

## LRRK2 (R1441C)

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

- **CAT#:** LRRK2 (R1441C)
- **Enzyme:** Human LRRK2 (R1441C)
- **Substrate:** Peptide substrate, [RLGRDKYKTLRQIRQ], 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	3.4
Ro 31-8220	230
GW5074	26
H-89	26,000
Rottlerin	>100,000

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

### Reference compound IC50 for LRRK2 (R1441C)

