

---

# 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

---