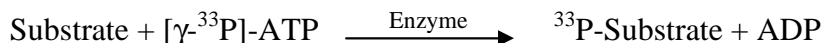


LRRK2

(Leucine-rich repeat serine/threonine-protein kinase 2, Dardarin, PARK8)

- **CAT#:** LRRK2
- **Enzyme:** Human LRRK2
- **Substrate:** Peptide substrate, [RLGRDKYKTLRQIRQ], 20 μ M
ATP 10 μ M
- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	6.8
Ro 31-8220	178
GW5074	50
H-89	23,000

- **Scientific Information:** Q5S007 (<http://www.expasy.org/uniprot/Q5S007>)

Reference compound IC50 for LRRK2

