

# PDE9A Phosphodiesterase Assay Service

PDE9A [Phosphodiesterase Assay](#) for compound screening and profiling via fluorescence polarization-based quantification of enzyme activity.

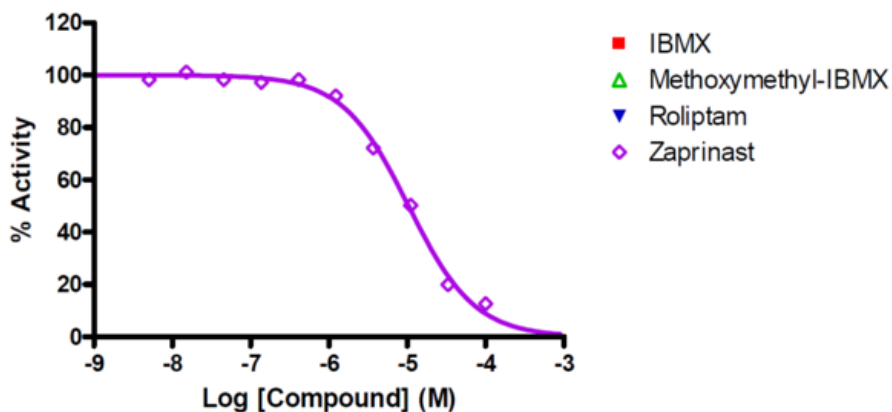
## Target Overview

CAT#	PDE9A
Enzyme	Human PDE9A
Synonym	High affinity cGMP-specific 3',5'-cyclic phosphodiesterase 9A
Substrate	0.2 $\mu$ M cGMP

## Assay Properties

Reaction	Nucleoside 3',5'-cyclic phosphate + H <sub>2</sub> O = nucleoside 5'-phosphate	
Readout	Enzymatic activity	
Reference Compound IC50s	<b>Compounds</b>	<b>IC50 (nM)</b>
	IBMX	>100,000
	Methoxymethyl-IBMX	>100,000
	Roliptam	>100,000
	Zaprinast	10,500
Scientific Information	<a href="http://www.uniprot.org/uniprot/O76083">O76083</a> ( <a href="http://www.uniprot.org/uniprot/O76083">http://www.uniprot.org/uniprot/O76083</a> )	
Screening Location	Malvern, USA	

## Reference Compound IC50 for PDE9A



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/pde9a\\_pde\\_malvern](https://www.reactionbiology.com/datasheet/pde9a_pde_malvern)

