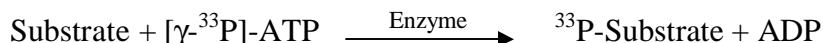


RON

(Macrophage-stimulating protein receptor [Precursor], MSP receptor, p185-Ron, MST1R)

- **CAT#:** RON
- **Enzyme:** Human RON
- **Substrate:** Peptide substrate, [KKS_RRGDYMTMQIG], 5 μM
ATP 1 μM

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	22
AG1478	94,000
PDGFR I-III	1,900
D-64406	17,200

- **Scientific Information:** Q04912 (<http://www.expasy.org/uniprot/Q04912>)

