

LabChip[®] EZ Reader Substrate Selection Guide

Substrate name	Cat. No.	Sequence	Recommended Enzymes
ProfilerPro Peptide 1	760345	5-FAM-AKRRRLSSLRA-COOH	IKK α , IKK β , IKK ϵ , LOK, PAK2, PAK3, PKG1 α , PKG1 β , ROCK2, ROCK1, DAPK1, ZIPK, PIM3, CLK3
ProfilerPro Peptide 2	760346	5-FAM-EAIYAAPFAKKK-CONH2	Abl, Abl(Q252H), Abl(T315I), Abl(H396P), Abl(E255K), Abl(Y253F), Abl(G250E), FGR, Arg, BTK, Flt-3, Flt-3(D835Y), MET, KIT, RET
ProfilerPro Peptide 4	760348	5-FAM-EGIYGVLFKKK-CONH2	FYN, LCK, LYN, SRC, Hck, BLK
ProfilerPro Peptide 5	760349	5-FAM-GQLIDSMANSFVGTR-CONH2	cRAF
ProfilerPro Peptide 6	760350	5-FAM-GRPRTSSFAEG-CONH2	AKT1, AKT2, AKT3, MSK1, MSK2, SGK1
ProfilerPro Peptide 7	760351	5-FAM-HMRSAMSGLHLVKRR-COOH	AMPK α 1
ProfilerPro Peptide 8	760352	5-FAM-IPTSPITTTYFFFKKK-COOH	Erk1, Erk2, CDK6, IRAK4, p38 α , p38 α (T106M), p38 β 2, p38 δ , p38 γ
ProfilerPro Peptide 10	760354	5-FAM-KKKVSRSGLYRSPSPENLNRPR-COOH	AMPK α 2/b1/1, BRSK1, BRSK2, CHK1, CHK2, MARK1, MARK2, MARK3, MARK4, MNK1, NuaK1, c-TAK1, TSSK1, TSSK2
ProfilerPro Peptide 11	760355	5-FAM-KKLNRTLVA-COOH	CAMK2 α , CAMK2 β , CAMK2 γ , CAMK4, RSK1, RSK2, RSK3, MAPKAPK3, NEK2, SGK2, SGK3, PhKg1, PhKg2, DCAML1, PKN1, PKN2, PASK
ProfilerPro Peptide 12	760356	5-FAM-KKLRRTLVA-COOH	Camk1a, Camk1d, MAPKAPK2, PKD1, PKD2, PDK3, MAPKAPK5 (PRAK), MELK
ProfilerPro Peptide 13	760357	5-FAM-KKSRGDYMTMQIG-CONH2	DDR2, INSR, IGF1R, MST2, FGFR1, ALK

Substrate name	Cat. No.	Sequence	Recommended Enzymes
ProfilerPro Peptide 14	760358	5-FAM-KRELVEPLTPSGEAPNQALLR-CONH2	JNK2
ProfilerPro Peptide 15	760359	5-FAM-KRREILSRRPpSYR-COOH	GSK3 β , GSK3a
ProfilerPro Peptide 16	760360	5-FAM-KRRRALpSVASLPGL-CONH2	CK1 δ , CK1 ϵ , CSNK1a, CK1g1, CK1g2, CK1g3
ProfilerPro Peptide 18	760362	5-FAM-QSPKKG-CONH2	CDK2
ProfilerPro Peptide 19	760363	5-FAM-RFARKGSLRQKNV-COOH	PKC α , PKC β 1, PKC β 2, PKC γ , PKC δ , PKC ϵ , PKC ξ , PKC θ , PKC ι
ProfilerPro Peptide 20	760364	5-FAM-RSRHSSYPAGT-CONH2	PIM1, PIM2
ProfilerPro Peptide 21	760365	5-FAM-LRRASLG-CONH2	PKAC α , AurA, AurB, AurC, PRKX
ProfilerPro Peptide 22	760366	5-FAM-EEPLYWSFPAKKK-CONH2	KDR (VEGFR2), EGFR, EGFR(T790M), EGFR(T790M,L856R), FER, BMX, FGFR1(V561M), FGFR2, FGFR2(N549H), FGFR3, FGFR3(K650E), FGFR4, FRK, HER4, JAK2, LTK, ROS, RET, RET(Y791F), RET(V804L), SYK,TEC, YES
ProfilerPro Peptide 24	760387	5-FAM-KKISGRLSPIMTEQ-CONH2	DYRK1 α , DYRK1b, DYRK3, DYRK4, HIPK1, HIPK2, HIPK3
ProfilerPro Peptide 25	760388	5-FAM-VDGKEIYNTIRRK-CONH2	HGK, GCK, MINK1, MST1, MST3, NEK1, CLK2
ProfilerPro Peptide 26	760389	5-FAM-ARKRERTYSFGHHA-COOH	P70S6K
ProfilerPro Peptide 27	760424	5-FAM-EFPIYDFLPAKKK-CONH2	FAK2, EphA1, EphA2, EphA3, EphA4, EphA5, EpHa8, EpHB1, EphB2, EphB3, EphB4, FGFR3, LynB, FES, PYK2, TXK, ITK, TYRO3, TRKB, TRKC, MER, SRM
ProfilerPro Peptide 29	760429	5-FAM-GGGPATPKKAKKL-CONH2	CDK2/-25, CDC2/cyclinB1, CDK3

Substrate name	Cat. No.	Sequence	Recommended Enzymes
ProfilerPro Peptide 30	760430	5-FAM-KKKKKEEYFFF-CONH2	FLT1, FLT4, FMS, AXL, PDGFRa, PDGFRa(D842V), PDGFRb, MSTR, KIT, KIT(T670I), MET(M1250T)
ProfilerPro Peptide 31	760480	5-FAM-RRRLSFAEPG-CONH2	PAK4, PAK5(PAK7)
ProfilerPro Peptide 32	760641	5-FAM-FLAKSFGSPNRAYKK-CONH2	NEK6 NEK7 NEK9
ProfilerPro Peptide 33	760642	5-FAM-C6-RLGRDKYKTLRQIRQ-CONH2	LRRK2 TAOK3 YSK SLK
ProfilerPro Peptide 34	760643	5-FAM-RRRFRPASPLRGPPK-COOH	DYRK2 MSSK1 CDK1 CDK4 CDK9 HIPK3 HIPK4 CLK1 CLK3 SRPK2
iFL-cAMP	760444	Proprietary	Phosphodiesterase Assays
iFL-cAMP	760445	Proprietary	Phosphodiesterase Assays
Histone 3 Acetylated Peptide	760511	FITC-TGGK(Ac)APR-COOH	Epigenetic Assays
p53 Acetylated Peptide	760512	FITC-TSPQPKK(Ac)-CONH2	Epigenetic Assays
Histone 4 Acetylated Peptide	760513	FITC-LGKGGAK(Ac)-CONH2	Epigenetic Assays

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2012, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

Version created 26 June 2012.