

---

---

## KIT[V654A]

Product code: 08-511

Human KIT, cytoplasmic domain [544-976(end) amino acids and V654A 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	: 220 / 250
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 4.5
IC50 at 1 mM ATP (nM)	: 8.2

