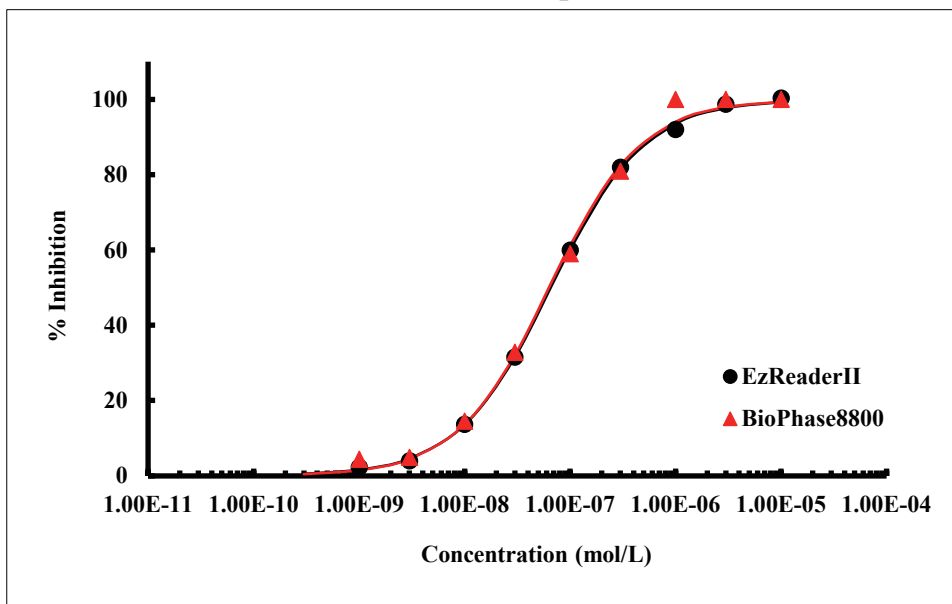


List of Substrates used in Mobility Shift Assay

A	ABLtide.....	01	K	Kemptide	13
	AMARA peptide	01	L	LIMKtide.....	13
B	Blk/Lyntide	02		LRRKtide	14
C	CDC25ctide	02	M	MCM2 peptide.....	14
	CDK7 peptide	03		MLCtide	15
	CDK9 substrate.....	03		Modified Erktide	15
	CHKtide	04		Modified Histone H1	16
	CK2tide	04		Modified IκBα-derived peptide	16
	CKtide.....	05		Moesin-derived peptide.....	17
	CREBtide-p.....	05	P	PDHKtide	17
	Crosstide	06		PKC peptide.....	18
	CSKtide	06	R	RS peptide	18
	CTD3 peptide.....	07	S	S6K peptide (N-FL)	19
D	DAPK1tide	07		S6K2 peptide	19
	DYRKtide-F.....	08		SAMS peptide	20
E	EEF2Ktide	08		SGKtide	20
	EGFR-derived peptide	09		SPAKtide.....	21
G	GS peptide.....	09		Srctide	21
H	H2A peptide.....	10		Synapsin peptide	22
	Histone H3 peptide	10	T	T308tide	22
I	IRAK1 peptide.....	11		TAOKtide.....	23
	IRS1	11	W	WASP peptide	23
	IκBα peptide	12			
J	JAK1 substrate peptide	12			

Substrate : **ABLtide**

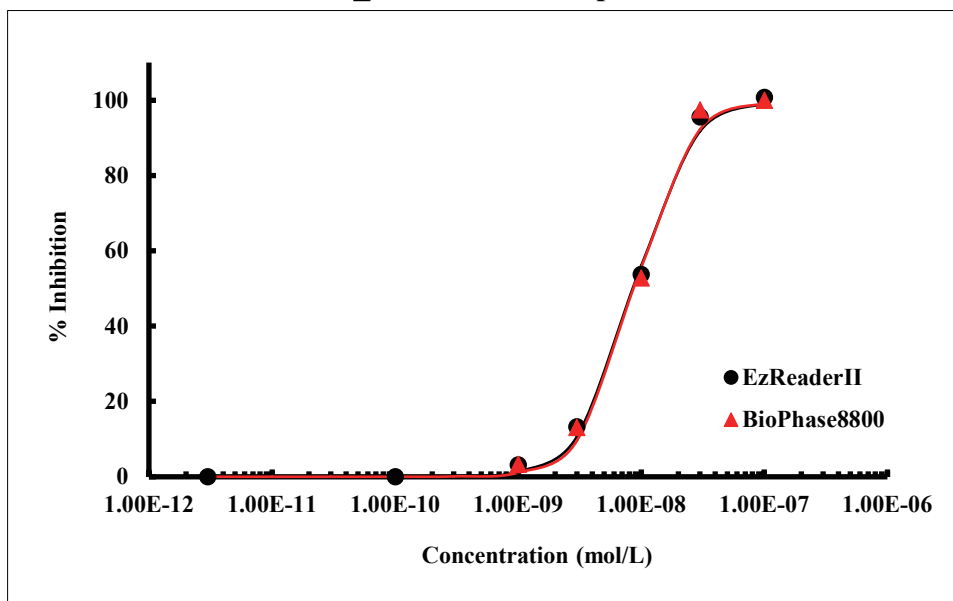
ABL vs. Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	66	63

Substrate : **AMARA peptide**

QIK_1mM vs Staurosporine

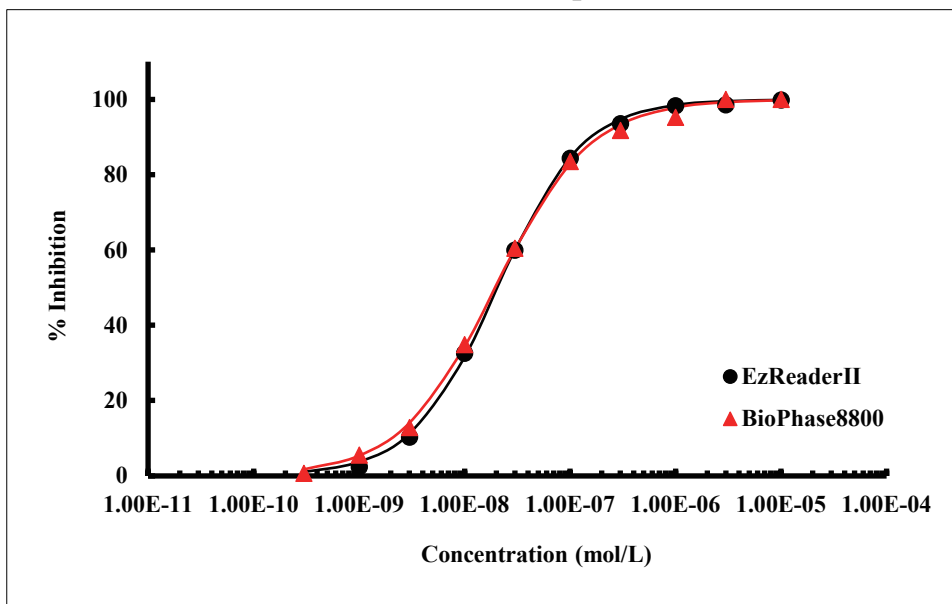


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	8.8	9.0



Substrate : **Blk/Lyntide**

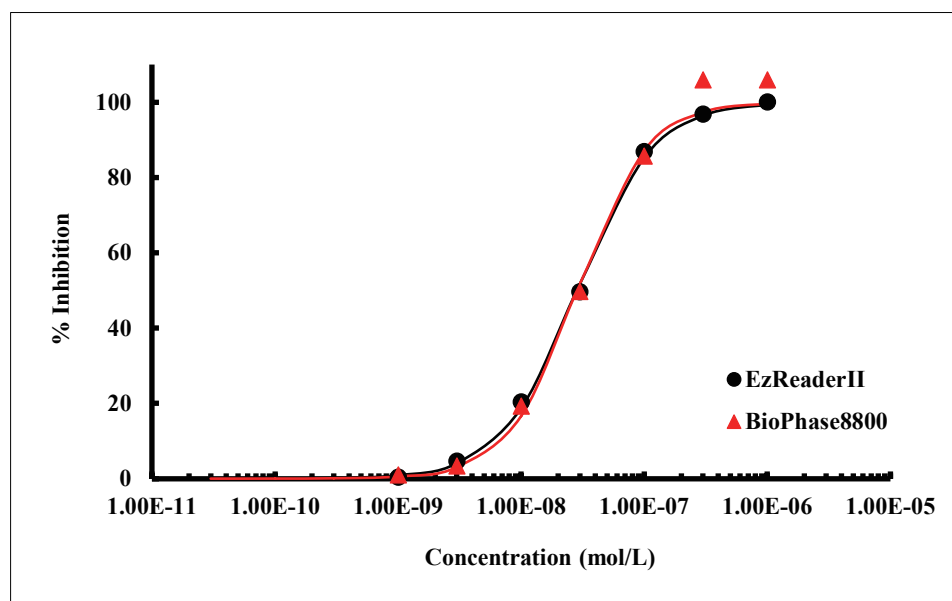
EPHA1 vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	21	20

Substrate : **CDC25ctide**

PLK3 vs GW843682X

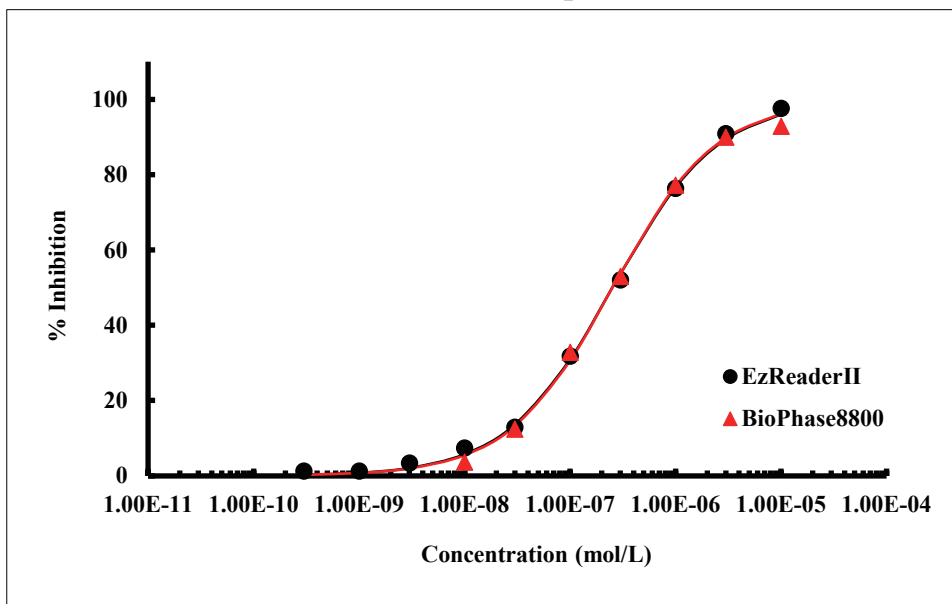


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	29	29



Substrate : CDK7 peptide

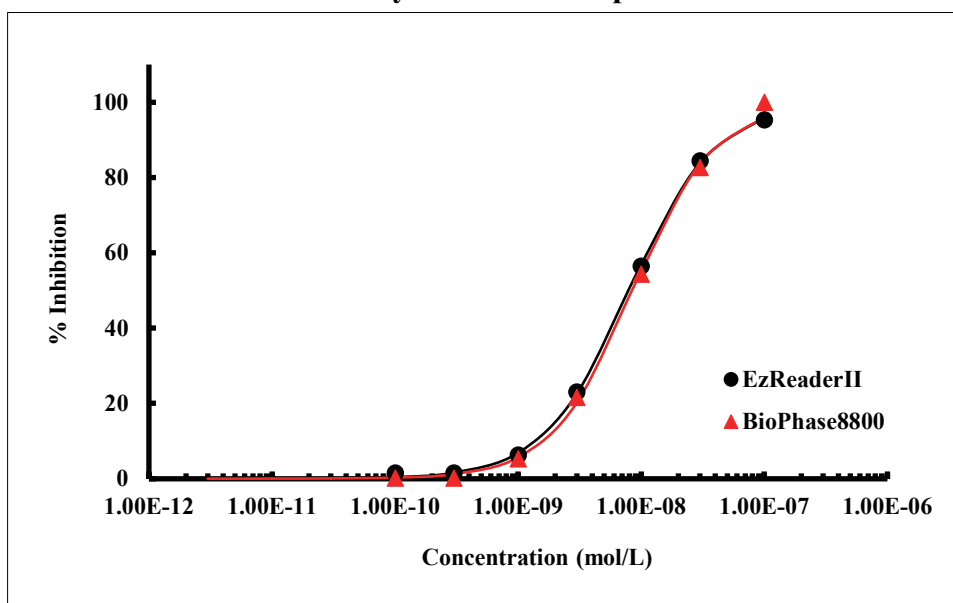
NEK9 vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	250	250

Substrate : CDK9 substrate

CDK9/CycT1 vs Staurosporine

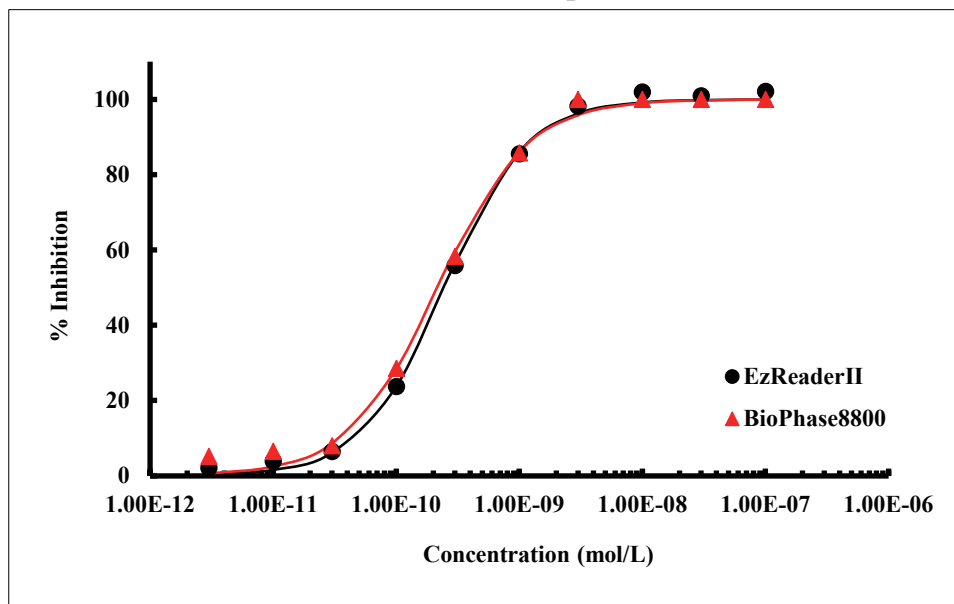


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	8.0	8.6



Substrate : **CHKtide**

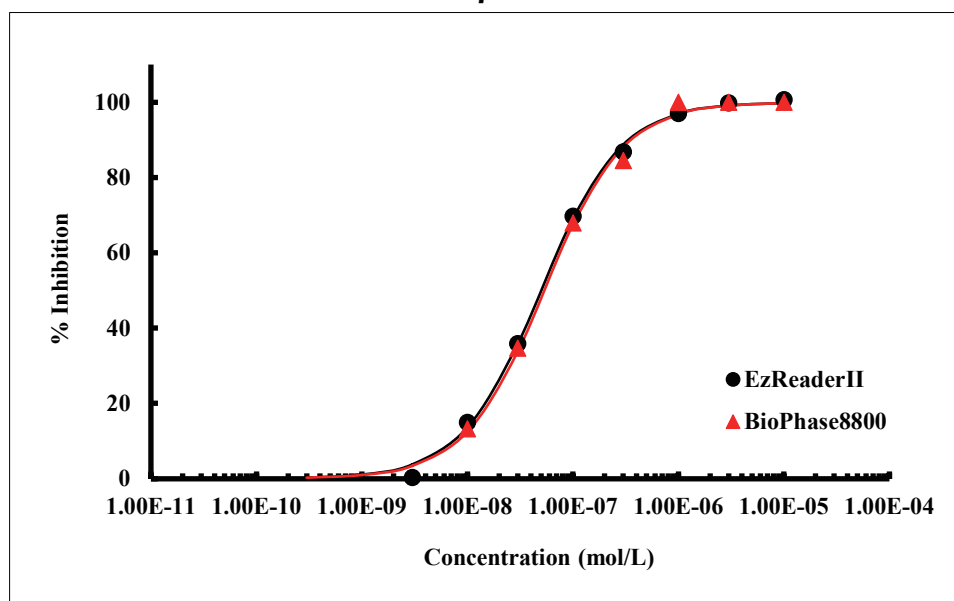
CHK1 vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	0.25	0.22

Substrate : **CK2tide**

CK2 α 1/ β vs TBB

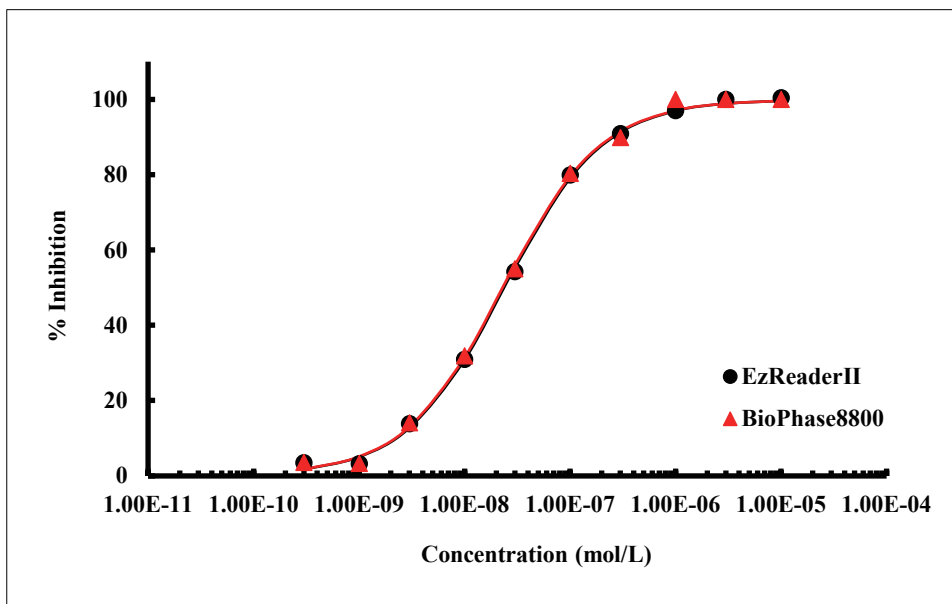


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	50	54



Substrate : CKtide

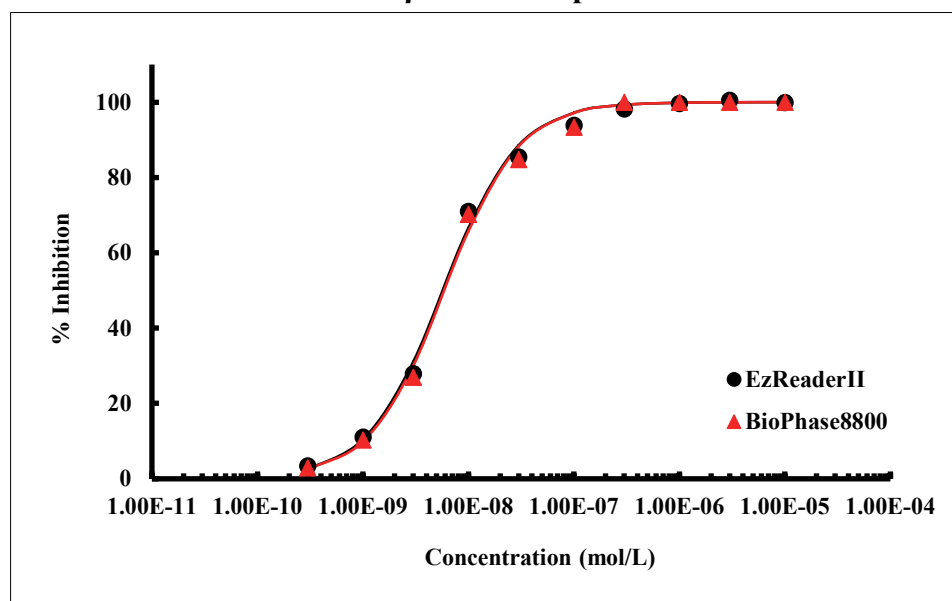
CK1δ vs 5-Iodotubercidin



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	24	23

Substrate : CREBtide-p

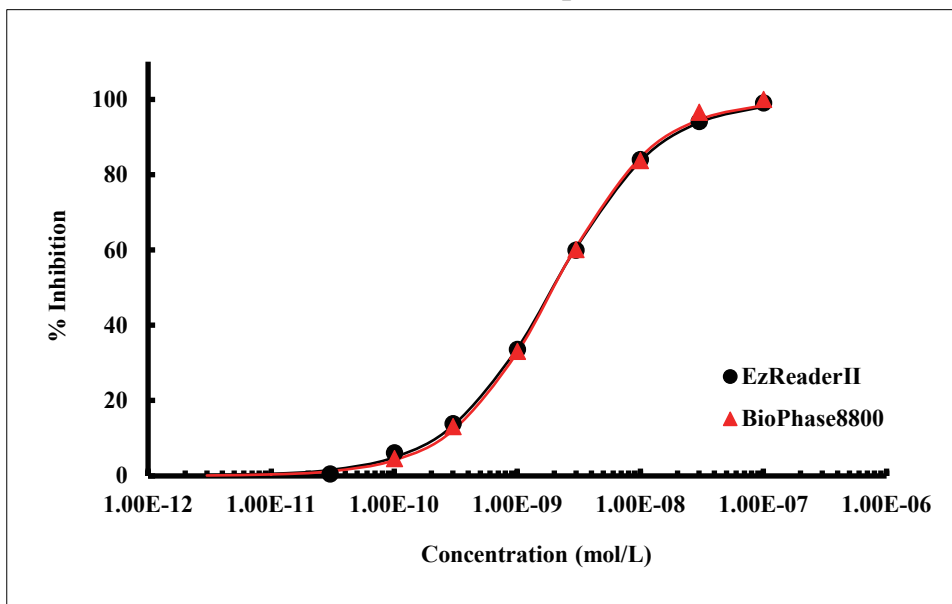
GSK3β vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	5.7	5.9

Substrate : Crosstide

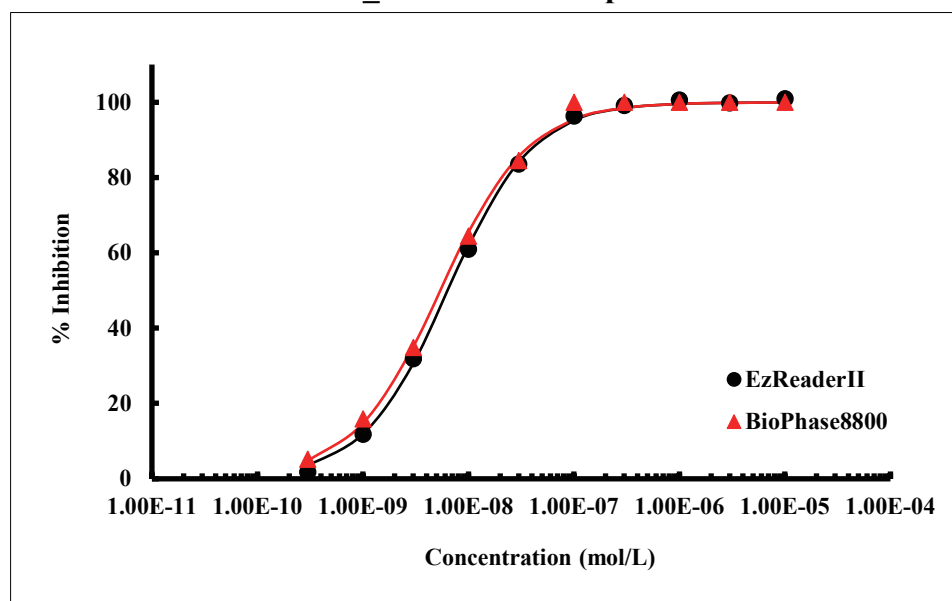
AKT1 vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	2.0	2.0

Substrate : CSKtide

FLT1_1mM vs Staurosporine

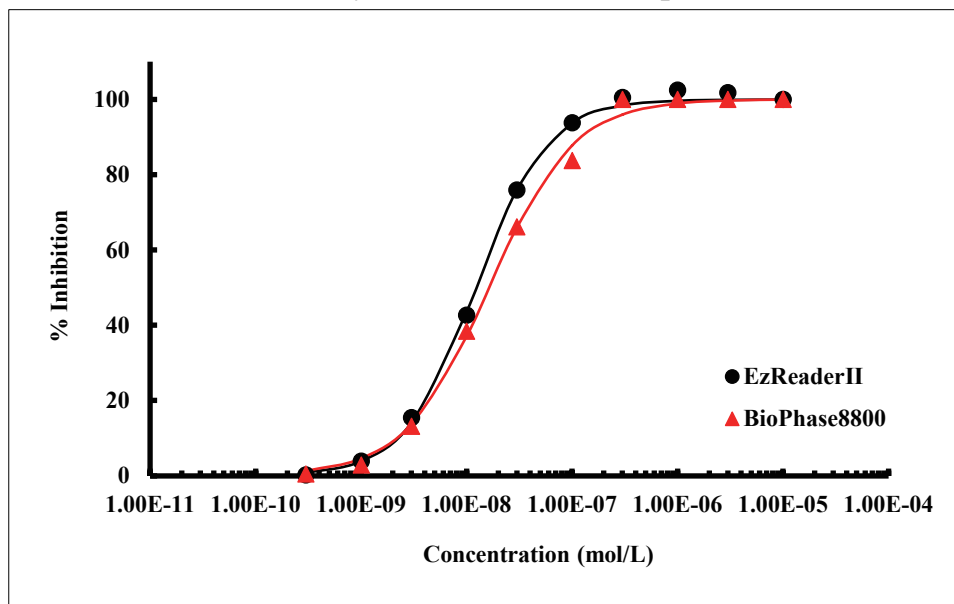


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	6.4	5.4



Substrate : CTD3 peptide

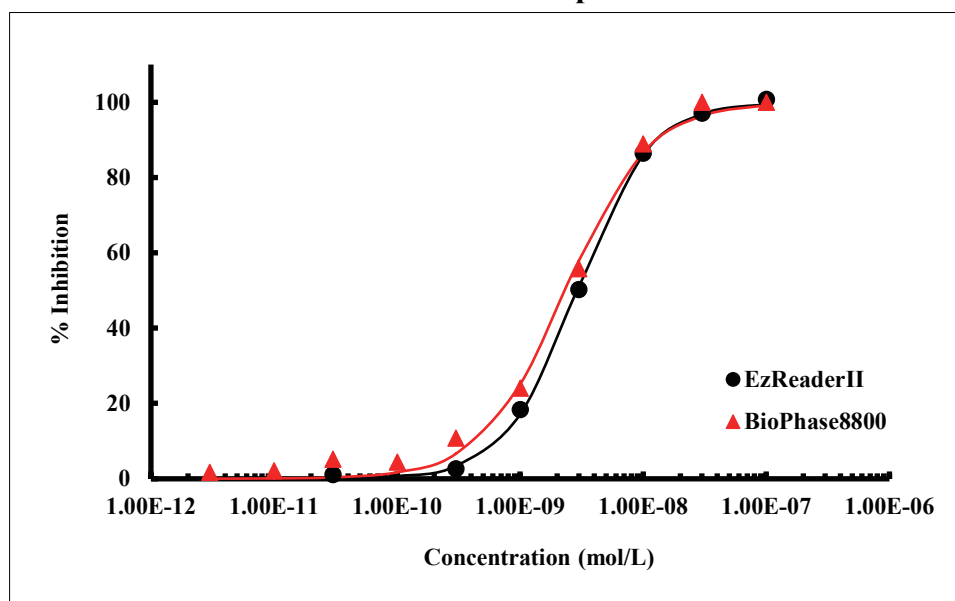
CDK7/CycH/MAT1 vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	12	16

Substrate : DAPK1tide

DAPK1 vs Staurosporine

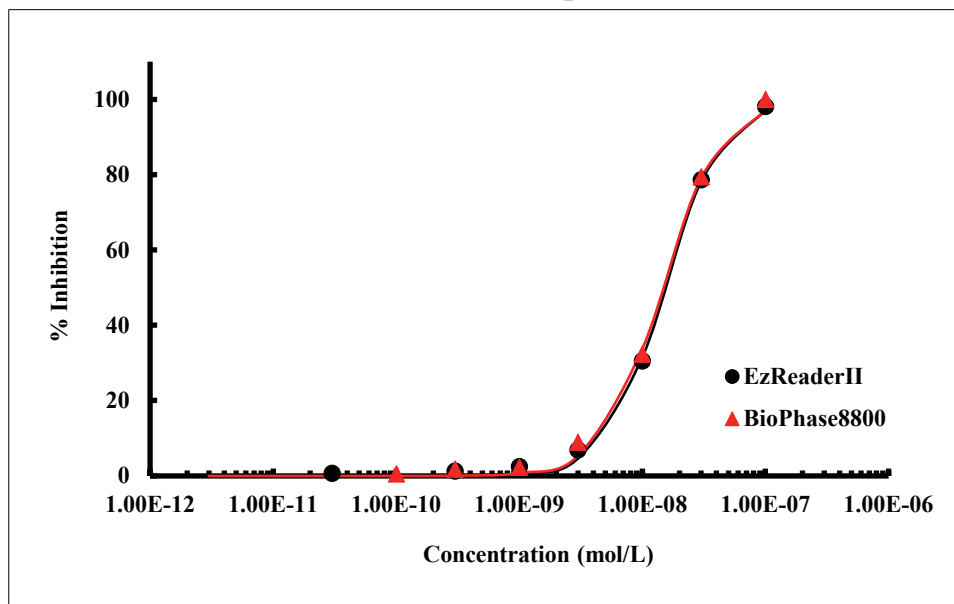


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	2.9	2.3



Substrate : **DYRKtide-F**

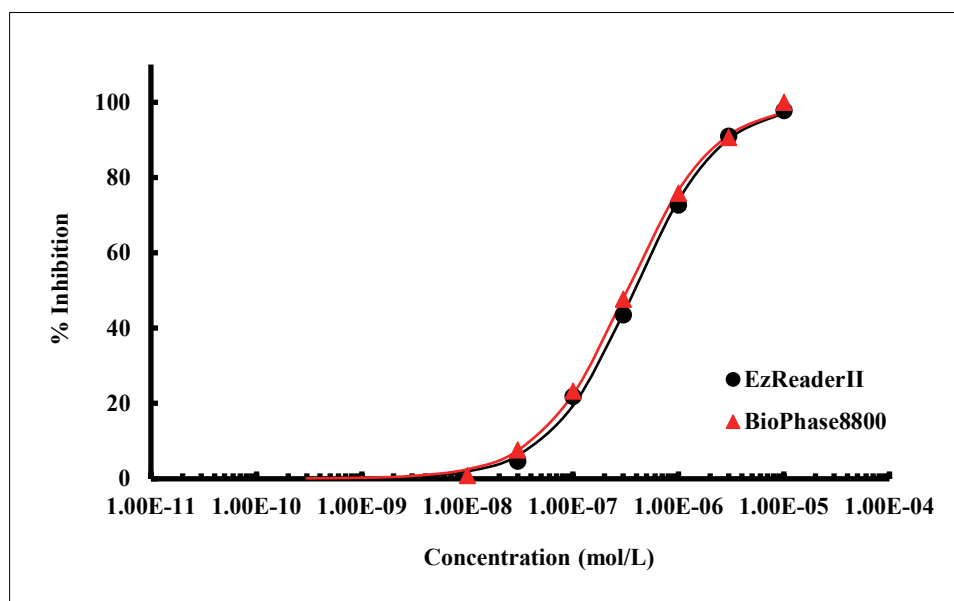
CLK1 vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	15	14

Substrate : **EEF2Ktide**

EEF2K vs A-484954

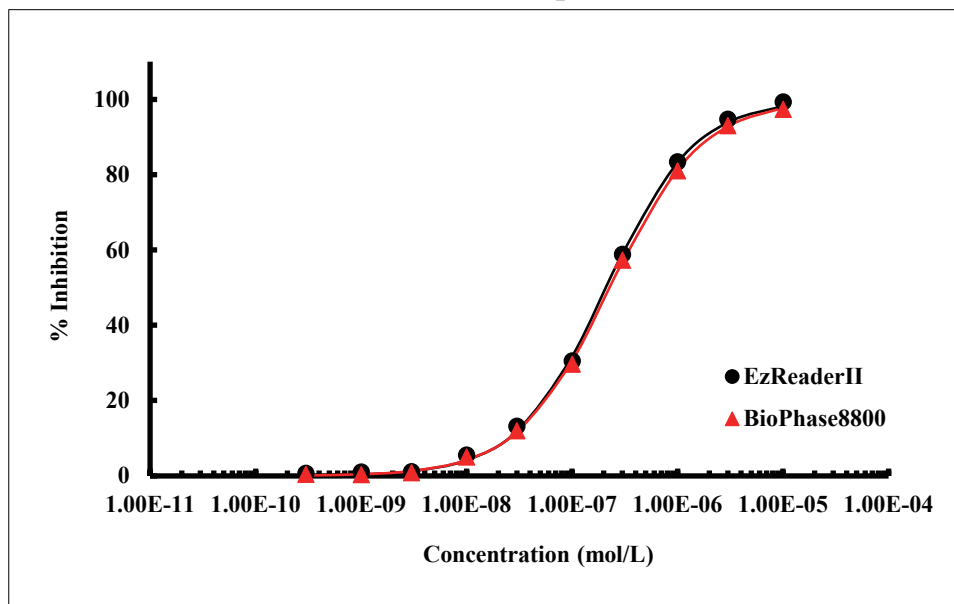


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	380	330



Substrate : EGFR-derived peptide

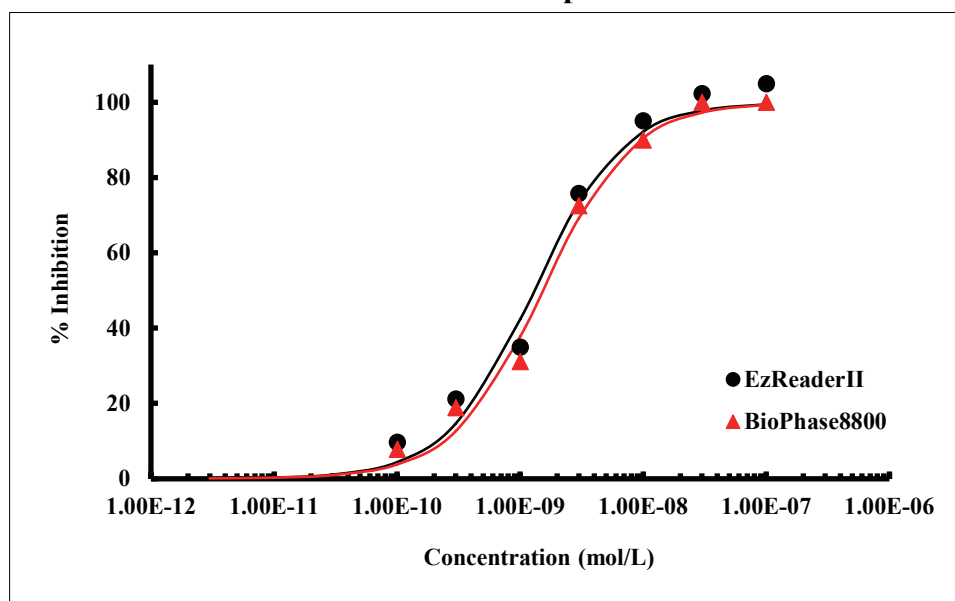
Erk5 vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	210	230

Substrate : GS peptide

PKD1 vs Staurosporine

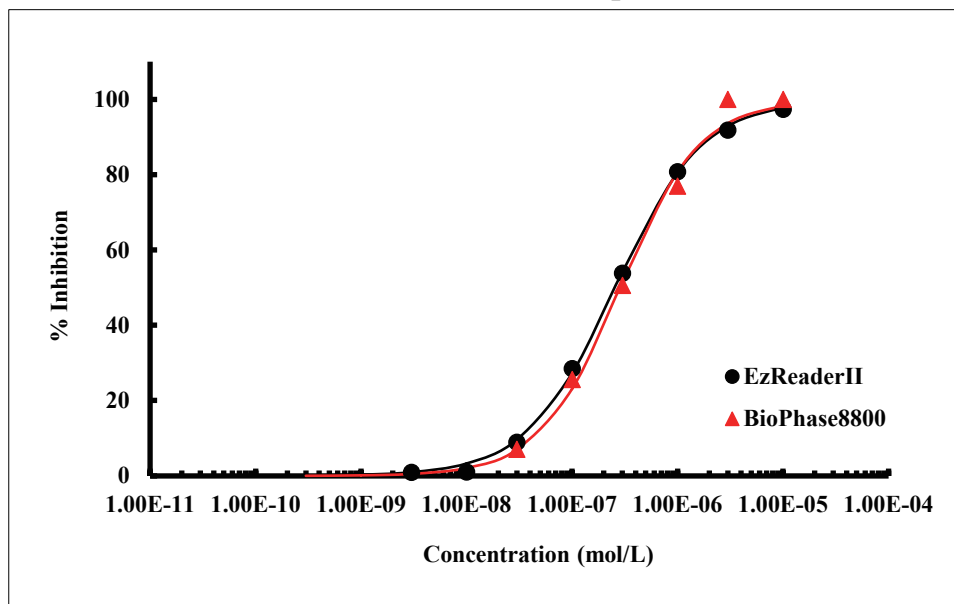


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	1.3	1.5



Substrate : H2A peptide

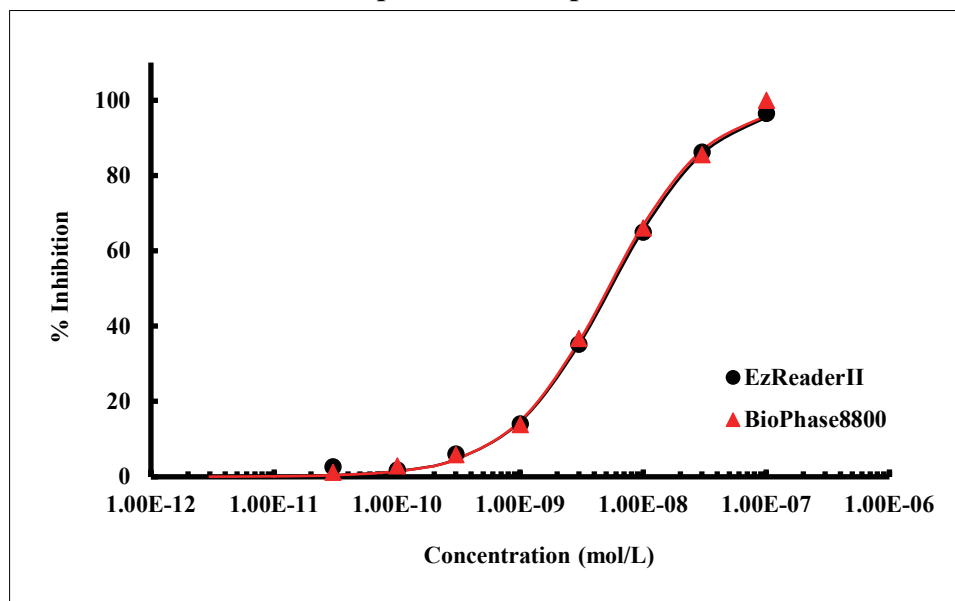
BUB1/BUB3 vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	260	290

Substrate : Histone H3 peptide

Haspin vs Staurosporine

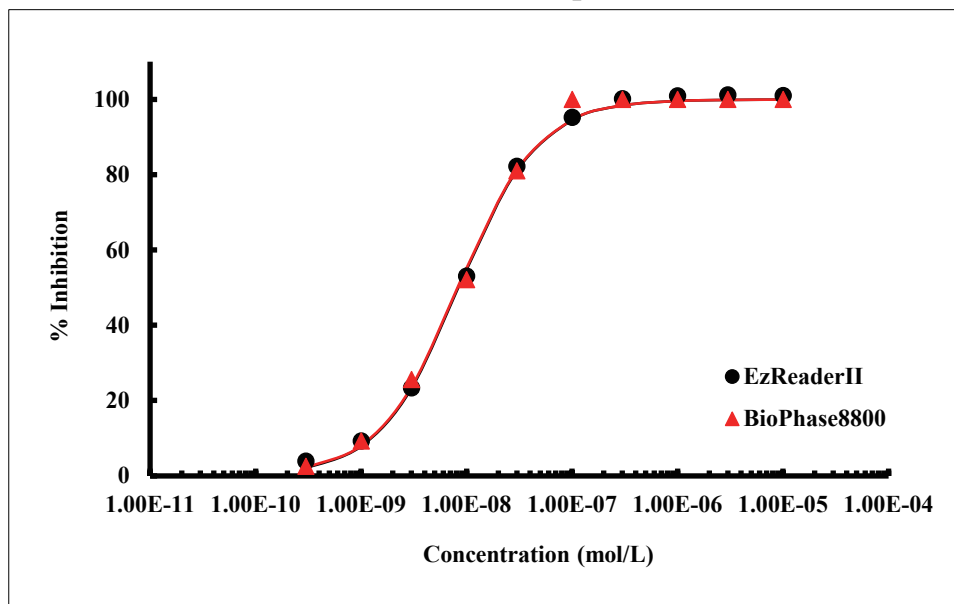


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	5.4	5.2



Substrate : IRAK1 peptide

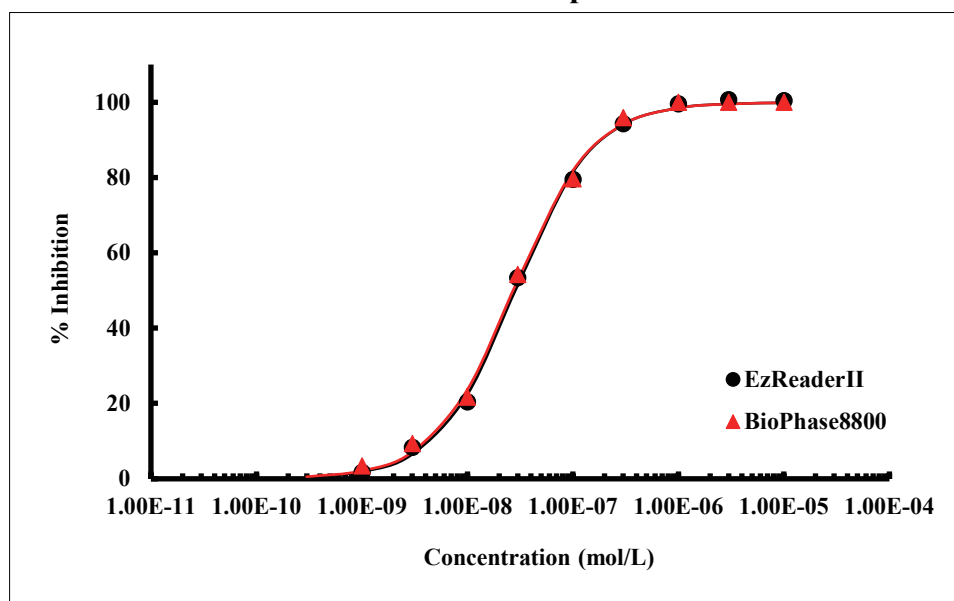
IRAK4 vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	8.5	8.3

Substrate : IRS1

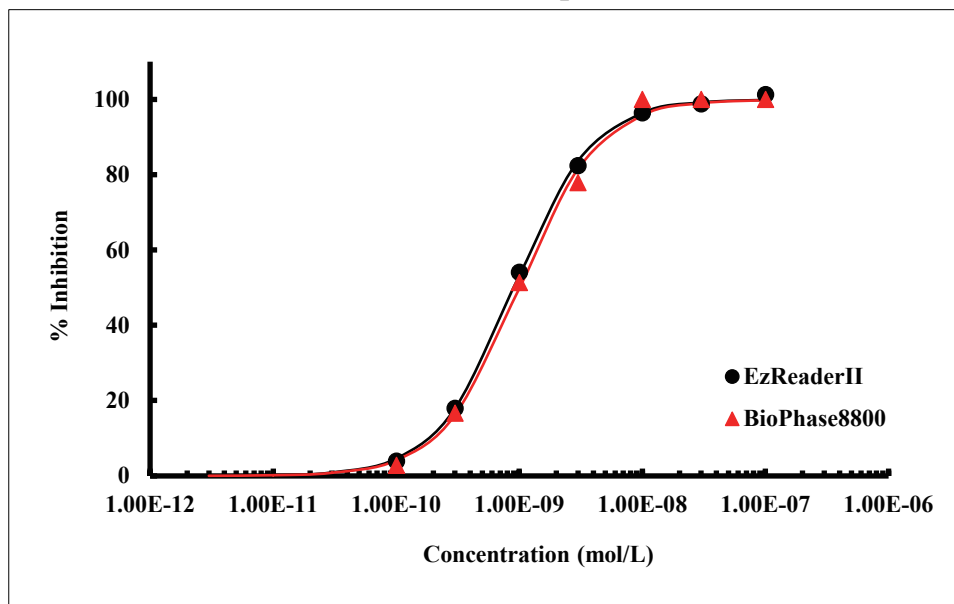
IGF1R vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	29	27

Substrate : **I κ B α peptide**

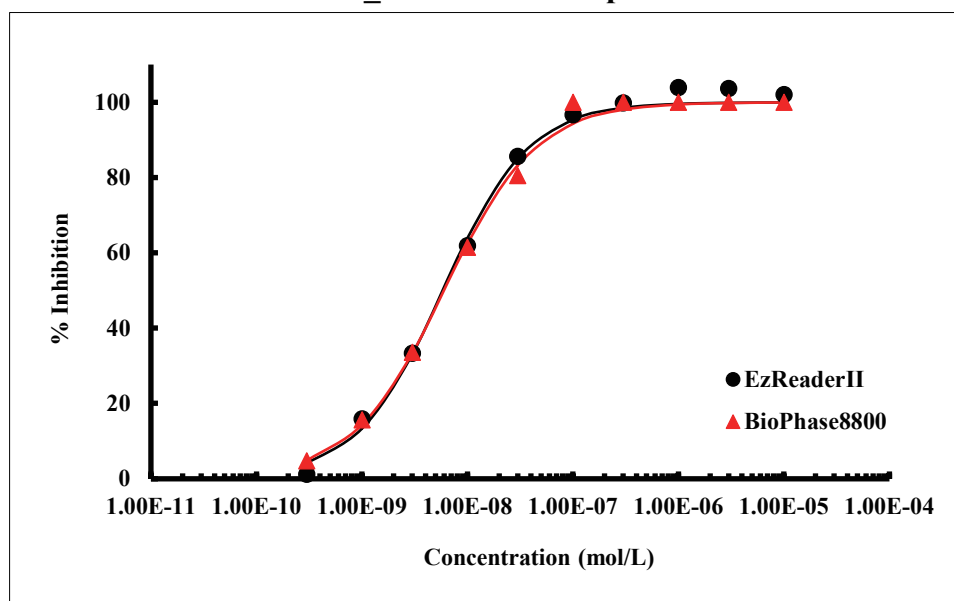
IKK ϵ vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	0.91	1.0

Substrate : **JAK1 substrate peptide**

JAK1_1mM vs Staurosporine

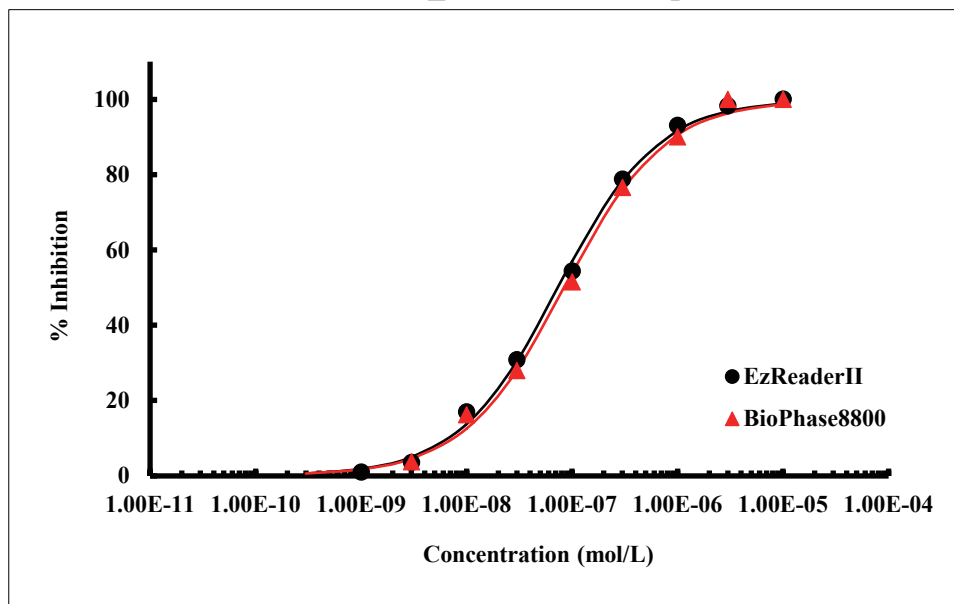


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	5.8	6.0



Substrate : **Kemptide**

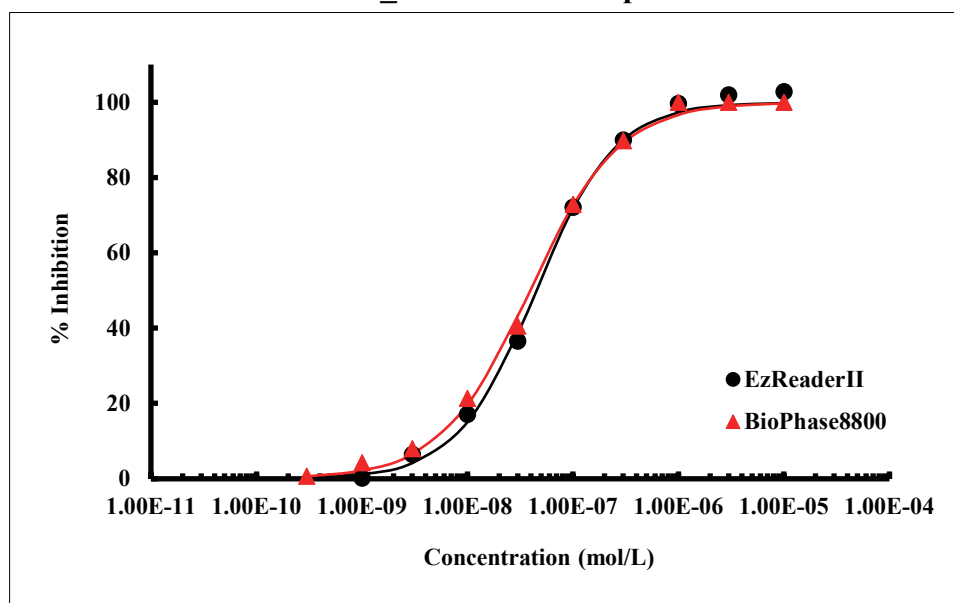
AurB/INCENP_1mM vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	74	85

Substrate : **LIMKtide**

ROCK1_1mM vs Staurosporine

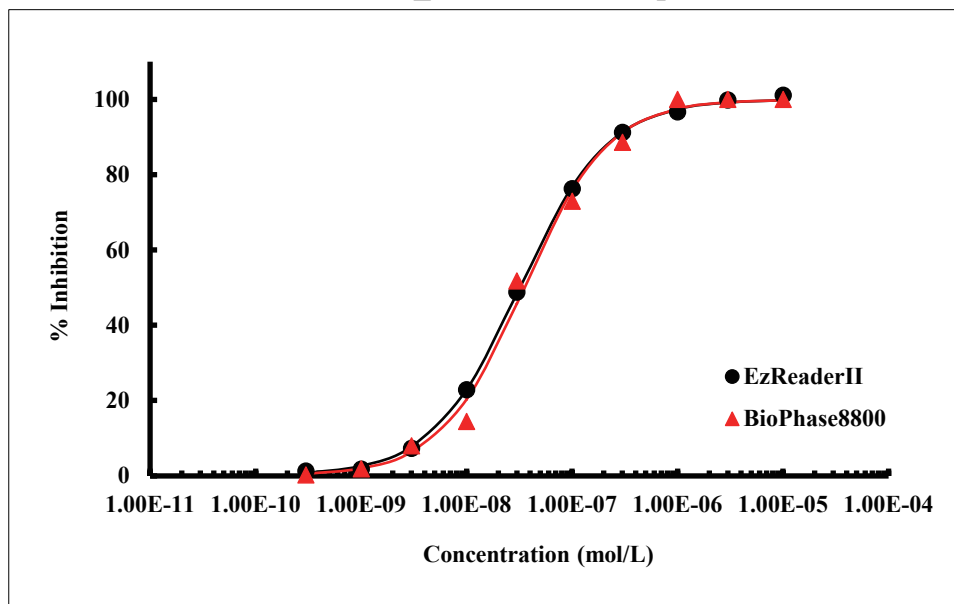


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	45	39



Substrate : LRRKtide

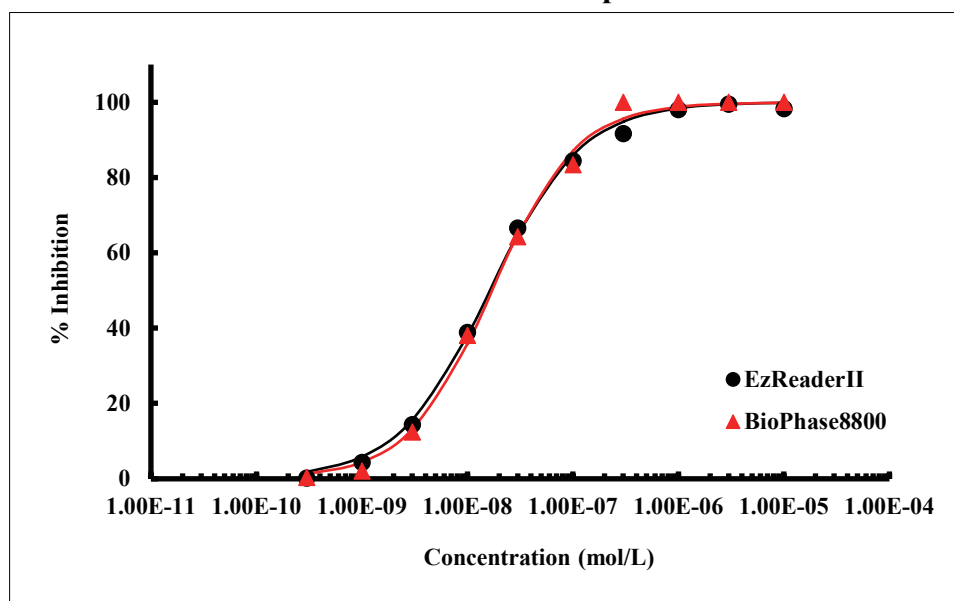
TAK1-TAB1_1mM vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	32	35

Substrate : MCM2 peptide

CDC7/ASK vs Staurosporine

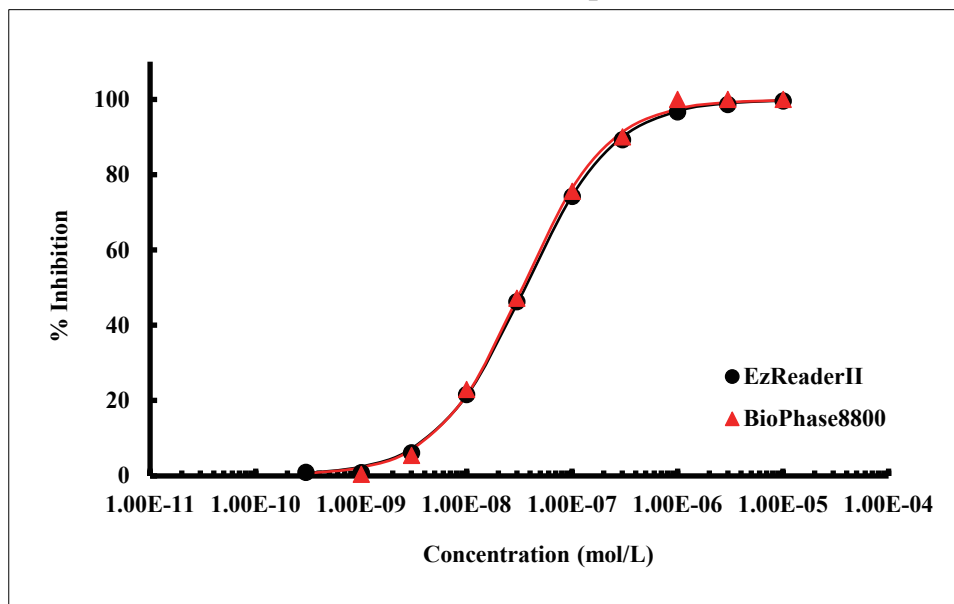


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	16	17



Substrate : **MLCtide**

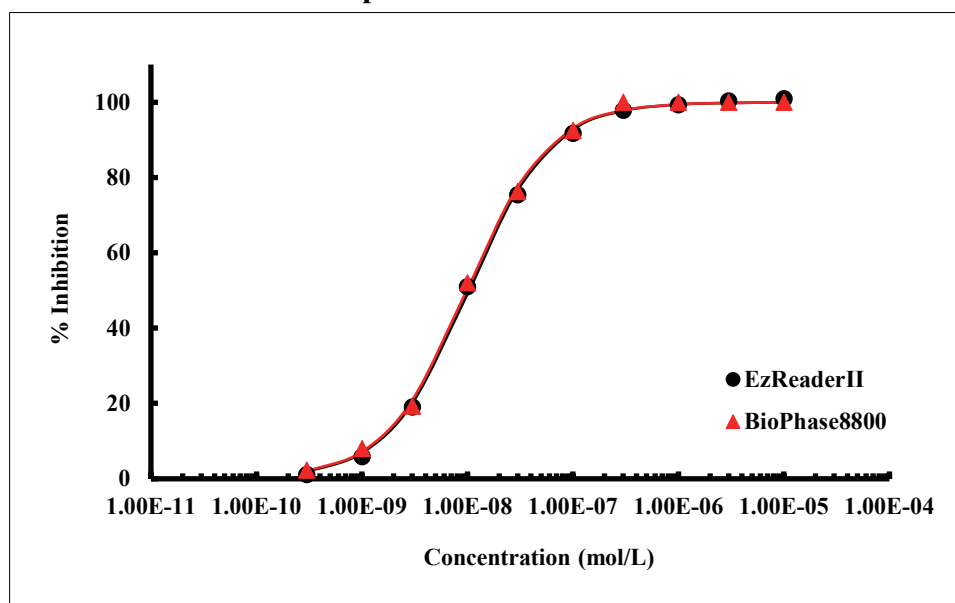
skMLCK vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	36	34

Substrate : **Modified Erktide**

p38α vs SB202190

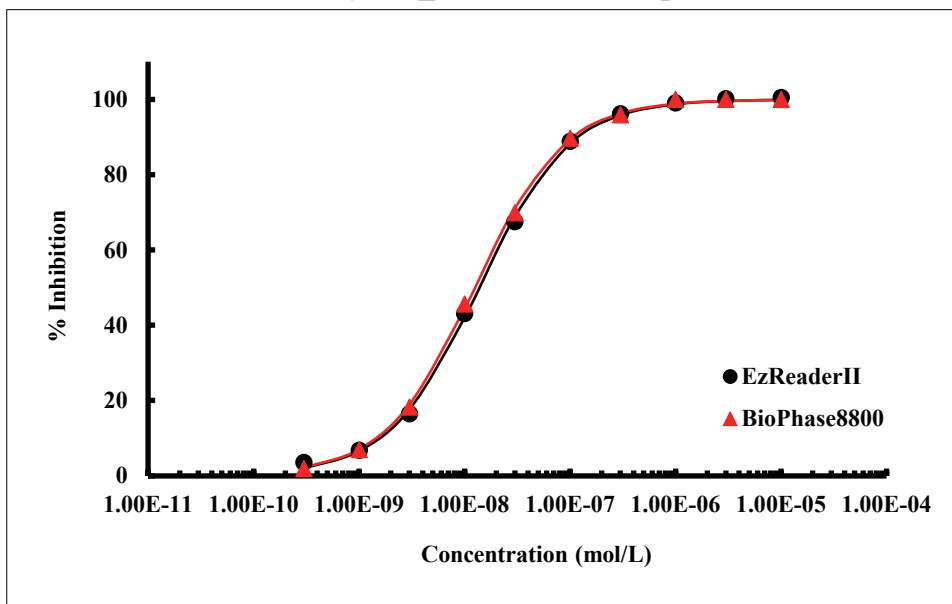


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	10	9.9



Substrate : **Modified Histone H1**

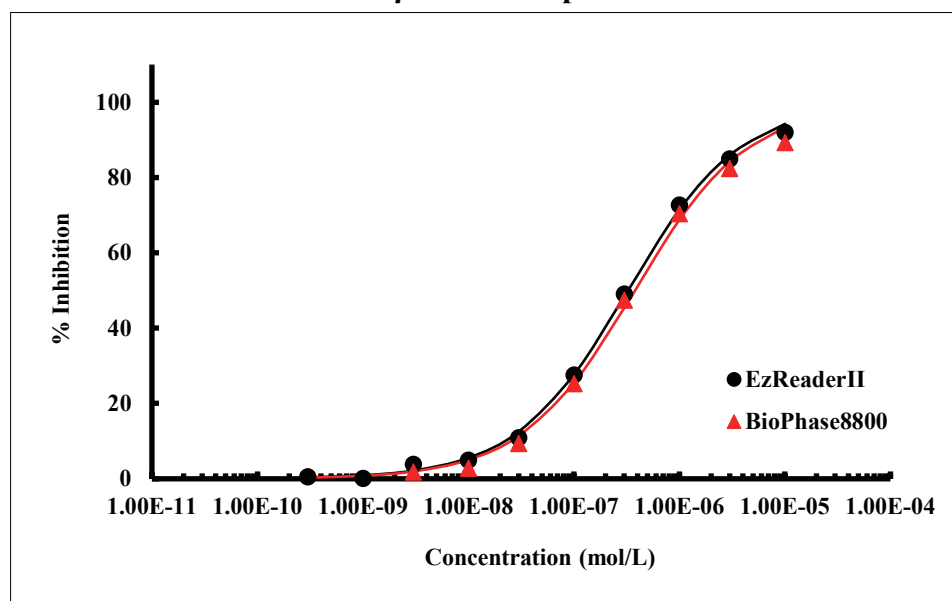
CDK2/CycA2_1mM vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	14	12

Substrate : **Modified IκBα-derived peptide**

IKKβ vs Staurosporine

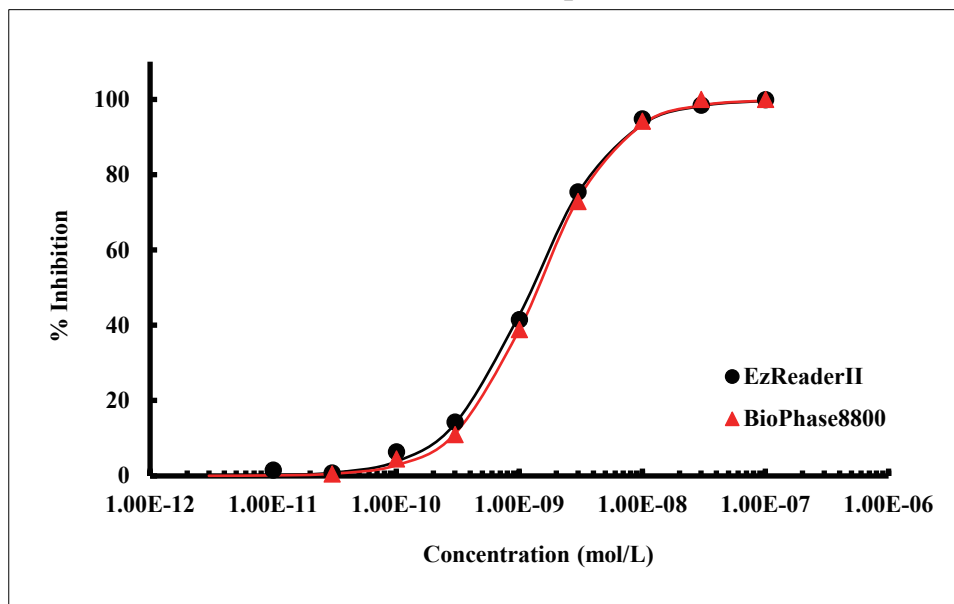


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	320	380



Substrate : Moesin-derived peptide

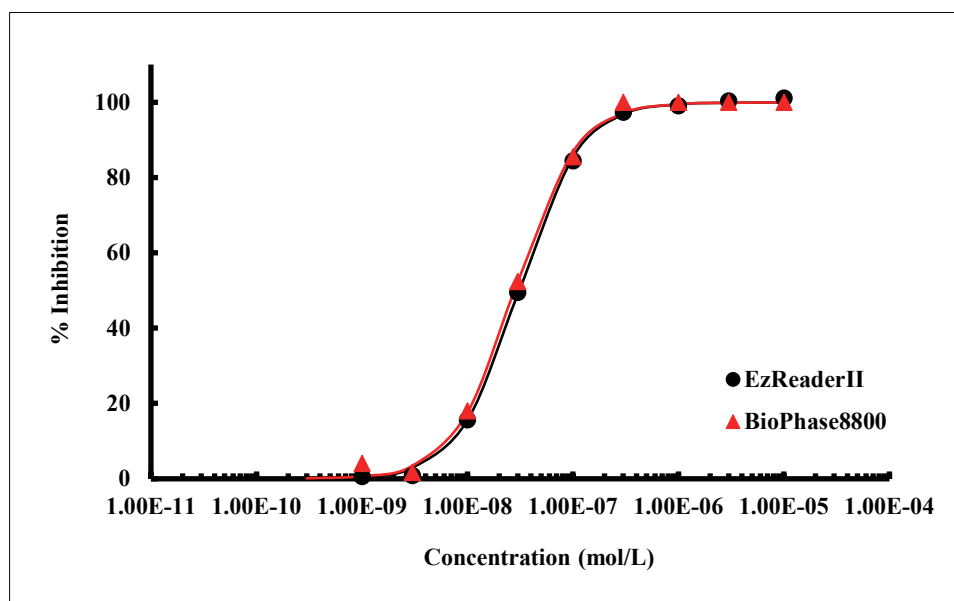
TNIK vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	1.3	1.4

Substrate : PDHKtide

PDHK2 vs VER-246608

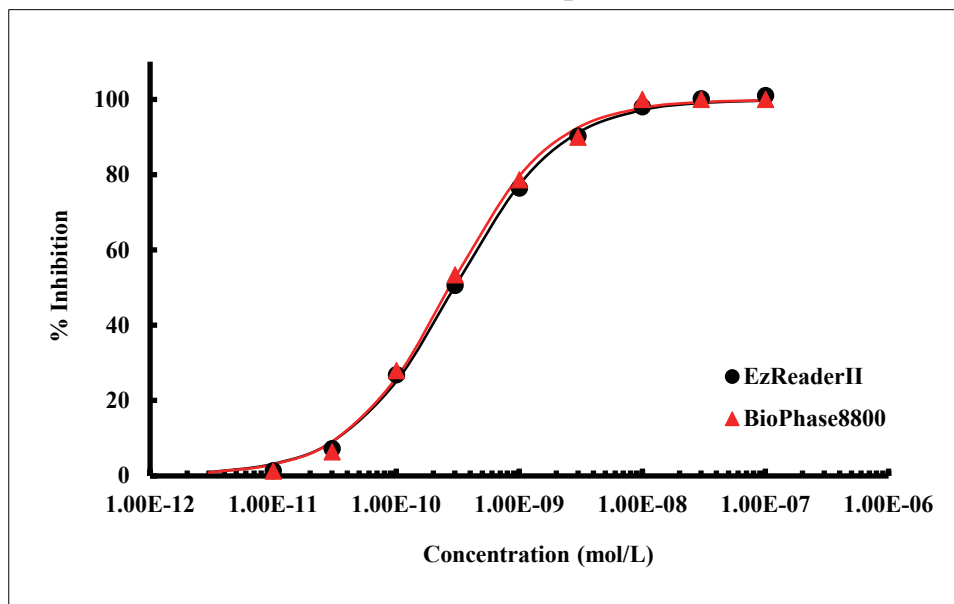


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	31	28



Substrate : PKC peptide

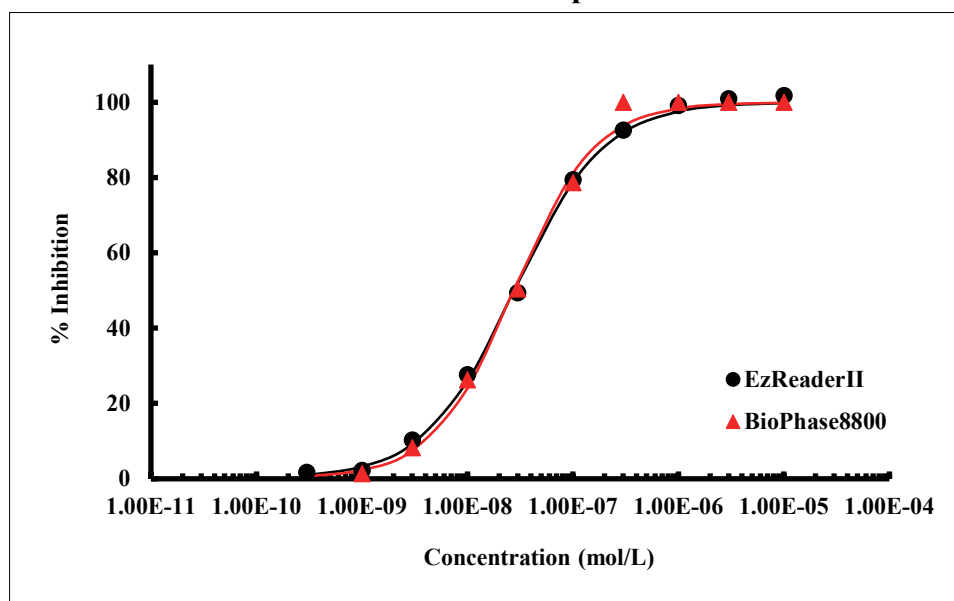
PKC α vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	0.30	0.27

Substrate : RS peptide

MNK1 vs Staurosporine

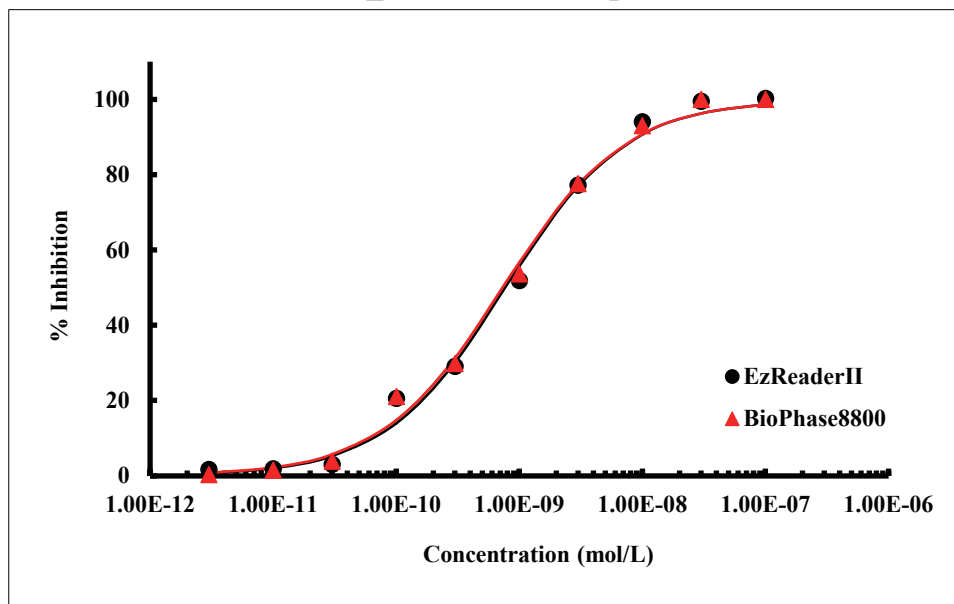


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	28	28



Substrate : S6K peptide(N-FL)

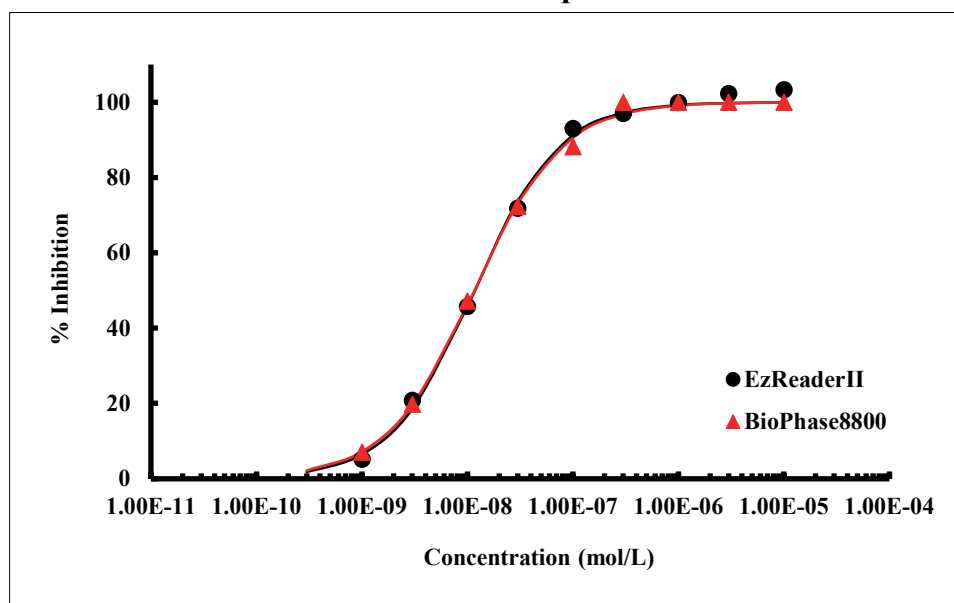
RSK4_1mM vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	0.77	0.73

Substrate : S6K2 peptide

PIM1 vs Staurosporine

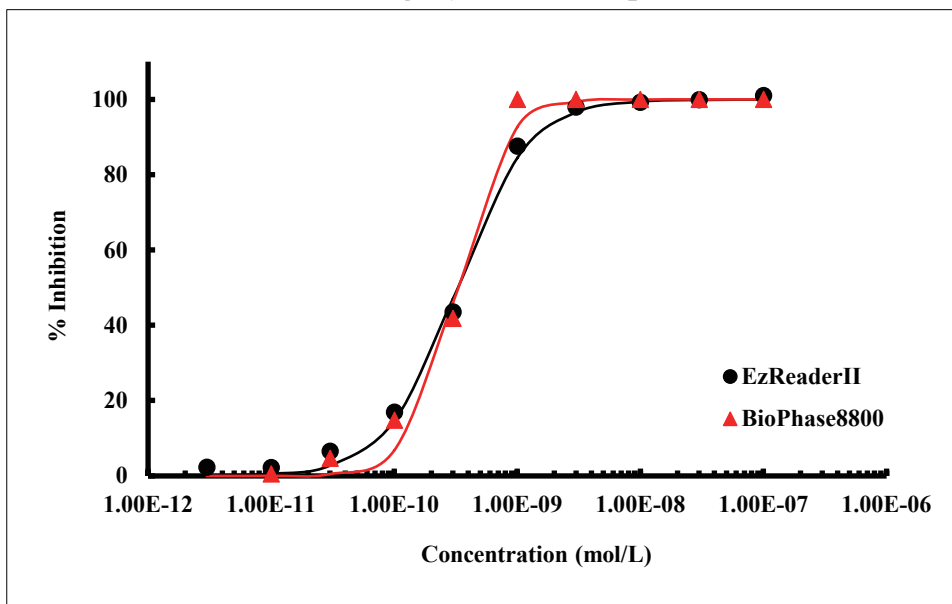


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	12	12



Substrate : SAMS peptide

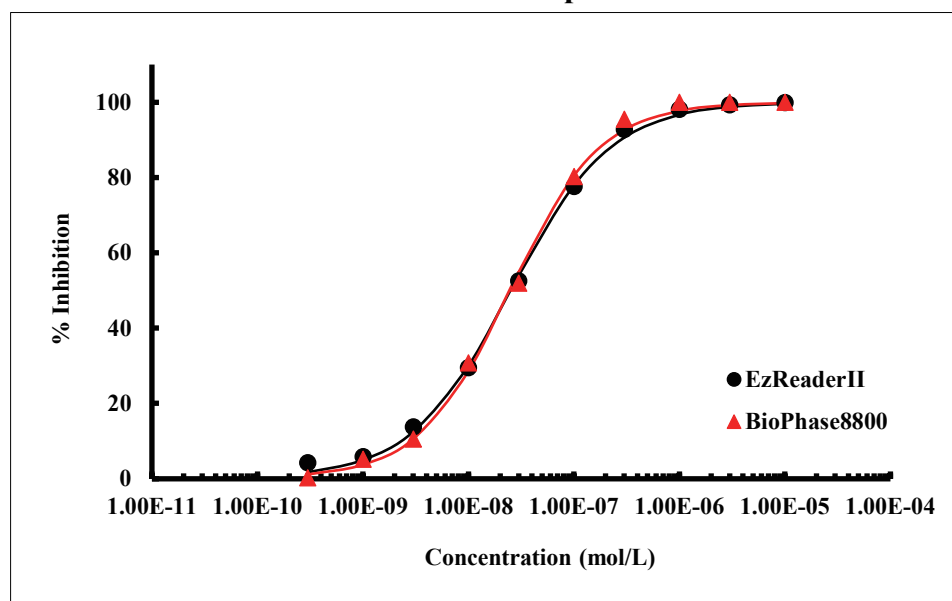
AMPK α 1/ β 1/ γ 1 vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	0.32	0.32

Substrate : SGKtide

SGK2 vs Staurosporine

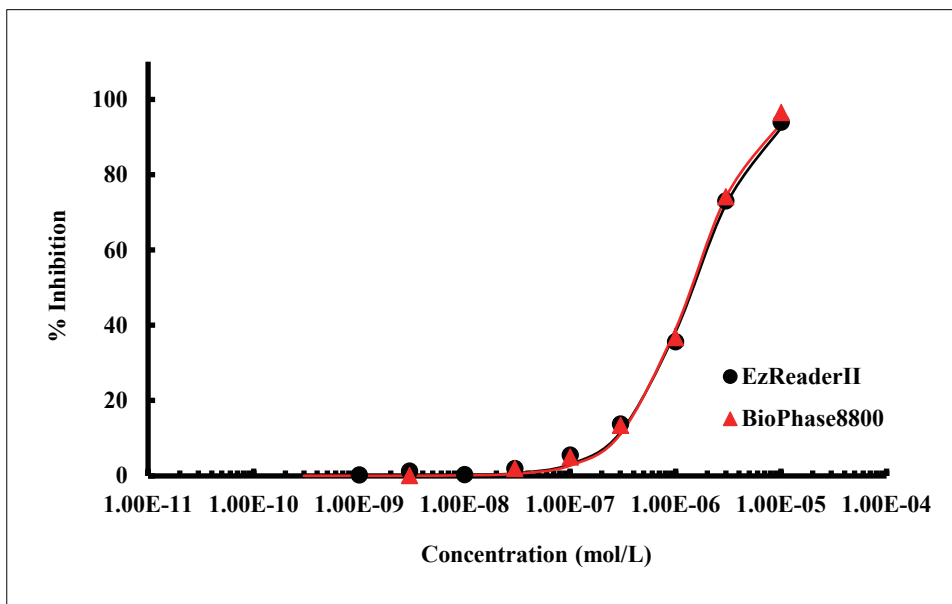


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	25	25



Substrate : **SPAKtide**

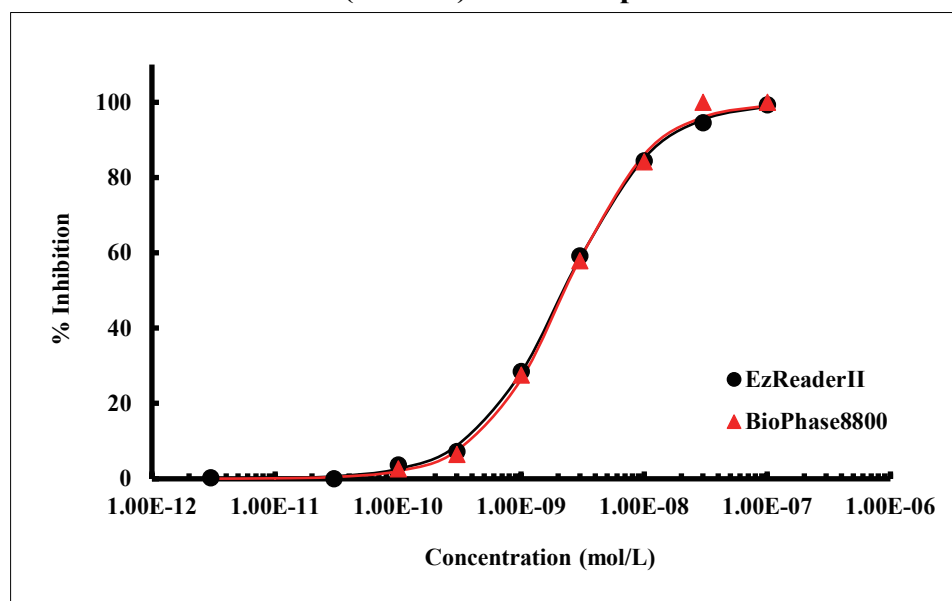
WNK3 vs K252a



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	1400	1400

Substrate : **Srctide**

ALK(F1174L) vs Staurosporine

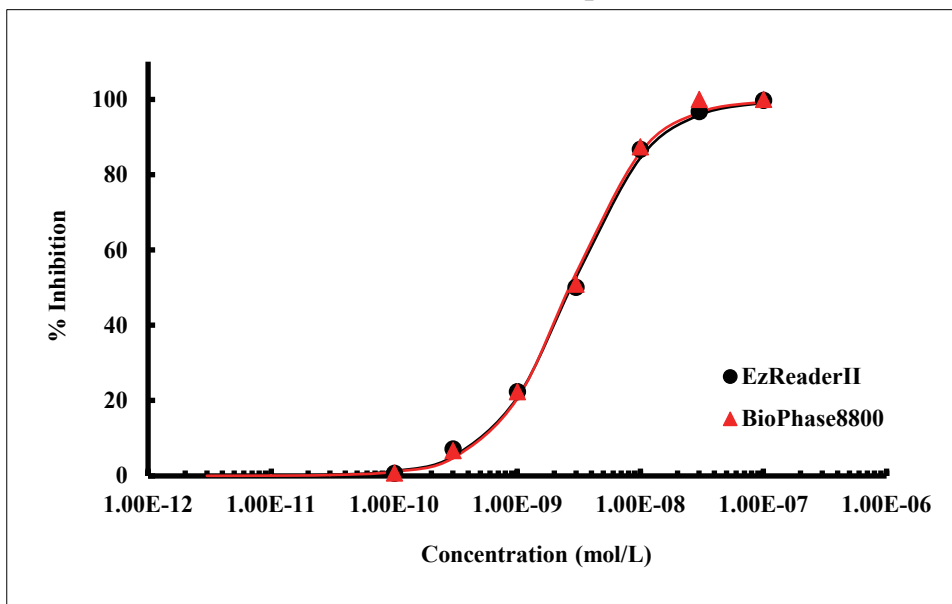


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	2.2	2.3



Substrate : Synapsin peptide

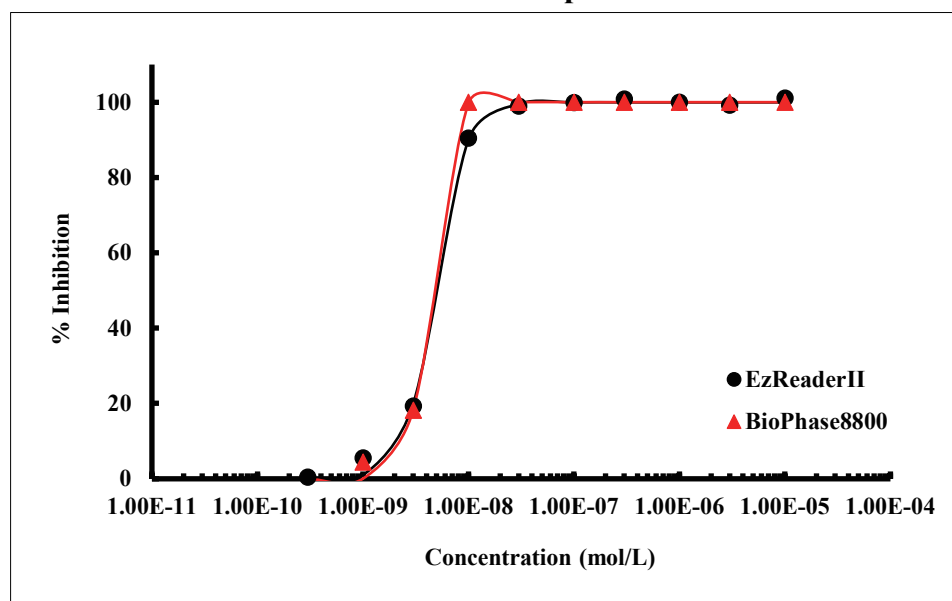
CaMK1δ vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	2.8	2.7

Substrate : T308tide

PDK1 vs Staurosporine

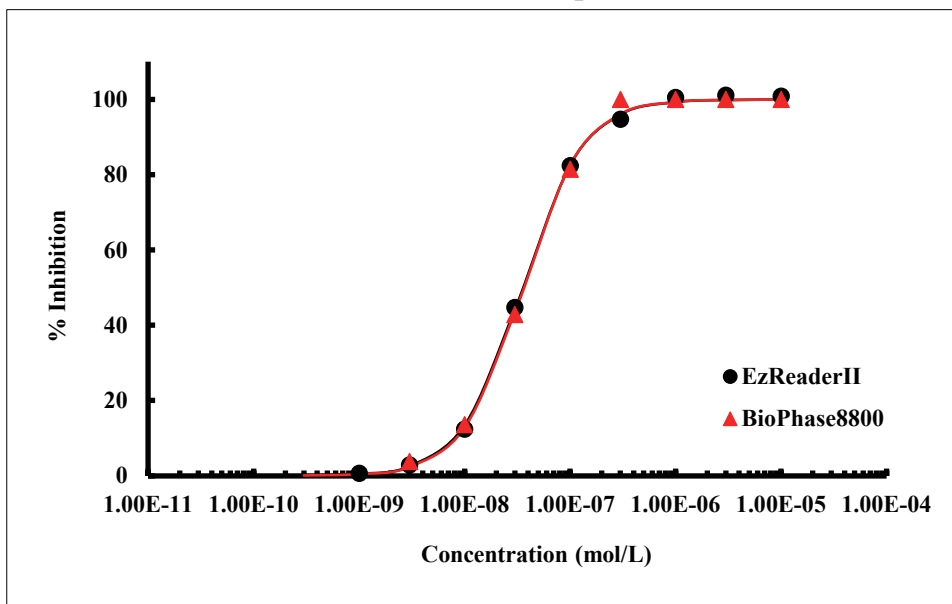


	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	4.8	3.9



Substrate : **TAOKtide**

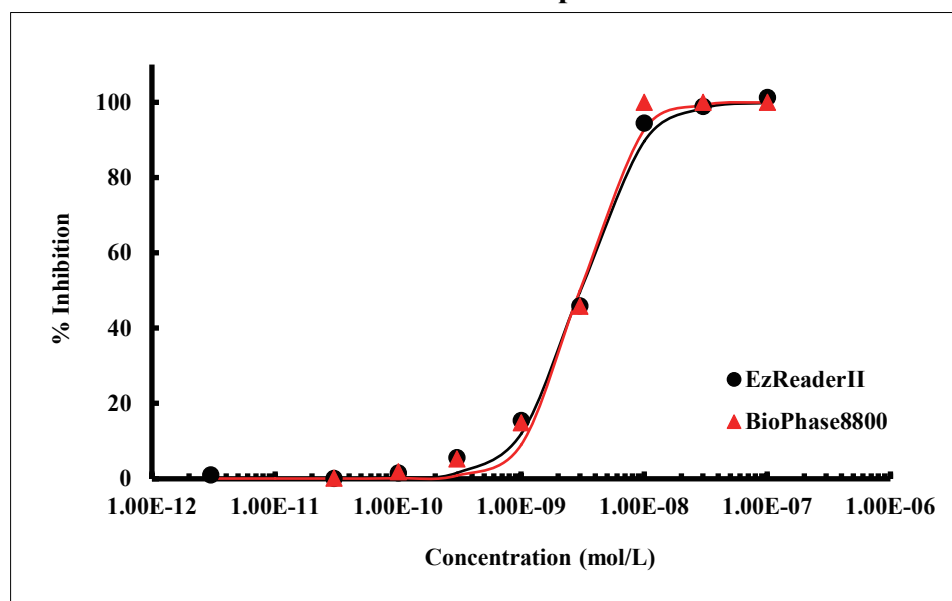
TAOK2 vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	35	36

Substrate : **WASP peptide**

ACK vs Staurosporine



	EzReaderII	BioPhase8800
IC ₅₀ (nmol/L)	3.0	3.0



Kinase-based Substrate List

弊社 Mobility Shift Assay で使用しているキナーゼと基質の対応表です。各基質名をクリックして頂くと新旧システム (BioPhase™ 8800 & EZReader) で取得された該当基質の IC₅₀ データへ移動します。

Kinase	Substrate	Kinase	Substrate
A	ABL	CDK5/p25	Modified Histone H1
	ABL [E255K]	CDK6/CycD3	DYRKtide-F
	ABL [T315I]	CDK7/CycH/MAT1	CTD3 peptide
	ACK	CDK9/CycT1	CDK9 substrate
	AKT1	CGK2	Kemptide
	AKT2	CHK1	CHKtide
	AKT3	CHK2	CHKtide
	ALK	CK1α	CKtide
	ALK [C1156Y]	CK1γ1	CKtide
	ALK [F1174L]	CK1γ2	CKtide
	ALK [G1202R]	CK1γ3	CKtide
	ALK [G1269A]	CK1δ	CKtide
	ALK [L1196M]	CK1ε	CKtide
	ALK [R1275Q]	CK2α1/β	CK2tide
	ALK [T1151_L1152insT]	CK2α2/β	CK2tide
	AMPKα1/β1/γ1	CLK1	DYRKtide-F
	AMPKα2/β1/γ1	CLK2	DYRKtide-F
	ARG	CLK3	DYRKtide-F
	AurA	COT	Modified Erktide
	AurA/TPX2	CRIK	Histone H3 peptide
	AurB/INCENP	CSK	Srctide
	AurC		
	AXL		
B	BLK	D	DAPK1tide
	BMX		GS peptide
	BRAF		IRS1
	BRAF [V600E]		IRS1
	BRK		Modified Erktide
	BRSK1		DYRKtide-F
	BRSK2		DYRKtide-F
	BTK		DYRKtide-F
	BTK [C481S]		DYRKtide-F
	BUB1/BUB3		
C	CaMK1α	E	EEF2Ktide
	CaMK1δ		Srctide
	CaMK2α		Srctide
	CaMK2β		Srctide
	CaMK2γ		Srctide
	CaMK2δ		Srctide
	CaMK4		Srctide
	CDC7/ASK		Srctide
	CDK1(CDC2)/CycB1		Srctide
	CDK2/CycA2		Srctide
	CDK2/CycE1		Srctide
	CDK3/CycE1		Srctide
	CDK4/CycD3		Srctide
			EML4-ALK
			EPHA1
			EPHA2

Kinase-based Substrate List

弊社 Mobility Shift Assay で使用しているキナーゼと基質の対応表です。各基質名をクリックして頂くと新旧システム (BioPhase™ 8800 & EZReader) で取得された該当基質の IC₅₀ データへ移動します。

Kinase	Substrate	Kinase	Substrate
E	EPHA3	H	HIPK2
	EPHA4		HIPK3
	EPHA5		HIPK4
	EPHA6		HPK1
	EPHA7	I	IGF1R
	EPHA8		IKKβ
	EPHB1		IKKε
	EPHB2		INSR
	EPHB3		IRAK4
	EPHB4		IRR
	Erk1		ITK
	Erk2	J	JAK1
	Erk5		JAK2
F	FAK		JAK3
	FER		JNK1
	FES		JNK2
	FGFR1		JNK3
	FGFR1 [V561M]	K	KDR
	FGFR2		KIT
	FGFR2 [V564I]		KIT [D816E]
	FGFR3		KIT [D816V]
	FGFR3 [K650E]		KIT [D816Y]
	FGFR3 [K650M]		KIT [T670I]
	FGFR3 [V555L]		KIT [V560G]
	FGFR3 [V555M]		KIT [V654A]
	FGFR4	L	LATS1/MOBKL1A
	FGFR4 [N535K]		LATS2/MOBKL1A
	FGFR4 [V550E]		LCK
	FGFR4 [V550L]		LOK
	FGR		LTK
	FLT1		LYNa
	FLT3		LYNb
	FLT4	M	MAP2K1
	FMS		MAP2K2
	FRK		MAP2K3
	FYN[isoform a]		MAP2K4
	FYN[isoform b]		MAP2K5
G	GSK3α		MAP2K6
	GSK3β		MAP2K7
H	Haspin		MAP3K1
	HCK		MAP3K2
	HER2		MAP3K3
	HER4		MAP3K4
	HGK		
	HIPK1		

Kinase-based Substrate List

弊社 Mobility Shift Assay で使用しているキナーゼと基質の対応表です。各基質名をクリックして頂くと新旧システム (BioPhase™ 8800 & EZReader) で取得された該当基質の IC₅₀ データへ移動します。

	Kinase	Substrate		Kinase	Substrate
M	MAP3K5	Modified Erktide	P	p38γ	Modified Erktide
	MAP4K2	S6K2 peptide		p38δ	Modified Erktide
	MAPKAPK2	GS peptide		p70S6K	S6K2 peptide
	MAPKAPK3	GS peptide		p70S6Kβ	S6K2 peptide
	MAPKAPK5	GS peptide		PAK1	LIMKtide
	MARK1	CHKtide		PAK2	DAPK1tide
	MARK2	CHKtide		PAK4	SGKtide
	MARK3	CHKtide		PAK5	DAPK1tide
	MARK4	CHKtide		PAK6	SGKtide
	MELK	GS peptide		PASK	GS peptide
	MER	CSKtide		PBK	Histone H3 peptide
	MET	Srctide		PDGFRα	CSKtide
	MET [D1228H]	Srctide		PDGFRα [D842V]	CSKtide
	MET [M1250T]	Srctide		PDGFRα [T674I]	CSKtide
	MET [Y1235D]	Srctide		PDGFRα [V561D]	CSKtide
	MINK	Modified Erktide		PDGFRβ	CSKtide
	MLK1	Modified Erktide		PDHK2	PDHKtide
	MLK2	Modified Erktide		PDHK4	PDHKtide
	MLK3	Modified Erktide		PDK1	T308tide
	MNK1	RS peptide		PGK	Kemptide
	MNK2	RS peptide		PHKG1	GS peptide
	MOS	Modified Erktide		PHKG2	GS peptide
	MRCKα	DAPK1tide		PIM1	S6K2 peptide
	MRCKβ	DAPK1tide		PIM2	S6K2 peptide
	MSK1	Crosstide		PIM3	S6K2 peptide
	MSK2	Crosstide		PKACα	Kemptide
	MSSK1	DYRKtide-F		PKACβ	Kemptide
	MST1	IRS1		PKACγ	Kemptide
	MST2	IRS1		PKCα	PKC peptide
	MST3	Moesin-derived peptide		PKCβ1	PKC peptide
	MST4	Moesin-derived peptide		PKCβ2	PKC peptide
	MUSK	CSKtide		PKCγ	PKC peptide
N	NDR1	SGKtide	PKCδ	PKC peptide	
	NDR2	SGKtide	PKCε	PKC peptide	
	NEK1	CDK7 peptide	PKCζ	PKC peptide	
	NEK2	CDK7 peptide	PKCη	PKC peptide	
	NEK4	GS peptide	PKCθ	PKC peptide	
	NEK6	CDK7 peptide	PKCι	PKC peptide	
	NEK7	CDK7 peptide	PKD1	GS peptide	
	NEK9	CDK7 peptide	PKD2	GS peptide	
	NIM1K	CHKtide	PKD3	GS peptide	
	NPM1-ALK	Srctide	PLK1	CDC25ctide	
	NuaK1	CHKtide	PLK3	CDC25ctide	
	NuaK2	CHKtide	PRKX	Kemptide	
P	p38α	Modified Erktide	Q	QIK	AMARA peptide
	p38β	Modified Erktide			

Kinase-based Substrate List

弊社 Mobility Shift Assay で使用しているキナーゼと基質の対応表です。各基質名をクリックして頂くと新旧システム (BioPhase™ 8800 & EZReader) で取得された該当基質の IC₅₀ データへ移動します。

	Kinase	Substrate		Kinase	Substrate
R	RAF1	Modified Erktide	Y	YES [T348I]	Srctide
	RET	CSKtide			
	RET [G691S]	CSKtide	Z	ZAP70	Blk/Lyntide
	RET [M918T]	CSKtide			
	RET [S891A]	CSKtide			
	RET [Y791F]	CSKtide			
	ROCK1	LIMKtide			
	ROCK2	LIMKtide			
	RON	Srctide			
	ROS	IRS1			
	RSK1	S6K peptide(N-FL)			
	RSK2	S6K peptide(N-FL)			
	RSK3	S6K peptide(N-FL)			
	RSK4	S6K peptide(N-FL)			
S	SGK	SGKtide			
	SGK2	SGKtide			
	SGK3	SGKtide			
	SIK	AMARA peptide			
	skMLCK	MLCtide			
	SLK	Moesin-derived peptide			
	SRC	Srctide			
	SRM	Blk/Lyntide			
	SRPK2	DYRKtide-F			
	SYK	Blk/Lyntide			
T	TAK1-TAB1	LRRKtide			
	TAOK2	TAOKtide			
	TBK1	CKtide			
	TEC	Srctide			
	TIE2	Blk/Lyntide			
	TNIK	Moesin-derived peptide			
	TNK1	CSKtide			
	TRKA	CSKtide			
	TRKB	Srctide			
	TRKC	Srctide			
	TSSK1	GS peptide			
	TSSK2	GS peptide			
	TSSK3	GS peptide			
	TXK	Srctide			
	TYK2	Srctide			
	TYRO3	CSKtide			
W	WNK1	SPAKtide			
	WNK2	SPAKtide			
	WNK3	SPAKtide			
Y	YES	Srctide			