

Info Sheet KinaseFinder

(Version 16)

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 in ProQinase's radiometric ³³PanQinase® Assay.

1. Submission of Test Samples

1.1 Protein Samples

Please provide the indicated volumes of a 1 mg/ml stock solution in H₂O or 50 mM HEPES, pH 7.5. Glycerol up to 10%, beta-mercaptoethanol/DTT up to 5 mM and NaCl up to 300 mM are tolerable. The buffer must not contain detergents or imidazole above 50 mM since these will interfere with the assay. If your test sample is not soluble in aqueous solution please contact us. We recommend to ship samples on dry ice to avoid damage of the proteins during transport.

Primary screening will be performed at 5 µg/50 µl sample assay concentration (100 µg/ml).

Hit confirmation will be performed at 2.5 µg/50 µl, 5 µg/50 µl and 10 µg/50 µl sample assay concentration.

Please note: The sample(s) may be returned at customer's expense. From the date of report, the sample(s) will be stored at -20° Celsius for 6 months and will then be discarded without further notice.

KinaseFinder with 1 assay plate (Tyr version):

- 1 ml ⇒ primary hit finding and hit confirmation with max. 3 hits
- 1.5 ml ⇒ primary hit finding and hit confirmation with max. 7 hits

KinaseFinder with 3 assay plates (Ser/Thr version):

- 2.40 ml ⇒ primary hit finding and hit confirmation with max. 3 hits
- 3.25 ml ⇒ primary hit finding and hit confirmation with max. 7 hits

KinaseFinder with 4 assay plates (complete version):

- 3.25 ml ⇒ primary hit finding and hit confirmation with max. 3 hits
- 3.75 ml ⇒ primary hit finding and hit confirmation with max. 7 hits

Protein stock solutions of other concentrations:

Primary screening will be performed at a sample assay concentration 10 fold lower than the provided sample stock solution concentration.

Hit confirmation will be performed at sample assay concentrations 20-fold, 10-fold and 5-fold lower than the sample stock solution concentration.

1.2 Biotinylated Peptide Samples

Please provide a 200 μM stock solution of the biotinylated peptide, depending on the assay setup, see section 2) in H_2O or 50 mM HEPES, pH 7.5. If your test sample is not soluble in aqueous solution please contact us. Volumes required are a tenth of the volumes indicated above for protein test samples. We recommend to ship samples on dry ice to avoid damage of the peptides during transport.

Primary screening will be performed at 1 μM sample assay concentration.

Hit confirmation will be performed at 1 μM , 0.5 μM and 0.25 μM sample assay concentration.

Please note: The sample(s) may be returned at customer's expense. From the date of report, the sample(s) will be stored at -20° Celsius for 6 months and will then be discarded without further notice.

Peptide stock solutions of other concentrations:

Primary screening will be performed at a sample assay concentration which is 200-fold lower than the sample stock solution concentration.

Hit confirmation will be performed at a sample assay concentration which is 200-fold, 400-fold and 800-fold lower than the sample stock solution concentration.

Please note: Testing of biotinylated peptides above 1 μM assay concentration is not possible, due to the limited binding capacity of the streptavidine-coated HTS Plus FlashPlates.

2. Assay Conditions

The KinaseFinder experiments are performed with a radiometric filter plate assay (protein samples) or a radiometric assay based on streptavidin-coated FlashPlate[®] HTS PLUS plates (biotinylated peptide samples).

2.1 Phosphorylation reaction

The samples are tested against each kinase in singlicate in a 96-well polypropylene microtiter plate in a 50 µl reaction volume. The reaction cocktail is pipetted in the following order:

- 10 µl of kinase
- 40 µl of buffer/ATP/test sample mixture

The reaction cocktail contains 60 mM HEPES-NaOH, pH 7.5, 3 mM MgCl₂, 3 mM MnCl₂, 3 µM Na-orthovanadate, 1.2 mM DTT, 1 µM ATP/[γ-³³P]-ATP (approx. 7.5 x 10⁵ cpm per well), protein kinase (1-300 ng/50µl, according to the ³³PanKinase[®] screening assay concentration of the respective kinase/kinase lot) and the protein/peptide of interest.

All PKC assays (except the PKC-mu and the PKC-nu assay) additionally contained 1 mM CaCl₂, 4 mM EDTA, 5 µg/ml Phosphatidylserine and 1 µg/ml 1,2-Dioleoyl-glycerol.

The CAMK1D, CAMK2A, CAMK2B, CAMK2D, CAMK4, CAMKK1, CAMKK2, DAPK2, EEF2K, MYLK, MYLK2 and MYLK3 assays additionally contained 1 µg/ml Calmodulin and 0.5 mM CaCl₂.

The PRKG1 and PRKG2 assays additionally contained 1 µM cGMP.

The DNA-PK assay additionally contained 2.5 µg/ml DNA.

The reaction cocktails are incubated at 30° C for 60 minutes.

Subsequently, depending on the type of test sample (protein or biotinylated peptide), the reaction cocktails are processed differently (see below).

2.2 Detection of phosphorylated proteins

Reaction cocktails with a protein or peptide as potential substrate are stopped with 20 µl of 10% (v/s) H₃PO₄. Subsequently, the reaction cocktails are transferred to pre-wetted 96-well filter plates (Millipore MSFC glass filter), followed by 10 min incubation at room temperature. Subsequently, the plates are washed three times with 250 µl of 150 mM H₃PO₄ and once with 20 µl of 100% ethanol. After drying the plates for 30 min at 40° C, 50 µl of scintillator are added and incorporation of ³³P_i ("counting of cpm") is determined with a microplate scintillation counter.

2.3 Detection of phosphorylated, biotinylated peptides

Reaction cocktails with a biotinylated peptide are stopped with 20 µl of 4.7 M NaCl/35 mM EDTA. Subsequently, the reaction cocktails are transferred to streptavidin-coated FlashPlate[®] HTS PLUS plates followed by 30 min incubation at room temperature on a shaker. Then, the plates are washed three times with 250 µl of 0.9% NaCl. Incorporation of ³³P_i ("counting of cpm") is determined with a microplate scintillation counter.

3. Evaluation of Raw Data

For evaluation of the results of the filter assay (testing of protein or peptide test samples), kinase autophosphorylation has previously been determined in three independent experiments and a mean autophosphorylation value has been calculated for each kinase. These values (normalized on the radioactivity input of the current experiment) are subtracted from the corresponding raw data obtained with each kinase. Moreover, the median of three background values (sample protein without enzyme) is subtracted from the raw data.

For evaluation of the results from the streptavidin-coated FlashPlate® HTS PLUS plate assay, the activity value (raw counts of the kinase assay plus biotinylated sample peptide as measured in the FlashPlate® HTS PLUS plate assay), the corresponding kinase background value (kinase w/o sample peptide) and the median of three background values of the sample peptide (sample peptide w/o enzyme) are measured. The median of three background values is subtracted from the raw data of the results of the testings of biotinylated sample peptide plus kinase.

4. Hit confirmation

Hits detected in a KinaseFinder experiment can be verified in a hit confirmation experiment, where the test sample will be tested at three different concentrations in triplicate (see section 1.1 and 1.2) at enzyme concentrations according to the standard ³³PanQinase® screening assay concentration of the respective kinase/kinase lot.

5. Protein Kinases

All kinases currently available for the KinaseFinder Service are listed on the following pages.

5.1 Ser/Thr Kinases

No.	Kinase Name	No.	Kinase Name	No.	Kinase Name	No.	Kinase Name
1	ACV-R1	63	COT	125	MEK2	187	PKMYT1
2	ACV-R1B	64	DAPK1	126	MEK5	188	PKN3
3	ACV-R2A	65	DAPK2	127	MEKK2	189	PLK1
4	ACV-R2B	66	DAPK3	128	MEKK3	190	PLK3
5	ACV-RL1	67	DCAMKL2	129	MELK	191	PRK1
6	AKT1	68	DMPK	130	MINK1	192	PRK2
7	AKT2	69	DNA-PK	131	MKK4	193	PRKD2
8	AKT3	70	DYRK1A	132	MKK6SDTD	194	PRKG1
9	AMPK-alpha1	71	DYRK1B	133	MKK7	195	PRKG2
10	ARK5	72	DYRK2	134	MKNK1	196	PRKX
11	ASK1	73	DYRK3	135	MKNK2	197	RAF1 YDYD
12	Aurora-A	74	DYRK4	136	MLK4	198	RIPK2
13	Aurora-B	75	EEF2K	137	MST1	199	RIPK5
14	Aurora-C	76	EIF2AK2	138	MST2	200	ROCK1
15	BMPR1A	77	EIF2AK3	139	MST3	201	ROCK2
16	B-RAF VE	78	ERK1	140	MST4	202	RPS6KA1
17	B-RAF wt	79	ERK2	141	mTOR	203	RPS6KA2
18	BRSK1	80	ERK7	142	MYLK	204	RPS6KA3
19	BRSK2	81	GRK2	143	MYLK2	205	RPS6KA4
20	BUB1B	82	GRK3	144	MYLK3	206	RPS6KA5
21	CAMK1D	83	GRK4	145	NEK1	207	RPS6KA6
22	CAMK2A	84	GRK5	146	NEK11	208	S6K
23	CAMK2B	85	GRK6	147	NEK2	209	S6K-beta
24	CAMK2D	86	GRK7	148	NEK3	210	SAK
25	CAMK2G	87	GSG2	149	NEK4	211	SGK1
26	CAMK4	88	GSK3-alpha	150	NEK6	212	SGK2
27	CAMKK1	89	GSK3-beta	151	NEK7	213	SGK3
28	CAMKK2	90	HIPK1	152	NEK9	214	SIK1 aa1-350
29	CDC42BPA	91	HIPK2	153	NIK	215	SIK2
30	CDC42BPB	92	HIPK3	154	NLK	216	SLK
31	CDC7/ASK	93	HIPK4	155	p38-alpha	217	SNARK
32	CDK1/CycA	94	HRI	156	p38-beta	218	SNK
33	CDK1/CycE	95	IKK-alpha	157	p38-delta	219	SRPK1
34	CDK16/CycK	96	IKK-beta	158	p38-gamma	220	SRPK2
35	CDK1CycB1	97	IKK-epsilon	159	PAK1	221	STK17A
36	CDK2/CycA	98	IRAK1	160	PAK2	222	STK23
37	CDK2/CycE	99	IRAK4	161	PAK3	223	STK25
38	CDK3/CycE	100	JNK1	162	PAK4	224	STK33
39	CDK4/CycD1	101	JNK2	163	PAK6	225	STK39
40	CDK4/CycD3	102	JNK3	164	PAK7	226	TAOK2
41	CDK5/p25NCK	103	LIMK1	165	PASK	227	TAOK3
42	CDK5/p35NCK	104	LIMK2	166	PBK	228	TBK1
43	CDK6/CycD1	105	LRRK G2019S	167	PDK1	229	TGFB-R1
44	CDK6/CycD3	106	LRRK2 I2020T	168	PHKG1	230	TGFB-R2
45	CDK7CycH	107	LRRK2 R1441C	169	PHKG2	231	TLK1
46	CDK8/CycC	108	LRRK2 wt	170	PIM1	232	TLK2
47	CDK9/CycK	109	MAP3K1	171	PIM2	233	TSF1
48	CDK9/CycT	110	MAP3K10	172	PIM3	234	TSK2
49	CHK1	111	MAP3K11	173	PKA	235	TSSK1
50	CHK2	112	MAP3K7/MAP3K7IP1	174	PKC-alpha	236	TTBK1
51	CK1-alpha1	113	MAP3K9	175	PKC-beta1	237	TTBK2
52	CK1-delta	114	MAP4K2	176	PKC-beta2	238	TTK
53	CK1-epsilon	115	MAP4K4	177	PKC-delta	239	VRK1
54	CK1-gamma1	116	MAP4K5	178	PKC-epsilon	240	VRK2
55	CK1-gamma2	117	MAPKAPK2	179	PKC-eta	241	WEE1
56	CK1-gamma3	118	MAPKAPK3	180	PKC-gamma	242	WNK1
57	CK2-alpha1	119	MAPKAPK5	181	PKC-iota	243	WNK2
58	CK2-alpha2	120	MARK1	182	PKC-mu	244	WNK3
59	CLK1	121	MARK2	183	PKC-nu	245	ZAK
60	CLK2	122	MARK3	184	PKC-theta		
61	CLK3	123	MARK4	185	PKC-zeta		
62	CLK4	124	MEK1 wt	186	PKC-zeta wt aa184-592		

5.2 Tyr Kinases

No.	Kinase Name	No.	Kinase Name
1	ABL1 T315I	63	MERTK
2	ABL1 wt	64	MET T1250M
3	ABL2	65	MET wt
4	ACK1	66	MET Y1235D
5	ALK F1174L (GST-HIS-tag)	67	MUSK
6	ALK L1196M (GST-HIS-tag)	68	PDGFR-alpha D842V
7	ALK wt	69	PDGFR-alpha wt
8	AXL	70	PDGFR-beta
9	BLK	71	PYK2
10	BMX	72	RET V804M
11	BRK	73	RET wt
12	BTK	74	RON
13	CSF1-R	75	ROS
14	CSK	76	SRC
15	DDR2 T654M	77	SRMS
16	DDR2 wt	78	SYK
17	EGF-R T790M	79	TEC
18	EGF-R wt	80	TIE2 wt
19	EPHA1	81	TIE2 Y897S
20	EPHA2	82	TNK1
21	EPHA3	83	TRK-A
22	EPHA4	84	TRK-B
23	EPHA5	85	TRK-C
24	EPHA6	86	TXK
25	EPHA7	87	TYK2
26	EPHA8	88	TYRO3
27	EPHB1	89	VEGF-R1
28	EPHB2	90	VEGF-R2
29	EPHB3	91	VEGF-R3
30	EPHB4	92	WEE1
31	ERBB2 wt	93	YES
32	ERBB4	94	ZAP70
33	FAK		
34	FER		
35	FES		
36	FGF-R1 V561M		
37	FGF-R1 wt		
38	FGF-R2		
39	FGF-R3 K650E		
40	FGF-R3 wt		
41	FGF-R4		
42	FGR		
43	FLT3 D835Y		
44	FLT3 ITD		
45	FLT3 wt		
46	FRK		
47	FYN		
48	HCK		
49	IGF1-R		
50	INS-R		
51	INSR-R		
52	ITK		
53	JAK1		
54	JAK2		
55	JAK3		
56	KIT D816V		
57	KIT T670I		
58	KIT wt		
59	LCK		
60	LTK		
61	LYN		
62	MATK		