






NON-KINASE ENZYME PROFILE

Ref. **P7**Suggested testing: 10 μ M (n=2)
 3 weeks

















The Non-kinase enzyme profile includes 42 robust and diverse enzymes broadly chosen among various enzyme families with a particular focus on specific enzymes involved in cell cycle regulation.

Due to its diversity, this profile allows an optimal assessment of enzymatic specificity of drug candidates.

PHOSPHATASES

Sub-family	Enzyme	Ref.
	phosphatase 1B (PTP1B)	 2593
	phosphatase 2A (PP2A)	 2584
	 phosphatase CDC25A	 2561
	phosphatase MKP1	 2652

PROTEASES

Sub-family	Enzyme	Ref.
SERINE PROTEASES	elastase	 0183
	trypsin	 0512
CYSTEINE PROTEASES	caspase-3	 0774
	caspase-8	 0668
	cathepsin L	 0802
ASPARTIC PROTEASES	BACE-1 (β -secretase)	 0701
	cathepsin D	 0650
	HIV-1 protease	 0346
METALLOPROTEASES	ACE	 0979
	ACE-2	 1954
	ECE-1	 0184
	MMP-1	 0510
	MMP-3	 0490
	MMP-9	 0491
	neutral endopeptidase	 0509
	TACE	 0702









PHOSPHODIESTERASES







Sub-family	Enzyme	Ref.
	PDE2A	 2426
	PDE3A	 2432
	PDE4D	 2434
	 PDE5 (non-selective)	 0204
	PDE6 (non-selective)	0478

NO SYNTHASES

Sub-family	Enzyme	Ref.
	inducible NOS	0196

ARACHIDONIC ACID METABOLISM

Sub-family	Enzyme	Ref.
PHOSPHOLIPASE	 PLA ₂	2999
CYCLOOXYGENASES	COX ₁	 0726
	COX ₂	 0727
LIPOXYGENASE	 5-Lipoxygenase 	 0772
	 12-lipoxygenase	 0883

 profile changed in 2010  Assay added in 2010  new assay  new protocol  human  standard turnaround time

MONOAMINE & NEUROTRANSMITTER SYNTHESIS & METABOLISM

Sub-family	Enzyme	Ref.
	acetylcholinesterase	0363
	MAO-A	0191
	MAO-B	0621

MISCELLANEOUS ENZYMES

Sub-family	Enzyme	Ref.
	CENP-E	2150
	Eg5	2151
	HDAC3	2083
	HDAC4	2493
	HDAC6	2495
	HDAC11	2663
	sirtuin 1	2581
	sirtuin 2	2582

SAMPLE SIZE (including IC₅₀ follow-up studies)

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

	SCREENING		SCREENING + FOLLOW UP with highest concentration at 10 μM ¹		SCREENING + FOLLOW UP with highest concentration at 100 μM ¹	
	WEIGHT (pre-weighed)	VOLUME (100% DMSO)	WEIGHT (pre-weighed)	VOLUME (100% DMSO)	WEIGHT (pre-weighed)	VOLUME (100% DMSO)
NON-KINASE ENZYME PROFILE						
Number of assays ▼	42	1 mg	185 μL@10 mM	1.5 mg	300 μL@10 mM	3 mg

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

ASSAY CATALOG REFERENCES

In 2009, Cerep finalized one of the milestones in the industrialization process: the implementation of new referencing and supply chain management systems. The references of each of the assays have thus been simplified: **each reference will now be displayed as 4 digits.**

A correlation table between old and new assay references is available at [www.cerep.com CATALOG ONLINE](http://www.cerep.com/CATALOG_ONLINE)
<<http://www.cerep.com/Cerep/Users/pages/Catalog/Assay/catalog.asp>>

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

(Headquarters)

155 boulevard Haussmann
75008 PARIS
tel. +33 (0)1 45 64 44 60

USA

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

JAPAN

Namiki Shoji Co., Ltd.
Kenseishinjuku Bldg. 5-5-3
Shinjuku, Shinjuku-ku
TOKYO, 160-0022
tel. +81 (0)3 3354 4026
fax +81 (0)3 3352 2196

CHINA

Ai Di Sheng (Edison) Road 326,
302-1 room
Zhangjiang High-Tech Park
SHANGHAI
tel. +86 18702160370