

# PDE2A Phosphodiesterase Assay Service

PDE2A [Phosphodiesterase Assay](#) for compound screening and profiling via fluorescence polarization-based quantification of enzyme activity.

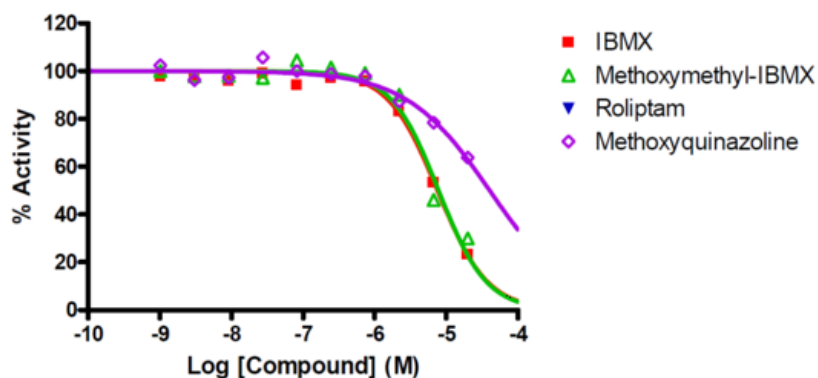
## Target Overview

<b>CAT#</b>	PDE2A
<b>Enzyme</b>	Human PDE2A
<b>Synonym</b>	cGMP-dependent 3',5'-cyclic phosphodiesterase, Cyclic GMP-stimulated phosphodiesterase, cGS-PDE)
<b>Substrate</b>	1 $\mu$ M cGMP

## Assay Properties

<b>Reaction</b>	Nucleoside 3',5'-cyclic phosphate + H <sub>2</sub> O = nucleoside 5'-phosphate	
<b>Readout</b>	Enzymatic activity	
<b>Reference Compound IC50s</b>	<b>Compounds</b>	<b>IC50 (nM)</b>
	IBMX	7,600
	Methoxymethyl-IBMX	7,900
	Roliptam	>50,000
	Methoxyquinazoline	>41,000
<b>Scientific Information</b>	<a href="http://www.uniprot.org/uniprot/O00408">O00408</a> ( <a href="http://www.uniprot.org/uniprot/O00408">http://www.uniprot.org/uniprot/O00408</a> )	
<b>Screening Location</b>	Malvern, USA	

## Reference Compound IC50 for PDE2A



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/pde2a\\_pde\\_malvern](https://www.reactionbiology.com/datasheet/pde2a_pde_malvern)

