

# PDE5A Phosphodiesterase Assay Service

PDE5A [Phosphodiesterase Assay](#) for compound screening and profiling via fluorescence polarization-based quantification of enzyme activity.

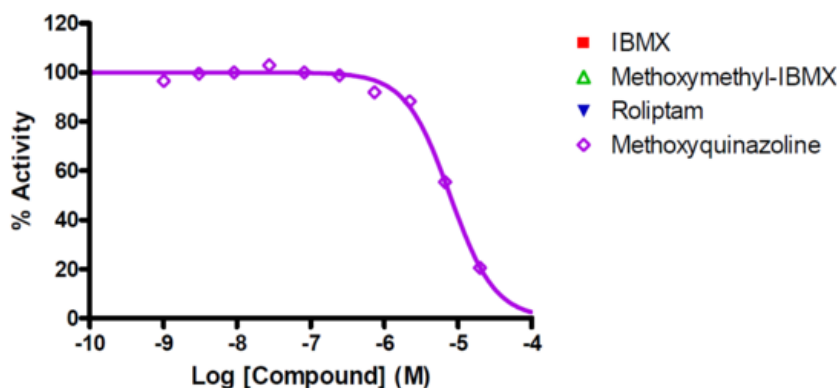
## Target Overview

CAT#	PDE5A
Enzyme	Human PDE5A
Synonym	cGMP-specific 3',5'-cyclic phosphodiesterase, CGB-PDE)
Substrate	1 $\mu$ 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	>50,000
	Methoxymethyl-IBMX	>50,000
	Roliptam	>50,000
	Methoxyquinazoline	7,900
Scientific Information	<a href="http://www.uniprot.org/uniprot/076074">076074</a> ( <a href="http://www.uniprot.org/uniprot/076074">http://www.uniprot.org/uniprot/076074</a> )	
Screening Location	Malvern, USA	

## Reference Compound IC50 for PDE5A



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/pde5a\\_pde\\_malvern](https://www.reactionbiology.com/datasheet/pde5a_pde_malvern)

