
BMX

Product code: 08-179

Full-length human BMX [1-675(end) amino acids of accession number NP_001712.1] was expressed as N-terminal GST-fusion protein (105 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP (μ M) Km app / Bin	: 75 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 10
IC50 at 1 mM ATP (nM)	: 45

