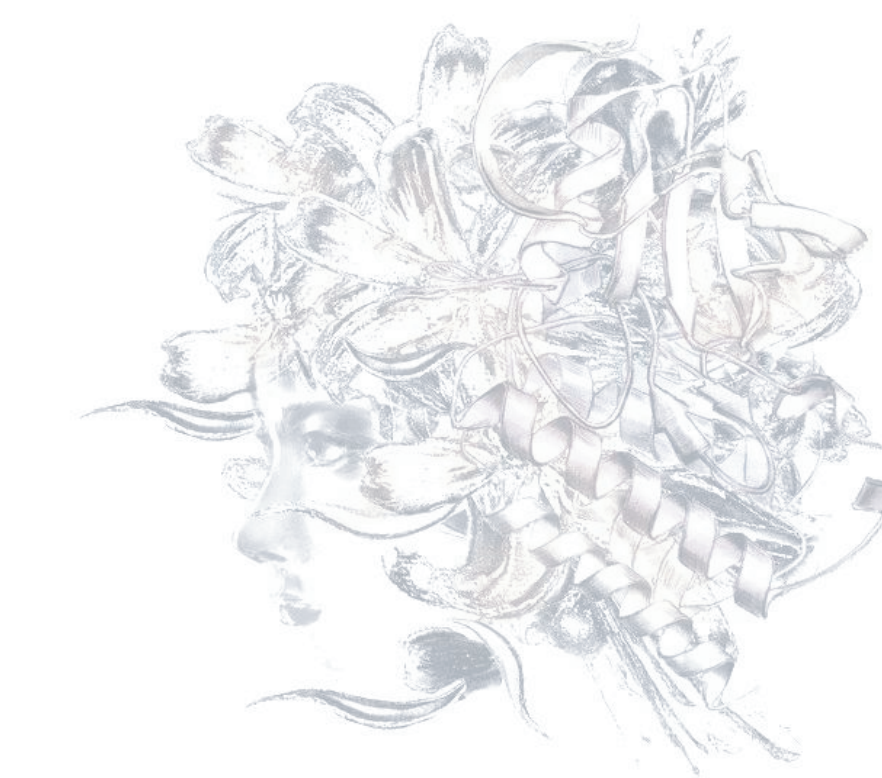


QuickScout™

A Comprehensive Tyrosine Kinase Profiling!



Kinase Inhibitor Profiling Service -- Comparative Inhibitory Activity

Our newest profiling services use pre-selected Tyrosine Kinase panels to permit the most rapid and thorough evaluation of potential kinase inhibitor selectivity.

QuickScout™ Comprehensive TK Panel ver 1.0

Carna Biosciences currently maintains stocks of, and has developed specific assays for 77 of the approximately 90 reported human tyrosine kinases, with more being developed continuously. QuickScout Comprehensive tests your activity modulator against each of these 77 enzymes, and offers the most complete specificity and selectivity assessment on the planet.

QuickScout™ TK Screen Panel ver 1.0

For the customer on a tighter budget, our QuickScout TK Screening Panel provides similar information, but on a smaller 20-enzyme panel of typical tyrosine kinases (contact us for detailed information).

-- Available Kinases for Screening --

ABL	JAK3	EGFR	FLT3	PDGFR α
ACK	SRC	EphA2	IGF1R	TIE2
CSK	SYK	EphB4	KDR	TRKA
FAK	TEC	FGFR1	MET	TYRO3

77 Tyrosine kinases are available NOW!

--- Also, we are providing custom profiling service such as IC₅₀ determinations. Please contact us for further information! ---

