
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. GST-BMX was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Srctide

ATP (μ M) K_{mapp} / B_{in} : 75 / 75

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 10

IC₅₀ at 1 mM ATP (nM) : 45
