
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 (μ 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 |

