

# 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	

Study Information	
Assay Type	<input checked="" type="radio"/> IC50 determination
Physical State	<input checked="" type="radio"/> Solution <input type="radio"/> Solid
Unit of Concentration	<input type="radio"/> $\mu$ M <input checked="" type="radio"/> $\mu$ g/mL
Residual materials	<input checked="" type="radio"/> Dispose after completion of study. <input type="radio"/> Return to the customer (at customer's cost)

Additional Information
<div style="border: 1px solid black; padding: 5px; text-align: center;"> For special request, please write here. </div>

Notice
<p>Detailed information about each kinase and assay condition can be checked on the latest Kinase Profiling Book.  <a href="http://www.carnabio.com/output/pdf/ProfilingProfilingBook_en.pdf">http://www.carnabio.com/output/pdf/ProfilingProfilingBook_en.pdf</a></p> <p><b>Sample Preparation</b>  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 : <math>\leq 100</math> target kinases = 500 <math>\mu</math>L , <math>&gt; 100</math> target kinases = 1000 <math>\mu</math>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.</p> <p>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 <math>&gt;50</math> <math>\mu</math>M test concentration, please contact us in advance)</p> <p>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: <a href="mailto:info@carnabio.com">info@carnabio.com</a></p> <p><b>Storage</b>  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.</p> <p><b>Disposal</b>  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.</p> <p>Please feel free to contact us if you have any questions or comments.</p>

## 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. (mg/mL)	Volume ( $\mu$ L)	Purity (%)	Storage Temp.	Start Concentration (ug/mL) 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																			

## Target Kinase Selection

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

The total number of assays you have chosen

**7 assays selected**

Check the box of kinase you would like to request.

☐ Hide All Unselected Kinases.

### Tyrosine Kinases

	Km	1mM		Km	1mM		Km	1mM		Km	1mM
ABL	<input type="checkbox"/>	<input type="checkbox"/>	ABL [E255K]	<input type="checkbox"/>	<input type="checkbox"/>	ABL [T315I]	<input type="checkbox"/>	<input type="checkbox"/>			
ALK	<input checked="" type="checkbox"/>	<input type="checkbox"/>	ALK [C1156Y]	<input checked="" type="checkbox"/>	<input type="checkbox"/>	ALK [F1174L]	<input checked="" type="checkbox"/>	<input type="checkbox"/>	ALK [G1202R]	<input checked="" type="checkbox"/>	<input type="checkbox"/>
ALK [G1269A]	<input checked="" type="checkbox"/>	<input type="checkbox"/>				ALK [L1196M]	<input checked="" type="checkbox"/>	<input type="checkbox"/>	ALK [R1275Q]	<input checked="" type="checkbox"/>	<input type="checkbox"/>
			NPM1-ALK	<input type="checkbox"/>	<input type="checkbox"/>	ARG	<input type="checkbox"/>	<input type="checkbox"/>			
			BMX	<input type="checkbox"/>	<input type="checkbox"/>						
EGFR	<input 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"/>
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"/>
EPHA2	<input 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"/>
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"/>
EPHB2	<input type="checkbox"/>	<input type="checkbox"/>	EPHB3	<input type="checkbox"/>	<input type="checkbox"/>	EPHB4	<input type="checkbox"/>	<input type="checkbox"/>			
FER	<input type="checkbox"/>	<input type="checkbox"/>	FES	<input type="checkbox"/>	<input type="checkbox"/>	FGFR1	<input type="checkbox"/>	<input type="checkbox"/>			
FGFR2	<input type="checkbox"/>	<input type="checkbox"/>									
FGFR4	<input type="checkbox"/>	<input type="checkbox"/>							FGFR4 [V550L]	<input type="checkbox"/>	<input type="checkbox"/>
FGR	<input type="checkbox"/>	<input type="checkbox"/>	FLT1	<input type="checkbox"/>	<input type="checkbox"/>				FLT4	<input type="checkbox"/>	<input type="checkbox"/>
			FRK	<input type="checkbox"/>	<input type="checkbox"/>						
						HER4	<input type="checkbox"/>	<input type="checkbox"/>	IGF1R	<input type="checkbox"/>	<input type="checkbox"/>
INSR	<input type="checkbox"/>	<input type="checkbox"/>									
JAK2	<input type="checkbox"/>	<input type="checkbox"/>	JAK3	<input type="checkbox"/>	<input type="checkbox"/>	KDR	<input type="checkbox"/>	<input type="checkbox"/>			
						LCK	<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 type="checkbox"/>	<input 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"/>			
PDGFRα	<input type="checkbox"/>	<input type="checkbox"/>	PDGFRα [D842V]	<input type="checkbox"/>	<input type="checkbox"/>				PDGFRα [V561D]	<input type="checkbox"/>	<input type="checkbox"/>
PDGFRβ	<input type="checkbox"/>	<input type="checkbox"/>	PYK2	<input type="checkbox"/>	<input type="checkbox"/>	RET	<input type="checkbox"/>	<input type="checkbox"/>	RET [G691S]	<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"/>
ROS	<input type="checkbox"/>	<input type="checkbox"/>	SRC	<input type="checkbox"/>	<input type="checkbox"/>	SRM	<input type="checkbox"/>	<input type="checkbox"/>			
TEC	<input type="checkbox"/>	<input type="checkbox"/>	TIE2	<input type="checkbox"/>	<input type="checkbox"/>						
TRKB	<input type="checkbox"/>	<input type="checkbox"/>	TRKC	<input type="checkbox"/>	<input type="checkbox"/>						
TYRO3	<input type="checkbox"/>	<input type="checkbox"/>	YES	<input type="checkbox"/>	<input type="checkbox"/>	YES [T348I]	<input type="checkbox"/>	<input type="checkbox"/>			

## Serine/Threonine Kinases

	Km	1mM		Km	1mM		Km	1mM		Km	1mM
AKT1	<input type="checkbox"/>	<input type="checkbox"/>	AKT2	<input type="checkbox"/>		AKT3	<input type="checkbox"/>		AMPK $\alpha$ 1/ $\beta$ 1/ $\gamma$ 1	<input type="checkbox"/>	<input type="checkbox"/>
AMPK $\alpha$ 2/ $\beta$ 1/ $\gamma$ 1	<input type="checkbox"/>		AurA	<input type="checkbox"/>	<input type="checkbox"/>	AurA/TPX2	<input type="checkbox"/>		AurB	<input type="checkbox"/>	<input type="checkbox"/>
			BRSK1	<input type="checkbox"/>	<input type="checkbox"/>	BRSK2	<input type="checkbox"/>		BUB1/BUB3	<input type="checkbox"/>	
						CaMK2 $\alpha$	<input type="checkbox"/>		CaMK2 $\beta$	<input type="checkbox"/>	
CaMK2 $\gamma$	<input type="checkbox"/>		CaMK2 $\delta$	<input type="checkbox"/>		CaMK4	<input type="checkbox"/>	<input type="checkbox"/>	CDC2/CycB1	<input type="checkbox"/>	<input type="checkbox"/>
			CDK2/CycA2	<input type="checkbox"/>	<input type="checkbox"/>	CDK2/CycE1	<input type="checkbox"/>	<input type="checkbox"/>	CDK3/CycE1	<input type="checkbox"/>	Km = 1mM
			CDK5/p25	<input type="checkbox"/>	<input type="checkbox"/>						
			CGK2	<input type="checkbox"/>		CHK1	<input type="checkbox"/>	<input type="checkbox"/>	CHK2	<input type="checkbox"/>	<input type="checkbox"/>
			CK1 $\gamma$ 1	<input type="checkbox"/>		CK1 $\gamma$ 2	<input type="checkbox"/>		CK1 $\gamma$ 3	<input type="checkbox"/>	
CK1 $\delta$	<input type="checkbox"/>					CK2 $\alpha$ 1/ $\beta$	<input type="checkbox"/>	<input type="checkbox"/>	CK2 $\alpha$ 2/ $\beta$	<input type="checkbox"/>	
CLK1	<input type="checkbox"/>	<input type="checkbox"/>	CLK2	<input type="checkbox"/>	<input type="checkbox"/>	CLK3	<input type="checkbox"/>				
DAPK1	<input type="checkbox"/>	<input type="checkbox"/>				DYRK1A	<input type="checkbox"/>	<input type="checkbox"/>	DYRK1B	<input type="checkbox"/>	<input type="checkbox"/>
DYRK2	<input type="checkbox"/>		DYRK3	<input type="checkbox"/>					Erk1	<input type="checkbox"/>	<input type="checkbox"/>
Erk2	<input type="checkbox"/>	<input type="checkbox"/>				GSK3 $\alpha$	<input type="checkbox"/>	<input type="checkbox"/>	GSK3 $\beta$	<input type="checkbox"/>	<input type="checkbox"/>
Haspin	<input type="checkbox"/>					HIPK1	<input type="checkbox"/>		HIPK2	<input type="checkbox"/>	
HIPK3	<input type="checkbox"/>		HIPK4	<input type="checkbox"/>	<input type="checkbox"/>				IKK $\beta$	<input type="checkbox"/>	<input type="checkbox"/>
									JNK1	<input type="checkbox"/>	<input type="checkbox"/>
JNK2	<input type="checkbox"/>	<input type="checkbox"/>	JNK3	<input type="checkbox"/>	<input type="checkbox"/>						
			MAPKAPK2	<input type="checkbox"/>	<input type="checkbox"/>	MAPKAPK3	<input type="checkbox"/>		MAPKAPK5	<input type="checkbox"/>	
MARK1	<input type="checkbox"/>		MARK2	<input type="checkbox"/>		MARK3	<input type="checkbox"/>		MARK4	<input type="checkbox"/>	
						MNK1	<input type="checkbox"/>		MNK2	<input type="checkbox"/>	
						MSK1	<input type="checkbox"/>			<input type="checkbox"/>	
NEK2	<input type="checkbox"/>	<input type="checkbox"/>	NEK4	<input type="checkbox"/>							
			NIM1K	<input type="checkbox"/>		NuaK1	<input type="checkbox"/>		NuaK2	<input type="checkbox"/>	
p38 $\alpha$	<input type="checkbox"/>	<input type="checkbox"/>	p38 $\beta$	<input type="checkbox"/>	<input type="checkbox"/>	p38 $\gamma$	<input type="checkbox"/>	<input type="checkbox"/>	p38 $\delta$	<input type="checkbox"/>	<input type="checkbox"/>
p70S6K	<input type="checkbox"/>	<input type="checkbox"/>	p70S6K $\beta$	<input type="checkbox"/>		PAK1	<input type="checkbox"/>		PAK2	<input type="checkbox"/>	<input type="checkbox"/>
			PAK5	<input type="checkbox"/>							
PIM1	<input type="checkbox"/>	<input type="checkbox"/>									
						PKC $\delta$	<input type="checkbox"/>		PKC $\epsilon$	<input type="checkbox"/>	<input type="checkbox"/>
PKC $\zeta$	<input type="checkbox"/>		PKC $\eta$	<input type="checkbox"/>							
PKD1	<input type="checkbox"/>		PKD2	<input type="checkbox"/>	<input type="checkbox"/>						

