

PDE7A Phosphodiesterase Assay Service

PDE7A [Phosphodiesterase Assay](#) for compound screening and profiling via fluorescence polarization-based quantification of enzyme activity.

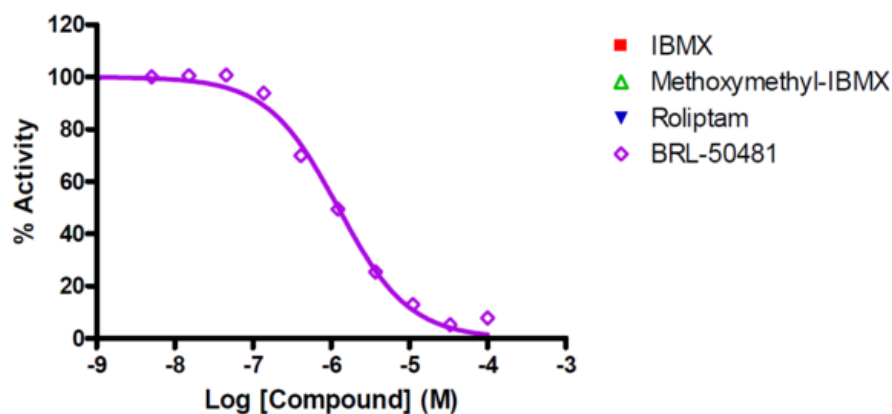
Target Overview

CAT#	PDE7A
Enzyme	Human PDE7A
Synonym	High affinity cAMP-specific 3',5'-cyclic phosphodiesterase 7A, HCP1, TM22
Substrate	0.2 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	>100,000
	Methoxymethyl-IBMX	>100,000
	Roliptam	>100,000
	BRL-50481	1,200
Scientific Information	Q13946 (http://www.uniprot.org/uniprot/Q13946)	
Screening Location	Malvern, USA	

Reference Compound IC50 for PDE7A



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/pde7a_pde_malvern

