

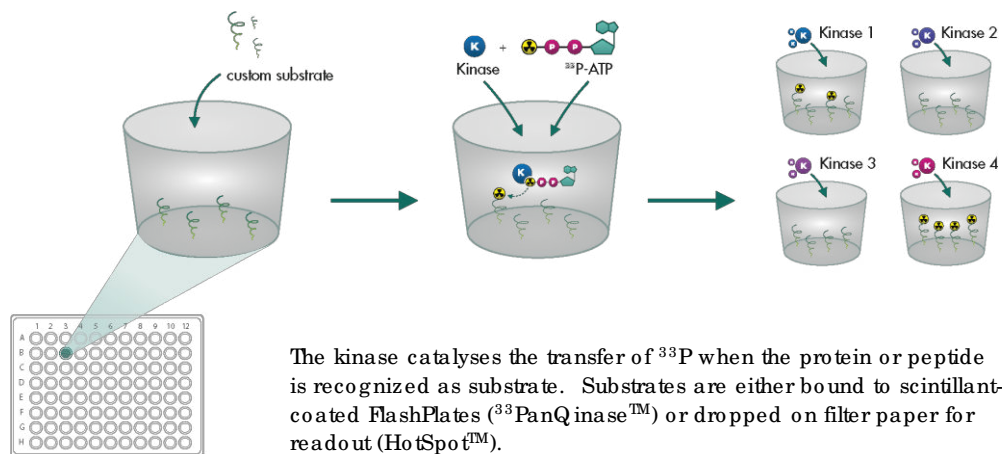
Purpose of the KinaseFinder Service

The KinaseFinder Service is intended to identify protein kinases, which accept a given test sample (protein or biotinylated peptide) as a substrate.

Type of KinaseFinder	Number of protein kinases	Assay Format
Tyr protein kinases	94	³³ PanQinase™
Ser/Thr protein kinases	245	³³ PanQinase™
Tyr and Ser/Thr protein kinases	339	³³ PanQinase™
Wild Type Kinase Panel	376	HotSpot™
Custom panel	Selected by customer	HotSpot™

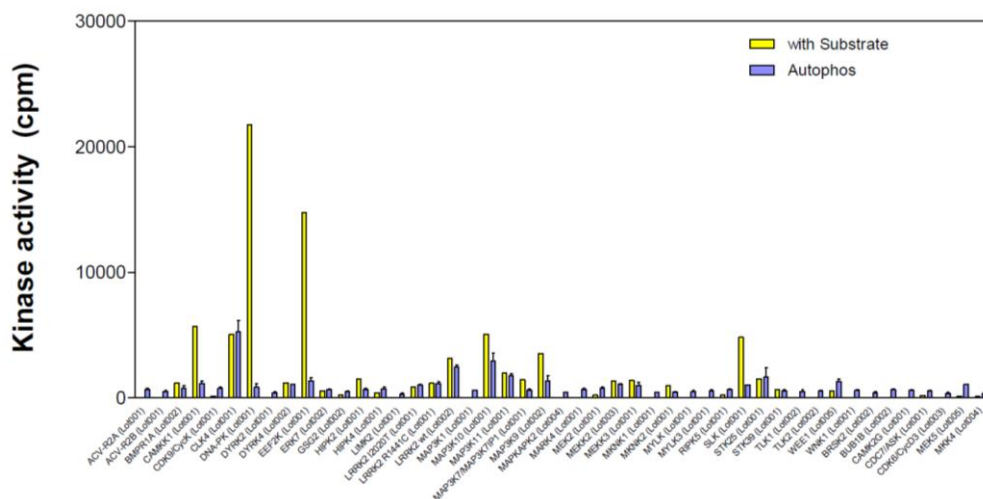
Assay principle

The target substrate is incubated with different kinases in every well of a multi-well plate together with ³³P-ATP that serves as phosphate donor.



The kinase catalyses the transfer of ³³P when the protein or peptide is recognized as substrate. Substrates are either bound to scintillant-coated FlashPlates (³³PanQinase™) or dropped on filter paper for readout (HotSpot™).

Example Data



Example data with 50 Ser/Thr kinases. Activity of kinases with example substrate protein are shown (corrected for substrate background activity). Substrate protein was tested at one concentration. Yellow: Kinase with substrate protein; blue: Kinase without substrate protein (autophosphorylation). A ratio value between substrate phosphorylation and apparent autophosphorylation > 3 may be considered as significant. Performed with ³³PanQinase™ assay. cpm = counts per minute