

PHOSPHODIESTERASES

Phosphodiesterases (PDEs) play a central role in the regulation of various signaling cascades by specifically hydrolyzing the cyclic nucleotides, cAMP and cGMP into their nucleoside 5' – monophosphate form i.e. AMP and GMP. Hydrolysis of these nucleotides results in the breakdown of these second messengers and the abrogation of signaling. To date, eleven members of the mammalian PDE superfamily have been identified, PDE 1 – 11 respectively. Given the importance of these enzymes in the regulation of cAMP/cGMP signaling, SignalChem is pleased to announce that various members of the PDE superfamily are now available.

Product Name	Cat #
PDE1A, Active	P89-30H
PDE1B, Active	P89-30BG
PDE1C, Active	P89-30CG
PDE2A, Active	P90-30G
PDE3A, Active	P91-31G
PDE3B, Active	P91-31BG
PDE4A, Active	P92-31G

Product Name	Cat #
PDE4B, Active	P92-31BG
PDE4C, Active	P92-31CG
PDE4D, Active	P92-31DG
PDE5A, Active	P93-31G
PDE6A, Active	P94-31G
PDE6B, Active	P94-30BG
PDE6C Protein	P94-34CG

Product Name	Cat #
PDE7A, Active	P95-31G
PDE7B, Active	P95-30BG
PDE8A, Active	P96-30G
PDE8B, Active	P96-31BG
PDE9A, Active	P97-30G
PDE10A, Active	P98-31G
PDE12 Protein	P100-30G