
KIT

Product code: 08-156

Human KIT, cytoplasmic domain [544-976(end) amino acids of accession number NP_000213.1] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP (μ M) Km app / Bin	: 370 / 400
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.2
IC50 at 1 mM ATP (nM)	: 2.0