

---

# KDR

Product code: 08-191

Human KDR, cytoplasmic domain [790-1356(end) amino acids of accession number NP\_002244.1] was expressed as N-terminal GST-fusion protein (90 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CSKtide
ATP ( $\mu$ M) Km app / Bin	: 74 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.1
IC50 at 1 mM ATP (nM)	: 13