
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) K_m app / B_{in} : 32 / 50
Metal : Mg
Reference compound : Staurosporine
IC₅₀ at ATP B_{in} (nM) : 1.1
IC₅₀ at 1 mM ATP (nM) : 7.9
