
PLK3

Product code: 05-159

Human PLK3, catalytic domain [58-340 amino acids of accession number NP_004064.2] was expressed as N-terminal GST-fusion protein (59 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : CDC25ctide

ATP (μ M) K_m app / B_{in} : 6.8 / 5

Metal : Mg

Reference compound : GW843682X

IC₅₀ at ATP B_{in} (nM) : 33

IC₅₀ at 1 mM ATP (nM) : 450
