

Target	Product Name	Partner Protein	Select for Profiling
Receptor Tyrosine Kinase Assays			
C-MET*	PathHunter™ C-MET Functional Assay	GrB2	
DDR2	PathHunter™ DDR2 Functional Assay	PLCG1	
EphB4*	PathHunter™ EphB4 Functional Assay	SHC1	
ErbB1*	PathHunter™ ErbB1 Functional Assay	PLCG1	
ErbB2/ErbB3*	PathHunter™ ErbB2/ErbB3 Functional Assay	GrB2	
ErbB4*	PathHunter™ ErbB4 Functional Assay	SHC1	
Flt3*	PathHunter™ Flt3 Functional Assay	P85	
IGFR1*	PathHunter™ IGFR1 Functional Assay	SHC1	
INSR*	PathHunter™ INSR Functional Assay	PLCG1	
PDGFRb*	PathHunter™ PDGFRb Functional Assay	PLCG1	
TrkA*	PathHunter™ TrkA Functional Assay	SHC1	
TrkA-P75*	PathHunter™ TrkA –P75Functional Assay	SHC1	
TrkB*	PathHunter™ TrkB Functional Assay	SHC1	
TrkB-P75*	PathHunter™ TrkB-P75 Functional Assay	SHC1	
TrkC*	PathHunter™ TrkC Functional Assay	SHC1	
TrkC-P75*	PathHunter™ TrkC-P75 Functional Assay	SHC1	
Cytosolic Tyrosine Kinase Assays			
PRLR-JAK1*	PathHunter™ PRLR-JAK1 Functional Assay	PLCG2	
PRLR-JAK2*	PathHunter™ PRLR-JAK2 Functional Assay	PLCG2	
CSF3R-JAK1	PathHunter™ PRLR-JAK1 Functional Assay	SHC1	
GHR-JAK1	PathHunter™ PRLR-JAK2 Functional Assay	PLCG2	

* ALSO AVAILABLE IN A SCHEDULED PANEL