

# PDE4A Phosphodiesterase Assay Service

PDE4A [Phosphodiesterase Assay](#) for compound screening and profiling via fluorescence polarization-based quantification of enzyme activity.

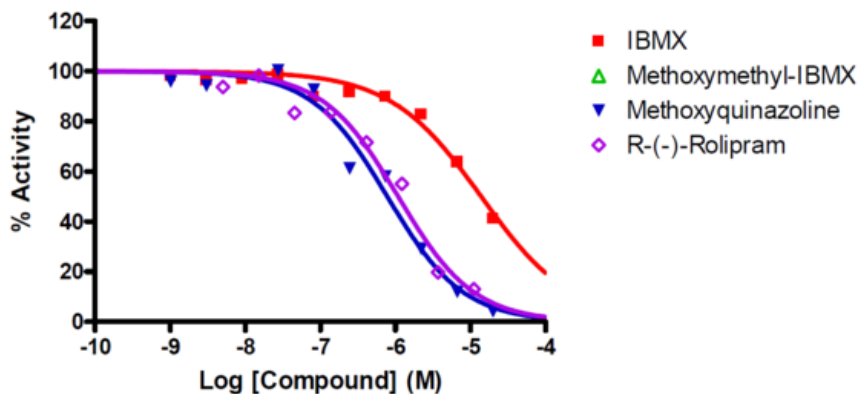
## Target Overview

<b>CAT#</b>	PDE4A
<b>Enzyme</b>	Human PDE4A
<b>Synonym</b>	cAMP-specific 3',5'-cyclic phosphodiesterase 4A, DPDE2, PDE46
<b>Substrate</b>	1 $\mu$ M cAMP

## 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	13,900
	Methoxymethyl-IBMX	>50,000
	Roliptam	>780
	Methoxyquinazoline	>1,100
<b>Scientific Information</b>	<a href="http://www.uniprot.org/uniprot/P27815">P27815</a> ( <a href="http://www.uniprot.org/uniprot/P27815">http://www.uniprot.org/uniprot/P27815</a> )	
<b>Screening Location</b>	Malvern, USA	

## Reference Compound IC50 for PDE4A



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/pde4a\\_pde\\_malvern](https://www.reactionbiology.com/datasheet/pde4a_pde_malvern)

