

PDE HIGH-THROUGHPUT PROFILE

▲ Ref. P23

Suggested testing: 10 μ M (n=2)

🕒 3 weeks

The PDE High-throughput profile is designed to determine the inhibitory activity and selectivity of compounds on the PDE superfamily. The robustness, reproducibility and relevance of this 13 enzyme assay profile were validated by determining the potency and selectivity of a broad panel of commercially available reference inhibitors and known clinical drugs in three independent experiments (poster available on the website).

→ For Cellular phosphodiesterase assays, please contact us at customresarch@cerep.com

■ ENZYMES

Family	Assay	Ref.
PHOSPHODIESTERASES	📄 PDE1B	🇫🇷 2431
	PDE2A ₁	🇫🇷 2426
	PDE3A	🇫🇷 2432
	PDE3B	🇫🇷 2705
	PDE4A _{1A}	🇫🇷 2342
	PDE4B ₁	🇫🇷 2413
	📄 PDE4D ₂	🇫🇷 2434
	PDE5 (non-selective)	🇫🇷 0204
	PDE6 (non-selective)	0478
	PDE7A	🇫🇷 2351
	PDE8A ₁	🇫🇷 2355
	PDE10A ₁	🇫🇷 2357
	PDE11A ₄	🇫🇷 2358

📄 new profile 📄 new assay conditions 🇫🇷 human 🕒 standard turnaround time

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)
PDE HIGH-THROUGHPUT PROFILE						
Number of assays ▼						
13	1 mg	200 μ L@10 mM	1.4 mg	275 μ L@10 mM	inquire	inquire

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.

Anticipating clinical effects from *in vitro* data

■ 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