
RET

Product code: 08-159

Human RET, cytoplasmic domain [658-1114(end) amino acids of accession number NP_066124.1] was expressed as N-terminal GST-fusion protein(79 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : CSKtide

ATP (μ M) Km app / Bin : 7.5 / 10
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 1.3
IC50 at 1 mM ATP (nM) : 20
