

EXPRESS DIVERSITY KINASE PROFILE

▲ Ref. P10

Suggested testing: 10 μ M (n=2)

📅 5 business days

(Starting on Monday: Compounds must be received at Cerep France site on Friday before noon.)

The **ExpresSDiversity kinase profile** is a fast turnaround profile containing **48 biochemical kinase assays**. The assays were selected taking into consideration the diversity of assay families and the automation level of each assay to maintain the one week turnaround time. As Kinase-inhibitory compounds often interact with a spectrum of kinases, identification of those with improved selectivity is essential. This need can be supported by rapid profiling of your libraries on ExpresSDiversity kinase profile.

→ All assays are part of the **Comprehensive kinase profile**

■ PROTEIN-TYROSINE KINASES

Family/Sub-family	Assay	Ref.
RTK	c-Met kinase	📅 2867
	📅 EGFR kinase	📅 2865
	📅 EphA2 kinase	📅 3055
	📅 EphA3 kinase	📅 3066
	EphB4 kinase	📅 3059
	📅 FGFR1 kinase	📅 2868
	📅 FGFR2 kinase	📅 2893
	FGFR3 kinase	📅 2894
	IRK (InsR)	📅 2898
	KDR kinase (VEGFR2)	📅 2864
	TRKA	📅 2901
CTK	Abl kinase	📅 3056
	📅 JAK3	📅 2905
	Lck kinase	📅 2906
	Src kinase	📅 2907

■ PROTEIN-SERINE/THREONINE KINASES

Family/Sub-family	Assay	Ref.
CMGC	📅 CDC2/CDK1 (cycB)	📅 2875
	📅 CDK2 (cycA)	📅 2908
	ERK ₂ (P42 ^{mapk})	📅 2878
	📅 GSK3 β	📅 2879
	JNK1	📅 2880
	p38 α kinase	📅 2881
CAMK	CaMK2 α	📅 3024
	CHK1	📅 2917
	📅 CHK2	📅 2882
	📅 MAPKAPK2	📅 3367
	📅 MARK1	📅 3362
	MNK2	📅 2918
	📅 Pim2 kinase	📅 2920
AGC	📅 SIK	📅 2921
	📅 Akt1/PKB α	📅 2923
	📅 PDK1	📅 2926
	PKA	📅 2927
	PKC β 2	📅 2750
	📅 ROCK1	📅 3364
📅 SGK1	📅 2929	

▲ Profile changed as compared to 2010 📅 assay developed in 2010 📅 new assay conditions 📅 human 📅 standard turnaround time

Anticipating clinical effects from *in vitro* data

PROTEIN-SERINE/THREONINE KINASES (cont'd)

Family/Sub-family	Assay	Ref.
CK1	CK1 α	1716
STE	HGK (MAP4K4)	3365
	MKK6	2931
	MST4 kinase	3366
	PAK2	2932
	PAK4	3363
TKL	TAOK2 (TAO1)	2586
	IRAK4	2933
	RAF-1 kinase	2936

OTHER KINASES

Family/Sub-family	Assay	Ref.
OTHER KINASES	AurA/Aur2 kinase	2885
	IKK α	2937
	NEK2	2939
	PLK1	2049

SAMPLE SIZE (including IC₅₀ follow-up studies)

Values are based on a molecular weight ≤ 500 g/mol and a typical testing concentration of 10 μ M in duplicate (including a possible retest).

	SCREENING		SCREENING + FOLLOW UP with highest concentration at 10 μ M ¹		SCREENING + FOLLOW UP with highest concentration at 100 μ M ¹	
	WEIGHT (pre-weighed)	VOLUME (100% DMSO)	WEIGHT (pre-weighed)	VOLUME (100% DMSO)	WEIGHT (pre-weighed)	VOLUME (100% DMSO)
EXPRESS DIVERSITY KINASE PROFILE						
Number of assays ▼	48	140 μ L@10 mM	1 mg	190 μ L@10 mM	2.7 mg	540 μ L@10 mM

¹ Assuming ~10% of test in IC₅₀. Usually, for 1 IC₅₀: 30 μ L@10 mM and + 25 μ L@10 mM by additional IC₅₀.

QUESTIONS OR CONCERNS?

Please contact us: sales@cerep.com



FRANCE
Le Bois l'Evêque
86600 CELLE L'EVESCAULT
tel. +33 (0)5 49 89 30 00

USA
15318 N.E. 95th Street
REDMOND, WA 98052
tel. +1 (425) 895 8666

CHINA
326 Aidisheng Road, B 302-1
Zhangjiang High-Tech Park
SHANGHAI 201203
tel. +86 21 5132 0568

sales@cerep.com
www.cerep.com