

---

---

# FMS(CSF1R)

Product code: 08-155

Human FMS, cytoplasmic domain [538-972(end) amino acids of accession number NP\_005202.2] 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	: 26 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.26
IC50 at 1 mM ATP (nM)	: 0.70

