

---

# FLT3

Product code: 08-154

Human FLT3, cytoplasmic domain [564-993(end) amino acids of accession number NP\_004110.2] was expressed as N-terminal GST-fusion protein (77 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP ( $\mu$ M) Km app / Bin	: 94 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.20
IC50 at 1 mM ATP (nM)	: 0.34

