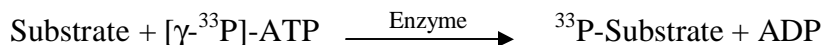


FMS

(Macrophage colony-stimulating factor 1 receptor [Precursor], CSF-1-R
Fms proto-oncogene, c-fms, CSF1R)

- **CAT#:** FMS
- **Enzyme:** Human FMS
- **Substrate:** Peptide substrate, [KKKSPGEYVNIEFG], 5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	15.3
PP2	4,900
AG1478	48,700
PDGFR I-III	12,000
D-64406	6,000

- **Scientific Information:** P07333 (<http://www.expasy.org/uniprot/P07333>)

Reference Compound IC50 for FMS

