
RET[M918T]

Product code: 08-508

Human RET, cytoplasmic domain [658-1114(end) amino acids and M918T 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	: 4.2 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.4
IC50 at 1 mM ATP (nM)	: 81