

---

---

## KIT[V560G]

Product code: 08-504

Human KIT, cytoplasmic domain [544-976(end) amino acids and V560G 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 ( $\mu$ M) Km app / Bin	: 110 / 250
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.2
IC50 at 1 mM ATP (nM)	: 1.6

---

---