
AXL

Product code: 08-107

Human AXL, cytoplasmic domain [464-885(end) amino acids of accession number NP_001690.2] was expressed as N-terminal GST-fusion protein (74 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : CSKtide

ATP (μ M) Km app / Bin : 32 / 50

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 1.1

IC50 at 1 mM ATP (nM) : 7.9
