

# PDE4C Phosphodiesterase Assay Service

PDE4C [Phosphodiesterase Assay](#) for compound screening and profiling via fluorescence polarization-based quantification of enzyme activity.

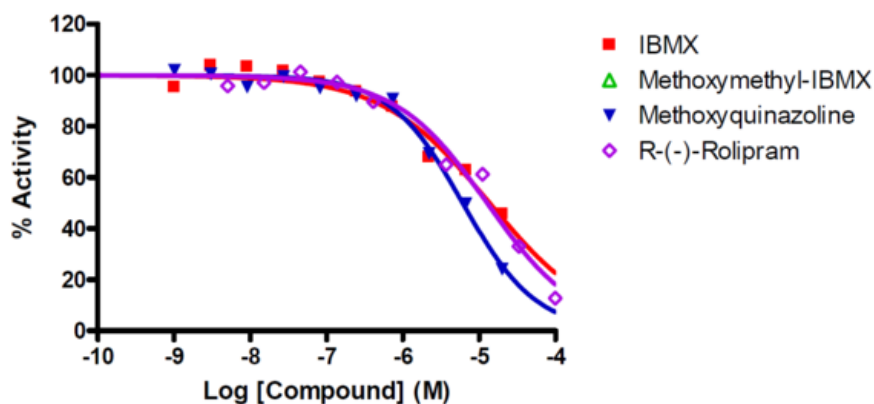
## Target Overview

CAT#	PDE4C
Enzyme	Human PDE4C
Synonym	cAMP-specific 3',5'-cyclic phosphodiesterase 4C, DPDE1, PDE21)
Substrate	1 $\mu$ M cAMP

## 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	13,700
	Methoxymethyl-IBMX	>50,000
	Roliptam	6,100
	Methoxyquinazoline	12,700
Scientific Information	<a href="http://www.uniprot.org/uniprot/Q08493">Q08493</a> ( <a href="http://www.uniprot.org/uniprot/Q08493">http://www.uniprot.org/uniprot/Q08493</a> )	
Screening Location	Malvern, USA	

## Reference Compound IC50 for PDE4C



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/pde4c\\_pde\\_malvern](https://www.reactionbiology.com/datasheet/pde4c_pde_malvern)

