox"/>	<input type="checkbox"/>	LTK	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
LYNa	<input type="checkbox"/>	<input type="checkbox"/>	LYNb	<input type="checkbox"/>	<input type="checkbox"/>	MER	<input type="checkbox"/>	<input type="checkbox"/>	MET	<input checked="" type="checkbox"/>	<input type="checkbox"/>		<input checked="" type="checkbox"/>	<input type="checkbox"/>		<input checked="" type="checkbox"/>	
MET [D1228H]	<input type="checkbox"/>	<input type="checkbox"/>	MET [M1250T]	<input type="checkbox"/>	<input type="checkbox"/>	MET [Y1235D]	<input type="checkbox"/>	<input type="checkbox"/>	MUSK	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
PDGFRα	<input checked="" type="checkbox"/>	<input type="checkbox"/>	PDGFRα [D842V]	<input type="checkbox"/>	<input type="checkbox"/>	PDGFRα [T674I]	<input type="checkbox"/>	<input type="checkbox"/>	PDGFRα [V561D]	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
PDGFRβ	<input type="checkbox"/>	<input type="checkbox"/>	PYK2	<input checked="" type="checkbox"/>	<input type="checkbox"/>	RET	<input type="checkbox"/>	<input type="checkbox"/>	RET [G691S]	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
RET [M918T]	<input type="checkbox"/>	<input type="checkbox"/>	RET [S891A]	<input type="checkbox"/>	<input type="checkbox"/>	RET [Y791F]	<input type="checkbox"/>	<input type="checkbox"/>	RON	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
ROS	<input type="checkbox"/>	<input type="checkbox"/>	SRC	<input checked="" type="checkbox"/>	<input type="checkbox"/>	SRM	<input type="checkbox"/>	<input type="checkbox"/>	SYK	<input checked="" type="checkbox"/>	<input type="checkbox"/>		<input checked="" type="checkbox"/>	<input type="checkbox"/>		<input checked="" type="checkbox"/>	
TEC	<input type="checkbox"/>	<input type="checkbox"/>	TIE2	<input checked="" type="checkbox"/>	<input type="checkbox"/>	TNK1	<input type="checkbox"/>	<input type="checkbox"/>	TRKA	<input checked="" type="checkbox"/>	<input type="checkbox"/>		<input checked="" type="checkbox"/>	<input type="checkbox"/>		<input checked="" type="checkbox"/>	
TRKB	<input type="checkbox"/>	<input type="checkbox"/>	TRKC	<input type="checkbox"/>	<input type="checkbox"/>	TXK	<input type="checkbox"/>	<input type="checkbox"/>	TYK2	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	
TYRO3	<input checked="" type="checkbox"/>	<input type="checkbox"/>	YES	<input type="checkbox"/>	<input type="checkbox"/>	YES [T348I]	<input type="checkbox"/>	<input type="checkbox"/>	ZAP70	<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	

Serine/Threonine Kinases

	Km	1mM		Km	1mM		Km	1mM		Km	1mM
AKT1			AKT2			AKT3			AMPKα1/β1/γ1		
AMPKα2/β1/γ1			AurA			AurA/TPX2			AurB		
AurC			BRSK1			BRSK2			BUB1/BUB3		
CaMK1α		Km = 1mM	CaMK1δ			CaMK2α			CaMK2β		
CaMK2γ			CaMK2δ			CaMK4			CDC2/CycB1		
CDC7/ASK			CDK2/CycA2			CDK2/CycE1			CDK3/CycE1		Km = 1mM
CDK4/CycD3			CDK5/p25			CDK6/CycD3			CDK7/CycH/MAT1		
CDK9/CycT1			CGK2			CHK1			CHK2		
CK1α			CK1γ1			CK1γ2			CK1γ3		

How to use "Kinase List" function

- Identify the Complete Kinase List (CKL) sheet to the right (the fourth sheet).
- Extract your request kinase names by using the alphabetized kinase list, and copy the target names from the CKL and paste them to the cells below. For Km testing, the target name is sufficient. For 1mM ATP testing, choose the target name followed by 1mM on the CKL.

All kinase names chosen are indicated in red and the total target number is shown on the application sheet.

Please be noted that any kinase names other than our official product names cannot be entered.

Kinase List

Complete Kinase List sheet (on the fourth page)

Kinase Name

ABL
ABL_1mM
ABL [E255K]
ABL [E255K]_1mM
ABL [T315I]
ABL [T315I]_1mM
ACK
ACK_1mM
AKT1
AKT1_1mM
AKT2
AKT3
ALK
ALK_1mM
ALK [C1156Y]
ALK [C1156Y]_1mM
ALK [F1174L]
ALK [F1174L]_1mM
ALK [G1202R]
ALK [G1202R]_1mM
ALK [G1269A]
ALK [G1269A]_1mM
ALK [L1152insT]
ALK [L1152insT]_1mM
ALK [L1196M]
ALK [L1196M]_1mM
ALK [R1275Q]
ALK [R1275Q]_1mM
AMPK α 1/ β 1/ γ 1
AMPK α 1/ β 1/ γ 1_1mM
AMPK α 2/ β 1/ γ 1
ARG
ARG_1mM
AurA
AurA_1mM
AurA/TPX2
AurB
AurB_1mM
AurC
AurC_1mM
AXL
AXL_1mM
BLK
BLK_1mM
BMX
BMX_1mM
BRAF [V600E]_Cascade
BRAF_Cascade
BRK
BRK_1mM
BRSK1
BRSK1_1mM
BRSK2
BTK
BTK_1mM
BTK [C481S]
BTK [C481S]_1mM
BUB1/BUB3
CaMK1 α
CaMK1 δ
CaMK2 α
CaMK2 β
CaMK2 γ
CaMK2 δ
CaMK4
CaMK4_1mM

Target for 1mM ATP testing

How to use the Complete Kinase List (CKL) sheet

- 1) Extract your request kinase names by filter function of the A row.
- 2) Copy the extracted kinase names and paste them onto the "Kinase List" on the first sheet

For 1mM ATP testing, choose the target name followed by _1mM [underscore1mM]