

---

# 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          |

