

PDE3A Phosphodiesterase Assay Service

PDE3A [Phosphodiesterase Assay](#) for compound screening and profiling via fluorescence polarization-based quantification of enzyme activity.

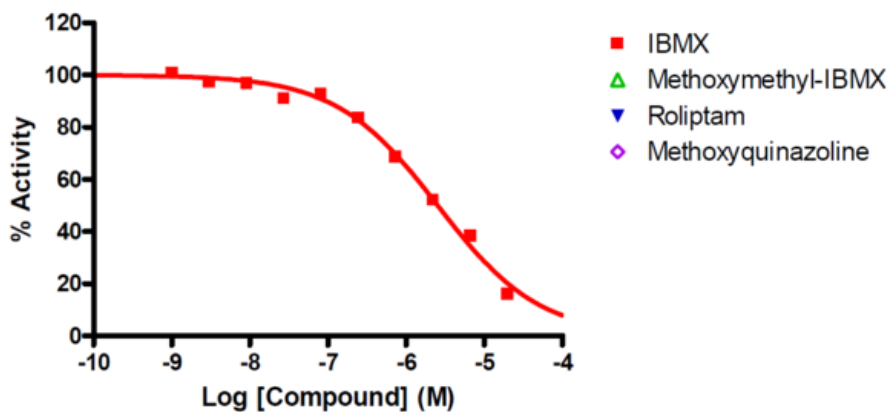
Target Overview

CAT#	PDE3A
Enzyme	Human PDE3A
Synonym	cGMP-inhibited 3',5'-cyclic phosphodiesterase A, CGI-PDE A
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	2,500
	Methoxymethyl-IBMX	>50,000
	Roliptam	>50,000
	Methoxyquinazoline	>50,000
Scientific Information	Q14432 (http://www.uniprot.org/uniprot/Q14432)	
Screening Location	Malvern, USA	

Reference Compound IC50 for PDE3A



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/pde3a_pde_malvern

