

PDE1A Phosphodiesterase Assay Service

PDE1A [Phosphodiesterase Assay](#) for compound screening and profiling via fluorescence polarization-based quantification of enzyme activity.

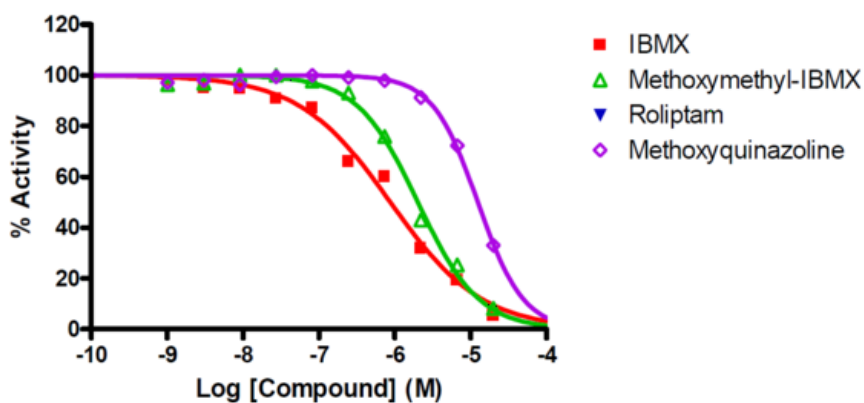
Target Overview

CAT#	PDE1A
Enzyme	Human PDE1A
Synonym	Calcium/calmodulin-dependent 3',5'-cyclic nucleotide phosphodiesterase 1A, 61 kDa Cam-PDE, hCam-1)
Substrate	1 μ M cGMP, Ca-CaM

Assay Properties

Reaction	Nucleoside 3',5'-cyclic phosphate + H ₂ O = nucleoside 5'-phosphate	
Readout	Enzymatic activity	
Reference Compound IC50s	Compounds	IC50 (nM)
	IBMX	880
	Methoxymethyl-IBMX	2,030
	Roliptam	>50,000
	Methoxyquinazoline	12,500
Scientific Information	P54750 (http://www.uniprot.org/uniprot/P54750)	
Screening Location	Malvern, USA	

Reference Compound IC50 for PDE1A



Let's Discover Together.

© Oct 2022 **Reaction Biology**. All rights reserved.

https://www.reactionbiology.com/datasheet/pde1a_pde_malvern

