

PDE4B Phosphodiesterase Assay Service

PDE4B [Phosphodiesterase Assay](#) for compound screening and profiling via fluorescence polarization-based quantification of enzyme activity.

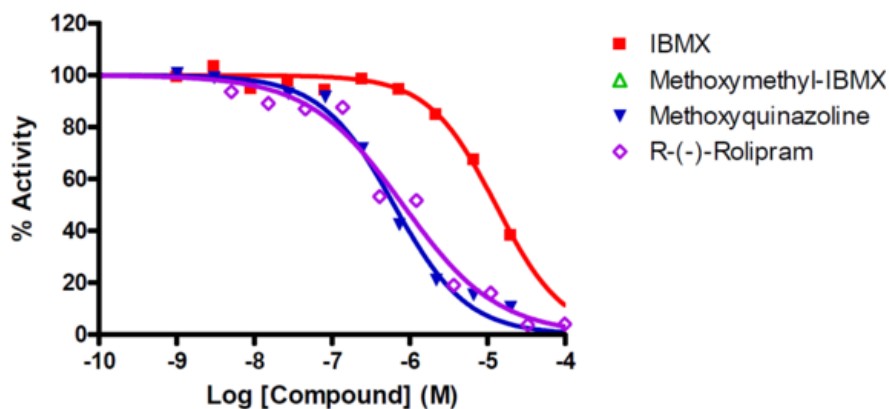
Target Overview

CAT#	PDE4B
Enzyme	Human PDE4B
Synonym	cAMP-specific 3',5'-cyclic phosphodiesterase 4B, DPDE4, PDE32
Substrate	1 μ M cAMP

Assay Properties

Reaction	Nucleoside 3',5'-cyclic phosphate + H ₂ O = nucleoside 5'-phosphate	
Readout	Enzymatic activity	
Reference Compound IC50s	Compounds	IC50 (nM)
	IBMX	12,900
	Methoxymethyl-IBMX	>50,000
	Roliptam	630
	Methoxyquinazoline	830
Scientific Information	Q07343 (http://www.uniprot.org/uniprot/Q07343)	
Screening Location	Malvern, USA	

Reference Compound IC50 for PDE4B



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/pde4b_pde_malvern

