

PDE3B Phosphodiesterase Assay Service

PDE3B [Phosphodiesterase Assay](#) for compound screening and profiling via fluorescence polarization-based quantification of enzyme activity.

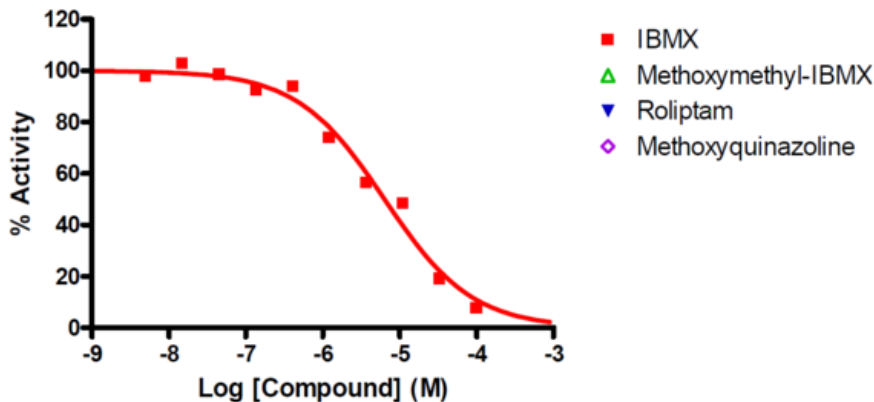
Target Overview

CAT#	PDE3B
Enzyme	Human PDE3B
Synonym	cGMP-inhibited 3',5'-cyclic phosphodiesterase B, CGI-PDE B
Substrate	1 μ M cAMP

Assay Properties

Reaction	Nucleoside 3',5'-cyclic phosphate + H ₂ O = nucleoside 5'-phosphate	
Readout	Enzymatic activity	
Reference Compound IC50s	Compounds	IC50 (nM)
	IBMX	6,400
	Methoxymethyl-IBMX	>100,000
	Roliptam	>100,000
	Methoxyquinazoline	>100,000
Scientific Information	Q13370 (http://www.uniprot.org/uniprot/Q13370)	
Screening Location	Malvern, USA	

Reference Compound IC50 for PDE3B



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/pde3b_pde_malvern

