

---

---

## RET[G691S]

Product code: 08-522

Human RET, cytoplasmic domain [658-1114(end) amino acids and G691S of accession number NP\_066124.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CSKtide
ATP ( $\mu$ M) Km app / Bin	: 13 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.1
IC50 at 1 mM ATP (nM)	: 24