

PDE1B Phosphodiesterase Assay Service

PDE1B [Phosphodiesterase Assay](#) for compound screening and profiling via fluorescence polarization-based quantification of enzyme activity.

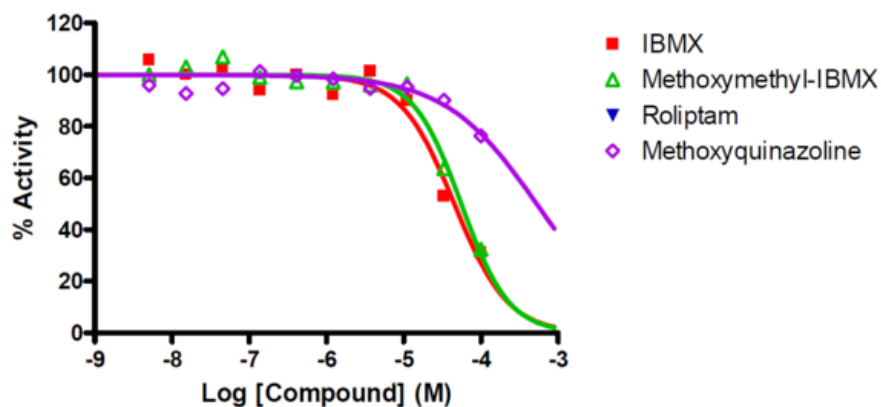
Target Overview

CAT#	PDE1B
Enzyme	Human PDE1B
Synonym	Calcium/calmodulin-dependent 3',5'-cyclic nucleotide phosphodiesterase 1B,63 kDa Cam-PDE, PDE1B1, PDES1B)
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	45,000
	Methoxymethyl-IBMX	56,000
	Roliptam	>100,000
	Methoxyquinazoline	>100,000
Scientific Information	Q01064 (http://www.uniprot.org/uniprot/Q01064)	
Screening Location	Malvern, USA	

Reference Compound IC50 for PDE1B



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/pde1b_pde_malvern

